

ICP-OES Agilent séries 5800 e 5900

Guia do usuário



Avisos

Nº de peça do manual

G8020-99002 1ª Edição, outubro de 2019

Direitos autorais

© Agilent Technologies, Inc. 2019

Nenhuma parte deste material pode ser reproduzida de qualquer forma ou por qualquer meio (incluindo armazenamento e recuperação eletrônica ou a tradução para outro idioma) sem autorização prévia por escrito da Agilent Technologies, Inc. De acordo com as leis de direitos autorais americanas e internacionais.

Impresso na Malásia

Agilent Technologies, Austrália (M) Pty Ltd 679 Springvale Road Mulgrave, Victoria, 3170

Austrália

www.agilent.com

Garantia

O material contido neste documento é fornecido "no estado em que se encontra" e está sujeito a alterações, sem aviso prévio em edicões futuras. Além disso. com o máximo rigor permitido pelas leis aplicáveis, a Agilent isenta-se de responsabilidade em relação a garantias, expressas ou implícitas, em relação a este manual e a qualquer informação contida nele, incluindo mas não limitado as garantias implícitas de adequação comercial e adequação a um propósito específico. A Agilent não será responsável por erros ou por danos incidentais ou consequenciais relacionados ao fornecimento, ao uso ou ao desempenho deste documento ou de qualquer informação nele contida. Se a Agilent e o usuário possuírem um acordo por escrito em separado com os termos de garantia cobrindo o material neste documento que entrem em conflito com esses termos, os termos de garantia do documento em separado prevalecerão.

Licenças de tecnologia

O hardware e/ou software descrito(s) neste documento é/são fornecido(s) sob licença, podendo ser usado(s) ou copiado(s) somente de acordo com os termos dessa licença.

Legenda de direitos restritos

Direitos restritos ao governo dos EUA. Os direitos de software e dados técnicos concedidos ao governo federal incluem apenas os direitos normalmente concedidos aos clientes usuários finais. A Agilent fornece esta licenca comercial habitual relativamente a software e dados técnicos de acordo com a FAR 12.211 (Dados técnicos) e 12.212 (Software de computador) e, para o Departamento de Defesa, segundo a DFARS 252.227-7015 (Dados técnicos – Itens comerciais) e DFARS 227.7202-3 (Direitos relativos a software de computador comercial ou documentação de software de computador).

Avisos de segurança

CUIDADO

Um aviso de **CUIDADO** representa um perigo. É utilizado para chamar a atenção para um procedimento de operação, prática ou ação similar, que pode resultar em danos ao produto ou perda de dados importantes se não for executado ou seguido corretamente. Não prossiga após uma indicação de **CUIDADO** até que as condições indicadas sejam completamente compreendidas e atendidas.

AVISO

Um AVISO representa um perigo. Ele chama a atenção para uma prática, um procedimento operacional ou similares que, se não forem seguidos corretamente poderão resultar em lesões pessoais ou fatais. Não prossiga após uma indicação de AVISO até que as condições indicadas tenham sido totalmente compreendidas e atendidas.

Índice

1	Riscos e práticas de segurança	7
	Geral	7
	Verificação do estado de segurança	8
	Plasma	8
	Calor, vapores e fumaça	9
	Riscos de gás comprimido	10
	Riscos elétricos	11
	Outras precauções	11
	Símbolos de advertência	13
2	Introdução	15
	Requisitos de preparo do local	15
	Documentação do usuário	15
	Convenções	16
	Notas e dicas	16
	Suprimento de gás argônio para a ICP-OES	16
	Sistema de exaustão	17
	Suprimento de ar de resfriamento do instrumento da ICP-OES	18
	Sistema de resfriamento de água	18
	Recipiente de drenagem	19
3	Instalação	21
	Visão geral do instrumento ICP-OES Agilent	22
	Codificação por cores do LED de status do instrumento	24
	Botão frontal Liga/Desliga	25
	Software ICP Expert	25
	Conectar o software ICP Expert ao ICP-OES	26

	Executar as calibrações do detector e do comprimento de onda	27
	Salvar e visualizar os dados da calibração	28
	Substituição de componentes de hardware	28
	Tochas do ICP-OES	29
	Montagem e desmontagem da tocha	30
	Montagem da tocha semidesmontável	32
	Montagem da tocha completamente desmontável	35
	Desmontagem da tocha	35
	Acessórios	39
	SPS 4	40
	Sistema avançado de válvulas (AVS) – AVS 4, AVS 6 e AVS 7	40
	Adaptador de duto de entrada externo	41
	Acessório de geração de vapor (VGA)	41
	Sistema de introdução de amostras multimodal (MSIS)	41
	Umidificador de argônio	41
	Câmara de nebulização de temperatura programável IsoMist	42
4	Operação	43
	Lista de verificação de análises	43
	Ligar o instrumento e o software	44
	Ligar o instrumento pela primeira vez (ou após desligamento prolongado)	44
	Operar o instrumento em estado de espera	45
	Preparar para análise	46
	Executar as calibrações do detector e do comprimento de onda	46

Criar/abrir uma worksheet	47
Criar uma nova worksheet	47
Abrir uma worksheet existente	47
Criar uma nova worksheet a partir de um modelo	47
Desenvolver um método	48
Executar amostras	50
Imprimir um relatório	50
Desligar o instrumento	51
Modo de espera do instrumento com argônio ativado	51
Modo de espera do instrumento com argônio desativado	51
Configurar o instrumento no modo de espera com argônio ativado ou desativado	52
Desligar o instrumento para desligamento prolongado	54
Manutenção e solução de problemas Rotina	55 56
Limpeza	58
Limpeza da tocha	59
Importante	60
Importante Procedimento de limpeza por imersão com ácido para a tocha em amostras de base aquosa ou ácida (tochas de peça única e semidesmontáveis)	60 60
Importante Procedimento de limpeza por imersão com ácido para a tocha em amostras de base aquosa ou ácida (tochas de peça única e semidesmontáveis) Lavagem da tocha	60 60 63
Importante Procedimento de limpeza por imersão com ácido para a tocha em amostras de base aquosa ou ácida (tochas de peça única e semidesmontáveis) Lavagem da tocha Secagem da tocha	60 60 63 65
Importante Procedimento de limpeza por imersão com ácido para a tocha em amostras de base aquosa ou ácida (tochas de peça única e semidesmontáveis) Lavagem da tocha Secagem da tocha Verificações adicionais após a limpeza	60 60 63 65 67
Importante Procedimento de limpeza por imersão com ácido para a tocha em amostras de base aquosa ou ácida (tochas de peça única e semidesmontáveis) Lavagem da tocha Secagem da tocha Verificações adicionais após a limpeza Solução de problemas	60 60 63 65 67 68
Importante Procedimento de limpeza por imersão com ácido para a tocha em amostras de base aquosa ou ácida (tochas de peça única e semidesmontáveis) Lavagem da tocha Secagem da tocha Verificações adicionais após a limpeza Solução de problemas Peças sobressalentes	60 60 63 65 67 68 68
 Importante Procedimento de limpeza por imersão com ácido para a tocha em amostras de base aquosa ou ácida (tochas de peça única e semidesmontáveis) Lavagem da tocha Secagem da tocha Verificações adicionais após a limpeza Solução de problemas Peças sobressalentes Suporte técnico 	60 63 65 67 68 68 68

5

Esta página foi deixada em branco intencionalmente.

1 Riscos e práticas de segurança

Geral	7
Verificação do estado de segurança	8
Plasma	8
Calor, vapores e fumaça	9
Riscos de gás comprimido	10
Riscos elétricos	11
Outras precauções	11
Símbolos de advertência	13

Geral

Salvo especificação em contrário, as declarações deste manual se aplicam aos instrumentos ICP-OES Agilent 5800 e 5900.

A operação de um ICP-OES Agilent envolve o uso de gases comprimidos, energia de radiofrequência de alta tensão e materiais perigosos, incluindo fluidos corrosivos e líquidos inflamáveis. O uso descuidado, indevido ou não qualificado deste espectrômetro, ou de produtos químicos utilizados com ele, pode causar morte ou ferimentos graves ao pessoal e/ou danos graves ao equipamento e à propriedade. Apenas pessoal treinado deve usar este instrumento.

O espectrômetro incorpora travas e tampas projetadas para impedir o contato descuidado com possíveis riscos. Se o instrumento for usado de alguma forma não especificada pela Agilent, essa proteção oferecida pelo equipamento poderá ser comprometida. É recomendável desenvolver hábitos seguros de trabalho que não dependam do bom funcionamento das travas para uma operação segura. É essencial que nenhuma trava ou tampa seja contornada, danificada ou removida.

As práticas de segurança descritas abaixo são fornecidas para ajudar o usuário a operar o instrumento com segurança. Leia todos os tópicos de segurança na íntegra antes de tentar operar o instrumento e sempre opere o espectrômetro de acordo com essas práticas de segurança.

Verificação do estado de segurança

As seguintes precauções gerais de segurança devem ser seguidas durante todas as fases de operação, manutenção e reparo deste instrumento.

Para garantir a segurança contínua do instrumento após procedimentos de manutenção e reparo, verifique se o instrumento voltou a um estado seguro para o usuário. Isso inclui a execução de verificações de desempenho para se certificar se os sistemas de segurança dos instrumentos estão funcionando corretamente. Verifique a condição geral do instrumento durante a operação quanto a desgastes ou sinais de corrosão que podem inibir a funcionalidade ou a segurança.

O descumprimento destas precauções ou advertências específicas em qualquer parte deste manual viola os padrões de segurança de projeto, fabricação e uso previsto do instrumento. A Agilent Technologies não assume nenhuma responsabilidade se o cliente não cumprir esses requisitos.

Plasma

O plasma é extremamente quente (cerca de 10.000°C) e irradia níveis perigosos de energia ultravioleta (UV) e de radiofrequência (RF). A bobina de indução opera a 1.500 V RMS e a cerca de 27 MHz. A exposição à energia UV e de RF pode causar danos graves à pele e cataratas nos olhos, enquanto o contato próximo com o plasma em operação pode resultar em queimaduras graves na pele e uma descarga elétrica que percorre distância considerável e pode causar morte, choque elétrico grave ou queimaduras na derme da pele.

O plasma nunca deve ser operado, exceto se:

- a porta do compartimento do plasma e sua alça estiverem completamente fechadas, e
- o espaço acima da chaminé e da entrada de ar estiver livre de objetos.

A proteção ao redor do compartimento da tocha foi projetada para reduzir as radiações UV visíveis e de RF a níveis seguros, permitindo ainda fácil acesso à tocha, à sua instalação e visualização. O espectrômetro possui um sistema de intertravamento projetado para extinguir o plasma se houver falha na rede elétrica, se a alça da porta do compartimento da tocha for aberta ou se a alça de carregamento da tocha estiver aberta. Não tente contornar o sistema de intertravamento.

Antes de abrir a porta do compartimento da tocha, *sempre* extinga o plasma pressionando SHIFT + F5 no teclado ou clicando no ícone "Plasma Off" (Plasma apagado) na barra de ferramentas do software ICP Expert.

A tocha e seus arredores permanecem quentes por até cinco minutos após a extinção do plasma. Entrar em contato com esta área antes que ela tenha esfriado o suficiente pode resultar em queimaduras. Permita que a tocha e seu compartimento esfriem antes de realizar outro trabalho nessa área ou use luvas resistentes ao calor.

O sistema de plasma foi cuidadosamente projetado para operar com segurança e eficácia durante o uso de tochas e componentes relacionados que estejam em conformidade com os critérios de projeto da Agilent. O uso de componentes não aprovados no compartimento do plasma pode tornar o sistema inoperante e/ou perigoso. Pode também invalidar a garantia do instrumento. Use apenas tochas e componentes relacionados fornecidos ou autorizados pela Agilent.

Calor, vapores e fumaça

O calor, o ozônio, vapores e a fumaça gerados pelo plasma podem ser perigosos e devem ser extraídos do instrumento por meio de um sistema de exaustão. Verifique se o instrumento está equipado com um sistema de exaustão apropriado (conforme especificado no Guia de preparo do local). O sistema deve ser ventilado para o exterior, de acordo com os regulamentos locais, e nunca para dentro do edifício. Verifique regularmente o sistema de exaustão através de testes de gases, para garantir que o mesmo esteja funcionando corretamente. O sistema de exaustão deve sempre ser ligado *antes* da ignição do plasma.

Riscos de gás comprimido

Todos os gases comprimidos (com exceção do ar) podem gerar riscos se vazarem para a atmosfera. Até mesmo pequenos vazamentos nos sistemas de suprimento de gás podem ser perigosos. Qualquer vazamento (com exceção do ar ou oxigênio) pode resultar em uma atmosfera deficiente em oxigênio, podendo causar asfixia. A área em que os cilindros estão armazenados e aquela ao redor do instrumento devem ser adequadamente ventiladas para evitar tais acúmulos de gás.

Os cilindros de gás devem ser armazenados e manipulados em estrita conformidade com os regulamentos e códigos de segurança locais. Os cilindros devem ser usados e armazenados apenas na posição vertical e presos a uma estrutura imóvel ou a um suporte para cilindro construído adequadamente. Movimente os cilindros apenas ao colocá-los em um carrinho construído adequadamente.

Use apenas conectores de mangueiras e reguladores aprovados (consulte as instruções do fornecedor de gás). Mantenha os cilindros de gás resfriados e devidamente rotulados. (Todos os cilindros são equipados com um dispositivo de alívio de pressão que se romperá e esvaziará o cilindro caso a pressão interna se eleve acima do limite de segurança devido a temperaturas excessivas). Certifiquese de que o gás correto esteja sendo usado antes de conectá-lo ao instrumento.

O gás primário usado com o espectrômetro é o argônio, que é o gás condutor para o plasma. É possível usar argônio ou nitrogênio como gás de purga do policromador. Podem ser necessários outros gases em futuras opções e acessórios. Use apenas gases com "grau de instrumento" com o seu espectrômetro.

Se estiver usando gases criogênicos (por exemplo, argônio líquido), evite queimaduras graves usando roupas protetoras e luvas adequadas.

Riscos elétricos

O sistema do espectrômetro e alguns acessórios contêm circuitos elétricos, dispositivos e componentes que funcionam com tensões perigosas. O contato com esses circuitos, dispositivos e componentes pode resultar em morte, ferimentos graves ou choque elétrico doloroso. Os painéis ou tampas que são fixados por parafusos no espectrômetro, bem como os acessórios, podem ser abertos *apenas* por engenheiros de serviços de campo treinados, qualificados ou autorizados pela Agilent (salvo indicação contrária). Consulte os manuais ou as etiquetas do produto fornecidos com seu computador pessoal (PC), monitor, impressora ou sistema de resfriamento por água para saber quais partes podem ser acessadas pelo operador.

Conectar o ICP-OES Agilent a uma fonte de energia não equipada com aterramento de proteção gera riscos de choque para o operador e pode danificar o instrumento. Da mesma forma, interromper o condutor de proteção na parte interna ou externa do ICP-OES Agilent ou anular o aterramento do cabo elétrico gera riscos de choque para o operador e pode danificar o instrumento.

Outras precauções

O uso do sistema do espectrômetro e dos acessórios pode envolver materiais, solventes e soluções inflamáveis, corrosivos, tóxicos ou outros produtos perigosos. O uso indevido, descuidado ou não qualificado desses materiais, solventes e soluções pode gerar riscos de explosão, queimadura química, incêndio, intoxicação e outros perigos que podem resultar em morte, ferimentos graves ou danos ao equipamento. Tome todas as precauções necessárias, incluindo o uso de jalecos, óculos protetores e outras formas apropriadas de proteção pessoal. Todos os resíduos deverão ser descartados de acordo com os requisitos regulatórios locais.

A operação de um ICP-OES envolve análises de soluções que foram preparadas ou digeridas com ácidos ou, em alguns casos, amostras que foram preparadas em solventes orgânicos.

Em caso de dúvida sobre um fluido específico, esse fluido não deve ser utilizado até que o fabricante confirme que ele não representa um perigo.

A concentração de ácido na amostra medida varia de acordo com as etapas de digestão e os tipos de ácido utilizados. Os usuários do instrumento devem estar cientes dos riscos associados ao uso dos ácidos utilizados no preparo de amostras e devem tomar todas as precauções necessárias, incluindo o uso de jalecos, óculos protetores e outras formas adequadas de proteção pessoal. Os resíduos ácidos deverão ser descartados de acordo com os requisitos regulatórios locais.

O tipo, a volatilidade e a concentração dos solventes orgânicos usados na amostra medida variam de acordo com o solvente selecionado e com o preparo de amostras envolvido. Os usuários do instrumento devem estar cientes dos riscos associados ao uso dos solventes orgânicos utilizados no preparo de amostras e devem tomar todas as precauções necessárias, incluindo assegurar uma ventilação adequada durante a utilização e o uso de jalecos, óculos protetores, luvas e outras formas adequadas de proteção pessoal. Todos os resíduos orgânicos deverão ser descartados de acordo com os requisitos regulatórios locais.

O fluxo de ar até a porta de entrada de ar do espectrômetro e seus acessórios deve estar desobstruído. Não obstrua as grades de ventilação no espectrômetro e seus acessórios. Consulte os manuais fornecidos com seu PC, monitor, impressora e sistema de resfriamento por água relacionados aos requisitos específicos de ventilação.

É necessário muito cuidado ao trabalhar com peças de vidro ou de quartzo para evitar quebras e cortes. Essa questão é particularmente importante na conexão do nebulizador à câmara de nebulização ou na remoção e substituição de peças de uma tocha quebrada.

O espectrômetro pesa aproximadamente 90 kg. Para evitar ferimentos ao pessoal ou danos ao instrumento ou à propriedade, sempre use um dispositivo de elevação mecânica adequado para mover o instrumento.

Use apenas peças sobressalentes autorizadas ou fornecidas pela Agilent com o seu instrumento. Somente operadores treinados devem usar o instrumento.

Símbolos de advertência

Consulte a seguir uma lista de símbolos que podem aparecer juntamente com as advertências neste manual ou no espectrômetro. Além disso, é mostrado o perigo que eles descrevem. O início do texto de advertência é indicado por um ícone de advertência:

ADVERTÊNCIA

Um símbolo triangular indica uma advertência. O significado dos símbolos que podem aparecer juntamente com as advertências na documentação ou no próprio instrumento são os seguintes:





Superfície quente



Risco químico



Risco de frio extremo



Pesado (perigo para os pés)



Gases nocivos



Desligar a energia



Risco para os olhos



Pesado (perigo para as mãos)



Radiacão de RF

O símbolo a seguir pode ser usado em etiquetas de advertência afixadas ao instrumento. Ao ver este símbolo, consulte o manual de operação ou de serviço em questão para saber qual é o procedimento correto indicado por essa etiqueta de advertência.

Os símbolos seguintes aparecem no instrumento para fins de informação.



2 Introdução

Requisitos de preparo do local	15
Documentação do usuário	15
Suprimento de gás argônio para a ICP-OES	16
Sistema de exaustão	17
Suprimento de ar de resfriamento do instrumento da ICP-OES	18
Sistema de resfriamento de água	18
Recipiente de drenagem	19

Requisitos de preparo do local

Antes de receber o instrumento, você receberá o Guia de preparo do local para o ICP-OES Agilent 5800 e 5900 descrevendo os requisitos operacionais e ambientais do sistema ICP-OES. Antes de instalar o ICP-OES, você deve preparar o laboratório de acordo com essas instruções. Guarde o Guia de preparo do local para consultas futuras, por exemplo, se você decidir mudar o instrumento de lugar. Se você não tiver mais sua cópia, é possível obtê-la gratuitamente no escritório Agilent local.

Posicione o equipamento de modo a ter fácil acesso ao botão Liga/Desliga no lado esquerdo do instrumento.

Documentação do usuário

Você recebeu a seguinte documentação para ajudá-lo a configurar e operar seu sistema ICP-OES Agilent:

 Este Guia do usuário, com práticas de segurança e informações acerca de riscos, instruções para encontrar informações sobre instalação e manutenção dos componentes do ICP-OES e uma breve visão geral da operação.

 Um abrangente Help and Learning Center (Centro de treinamento e ajuda) contendo ajuda sensível ao contexto, com informações sobre como configurar métodos e executar testes de instrumentos, noções básicas sobre hardware e vídeos de manutenção, instruções passo a passo para as operações executadas com frequência e instruções para o uso de qualquer acessório solicitado.

Convenções

As seguintes convenções foram utilizadas em toda a documentação:

- Itens de menu, opções de menu e nomes de campos (por exemplo, clique em Copiar no menu Editar) foram digitados em negrito. O negrito também é usado para indicar os botões que aparecem no software (por exemplo, clique em OK).
- As LETRAS MAIÚSCULAS indicam comandos de teclado (por exemplo, pressione a tecla F2) e o texto que você deve digitar usando o teclado (por exemplo, digite SETUP quando solicitado).

Notas e dicas

A Nota é usada para fazer recomendações ou dar informações.

A Dica é usada para dar sugestões práticas que o ajudarão a alcançar o melhor desempenho possível de seu ICP-OES.

Suprimento de gás argônio para a ICP-OES

Tabela 1. Fluxo típico para os instrumentos da ICP-OES Agilent

Vazões típicas	Argônio (com gás de purga de nitrogênio)
Medição de comprimentos de onda > 189 nm (multiplicador de pressão do policromador desativado)	Fluxo de argônio 16,7 L/min Fluxo de nitrogênio 0,8 L/min
Medição de comprimentos de onda < 189 nm (multiplicador de pressão do policromador ativado)	Fluxo de argônio 16,7 L/min Fluxo de nitrogênio 4,4 L/min

O argônio pode ser usado como gás de plasma e de purga, por isso o fluxo total de argônio é maior.

Os usuários (ou outras pessoas autorizadas) devem realizar testes de vazamento adequados, necessários para garantir a segurança das conexões de gases e líquidos que o operador foi orientado a montar durante a instalação, uso normal ou manutenção.

Sistema de exaustão

Superfície quente

O plasma opera a temperaturas extremamente altas. A fumaça de exaustão do instrumento ICP-OES pode ser nociva ou corrosiva.

O sistema de exaustão deve ser inspecionado regularmente a fim de assegurar um fluxo adequado. Consulte o Guia de preparo do local do ICP-OES para conhecer as vazões apropriadas.

ADVERTÊNCIA



O respiradouro externo pode ficar quente durante a operação do instrumento da ICP-OES e permanecer quente por algum tempo após o desligamento do instrumento. Deixe o respiradouro externo esfriar por pelo menos cinco minutos antes de tentar remover a mangueira de exaustão. Use luvas resistentes ao calor.

Suprimento de ar de resfriamento do instrumento da ICP-OES

O instrumento da ICP-OES requer *ar limpo, seco e não corrosivo para fins de resfriamento.* Ele é fornecido ao instrumento através de um respiradouro de suprimento de ar localizado na parte superior esquerda do instrumento. O respiradouro contém um filtro de poeira para filtrar matéria particulada.

O suprimento de ar é usado para resfriar os componentes mecânicos e eletrônicos internos do instrumento. Vários desses conjuntos contêm peças propensas à corrosão. A introdução de ar de resfriamento contaminado, com altas concentrações de vapores ácidos ou outras substâncias corrosivas, pode causar danos ao instrumento.

Devido à natureza corrosiva de alguns trabalhos de análise, recomenda-se que as aplicações que exigem uso de grande quantidade de materiais corrosivos tenham um sistema de suprimento de ar de resfriamento externo. É *necessário* que o ar de resfriamento seja proveniente de uma área ambientalmente controlada que esteja longe da exaustão do instrumento e de qualquer outra área onde materiais corrosivos sejam armazenados ou utilizados. Não conduza ar quente e úmido para dentro de um instrumento que esteja em um ambiente de laboratório resfriado.

O sistema de ar de resfriamento, que possui conduto de gases, ventilador, dutos e coifa, deve fornecer na entrada do instrumento uma pressão positiva de ar de 4 m³/min quando se utiliza o conjunto adaptador de duto de entrada externo. Os dutos devem ser resistentes à corrosão e à prova de fogo.

Sistema de resfriamento de água

Os instrumentos ICP-OES Agilent requerem uma fonte de água para resfriamento. Consulte o Guia de preparo do local, para ver quais são os sistemas de resfriamento de água compatíveis, e a documentação fornecida com o sistema de resfriamento de água para obter informações sobre a instalação e as instruções de utilização.

NOTA

Recomenda-se a regulação da pressão dos fornecimentos em que a pressão da água para resfriamento possa estar sujeita a flutuações. A regulação da pressão é necessária em fornecimentos que possam exceder a pressão máxima admissível de 400 kPa.

O instrumento é equipado com um sensor de fluxo de água, que interromperá a operação do plasma caso o fluxo da água para resfriamento que passa pelo instrumento caia para menos de 1,7 L/min. Um segundo sensor de fluxo de água interromperá o conjunto de resfriamento Peltier da câmera caso o fluxo de água que passa pelo instrumento caia para menos de 0,2 L/min.

CUIDADO

Certifique-se sempre de que o sistema de resfriamento de água esteja ligado antes de acender o plasma.

Recipiente de drenagem

O sistema ICP-OES Agilent precisa de um recipiente de drenagem para despejar fluidos em excesso da câmara de nebulização e do amostrador automático. A tubulação adequada é fornecida com o espectrômetro para o uso com solventes inorgânicos. Ao utilizar solventes orgânicos, será necessário usar uma tubulação de drenagem diferente, que seja adequada para o solvente em uso.

Deverá ser fornecido pelo usuário um recipiente quimicamente inerte, que não seja de vidro ou tenha bocal estreito, para que se mantenha um mínimo de 2 litros de resíduos. O recipiente deve ficar embaixo do compartimento de amostras (ou no lado direito do instrumento), onde ficará protegido pela bancada e à vista do operador.

Esta página foi deixada em branco intencionalmente.

Visão geral do instrumento ICP-OES Agilent	22
Codificação por cores do LED de status do instrumento	24
Botão frontal Liga/Desliga	25
Software ICP Expert	25
Conectar o software ICP Expert ao ICP-OES	26
Executar as calibrações do detector e do comprimento de onda	27
Substituição de componentes de hardware	28
Montagem e desmontagem da tocha	30
Acessórios	39

O ICP-OES Agilent deve ser instalado por um engenheiro de serviços de campo treinado, qualificado ou autorizado pela Agilent.

Você deve preencher e devolver o formulário do Guia de preparo do local declarando que preparou o laboratório de acordo com os requisitos detalhados nesse manual. Um dos representantes da Agilent marcará com você uma data adequada para a instalação.

Os detalhes sobre como desembalar o instrumento e o que fazer em caso de danos durante o transporte também estão descritos no Guia de preparo do local.

Visão geral do instrumento ICP-OES Agilent



Figura 1. Visão frontal e lateral do instrumento ICP-OES

1. Exaustão	8. Alça do suporte da tocha	15. Alça do compartimento da tocha
2. Filtro de entrada de ar	9. Câmara de nebulização	16. Conjunto de entrada de água
3. Janela do cone pré-óptico e de visão axial (não exibida)	10. Nebulizador	17. Filtro óptico de purga para argônio ou nitrogênio
4. Janela de visualização radial e da ponta cônica (não exibida)	11. Bomba peristáltica	18. Conjunto de suprimento de gás
5. Bobina de indução	12. Botão de alimentação e conexão do cabo de energia	19. Localização do acessório Válvula seletora AVS 4, AVS 6 ou AVS 7 opcional
6. Tocha	13. Botão Liga/Desliga do painel frontal	20. Dreno de transbordamento de líquido
7. Conexões do nebulizador e do gás make up	14. LED indicador de status do instrumento	

Todas as conexões de serviços do instrumento ICP-OES são feitas no lado direito do instrumento, exceto a de energia elétrica, que é feita no lado esquerdo. Remova manualmente a tampa do lado direito para ter acesso às conexões da rede Ethernet, de acessórios, de água e de gás. O botão Liga/Desliga está localizado na parte frontal do instrumento, do lado inferior esquerdo, e o botão de alimentação elétrica está localizado na parte de trás, do lado esquerdo.



Figura 2. Conexões de entrada e saída na lateral do instrumento ICP-OES

	Item	Descrição
1	Conexões de	Conexões de acessórios Agilent e do cabo Ethernet
	acessórios e do cabo	blindado para comunicação entre o computador PC e o
	de rede LAN	instrumento
2	Entrada de gás argônio	Entrada de gás argônio, padrão
3	Entrada de gás	Entrada de gás opcional (mistura 80% Ar/20% O2). Entrada
	opcional	de gás nitrogênio opcional (não exibida)
4	Filtro óptico de purga	Filtro de gás argônio ou nitrogênio
5	Saída de água	Conecta-se para retorno do resfriador de água
6	Filtro de água	Filtro de água com partículas grossas
7	Entrada de água	Conecta-se para fornecimento do resfriador de água

Codificação por cores do LED de status do instrumento

A luz indicadora de status do instrumento, na parte frontal superior direita do ICP-OES Agilent, exibe diferentes cores que representam o estado do instrumento:

- A luz verde indica o seguinte:
 - O instrumento e o software estão conectados e prontos para tentar uma sequência de ignição do plasma; ou
 - O instrumento e o software estão conectados, e a sequência de ignição do plasma está em andamento; ou
 - O instrumento e o software estão conectados, o plasma está aceso e operante, sendo possível realizar a análise de amostras.
- Uma luz amarela piscante indica que o instrumento e o software não estão conectados (o instrumento está desativado).
- A luz amarela indica que o instrumento e o software estão conectados, mas o instrumento não está pronto para dar início a uma sequência de ignição do plasma devido a uma porta bloqueada ou ao travamento da alça da tocha, baixa pressão do gás, baixo fluxo de água ou de ar de resfriamento. É necessária a intervenção do usuário para resolver o problema.
- Uma luz laranja pulsante (intensa ou diminuindo) indica que o instrumento está inicializando ou realizando uma atualização de firmware.
- Uma luz vermelha indica que um erro no firmware do instrumento está causando uma interrupção na análise ou restringindo a operação normal do instrumento. É necessária a intervenção do usuário para resolver o problema.

Botão frontal Liga/Desliga

O LED do botão Liga/Desliga frontal indica o estado ativo ou não ativo básico do instrumento.

- Desligado indica que o instrumento está completamente sem energia quando o botão de alimentação elétrica do lado esquerdo do instrumento também está desligado.
- Uma luz verde que pisca a cada 10 segundos indica que a maior parte do instrumento está sem energia, exceto o circuito sensor do botão Liga/Desliga no painel frontal, que fica dentro do módulo de alimentação elétrica. Pressione o botão Liga/Desliga no painel frontal para ligar o instrumento por completo.
- Luz verde piscando devagar (1 Hz). O módulo de alimentação elétrica está sendo inicializado e executado através da verificação automática de inicialização ou está desligando o instrumento.
- Luz verde piscando rápido (2 Hz). Houve falha na verificação automática de inicialização do módulo de alimentação elétrica, e o instrumento não pode ser utilizado. A fonte de alimentação elétrica não é adequada ou o módulo de alimentação elétrica interno precisa ser substituído.
- Uma luz verde constantemente acesa indica que o instrumento está ligado.

Software ICP Expert

O engenheiro de serviços de campo treinado, qualificado ou autorizado pela Agilent instalará o software ICP Expert para você durante o processo de instalação. Contudo, você terá de instalar o software por si só posteriormente, por exemplo, se você trocar de computador. As instruções são fornecidas.

Existem diversas situações de instalação:

- Arquivo plano independente ou de rede
- Banco de dados independente ou de rede para instalações conforme a 21 CFR Parte 11
- Upgrade de banco de dados independente ou de rede para instalações do ICP Expert

A instalação inclui:

- Instalação do software ICP Expert
- Instalação da Ajuda do ICP Expert
- Instalação do cabo de rede LAN
- Configuração do endereço IP do instrumento
- Instalação e configuração da SDA para conformidade com a 21 CFR Parte 11
- Instalação e configuração da SCM para conformidade com a 21 CFR Parte 11
- Conclusão de uma varredura de corrente escura e de uma calibração de comprimento de onda

Para obter instruções sobre como instalar o seu software ICP Expert, consulte a documentação fornecida com o software:

- Instruções de instalação do software ICP Expert
- Instruções de instalação do software ICP Expert para ambientes conforme a 21 CFR Parte 11

Após instalar o software apropriado, conecte o cabo Ethernet, configure o endereço IP do instrumento, se necessário, e realize a calibração do detector e do comprimento de onda.

Conectar o software ICP Expert ao ICP-OES

Para conectar o software e o instrumento:

- 1 Execute o software ICP Expert clicando duas vezes no ícone da área de trabalho.
- 2 Clique em Instrument (Instrumento).
- 3 Clique em Connect (Conectar).
- 4 Insira o endereço IP do instrumento ou selecione um instrumento existente.
- 5 Clique em Connect (Conectar).
- 6 Clique em Close (Fechar).

Executar as calibrações do detector e do comprimento de onda

Para realizar a calibração do detector e do comprimento de onda:

- 1 Ligue a exaustão externa.
- 2 Execute o software ICP Expert e abra a janela de configuração do instrumento.
- 3 Conecte o software ao instrumento, se necessário.
 - **a** Clique em **Instrument** (Instrumento) na barra de ferramentas do ICP Expert.
 - **b** Clique em **Connect** (Conectar).
 - c Selecione o instrumento na lista e clique em Connect (Conectar).
- 4 Clique em Calibrate (Calibrar) na seção Detector (Detector).

NOTA Assim que a calibração do detector terminar, a data e a hora da última calibração realizada com êxito serão exibidas, confirmando que a medição do detector foi concluída.

- 5 Certifique-se de que um nebulizador concêntrico de vidro convencional, uma câmara de nebulização de passagem dupla e a tocha apropriada do plasma (para exibição dupla ou radial) estejam instalados. Use a amostragem manual. A tubulação da bomba peristáltica recomendada é branca/branca para a bomba e azul/azul para o dreno.
- 6 Acenda o plasma.
- 7 Verifique se a purga com o multiplicador de pressão do policromador está ativada.
 - Na janela "Instrument" (Instrumento) do software ICP Expert, clique na guia Status e verifique, na seção Polychromator (Policromador), se o Boost (Multiplicador de pressão) está selecionado.
 - b Se o instrumento estiver em estado de espera e for ligado, a calibração de comprimento de onda nas etapas 7-10 poderá ser realizada imediatamente. Se o instrumento for ligado após ter ficado algumas horas sem alimentação elétrica, o policromador poderá levar várias horas para estabilizar a temperatura e fazer a purga.

NOTA

NOTA

- 8 Aspire a solução da calibração do comprimento de onda e selecione a guia **Calibration** (Calibração).
- 9 Clique em **Calibrate** (Calibrar) na seção "Instrument" (Instrumento) para realizar a calibração do comprimento de onda.

Assim que finalizada, a data e a hora da última calibração realizada com êxito será exibida, e uma marca de aprovação ou falha aparecerá juntamente com o erro do comprimento de onda, indicando se a calibração do instrumento foi concluída.

Salvar e visualizar os dados da calibração

Os dados da calibração estão armazenados em Logs (Registros). Para ver os dados, exporte os registros.

- Na janela principal do ICP Expert, clique em File > Logs > Export Logs (Arquivo > Registros > Exportar registros).
- 2 Salve o arquivo de registro compactado.
- **3** Abra o arquivo .zip para ver os arquivos do registro.

O local de salvamento padrão é \Users\'Nome do usuário'\Documents\Agilent\ICP Expert \Exported Results\ na unidade de disco em que o software ICP Expert está instalado.

Substituição de componentes de hardware

O ICP-OES ficará pronto para ser operado assim que o engenheiro de serviços de campo Agilent terminar de instalá-lo. Contudo, você terá de configurar os seguintes itens: bomba peristáltica, nebulizador, tocha ou câmara de nebulização. Para obter instruções para esse procedimento, consulte a seção "How to" (Como) no Help and Learning Center (Centro de treinamento e ajuda) do ICP Expert.

Tochas do ICP-OES

Muitas tochas do ICP-OES estão listadas abaixo, cada uma com propriedades únicas e adequadas para uma ampla gama de aplicações. Tochas adicionais podem estar disponíveis. Consulte o site da Agilent para obter mais informações.

	Tochas de exibição dupla			
Parâmetro	Tocha padrão	Tocha semidesmontável	Tocha inerte semidesmontável	Tocha semidesmontável de alto teor de sólidos
Construção	Uma peça	Duas peças desmontáveis	Duas peças desmontáveis	Duas peças desmontáveis
Diâmetro interno do injetor (mm)	1,8	1,8	1,8	2,4
Material do injetor	Quartzo	Quartzo	Cerâmica	Quartzo
Comprimento do tubo externo	Médio com fenda	Médio com fenda	Médio com fenda	Médio com fenda
Material do tubo externo	Quartzo	Quartzo	Quartzo	Quartzo
Tubo intermediário	Tulipa de quartzo	Tulipa de quartzo	Tulipa de quartzo	Tulipa de quartzo
Aplicação recomendada	Maioria das amostras	Maioria das amostras	Digestões com ácido fluorídrico (HF)	Amostras com alta concentração total de sólidos dissolvidos

Derâmetre	Tochas de visualização radial				
Parametro	Tocha padrão Tocha semidesmontável		Tocha inerte semidesmontável		
Construção	Uma peça	Duas peças desmontáveis	Duas peças desmontáveis		
Diâmetro interno do injetor (mm)	1,4	1,4	1,8		
Material do injetor	Quartzo	Quartzo	Cerâmica		
Comprimento do tubo externo	Curto (sem fenda)	Curto (sem fenda)	Curto (sem fenda)		
Material do tubo externo	Quartzo	Quartzo	Quartzo		
Tubo intermediário	Tulipa de quartzo	Tulipa de quartzo	Tulipa de quartzo		
Aplicação recomendada	Maioria das amostras	Maioria das amostras	Digestões com ácido fluorídrico (HF)		

			Tochas completame	ente desmontáveis	
Parâmetro	Padrão	Inerte	Alta concentração de sólidos	Orgânicos semivoláteis	Orgânicos voláteis
Construção	Completamen te desmontável	Completame nte desmontável	Completamente desmontável	Completamente desmontável	Completamente desmontável
Diâmetro interno do injetor (mm)	1,8	1,8	2,4	1,4	0,8
Material do injetor	Quartzo	Cerâmica	Quartzo	Quartzo	Quartzo
Comprimento do tubo externo	Médio com fenda	Médio com fenda	Médio com fenda	Médio com fenda	Curto (sem fenda)
Material do tubo externo	Quartzo	Quartzo	Quartzo	Quartzo de alta pureza (para solventes orgânicos)	Quartzo de alta pureza (para solventes orgânicos)
Tubo intermediário	Tulipa de quartzo	Tulipa de quartzo	Tulipa de quartzo	Tulipa de quartzo	Tulipa de quartzo
Aplicação recomendada	Maioria das amostras	Digestões com ácido fluorídrico (HF)	Amostras com alto teor total de sólidos dissolvidos	Solventes semivoláteis orgânicos, por exemplo, cetonas, querosene, Jet-A1	Solventes orgânicos voláteis, por exemplo, gasolina

Uma tocha completamente desmontável pode ser adaptada para outra aplicação através da substituição do injetor e/ou o tubo externo pela configuração recomendada para esse tipo de amostra. Consulte o site da Agilent Technologies para obter informações sobre pedidos.

Montagem e desmontagem da tocha

Ambos os modelos de tochas semi e completamente desmontáveis permitem que o conjunto de tubos externos/intermediários fique separado da base/injetor para facilitar a limpeza ou substituição. Além disso, as tochas completamente desmontáveis permitem que o injetor seja removido da base para facilitar a limpeza ou substituição.

Risco químico

AVISO



Superfície quente A tocha e seu compartimento podem ficar extremamente quentes durante a operação do instrumento e permanecem assim por algum tempo após o plasma ser apagado. Aguarde o compartimento do plasma esfriar por, no mínimo, cinco minutos, antes de acessar o compartimento do plasma. Aguarde a tocha esfriar por dois minutos antes de tocar o tubo externo ou tentar desmontar a tocha.

AVISO



Os ácidos nítricos e clorídricos e soluções alcalinas fortes são muito corrosivos e podem causar queimaduras graves quando entram em contato com a pele. O uso de roupas protetoras adequadas é sempre essencial durante o manuseio desses ácidos. Se o ácido entrar em contato com a pele, lave a pele imediatamente com água em abundância e procure um médico.

CUIDADO

Para evitar danos à tocha, sempre tome cuidado ao manuseá-la ou armazená-la. Para evitar danos ao instrumento, não use a tocha se esta estiver danificada ou úmida.



Figura 3. Tocha desmontável, onde: 1. Tocha, 2. Vedante superior e 3. Corpo da tocha

Montagem da tocha semidesmontável

Para montar a tocha:

1 Coloque o vedante superior de volta no conjunto de tubos de quartzo, deslize o vedante limpo de volta no conjunto de tubos a partir da extremidade inferior, certificando-se de que a face plana superior do vedante esteja posicionada voltada para a parte superior do conjunto de tubos. Consulte a Figura 4.



Figura 4. O vedante superior sendo recolocado no conjunto de tubos externos de quartzo

2 Deslize o vedante superior cerca de 1/3 acima da extensão do conjunto de tubos de quartzo e posicione-o de forma que a face superior do vedante fique alinhada com a parte inferior da marca de alinhamento "T", conforme mostrado na Figura 5.



Figura 5. O vedante inferior alinhado com a marca de alinhamento no tubo externo de quartzo

3 Deslize cuidadosamente o conjunto de tubos de quartzo no corpo da tocha, evitando o contato com o injetor.



Figura 6. Tocha inserida no corpo da tocha

4 Posicione o conjunto de tubos de quartzo de forma que a canaleta no corpo da tocha se alinhe com a marca de alinhamento "T" no conjunto de tubos de quartzo, conforme mostrado na Figura 7.



Figura 7. A marca de alinhamento no conjunto de tubos externos de quartzo deve estar alinhada com a canaleta no corpo da tocha

5 Pressione o conjunto de tubos externos de quartzo totalmente no corpo da tocha e pressione o vedante superior para fixar o conjunto de tubos no lugar. Consulte a Figura 8.



Figura 8. Certifique-se de que o vedante esteja encaixado corretamente pressionando firmemente as bordas superiores

Montagem da tocha completamente desmontável

Para montar a tocha:

- 1 Desaparafuse a porca de fixação na base da tocha.
- 2 Coloque o vedante superior de volta na parte superior da base da tocha e verifique se está encaixado corretamente pressionando firmemente as bordas superiores.
- 3 Insira o conjunto de tubos no corpo da tocha e posicione-o de forma que a canaleta no corpo da tocha fique alinhada com a marca de alinhamento "T" no conjunto de tubos. Pressione totalmente o conjunto de tubos externos no corpo da tocha. Consulte a Figura 7 e a Figura 8.
- 4 Insira totalmente o injetor na base do corpo da tocha até que a saliência entre em contato com o braço de retenção e "clique" na posição.



5 Substitua a porca de fixação e aperte para prender o injetor.

CUIDADO Aperte apenas manualmente. Não aperte demais nem use uma chave como alavanca.

Desmontagem da tocha

Para desmontar uma tocha semidesmontável:

Poderá ser necessária uma chave de fenda de ponta chata para soltar o anel de proteção, caso não seja possível desencaixar manualmente a tocha do corpo.

CUIDADO Use a chave de fenda para soltar o anel de proteção da tocha do corpo da tocha somente se não for possível fazê-lo manualmente.

Solte cuidadosamente o vedante superior pressionando-o com a unha do polegar (consulte a Figura 9) ou use uma chave de fenda de ponta chata para levantá-lo. Ao usar uma chave de fenda, insira a chave de fenda no local indicado, como mostrado na Figura 10, mantendo uma ação de cima para baixo para soltar o vedante. Repita a ação no lado oposto. O vedante deve ser separado do corpo da tocha.



Figura 9. Remoção do vedante superior de uma tocha semidesmontável pressionando-o com uma unha de polegar


Figura 10. Remoção do vedante superior com a ajuda de uma chave de fenda de ponta chata

2 Mantendo um movimento paralelo reto, deslize com cuidado o conjunto de tubos externos de quartzo para fora do corpo da tocha. Consulte a Figura 11.



Figura 11. Remoção do conjunto de tubos de quartzo do corpo da tocha

3 Para remover o vedante superior do conjunto de tubos de quartzo, deslize o vedante em direção à extremidade superior do conjunto de tubos. Consulte a Figura 12.



Figura 12. Remoção do vedante superior do conjunto de tubos de quartzo

Para desmontar a tocha completamente desmontável:

1 Desaparafuse a porca de fixação na base da tocha e remova o injetor puxando cuidadosamente a junta esférica.



2 Remova o conjunto de tubos de quartzo da parte superior da base da tocha. Se o vedante superior for removido com o conjunto de tubos externos de quartzo, deslize o vedante superior do conjunto de tubos em direção à extremidade superior.



Veja na página 59 o procedimento de limpeza da tocha.

Acessórios

Os acessórios a seguir podem estar disponíveis para uso com seu instrumento ICP-OES:

- Amostrador automático SPS 4
- Sistema avançado de válvulas (AVS) válvula de alternância AVS 4, AVS 6 e AVS 7
- Adaptador de duto de entrada externo
- Acessório de geração de vapor (VGA)
- Sistema de introdução de amostras multimodal (MSIS)
- Umidificador de argônio
- Câmara de nebulização de temperatura programável IsoMist

Para obter informações sobre instalação e segurança, consulte a documentação do acessório ou o Help and Learning Center (Centro de treinamento e ajuda) do ICP Expert.

SPS 4

Consulte as instruções fornecidas com o acessório para obter informações de segurança e para preparar a instalação do SPS 4.

O SPS 4 é compatível com uma ampla linha de racks de amostras autoclaváveis de baixo custo e comercialmente disponíveis.

A contaminação da amostra a partir de partículas aéreas é eliminada, e vapores corrosivos ou tóxicos são removidos durante a amostragem com o gabinete opcional para o ambiente.

Sistema avançado de válvulas (AVS) - AVS 4, AVS 6 e AVS 7

Para obter informações sobre segurança e como preparar o acessório para a instalação, consulte as instruções fornecidas com o acessório em questão.

A AVS 4 integrada (válvula seletora de 4 portas) enxágua o sistema de introdução de amostra enquanto a próxima amostra está sendo apresentada ao instrumento antes da medição. O excesso de amostra é desviado da câmara de nebulização e o enxágue é introduzido imediatamente depois da medição. Isso reduzirá o carryover da amostra e a frequência de limpeza necessária dos componentes de introdução de amostra.

A AVS 6 (válvula seletora de 6 portas) integrada e a AVS 7 (válvula seletora de 7 portas) aumentam a produção analítica, diminuem o tempo de espera e os custos operacionais. A válvula seletora está localizada entre o nebulizador e a bomba peristáltica do espectrômetro. As amostras são rapidamente carregadas no loop de amostra e ficam prontas para análise imediata pelo ICP-OES, reduzindo consideravelmente os atrasos na coleta de amostras. O enxágue preventivo da linha de amostragem indica que os tempos de análise de uma amostra para a seguinte foram reduzidos.

O AVS 6 e o AVS 7 contêm um injetor de bolhas que injeta bolhas automaticamente assim que a amostra é carregada no ciclo, isolando a amostra da solução de enxágue/portadora. Esse processo reduz o volume de amostra necessário para a medição (reduzindo, portanto, o tempo de análise), pois os efeitos de restos (ou de diluição) são minimizados.

Adaptador de duto de entrada externo

O adaptador de duto de entrada externo oferece um acoplamento para conduzir ar para a porta de entrada de ar, para uso em laboratórios com ambientes mais hostis.

Acessório de geração de vapor (VGA)

O VGA 77P é um acessório de geração de vapor de fluxo contínuo para instrumentos de espectroscopia de emissão óptica com plasma acoplado indutivamente (ICP-OES) que detecta o Hg e os elementos de formação de hidretos em concentrações de partes por bilhão (ppb). Todo o conjunto de encanamento está integrado a um módulo separado e pode ser simplesmente trocado durante a substituição dos elementos.

Sistema de introdução de amostras multimodal (MSIS)

Para obter informações sobre segurança e como preparar o MSIS para instalação, consulte as instruções fornecidas com este acessório e a Ajuda do ICP Expert.

O MSIS é usado com o instrumento ICP-OES para fornecer geração simultânea de vapor de vários elementos de formação de hidretos, permitindo a determinação com limites baixos de detecção de ppb.

O MSIS consiste em um nebulizador e uma câmara de nebulização ciclônica de vidro modificado que contém dois tubos cônicos verticais no centro da câmara. Isso permite que o redutor e a amostra se misturem de modo rápido e completo na câmara por meio da tecnologia de hidretos de película fina para formar os hidretos.

O MSIS pode ser operado de três modos: somente hidreto, hidreto simultâneo e nebulização convencional ou somente nebulização convencional.

Umidificador de argônio

O umidificador de argônio costuma ser usado na execução de amostras aquosas cujo conteúdo é composto por elementos sólidos ou sais altamente dissolvidos. Ao utilizar o acessório, o fluxo de gás de nebulização passa pelo umidificador para aumentar o vapor de água no gás. Descobriu-se que esse procedimento é benéfico ao reduzir o acúmulo de sal e outros elementos sólidos dissolvidos da ponteira do nebulizador no sistema de introdução de amostra. O umidificador de argônio, ao reduzir as obstruções, ajuda a garantir uma operação ininterrupta e livre de manutenções.

Câmara de nebulização de temperatura programável IsoMist

A IsoMist utiliza o efeito Peltier para fornecer resfriamento ou aquecimento elétrico à câmara de nebulização.

A "temperatura definida" é monitorada e controlada pelo software ICP Expert. O desempenho da IsoMist pode ser monitorado e controlado por meio de um computador de laboratório por USB ou tecnologia sem fio Bluetooth.

Lista de verificação de análises	43
Ligar o instrumento e o software	44
Preparar para análise	46
Executar as calibrações do detector e do comprimento de onda	46
Criar/abrir uma worksheet	47
Desenvolver um método	48
Executar amostras	50
Imprimir um relatório	50
Desligar o instrumento	51

Este capítulo oferece um breve guia para configurar o instrumento e executar amostras.

Você encontrará instruções passo a passo para operações comuns na Ajuda do ICP Expert. Para acessar essas informações:

- 1 Clique duas vezes no botão Help and Learning Center (Centro de treinamento e ajuda) do ICP Expert na área de trabalho.
- 2 Quando a Ajuda do ICP Expert for exibida, clique em **How to** (Como fazer) para ver as instruções passo a passo disponíveis.

Lista de verificação de análises

Você precisa realizar as seguintes etapas para medir uma ou mais amostras. Você encontrará informações sobre cada etapa neste capítulo.

- Ligar o computador, o instrumento e o software
- Conectar o software ao instrumento
- Preparar para análise

- Realizar uma varredura de corrente escura e uma calibração de comprimento de onda
- Criar/abrir uma worksheet
- Desenvolver um método
- Executar uma amostragem
- Imprimir um relatório

Ligar o instrumento e o software

Antes de iniciar o sistema, leia atentamente a seção Riscos e práticas de segurança no início deste manual para garantir que seu laboratório esteja configurado de acordo com os detalhes especificados no Guia de preparo do local.

Ligar o instrumento pela primeira vez (ou após desligamento prolongado)

Como ligar o instrumento pela primeira vez ou após o desligamento:

- 1 Verifique se as linhas de exaustão e a do adaptador de duto de entrada externo (se usado) estão devidamente conectadas ao instrumento ICP-OES e se o filtro de ar não está obstruído.
- 2 Ligue o sistema de exaustão do laboratório.
- **3** Certifique-se de que as linhas de gás e de água estejam conectadas ao instrumento ICP-OES.
- 4 Ligue o suprimento de gás.
- 5 Ligue o resfriador de água.
- 6 Verifique se os fornecimentos de gás e de água estão ligados e configurados com as pressões certas, e se o resfriador de água foi configurado com a temperatura correta.
- 7 Verifique se a tocha está limpa, em boas condições e instalada com a respectiva alça completamente fechada.
- 8 Verifique se todos os tubos na câmara de nebulização, nebulizador e bomba peristáltica estão instalados e conectados corretamente.

- 9 Verifique se a porta do compartimento do plasma está totalmente fechada.
- 10 Verifique se o cabo de rede LAN Ethernet está conectado ao computador ou à rede de área local.
- **11** Ligue o computador, o monitor e a impressora.
- 12 Conecte o cabo do ICP-OES à tomada na parede e coloque o botão de alimentação elétrica no lado esquerdo do instrumento em "On" (Ligado).
- 13 Pressione o botão Liga/Desliga na parte frontal do instrumento. O LED Liga/Desliga acenderá em verde quando tudo estiver ligado. O ICP-OES estará em estado de espera, realizando purga contínua e controlando o policromador pelo termostato.
- NOTA Para que o instrumento funcione, tanto o botão de alimentação elétrica no lado esquerdo do instrumento quanto o botão Liga/Desliga no painel frontal devem estar ligados.
 - 14 Clique duas vezes no ícone do ICP Expert na área de trabalho para iniciar o software ICP Expert. A página Start (Início) será exibida.

Operar o instrumento em estado de espera

Durante o estado de espera (com ambos os botões de energia ligados), a temperatura do policromador é regulada e a purga de gás do policromador é realizada para garantir que o instrumento fique pronto para análise 20 minutos depois que o plasma for aceso.

Como começar a usar o instrumento a partir do estado de espera:

- 1 Ligue o sistema de exaustão do laboratório.
- 2 Verifique se a tocha está limpa, em boas condições e instalada com a respectiva alça completamente fechada.
- **3** Verifique se todos os tubos na câmara de nebulização, nebulizador e bomba peristáltica estão conectados corretamente.
- 4 Verifique se a porta do compartimento da tocha está totalmente fechada.
- **5** Ligue o monitor e a impressora (se estiverem desligados).
- 6 Ligue o resfriador de água (se estiver desligado).
- 7 Se houver acessórios conectados, ligue-os.

Preparar para análise

Como preparar para análise:

1 Clique no botão **Plasma** no software ICP Expert. Como alternativa, pressione F5 ou selecione **Plasma on** (Plasma aceso) na seta do botão Plasma.

NOTA A sequência de ignição do plasma levará até 60 segundos para ser concluída. Se a ignição do plasma falhar, consulte a seção Troubleshooting (Solução de problemas) no Help and Learning Center (Centro de treinamento e ajuda) para obter mais informações.

NOTA

Para obter estabilidade e desempenho ideais, após a ignição do plasma, é recomendável aguardar um período de 20 minutos para que o instrumento ICP-OES aqueça.

Caso estejam sendo usados comprimentos de onda abaixo de 189 nm: a purga do policromador pode durar muitas horas. A estabilização térmica do policromador pode durar muitas horas caso o instrumento tenha ficado desligado por muito tempo.

- 2 Confira se a bomba peristáltica está configurada corretamente (consulte a seção Peristaltic Pump [Bomba Peristáltica] no Help and Learning Center (Centro de treinamento e ajuda) do ICP Expert). Caso ainda não tenha feito, ajuste as barras de pressão na bomba peristáltica para igualar o fluxo de amostra e a drenagem.
- **3** Coloque a tubulação de amostra da bomba peristáltica na solução de enxágue e a tubulação de drenagem no recipiente de drenagem.
- 4 Clique no botão Pump (Bomba) no software ICP Expert e escolha Normal (15 rpm) na seta abaixo do botão Pump (Bomba). A bomba será inicializada e a solução começará a ser aspirada.

Executar as calibrações do detector e do comprimento de onda

Consulte a página 27 para obter instruções.

Criar/abrir uma worksheet

Criar uma nova worksheet

Para criar uma nova worksheet, clique em New Quantitative (Novo quantitativo) ou Intelliquant Screening (Varredura Intelliquant) na página Start (Início) ou no menu File (Arquivo).

Será apresentada uma lista de arquivos usados recentemente quando for criada uma nova worksheet a partir de um modelo; caso contrário, é possível Browse (Procurar) mais arquivos. A caixa de diálogo "New From Template" (Novo a partir do modelo) será exibida neste caso.

Abrir uma worksheet existente

Como abrir uma worksheet existente:

- 1 Clique em Open (Abrir) na página Start (Início) ou no menu File (Arquivo).
- 2 Será exibida uma lista de arquivos usados recentemente. Caso contrário, é possível **Browse** (Procurar) mais arquivos. A caixa de diálogo **Open** (Abrir) será exibida neste caso.

Criar uma nova worksheet a partir de um modelo

Para criar uma nova worksheet a partir de um modelo, clique em **New From** (Novo a partir de) na página Start (Início) ou **New From Template** (Novo a partir de um modelo) no menu File (Arquivo).

Será apresentada uma lista de arquivos usados recentemente; caso contrário, é possível **Browse** (Procurar) mais arquivos. A caixa de diálogo New From Template (Novo a partir do modelo) será exibida neste caso.

A janela Worksheet será exibida com a nova worksheet carregada.

Desenvolver um método

NOTA

Para obter instruções sobre como configurar e usar acessórios, o IntelliQuant e as opções Extend QC, consulte o Help and Learning Center (Centro de treinamento e ajuda) do ICP Expert.

Como desenvolver um método:

- 1 Abra uma nova planilha ou use um modelo.
- 2 Selecione os recursos e acessórios na página Configuration (Configuração) a serem usados durante a análise. (Alguns recursos estão disponíveis apenas no software ICP Expert Pro).
- 3 Na página "Elements" (Elementos), selecione os elementos do menu suspenso "Element" (Elemento) ou digite o nome do elemento ou o símbolo e, em seguida, execute um dos métodos abaixo:
 - Clique em Add Analyte (Adicionar analito) para adicionar o comprimento de onda primário ao elemento selecionado.
 - Destaque o comprimento de onda que você deseja usar na lista de comprimentos de onda disponíveis. Clique em Add (Adicionar).
- NOTA Como alternativa, pressione CTRL e selecione o elemento da tabela periódica para adicionar o comprimento de onda primário ao elemento selecionado para o seu método.

O elemento será exibido na tabela com o comprimento de onda e as configurações padrão selecionadas.

- NOTA É recomendável adicionar vários comprimentos de onda a cada elemento. Resultados semelhantes para cada comprimento de onda indicam que não há interferências presentes. Se houver resultados diferentes, verifique o IntelliQuant para informar o comprimento de onda correto.
 - 4 Verifique se não há interferentes conhecidos ou outras linhas analíticas próximo à linha analítica selecionada. Sua intensidade relativa regulará a proximidade entre as linhas. Por exemplo, se sua matriz contém um elemento que não é de interesse analítico, mas é um interferente potencial e possui uma linha próxima a uma de suas linhas analíticas, a concentração desse elemento na sua matriz determinará se você precisa ou não escolher uma outra linha analítica.

- 5 Faça os ajustes necessários em cada elemento incluindo a seleção de um comprimento de onda diferente, inserindo informações adicionais na coluna "Label" (Legenda) e selecionando o tipo (analito, padrão interno ou interferente).
- 6 Clique **Conditions** (Condições) para modificar ambas as definições comuns de corrida e configurações para cada elemento. É possível usar até oito condições de medidas diferentes.
- 7 Clique em QC (CQ Controle de Qualidade) para inserir os limites de detecção do método, selecione os testes de CQ a serem usados e quais ações devem ser tomadas em caso de erro.
- 8 Clique em IEC para inserir a concentração dos elementos em seus padrões de analitos e interferentes.
- 9 Clique em Standards (Padrões) para inserir a concentração dos elementos em seus padrões e selecionar se serão usadas outras opções, como Standard Additions (Adições de padrão) ou MultiCal. Além disso, selecione também se será usada a solução em branco nas calibrações e se o reslope (cálculo da inclinação da curva da calibração) deverá ser habilitado.
- 10 Clique em Sequence (Sequência) para especificar o número de amostras, inserir os testes de CQ, selecionar o tipo de solução e editar as etiquetas das amostras e ações no final da corrida.
- **11** Clique em **FACT** para criar modelos FACT manualmente ou configurar o método para que eles sejam criados automaticamente durante a análise.
- 12 Se você estiver usando um amostrador automático, clique na guia "Autosampler" (Amostrador automático) para selecionar os racks e a profundidade da sonda (se necessário). As opções variam de acordo com o amostrador automático selecionado.
- NOTA Consulte o Help and Learning Center (Centro de treinamento e ajuda) do ICP Expert para obter uma descrição mais detalhada sobre a configuração de um método.

Executar amostras

Como executar amostras:

- 1 Configure o amostrador automático com todas as soluções, se necessário.
- 2 Clique na guia Analysis (Análise) e faça o seguinte:
 - a Confira se suas amostras estão selecionadas. A seleção é indicada por uma marca de seleção ao lado da coluna Rack:Tube (Rack:Tubo). Para selecionar todas as soluções, marque a caixa de seleção ao lado do título Rack:Tube (Rack:Tubo).

\bigcirc	Rack: Tube	Sample Label
~	S1:1	Blank
~	S1:2	Standard 1
~	S1:3	Standard 2

b Clique no ícone **Run** (Executar) na barra de ferramentas (ou pressione as teclas SHIFT+F8) para iniciar a análise e siga as instruções.

Para obter mais informações sobre como executar uma análise, consulte o Help and Learning Center (Centro de treinamento e ajuda) do ICP Expert.

Imprimir um relatório

Como imprimir um relatório:

- 1 Clique em **Report** (Relatório) na barra de ferramentas ou em **File > Report** (Arquivo > Relatório).
- 2 Escolha se deseja imprimir, visualizar ou salvar o relatório como arquivo PDF.

RECOMENDAÇÃO A visualização do relatório assegura-lhe que você incluiu todos os dados que você necessita.

- 3 Selecione um modelo de relatório e clique em Open (Abrir).
- 4 Clique no botão **Print** (Imprimir) para gerar um relatório, conforme especificado. Será exibida uma mensagem indicando brevemente o status do andamento do relatório em criação.
- 5 Será aberta a caixa de diálogo Print (Imprimir). Você pode optar por imprimir todo o relatório ou um intervalo de páginas. Os relatórios são impressos em sua impressora padrão, exceto se especificado o contrário. Você pode especificar as opções da impressora na caixa de diálogo Print Setup (Configurar impressão), acessada no menu File (Arquivo).

Desligar o instrumento

Existem três modos de espera e desligamento:

- Modo de espera com argônio ativado para uso rotineiro (Página 52)
- Modo de espera com argônio desativado para desligamento durante a noite, fins de semana e prolongado (Página 52)
- Desligamento total para um desligamento completo (Página 54)

Modo de espera do instrumento com argônio ativado

Esta opção do modo de espera deixa o ICP-OES 5800/5900 em estado de prontidão, consumindo apenas uma pequena quantidade de insumos (gás argônio e eletricidade). O benefício de estar no modo de espera é que o 5800/5900 pode estar operacional novamente em <20 minutos após a ignição do plasma. O modo de espera é ativado automaticamente assim que a análise anterior é concluída e o plasma é desligado. Siga as instruções "Configurar o instrumento no modo de espera com argônio ativado ou desativado" na página 52 para garantir que o instrumento esteja totalmente preparado para o modo de espera.

A recomendação típica para o modo de espera é deixar a energia do ICP-OES e o suprimento de gás argônio pressurizados, permitindo uma purga de argônio por gotejamento ou filete através do policromador, sem coletar dados analíticos.

Para voltar ao status operacional a partir deste modo, ligue o plasma e estabilize por 20 minutos.

Modo de espera do instrumento com argônio desativado

Para desligamento durante a noite, fins de semana e prolongado, recomenda-se que o instrumento esteja totalmente energizado, mas que o plasma esteja apagado. Deste modo, os sistemas de purga também são desativados. O sistema interno de resfriamento a ar opera com capacidade reduzida.

Esta opção do modo de espera reduzirá ainda mais o consumo de energia. O ICP-OES pode ficar com a energia ligada e o gás argônio pode ser completamente desligado. Siga as instruções "Configurar o instrumento no modo de espera com argônio ativado ou desativado" na página 52 para garantir que o instrumento esteja totalmente preparado para o modo de espera.

Para voltar ao status operacional a partir do gás desligado, ative o software e ligue o gás com o ICP-OES ligado e deixe purgar por 30 minutos antes da ativação do resfriador de água ou plasma. Essa purga de 30 minutos é necessária para remover qualquer vapor de água atmosférico da ótica.

NOTA Não é recomendável desativar o argônio, especialmente se você estiver analisando rotineiramente linhas abaixo de 190 nm. Desativar o gás argônio reduzirá o consumo de energia; no entanto, o ICP-OES levará mais tempo para retornar ao status operacional para linhas de UV profundas quando o gás argônio for ativado novamente.

NOTA Se você desativar o gás argônio, também desative o resfriador de água.

Configurar o instrumento no modo de espera com argônio ativado ou desativado

Como colocar o instrumento no modo de espera:

- 1 Enxágue a câmara de nebulização aspirando água por alguns minutos.
- NOTA Para a execução de amostras orgânicas recomenda-se que a câmara de nebulização esteja limpa e completamente seca entre as análises.
 - 2 Apague o plasma clicando no ícone Plasma Off (Plasma apagado), pressionando SHIFT + F5 no teclado ou escolhendo Plasma Off (Plasma apagado) no menu Analyze (Analisar). A bomba peristáltica para automaticamente quando o plasma é apagado.
 - 3 Desative o multiplicador de pressão do policromador durante a noite, desmarcando Polychromator > Boost (Policromador > Multiplicador de pressão) na aba Status da janela Instrument (Instrumento).
 - 4 Para o modo de espera apenas com argônio desativado: Desligue o suprimento de gás argônio.

- **5** Para aumentar a vida útil da tubulação da bomba, afrouxe os tubos da bomba peristáltica liberando as barras de pressão e levante os tubos para fora das canaletas. Para fazer isso:
 - a Levante os parafusos da barra de pressão. Isso os libera da barra de pressão (consulte a Figura 13).
 - **b** Deixe a barra de pressão solta.
 - c Levante os tubos para fora das canaletas.



Figura 13. Bomba peristáltica com as barras de pressão levantadas e a barra de pressão solta, permitindo que a tubulação seja liberada

- 6 Desligue o resfriador de água.
- 7 Feche a planilha clicando em "Close" (Fechar) no menu "File" (Arquivo), mas deixe o software ICP Expert em execução. Você pode desligar a impressora, o monitor e quaisquer acessórios, se desejar.

Verifique se a opção de economia de energia no seu PC está desativada (isso impedirá o desligamento do seu disco rígido). Se essa opção não estiver desativada, você poderá perder dados durante um desligamento inesperado.



Fumaça nociva

Perigo de inalação. O sistema de exaustão DEVERÁ permanecer ligado se o suprimento de gás estiver ligado.

Desligar o instrumento para desligamento prolongado

Para longos períodos de desuso, desligue completamente o instrumento. (Isso desativa toda a purga e o sistema de termostato do policromador).

Para desligar completamente seu instrumento:

CUIDADO

A unidade de controle de gás ICP-OES fornece purga ininterrupta de gás ao conjunto do policromador apenas durante os estados de funcionamento e inatividade somente para minimizar a entrada de umidade. Recomenda-se que os suprimentos de instrumentos e de gás permaneçam sempre ligados, exceto durante longos períodos sem uso.

- **1** Siga as etapas 1–7 do procedimento do modo de espera.
- 2 Desligue todos os acessórios (quando aplicável) e, em seguida, desligue o suprimento de gás argônio no cilindro.
- **3** Desligue a energia na parte frontal inferior esquerda do instrumento.
- 4 Aguarde o LED verde do botão frontal Liga/Desliga parar de piscar e, em seguida, desligue o botão de alimentação elétrica no lado esquerdo do instrumento. Isso desligará completamente o instrumento, inclusive o sistema de termostato do policromador.
- 5 Desligue o sistema de exaustão do laboratório.
- 6 Saia do software ICP Expert, se não for mais necessário, escolhendo Exit (Sair) no menu File (Arquivo). Desligue a impressora e o monitor.

Reiniciar o instrumento para estar pronto para análise novamente levará várias horas devido ao tempo de purga de gás e de estabilização térmica do policromador. Consulte a página 44 para obter instruções completas de inicialização.

NOTA Se o instrumento não for utilizado durante algum tempo, a tocha, o cone, a ponta cônica e o compartimento da tocha devem ser limpos de quaisquer depósitos, sujeira ou resíduo. Verifique a limpeza da janela axial removível quando o cone for removido.

Rotina	56
Limpeza	58
Limpeza da tocha	59
Solução de problemas	68
Peças sobressalentes	68
Suporte técnico	69

Este capítulo inclui os requisitos de manutenção do ICP-OES Agilent que podem ser realizados por um operador. Todos os procedimentos de manutenção não especificamente mencionados no presente capítulo devem ser efetuados apenas por engenheiros de serviço de campo treinados, qualificados ou autorizados pela Agilent.

AVISO



Risco para os olhos

Risco de choque

Sempre use óculos de segurança apropriados ao manusear soluções de amostra e outras substâncias químicas, ou quando o plasma estiver ativo, para minimizar o risco de ferimento aos olhos causados por líquidos perigosos e exposição aos raios ultravioleta.

AVISO



Este instrumento contém circuitos elétricos, dispositivos e componentes que funcionam com tensões perigosas. O contato com esses circuitos, dispositivos e componentes pode resultar em morte, provocar ferimentos graves ou choque elétrico doloroso.



Risco de RF e superfícies quentes

O plasma irradia níveis perigosos de energia de radiofrequência (RF). A exposição à energia de RF pode causar danos graves à pele e cataratas nos olhos, enquanto o contato próximo com o plasma em operação pode resultar em queimaduras graves na pele e em descarga elétrica que percorre distância considerável e é capaz de causar morte, choque elétrico grave ou queimaduras na derme da pele.

NOTA

Esta seção contém procedimentos de manutenção do instrumento ICP-OES. Você deve consultar os manuais dos acessórios, da impressora e do computador, para saber quais são seus procedimentos de manutenção, e a Ajuda do ICP Expert, para saber quais são os procedimentos de manutenção de todos os acessórios solicitados por você.

Rotina

As seguintes peças do ICP-OES, consumíveis e acessórios, necessitam de manutenção de rotina. As instruções de manutenção estão incluídas no Help and Learning Center (Centro de treinamento e ajuda) do ICP Expert. Para acessar essas instruções, clique duas vezes no ícone do Help and Learning Center (Centro de treinamento e ajuda) do ICP Expert na área de trabalho. Clicar ou pesquisar **Maintenance** (Manutenção).

A frequência de execução dessas tarefas de manutenção de rotina depende da frequência de uso e dos tipos de amostras analisadas. Com matrizes de amostra desafiadoras ou para sistemas operando continuamente, é necessária uma manutenção mais frequênte.

De hora em hora

□ Verifique e, se necessário, esvazie o recipiente de drenagem.

Diariamente

- Verifique o sistema de exaustão e a pressão do gás argônio.
- Verifique o nível de água do umidificador de argônio antes de cada uso (se aplicável).
- □ Limpe a superfície do ICP-OES (derramamentos devem ser limpos imediatamente).
- Inspecione a tocha quanto a obstruções do injetor ou outros danos.
- Verifique o nebulizador quanto a obstruções ou pulsações durante a operação.
- Inspecione a tubulação da bomba e substitua se estiver achatada ou se tiver perdido a elasticidade. Solte os tubos da bomba quando ela não estiver em uso.
- Faça uma verificação visual da câmara de nebulização para confirmar se todas as conexões estão seguras e sem vazamento e se os resíduos estão sendo drenados corretamente.

Semanalmente

- Limpe a tocha.
- □ Inspecione o cone (axial) ou a ponta cônica (radial); limpe se necessário.
- □ Verifique os outros tubos de introdução de amostras e O-rings.
- Substitua a tubulação da bomba peristáltica.
- Limpe o nebulizador.

Mensalmente

- Limpe a câmara de nebulização.
- Limpe o nebulizador.
- Verifique a outra introdução de amostras, tubos de transferência e O-rings (procure desgaste excessivo, má vedação ou dobras e substitua conforme necessário).
- Inspecione as janelas de pré-óptica axial e radial removíveis quanto à limpeza. Limpe ou substitua conforme necessário.

- Verifique as condições do filtro na entrada de ar na parte superior do seu instrumento. Remova o conjunto do filtro de entrada de ar e substitua o elemento do filtro se houver acúmulo de poeira e sujeira.
- Verifique o nível de água no resfriador de água (consulte o manual fornecido com o resfriador de água para obter detalhes).
- Verifique/limpe o trocador de calor (radiador) no resfriador para remover acúmulos de poeira e sujeira.
- **D** Realize uma calibração de comprimento de onda.
- Inspecione o sistema de suprimento de gás externo quanto a vazamentos, incluindo a tubulação conectada ao instrumento, e fissuras por tensão. Substitua todos os componentes danificados, desgastados ou com vazamento.

Anualmente

- Drene e substitua o líquido refrigerador.
- Substitua o filtro de gás da purga de argônio.
- Remova e limpe o filtro de água do lado direito do instrumento.
- Periodicamente, drene o líquido refrigerador do sistema de refrigeração e, em seguida, encha/trate com um algicida apropriado (conforme recomendação do fabricante).

Limpeza

Todo derramamento no compartimento de amostra deverá ser limpo imediatamente.

Os usuários (ou outras pessoas autorizadas) deverão realizar o procedimento de descontaminação adequado se for derramado material perigoso sobre ou dentro do ICP-OES.

As superfícies exteriores do ICP-OES devem ser mantidas limpas. Toda a limpeza deve ser feita com um pano macio. Se necessário, esse pano pode ser umedecido com água ou detergente suave. Não use solventes orgânicos ou agentes de limpeza abrasivos.

Antes de usar qualquer produto de limpeza, processo ou método de descontaminação, exceto aqueles especificados pela Agilent, os usuários (ou outras pessoas autorizadas) devem consultar o representante ou o engenheiro de serviços de campo local autorizado pela Agilent para confirmar que o método proposto não danificará o equipamento.

Limpeza da tocha

Para ajudá-lo a alcançar a vida útil máxima de todos os modelos da tocha de encaixe simples do ICP-OES 5800/5900, recomenda-se que os procedimentos de limpeza sejam seguidos assim que surgir qualquer descoloração no tubo externo da tocha.

Para maximizar a vida útil da tocha e evitar a contaminação, é necessário inspecionar a condição e a limpeza da tocha diariamente.

Consulte a Ajuda do ICP Expert para obter informações sobre como limpar a tocha de amostras a base de matéria orgânica e ter a acumulação de carbono no injetor.

Para acessar a Ajuda, clique duas vezes no ícone Ajuda do ICP Expert na área de trabalho do computador.

CUIDADO Não use fio para limpeza ou abrasivos, como escovas ou lã de aço, para limpar a tocha.

AVISO



Superfície quente e risco químico

A tocha e seu compartimento ficam extremamente quentes durante a operação do instrumento e permanecem assim por algum tempo após o instrumento ser desligado. Aguarde a tocha e seu compartimento esfriarem por pelo menos cinco minutos antes de tentar removê-la. Use luvas resistentes ao calor.

Os ácidos nítricos e clorídricos e soluções alcalinas fortes são muito corrosivos e podem causar queimaduras graves quando entram em contato com a pele. A preparação de soluções ácidas de imersão e qualquer imersão devem ser feitas sob um exaustor para a fumaça. O uso de roupas protetoras adequadas é sempre essencial durante o manuseio desses ácidos. Se o ácido entrar em contato com a pele, lave a pele imediatamente com água em abundância e procure um médico.

Importante

Use um béquer limpo, de diâmetro largo e topo aberto (de preferência com 100 mL de forma alta), ou um recipiente semelhante, para as soluções de imersão. Use o suporte de limpeza da tocha recomendado (consulte a Figura 15A) para as tochas de encaixe simples do ICP-OES 5800/5900 (número de peça G8010-68021) para manter a tocha invertida durante a imersão no processo de limpeza. O suporte de limpeza da tocha ajuda na limpeza da tocha, evitando danos à extremidade frágil do tubo externo de quartzo e limitando a exposição dos materiais de elastômero no corpo da tocha à fumaça ácida, reduzindo a degradação prematura.

- Use detergentes ou ácidos limpos/sem partículas para realizar a imersão.
- Certifique-se de que o ácido não entre em contato com o vedante entre o tubo externo de quartzo e a base plástica.
- Não molhe o vedante superior removível da tocha semidesmontável.
- Mantenha a tocha na vertical, com a junta esférica na parte superior, durante todas as etapas de limpeza e enxágue, a menos que seja indicado de outra forma.
- Não remova a tampa adesiva da tocha que cobre o vedante entre o quartzo e o plástico na tocha de peça única.
- Para limpar melhor, sempre desmonte a tocha semi ou completamente desmontável durante os procedimentos de enxágue e secagem.

Procedimento de limpeza por imersão com ácido para a tocha em amostras de base aquosa ou ácida (tochas de peça única e semidesmontáveis)



Consulte a página 30 para obter instruções de desmontagem da tocha desmontável.

Figura 14. Componentes da tocha, onde 1. Junta esférica, 2. Portas de gás, 3. Anel de proteção da tocha (apenas tocha desmontável) e 4. Tubo externo da tocha

Para limpar a tocha:

- 1 Prepare uma solução de 50% de água régia (1 parte de água deionizada para 1 parte de água régia) em um béquer aberto de grande diâmetro. Para fazer a água régia, combine 1:3 de ácido nítrico concentrado:ácido clorídrico.
- 2 Coloque o béquer que contém a solução ácida sob o suporte de limpeza da tocha.
- 3 Mergulhe as peças de quartzo da tocha em água régia a 50% por, no mínimo, 1 hora. O período de tempo necessário no procedimento de limpeza dependerá da extensão da contaminação. Não deixe a tocha no ácido por mais de 4 horas. Se os depósitos permanecerem após o uso de 50% de água régia, repita o processo de limpeza usando uma concentração mais alta de água régia.
- 4 Coloque a tocha na solução de água régia através do orifício fornecido no suporte de limpeza da tocha, como mostrado na Figura 15A, B, C e D. Verifique se o tubo externo de quartzo junto com o injetor está imerso na solução.

Para mergulhar o injetor sem o conjunto de tubos externos de quartzo, a placa do tubo injetor da tocha deve ser instalada no suporte de limpeza da tocha para minimizar a exposição do corpo da tocha à fumaça ácida.

Para tochas semi ou completamente desmontáveis, o conjunto de tubos externos de quartzo pode ser removido e embebido separadamente em um béquer de água régia (Figura 15D).

5 Pipete parte do ácido através da junta esférica do injetor para remover o acúmulo da parte inferior do injetor. Consulte a Figura 16A.

CUIDADO Evite o contato ácido com o vedante onde o quartzo encontra a base de plástico. Podem ocorrer danos aos vedantes e ao corpo da tocha.

- NOTA A Agilent possui um suporte para limpeza da tocha, que ajuda a manter a tocha de pé e estável durante a limpeza. Consulte o website de Agilent www.agilent.com para maiores detalhes.
 - 6 Mantenha a tocha de peça única invertida em todas as etapas de limpeza e enxágue, a menos que seja indicado de outra forma.



Figura 15A. Mergulhe a tocha na solução de água régia. Verifique se o injetor está imerso na solução.



Figura 15B e C. Suporte de limpeza da tocha montado com placas substituíveis para a tocha. Imersão do injetor em água régia.



Figura 15D. Imersão do conjunto de tubos externos de quartzo em água régia.

- 7 Enxágue a tocha e o conjunto de tubos externos e seque-os completamente antes de usar a tocha no instrumento.
- 8 Siga as etapas descritas abaixo "Lavagem da tocha" e "Secagem da tocha".

Lavagem da tocha

Para enxaguar a tocha (tochas de peça única e semidesmontáveis):

- 1 Segure a tocha com o conector da junta esférica para cima.
- 2 Lave minuciosamente a parte interna e externa da tocha com água deionizada (18 MΩ-cm) usando uma frasco de lavagem para direcionar o fluxo de água. Consulte a Figura 16A.
- 3 Inverta a tocha (consulte a Figura 16B), de modo que o tubo de quartzo ou o injetor fiquem para cima e o conector da junta esférica fique para baixo. Introduza a água de enxágue nos tubos de quartzo, para que a água flua através das portas de entrada de gás e do conector da junta esférica por, no mínimo, 1 minuto.



Figura 16A. Enxágue o injetor com água deionizada através do conector da junta esférica.



Figura 16B. Lave as portas de gás e o conector da junta esférica com água deionizada.

4 Se a tocha foi desmontada para limpeza, lave bem o conjunto de tubos externos de quartzo e o vedante superior removível. Consulte a Figura 17A e a Figura 17B.



Figura 17A e B. Lave o conjunto de tubos externos de quartzo e o vedante superior com água deionizada.

- **5** Seque a tocha completamente e monte novamente a tocha semidesmontável, se necessário, antes de usar a tocha no instrumento.
- NOTA Para amostras com alto teor de sólidos totais dissolvidos (TDS), como digestões de solo, mergulhe as partes de quartzo da tocha em uma solução de detergente a 5% v/v por 10 minutos após a etapa inicial de limpeza ácida. Sempre enxágue a tocha e seque-a completamente antes de usá-la no instrumento. Siga as etapas descritas acima em "Lavagem da tocha" e abaixo em "Secagem da tocha". Esta etapa adicional pode ajudar a reduzir a deposição no tubo externo da tocha.

CUIDADO

Não coloque as peças plásticas da tocha em um forno para secagem. Não é tão eficaz na remoção de umidade quanto o uso de ar comprimido, argônio ou nitrogênio e pode danificar a tocha.

Secagem da tocha

- 1 Segure a tocha de modo invertido (com o conector da junta esférica para cima).
- 2 Injete ar comprimido limpo, argônio ou nitrogênio nas portas de suprimento de gás na base e na abertura do conector da junta esférica para remover a umidade. Consulte a Figura 18A e a Figura 18B.



Figura 18A e B. Use ar comprimido, argônio ou nitrogênio para secar a tocha.

3 Injete ar comprimido, argônio ou nitrogênio limpo na extremidade aberta do conjunto de tubos externos de quartzo. Não insira o bico de ar dentro do conjunto de tubos. Consulte a Figura 19.



Figura 19. Sopre ar comprimido, argônio ou nitrogênio limpo na extremidade aberta do conjunto de tubos.

- 4 Se a tocha foi desmontada para limpeza, sopre ar comprimido, argônio ou nitrogênio limpo no conjunto de tubos externos de quartzo e no vedante superior para remover toda a umidade.
- 5 Retire com sopro, cuidadosamente, toda umidade deixada no corpo da tocha e nas peças de quartzo.
- 6 Verifique se toda a umidade foi removida antes de recolocar a tocha no instrumento.
- 7 Para a tocha semi e completamente desmontável, é aconselhável desmontála, facilitando o acesso ao corpo da tocha para uma secagem completa.

Verificações adicionais após a limpeza

Execute as seguintes verificações após a limpeza:

- 1 Inspecione a tocha quanto a danos, como conexões soltas dos tubos de quartzo na base de plástico, orifícios ou rachaduras significativas. Se forem encontrados danos, substitua a tocha imediatamente.
- 2 Verifique a ocorrência de excedentes após a reinstalação da tocha no instrumento, para determinar se o procedimento de limpeza foi suficiente. Se encontrar excedentes, repita o processo de limpeza.
- **3** Substitua a tocha quando a superfície externa do tubo externo de quartzo estiver áspera para a tocha (que são sinais de desgaste) ou se houver rachaduras visíveis.
- NOTA A exposição prolongada ao ácido durante a limpeza pode levar à descoloração da base de plástico. Essa alteração é apenas cosmética, e o desempenho não será impactado se a tocha estiver limpa e se os resultados das outras verificações da tocha estiverem satisfatórios.

RECOMENDAÇÃO Quando a tocha não estiver em uso, armazene-a na caixa original ou em um saco plástico, ou então use o rack de armazenamento da tocha de encaixe simples Agilent ICP-OES (nº de peça G8010-67000).

Solução de problemas

Para obter informações de soluções de problemas, consulte o Help and Learning Center (Centro de treinamento e ajuda) do ICP Expert:

- 1 Clique duas vezes no Help and Learning Center (Centro de treinamento e ajuda) do ICP Expert na área de trabalho do computador.
- 2 Clicar em **Troubleshooting** (Solução de problemas) para ver instruções de solução da problemas.

Peças sobressalentes

Para obter informações sobre peças sobressalentes e itens consumíveis, consulte o site da Agilent Technologies:

www.agilent.com

Para substituir os itens listados abaixo, você deve usar as peças fabricadas pela Agilent, que podem ser adquiridas *on-line* no site da Agilent ou através do seu representante de vendas local.

Ao substituir o cabo de comunicações, sempre use um cabo Ethernetcom classificação "Categoria 5E Patch Cord (cabo de interconexão curto), STP/FTP, blindado, trançado, branco, 14 pés, booted (inicializável)" para manter a compatibilidade eletromagnética (EMC).

Abaixo, há uma lista de peças sobressalentes que se deve ter à mão para reduzir o tempo de inatividade durante a manutenção e reparos:

- Tocha (uma peça) ou componentes sobressalentes, como o conjunto de tubos externos e o injetor, se estiver usando uma tocha desmontável
- Filtro de entrada de ar (filtro de poeira)
- Janela de pré-óptica axial
- Janela de pré-óptica radial
- Câmara de nebulização
- Nebulizador
- Tubulação da bomba peristáltica
- Tubulação de drenagem

Suporte técnico

Para obter informações de contato do suporte técnico, consulte o site da Agilent Technologies:

www.agilent.com

Esta página foi deixada em branco intencionalmente.

Neste manual

O manual descreve o seguinte:

- Riscos e práticas de segurança
- Introdução
- Instalação
- Operação
- Manutenção e solução de problemas

www.agilent.com

© Agilent Technologies, Inc. © Agilent Technologies 2019

1ª edição, 10/2019



