

Learn more about maximizing ICP-OES performance:

agilent.com/chem/ICP-OESSupplies

See our OneNeb Universal Nebulizer in action at:

agilent.com/chem/OneNebVideo

了解更多关于如何最大限度提升 ICP-OES 性能的信息。请访问: agilent.com/chem/ICP-OESSupplies

如需查看运行中的 OneNeb 通用雾化器。请访问:

agilent.com/chem/OneNebVideo

Erfahren Sie mehr darüber, wie Sie die Leistung Ihres ICP-OES maximieren können: agilent.com/chem/ICP-OESSupplies

Sehen Sie unseren OneNeb Universalzerstäuber in Aktion:

agilent.com/chem/OneNebVideo

Obtenga más información sobre cómo maximizar el rendimiento del ICP-OES: agilent.com/chem/ICP-OESSupplies

Vea el nebulizador universal OneNeb en acción en:

agilent.com/chem/OneNebVideo

Pour en savoir davantage sur l'optimisation des performances des systèmes ICP-OES: agilent.com/chem/ICP-OESSupplies

Regardez une démonstration de notre nébuliseur universel

OneNeb sur: agilent.com/chem/OneNebVideo

Maggiori informazioni sull'ottimizzazione delle prestazioni dei sistemi ICP-OES su: agilent.com/chem/ICP-OESSupplies

Guarda il nebulizzatore universale OneNeb in azione su:

agilent.com/chem/OneNebVideo

ICP-OES の性能を最大化する消耗品のご紹介:

agilent.com/chem/jp

OneNeb Universal Nebulizer 紹介ビデオ:

agilent.com/chem/OneNebVideo

Obtenha mais informações sobre como maximizar o desempenho do sistema ICP-OES: agilent.com/chem/ICP-OESSupplies

Veja nosso nebulizador universal OneNeb em ação em:

agilent.com/chem/OneNebVideo

Подробнее о повышении производительности ИСП-ОЭС:

agilent.com/chem/ICP-OESSupplies

Универсальный распылитель OneNeb в действии:

agilent.com/chem/OneNebVideo



ESSENTIAL SUPPLIES FOR ICP-OES

ICP-OES 的
必备备件

UNVERZICHTBARE VERBRAUCHSMATERIALIEN
FÜR ICP-OES

CONSUMIBLES ESENCIALES
PARA ICP-OES

CONSUMMABLES INDISPENSABLES
POUR L'ICP-OES

PRODOTTI DI CONSUMO
PER ICP-OES

ICP-OES
関連消耗品

CONSUMÍVEIS ESSENCIAIS
PARA ICP-OES

ОСНОВНЫЕ РАСХОДНЫЕ МАТЕРИАЛЫ
ДЛЯ ОПТИКО-ЭМИССИОННОГО
СПЕКТРОМЕТРА С ИНДУКТИВНО-
СВЯЗАННОЙ ПЛАЗМОЙ, ИСП-ОЭС

The Measure of Confidence



Agilent ICP-OES Systems

Agilent ICP-OES 系统 • Agilent ICP-OES-Systeme • Sistemas ICP-OES Agilent • Systèmes ICP-OES Agilent
Sistemi ICP-OES Agilent • Agilent ICP-OES システム • Sistemas ICP-OES Agilent • Системы ИСП-ОЭС Agilent

1. Nebulizer

雾化器
Zerstäuber
Nebulizador
Nébuliseur
Nebulizzatore
ネブライザ
Nebulizador
Распылитель

2. Spray Chamber

雾化室
Sprühkammer
Cámara de nebulización
Chambre de nébulisation
Camera di nebulizzazione
スプレーチャンバ
Câmara de nebulização
Распылительная камера

3. Torch

炬管
Fackel
Antorcha
Torche
Torchia
トーチ
Tocha
Горелка

4. Pump Tubing

泵管
Pumpenschlauch
Tubos para bomba
Tuyaux de pompe
Tubi pompa
ポンプチューブ
Tubulação da bomba
Трубки к насосу



5. Miscellaneous Supplies

其他备件
Diverses Zubehör
Consumibles varios
Consommables divers
Prodotti di consumo vari
その他の消耗品
Consumíveis diversos
Различные комплектующие и
расходные материалы

6. Standards

标样
Standards
Estándares
Étalons
Standard
標準試料
Padrões
Стандарты

7. SPS 3 Autosampler Supplies

SPS 3 自动进样器备件
Zubehör für automatischen
Probengeber SPS 3
Consumibles para el muestreador
automático SPS 3
Consommables pour passeur
automatique d'échantillons SPS 3
Prodotti di consumo per
autocampionatore SPS 3
SPS 3 オートサンプラ消耗品
Consumíveis para o amostrador
automático SPS 3
Комплектующие к
автосамплеру SPS 3