

# NexION ICP-MS CONSUMABLES AND SUPPLIES



## Introduction and Overview

NexION® ICP-MS consumables and supplies are designed with your instrument in mind. Each part is fit to perform and manufactured with the highest quality materials available. Whether you are looking for sample introduction components or standards, we have the consumables you need to keep your NexION up and running smoothly and efficiently.



NexION® 5000  
ICP-MS



NexION®  
1100/2200 ICP-MS



NexION®  
1000/2000 ICP-MS

**TRUST THE CONSUMABLES  
ENGINEERED FOR  
YOUR INSTRUMENTS**

## Table of Contents

Introduction and Overview.....	2
Sampler and Skimmer Cones .....	3
Quartz Torches .....	4
Demountable Injectors .....	4
Concentric Plus and PFA Nebulizers .....	5
DuraMist Nebulizers .....	5
MicroMist Nebulizers.....	5
SeaSpray Nebulizers .....	5
Direct Connection Nebulizer Gas Fitting .....	6
<i>pergo</i> Argon Humidifiers.....	6
<i>Elegra</i> Argon Humidifiers.....	6
Spray Chambers .....	7
Spray Chamber Components.....	7
Nebulizer Fittings and Capillaries.....	8
Dual-Stage Detector.....	8
Autosampler Probes .....	9-10
Autosampler Tubes .....	11
Pump Tubing .....	12-13
Additional Tubing.....	13
Internal Standard Components.....	13
Temperature Controlled Spray Chambers .....	14
Sample Introduction Kits .....	14-16
SMARTintro Sample Intro Modules and Cassettes ..	17
Reference Standards .....	18-20
Environmental Analysis Standards .....	21
Pharmaceutical Analysis Standards .....	22-23
Gold Nanoparticle Standards.....	23
Other Replacement Parts .....	24
NexION 1100/2200/5000 ICP-MS Consumable Kits .....	24
NexION 1000/2000 ICP-MS Consumable Kits .....	25
Single Cell Sample Introduction Kit .....	26

## Sampler and Skimmer Cones

Precision-designed and manufactured, large-orifice sampler and skimmer cones provide superior long-term stability and resist clogging, allowing analysis under both high and low sample-uptake conditions. Nickel is a very rugged, long-lasting material for the majority of sample types, while platinum is the material of choice for more corrosive samples. They have been designed to maximize signal stability and to minimize clogging during extended runs of samples containing high dissolved solids. Platinum cones are boron-free and have ultra-low levels of impurities.

### Nickel Sampler and Skimmer Cones

Type	Material	ICP-MS Model	Part No.
Sampler	Nickel	All NexION Models	W1033612
Skimmer	Nickel	NexION 1000/1100/2000	W1026356
Skimmer	Nickel	NexION 2200	N8171142
Skimmer	Nickel	NexION 5000	N8161117

### Platinum Sampler and Skimmer Cones

Type	Material	ICP-MS Model	Part No.
Sampler	Platinum	All NexION Models	N8161140
Sampler for HF and Organic Solvents	Platinum	All NexION Models	W1033614
Sampler for Sulfur Based Acids/Solvents	Platinum	All NexION Models	N8145028
Skimmer	Platinum	NexION 1000/1100/2000	W1026907
Skimmer	Platinum	NexION 2200	N8171143
Skimmer	Platinum	NexION 5000	N8161041

### Hyper-Skimmer Cones

Type	Material	ICP-MS Model	Part No.
Hyper-Skimmer Cone	Aluminum	NexION 1000/1100/2000	W1033995
Hyper-Skimmer Cone	Nickel	NexION 2200/5000	N8160120
Hyper-Skimmer for Geological Applications	Nickel	NexION 2200/5000	N8170140
Hyper-Skimmer Kit	Aluminum	NexION 1000/1100/2000	N8140533

## Replacement Parts

Description	ICP-MS Model	Part No.
Sampler Gasket	All NexION Models	W1040148
Hyper-Skimmer O-Ring	NexION 1000/1100/2000	09902123
Hyper-Skimmer O-Ring	NexION 2200/5000	09211389
Hyper-Skimmer Screw	NexION 1000/1100/2000	WE027484
Cone Removal Tool	NexION 1000/1100/2000	W1034694
Sampler and Skimmer Removal Tool	NexION 2200/5000*	N8160118
Hyper-Skimmer Removal Tool	NexION 2200/5000	N8160119
Torch Alignment Tool	NexION 1000/1100/2000/2200/5000	N8157042

\*Hyper-Skimmer Removal Tool (N8160119) also needed for NexION 5000.



Sampler Cone for NexION 1000/1100/2000/2200/5000 and Skimmer Cone for NexION 1000/1100/2000



Hyper-Skimmer Cone for NexION 1000/1100/2000



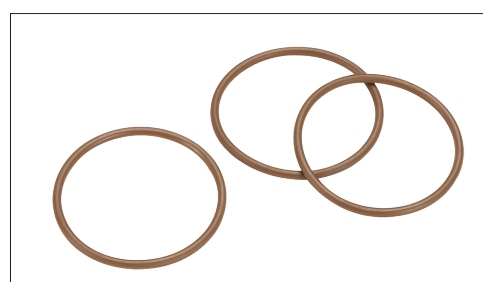
Hyper-Skimmer Cone for NexION 2200/5000



Cone Removal Tool for NexION 1000/1100/2000



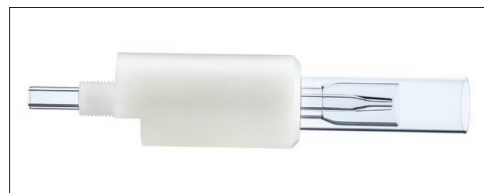
Hyper-Skimmer Removal Tool for NexION 2200/5000



Hyper-Skimmer O-Ring for NexION 1000/1100/2000

## Quartz Torches

Torch	Description	Part No.
Demountable Quartz Torch (Black)	Demountable Quartz Torch, PVDF O-Ring Free Holder for Demountable Injectors	N8152379
Demountable UHP Quartz Torch (Gray)	Demountable Ultra Pure Quartz Torch, PVDF O-Ring Free Holder for Demountable Injectors	N8152448
Demountable Quartz Torch (Organics)	Demountable Quartz Torch for Organic Analysis with PVDF O-Ring Free Holder	N8152596
Fixed 2.0 mm Injector UHP Quartz Torch (White)	One Piece Ultra Pure Quartz Torch with 2.0 mm Injector, PVDF O-Ring Free Holder for NexION 2000S ICP-MS	N8152428
Fixed 2.0 mm Injector Quartz Torch (Blue/Purple)	One Piece Quartz Torch with 2.0 mm Injector, PVDF O-Ring Free Holder for NexION 2000B ICP-MS	N8152472
Fixed 2.5 mm Injector Quartz Torch (Green)	One Piece Quartz Torch with 2.5 mm injector, PVDF O-Ring Free Holder for NexION 2000C ICP-MS	N8152473
Fixed 1.5 mm Injector Quartz Torch (Blue/Blue)	One Piece Quartz Torch with 1.5 mm Injector, PVDF O-Ring Free Holder	N8152427
Fixed 1.5 mm Injector Quartz Torch (Orange)	One Piece Quartz Torch with 1.5 mm Injector, PVDF O-Ring Free Holder for Organic Solvents	N8152616
Fixed 1.5 mm Injector UHP Quartz Torch (White/White)	One Piece Ultra Pure Quartz Torch with 1.0 mm Injector, PVDF O-Ring Free Holder	N8152601
Fixed 1.0 mm Injector Quartz Torch (Orange/Orange)	One Piece Quartz Torch with 1.0 mm Injector, PVDF O-Ring Free Holder for High Volatility Organic Solvents	N8152474
Fixed 1.0 mm Injector Quartz Torch (Blue/Blue/Blue)	One Piece Quartz Torch with 1.0 mm Injector, PVDF O-Ring Free Holder	N8152699
Fixed 1.0 mm Injector UHP Quartz Torch (White/White/White)	One Piece Ultra Pure Quartz Torch with 1.0 mm Injector, PVDF O-Ring Free Holder	N8152600
Torch Injector Nut		N8152451
Torch Injector Nut Removal Tool		N8152443
Torch O-Ring Kit	Viton O-Ring Kit Includes: 2.8 mm i.d. (Qty. 4), 14 mm ID (Qty. 1) and 16 mm i.d. (Qty. 1)	N8152476
PTFE Torch Light Shield Disk		N8152377



Fixed SiIQ 2 mm Torch (N8152428)



Fixed 1.5 mm Torch (N8152427)



One Piece Quartz Torch with 2.5 mm Injector (N8152473)



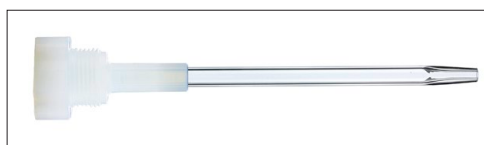
Torch Injector Nut (N8152451)



PTFE Torch Light Shield Disk (N8152377)

## Demountable Injectors

Description	Part No.
2.0 mm Demountable Quartz Injector (Black)	N8152376
1.5 mm Demountable Quartz Injector (Black/Orange)	N8152445
1.0 mm Demountable Quartz Injector (Black/Orange/Orange)	N8152444
1.0 mm Demountable Platinum Injector (Gray/Gray/Gray)	N8152603
2.0 mm Demountable Platinum Injector (Gray)	N8152429
1.5 mm Demountable Sapphire Injector (Black/Gold/Gold)	N8152606
2.0 mm Demountable Sapphire Injector (Black/Gold)	N8152447
Platinum Shield Disk for NexION Platinum and Sapphire Injectors	N8151106



2 mm Demountable Quartz Injector (N8152376)



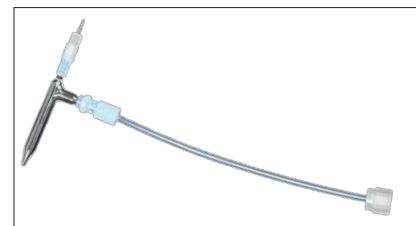
2 mm Demountable Platinum Injector (N8152429)

## Concentric Plus and PFA Nebulizers

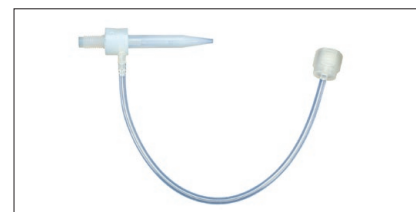
Concentric Plus nebulizers provide low dead volume and high sensitivity with excellent detection limits and washout. Each Plus nebulizer includes the FitKit+ gas and sample quick connect fittings. The connections are made with high-purity PFA, minimizing the chance of contamination during backflow events. The gas connection fits snugly over the maria on the sidearm of the nebulizer.

Material	Product Name	Description	Part No.
Glass	Concentric Type CT R +	High performance glass	N8152373
Quartz	Concentric Type CT Q +	High performance quartz	N8152372
UHP Quartz	Concentric Type CT SilQ +	High performance ultra-high purity quartz	N8152384
PFA	PFA ST3 MicroFlow	High performance, HF resistant	N8152378
PFA	PFA LC MicroFlow	For coupling HPLC-ICP-MS	N8152477

\*See page 8 for Nebulizer Fittings and Capillaries



Glass Low Volume Concentric Plus Nebulizer (N8152373)



PFA ST3 Nebulizer Assembly (N8152378)

## DuraMist Nebulizers

The DuraMist concentric nebulizer is an economical nebulizer for high precision analyses involving samples digested in HF.

Material	Uptake Rate	Part No.
PEEK	1.0 mL/min	N8151078

## MicroMist Nebulizers

The MicroMist micro-uptake glass concentric nebulizer offers high performance for limited sample volumes. Direct Connection (DC) nebulizers are fitted with a UniFit connector and an Argon connector. MicroMist nebulizers include a gas fitting with ratchet connector (N8151079).

Material	Uptake Rate	Part No.
Glass	0.2 mL/min	N8151074

## SeaSpray Nebulizers

The SeaSpray concentric nebulizer is ideal for samples with high dissolved solids and offers outstanding nebulization efficiency for trace level analysis. Direct Connection (DC) nebulizers are fitted with a UniFit connector and an Argon connector. SeaSpray nebulizers include a gas fitting with ratchet connector (N8151079).

Material	Uptake Rate	Part No.
Glass	0.4 mL/min	N8151070
Quartz	1.0 mL/min	N8151072
Glass	2.0 mL/min	N8151071



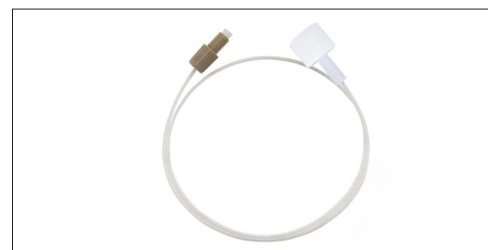
MicroMist Nebulizers



SeaSpray Nebulizers

## Direct Connection Nebulizer Gas Fitting

Description	Part No.
1/8 Gas Fitting with Ratchet Nebulizer Connector (DC)	N8151079



Gas Fitting with Ratchet Nebulizer Connector

## pergo Argon Humidifiers

The *pergo* improves performance of all concentric nebulizers for the routine analysis of brines, fusion digests and other high-TDS samples. *pergo 2* for ICP-MS humidifies both the nebulizer gas flow and the AMS gas flow for optimum performance.

Description	Part No.
<i>pergo 1</i> without Nebulizer	N8145315
<i>pergo 1</i> with Type C Plus Nebulizer (N8152373)	N8145316
<i>pergo 1</i> with PFA-ST Nebulizer	N8150427
<i>pergo 2</i> without Nebulizer	N8140743
<i>pergo 2</i> with Type C Plus Nebulizer (N8152373)	N8150428
<i>pergo 1</i> NexION 1000/2000 Gas Line Conversion Kit	N8152683
<i>pergo 2000</i> Nebulizer Gas and AMS Humidifier	N8150498
<i>pergo 2000</i> Nebulizer Gas Humidifier	N8150499

### FEATURES AND BENEFITS:

- Improved RSDs
- Improved short- and long-term stability
- Improved detection limits
- Faster washout
- Ability to use high-sensitivity nebulizers for high-TDS samples
- Reduce long-term drift
- Extends the length of analytical runs



*pergo 2000* Nebulizer Gas and AMS Humidifier

## Elegra Argon Humidifiers

The *Elegra* argon humidifier is used in ICP analyses involving samples with high concentrations of dissolved solids. It helps to alleviate salt deposits in the nebulizer and torch injector, allowing uninterrupted and maintenance-free operation.

- Highly efficient membrane humidification technology
- Improved signal stability for samples with high TDS
- *Elegra* Dual two-channel configuration available for ICP-MS instruments using auxiliary argon

Description	Part No.
<i>Elegra</i> Argon Humidifier	N8150493
<i>Elegra</i> Dual Two-Channel Argon Humidifier	N8150494



*Elegra* Argon Humidifier

## Spray Chambers

The cyclonic spray chamber relies on momentum to efficiently and quickly remove the larger droplets from the sample mist, decreasing any opportunity for occlusion or agglomeration of the smaller droplets. The result is higher efficiency, better precision and faster washouts than other spray chamber designs.

- Decreased sample introduction system installation and optimization time due to the preset nature of the spray chamber/nebulizer combination
- More samples per hour due to minimal dead volume and efficient draining

Type	Description	Part No.
Glass Cyclonic High Sensitivity Spray Chamber	For NexION 2000B and NexION 2200 ICP-MS	N8152375
Glass Cyclonic High Sensitivity Spray Chamber with Matrix Gas Port	For NexION 2000C/P and NexION 2200 ICP-MS	N8152389
SiIQ Cyclonic High Sensitivity Spray Chamber	For NexION 2000S and NexION 2200 ICP-MS	N8152424
SiIQ Cyclonic Spray Chamber with Matrix Gas Port	For NexION 5000 ICP-MS	N8152539
Quartz Cyclonic High Sensitivity Spray Chamber with Matrix Gas Port	For applications requiring low levels of Boron and Silicon detection	N8152383
PFA Barrel Spray Chamber	PFA Double Pass Spray Chamber for use with HF acid	N8152390
PFA Barrel Spray Chamber with Additional Gas Port	PFA Double Pass Spray Chamber for use with HF acid	N8150450

## Spray Chamber Components

Description	Part No.
UV Light Shield for Cyclonic Spray Chamber for NexION 2000B and NexION 2200 ICP-MS	N8152425
UV Light Shield for UV Shield for PC <sup>3x</sup> Peltier Cooler/Heater for NexION 2000C/P/S (for use with PC <sup>3</sup> ) and NexION 2200 ICP-MS	N8152475
Threaded Spray Chamber Drain Line 1/4 -28 Fitting	N8152456



Glass Cyclonic Chamber (N8152375)



Glass Cyclonic High Sensitivity Spray Chamber with Matrix Gas Port (N8152389)



PFA Barrel Spray Chamber with Additional Gas Port (N8150450)

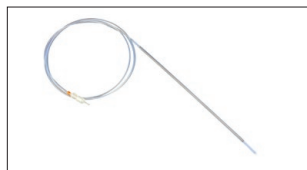


UV PC<sup>3</sup> Light Shield (N8152475)

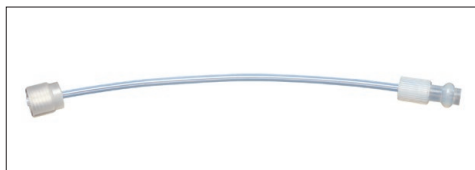


## Nebulizer Fittings and Capillaries

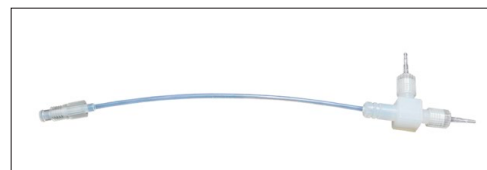
Description	Part No.
Internal Standard Tee for Concentric Nebulizers	N8152386
PFA-ST Nebulizer Internal Standard Addition Tee	N8152423
Nebulizer Fitting for Concentric Nebulizer (N8152378)	N8152387
Peripump Union - Male Barb CTFE Fitting	N8152453
Peripump Union - Female Barb CTFE Fitting	N8152388
Sample Introduction Line for PFA-ST Nebulizer	N8152455
Sample Introduction Line for Concentric Nebulizers	N8145503
Concentric Nebulizer 6" Gas Line	N8152385
Matrix Gas Line for Oxygen Addition	N8152426
AMS Gas Line	N8152374
Internal Standard Probe	N8152452
PTFE Torch Light Shield Disk	N8152377
Union Fittings for LC to PFA Nebulizer	N8152478
1/8 Gas Fitting with Ratchet Nebulizer Connector	N8151079



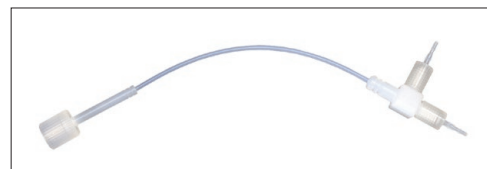
Internal Standard Probe (N8152452)



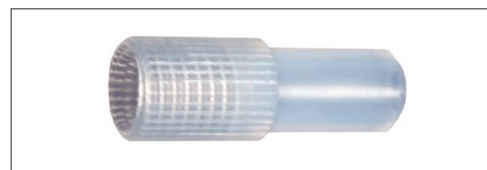
Nebulizer Gas Line (N8152385)



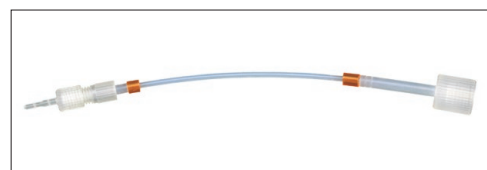
Internal Standard Tee for Concentric Nebulizer (N8152386)



Internal Standard Addition Tee (N8152423)



10-32 Union for Concentric Nebulizer (N8152373)

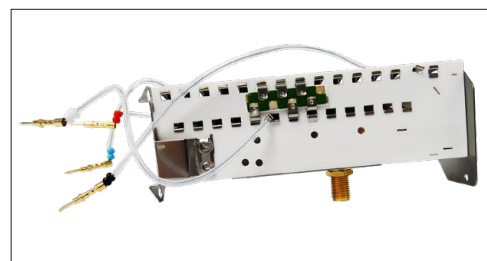


Sample Introduction Line for PFA-ST Nebulizer (N8152455)

## Dual-Stage Detector

The dual-stage detector is an integral part of the NexION ICP-MS SimulScan™ system. This discrete dynode electron multiplier measures both high- and low-level analytes simultaneously. This conserves valuable or limited samples, eliminates the need to perform time-consuming sample dilutions and allows you to quickly characterize unknown samples.

Description	Part No.
Dual-Stage Detector	N8151011



ICP-MS Detector (N8151011)



## Autosampler Probes

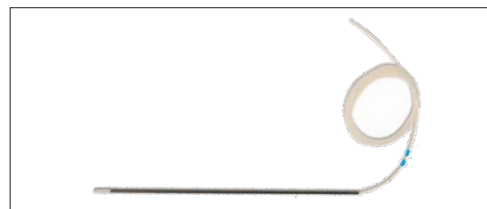
### Probes for S20 Series Autosamplers

Description	Part No.
1.0 mm i.d. Corrosion Resistant, FEP Covered Carbon-Fiber Probe with Attached Capillary Tubing (2 blue bands)	N0811956*
0.8 mm i.d. Corrosion Resistant, FEP Covered Carbon-Fiber Probe with Attached Capillary Tubing (red band)	N0811953*
0.5 mm i.d. Corrosion Resistant, FEP Covered Carbon-Fiber Probe with Attached Capillary Tubing (blue band)	N0777547*
Filtered Stainless Steel Autosampler Probe for Oils Applications	N0771529
Probe Fitting Kit - Includes one barb fitting (tan), one 1/16" flangeless fitting (blue) and two 1/16" CTFE ferrules (clear). To be purchased with autosampler probes N0811956, N0811953 and N0777547	N0830022

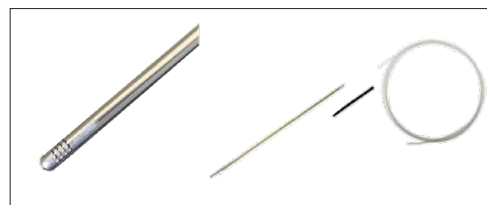
\*This probe must be purchased with the probe fitting kit N0830022.

### Probes for SC-DX, DX/DXCI, SampleSense FAST, prepFAST Autosamplers

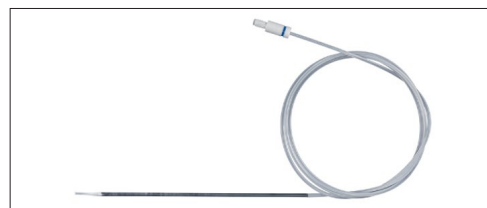
Description	Part No.
1.0 mm i.d., Corrosion-Resistant, Heavy-Walled, Self-Supporting Carbon-Fiber Probe with 130 cm Anti-Kinking Sample Transfer Line. Compatible with DX and DXCI Autosamplers	N8151113
0.25 mm i.d., Self-Aspirating, Carbon-Fiber Sampling Probe with 80 cm Capillary for ST Nebulizers. Compatible with SC-series FAST Autosamplers	N0777223
0.50 mm i.d., Self-Aspirating, Carbon-Fiber Sampling Probe with 80 cm Capillary for ST Nebulizers. Compatible with SC-series FAST Autosamplers	N0777225
0.25 mm i.d., Self-Aspirating, Carbon-Fiber Sampling Probe with 150 cm Capillary for ST Nebulizers. Compatible with SC-DX Autosamplers	N0777826
0.50 mm i.d., Self-Aspirating, Carbon-Fiber Sampling Probe with 150 cm Capillary for ST Nebulizers. Compatible with SC-DX Autosamplers	N0777827
Internal Standard Probe with Carbon Fiber - 0.5 mm i.d. (Orange), 150 cm Capillary. Includes Fluoropolymer Micro Fitting and can be used with Concentric and PFA Nebulizers	N8152452



Carbon Fiber Autosampler Probe for S20 Series Autosamplers (N0811956)



Filtered Stainless Steel Autosampler Probe for Oils Applications (N0771529)



Heavy-Walled, Anti-Kinking Probe for DXCI Autosamplers (N8151113)



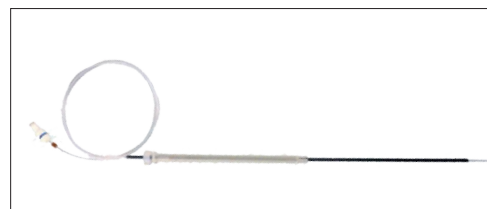
0.25 mm i.d., Self-Aspirating, Carbon-Fiber Sampling Probe for ST Nebulizers. (N0777223)

## Autosampler Probes *continued...*

### Probes for S10 Autosamplers

Description	Part No.
Sampling Probe, 0.6 mm i.d. - Includes PTFE encased stainless steel sampling probe and screw fitting (B3000152) and capillary tubing assembly (B3000157)	B3000159
FIAS Sampling Probe, 0.6 mm i.d. - Includes Stainless Steel Sampling Probe and Screw Fitting (B3000152) and FIAS Capillary Tubing Assembly (B3000158)	B3000160
FIAS Standard Sampling Probe, 1.0 mm i.d. - Includes PTFE encased stainless steel sampling probe and screw fitting (B3000152) and capillary tubing assembly (B0191060)	B3000161
Sampling Probe, 0.6 mm i.d. - Corrosion-resistant probe (black epoxy), suitable for inorganic acids and most organic solvents, except for NMP. Includes sampling probe with FEP tube and screw fitting (B3000055) and capillary tubing assembly (B3000157)	B3001770
FIAS Sampling Probe, 0.6 mm i.d. - Includes Sampling Probe with FEP Tube and Screw Fitting (B3000055) and FIAS Capillary Tubing Assembly (B3000158)	B3001771
FIAS Sampling Probe, 1.0 mm i.d. - Corrosion-resistant probe (black epoxy), suitable for inorganic acids and most organic solvents, except for NMP. Includes sampling probe with FEP tube and screw fitting (B3001769) and capillary tubing assembly (B0191060)	B3001772
Ultraclean Sampling Probe, 0.5 mm i.d. - For ultra-trace analysis. Operates at 400 µL/min. Requires CTFE union (N8122355) to connect to PFA-ST nebulizer	N8122361
Ultraclean Sampling Probe, 0.25 mm i.d.	N8122360
CTFE Union. Necessary to connect ultraclean autosampler probe (N8122360 & N8122361) with PFA-ST concentric nebulizer (N8122192)	N8122355
PTFE Sampling Probe Assembly with Carbon Fiber Support and Female Fitting, 0.25 mm i.d. for ICP-MS. Requires but does not include N8145017	N8145139
PTFE Sampling Probe Assembly with Carbon Fiber Support and Female Fitting, 0.5 mm i.d. Requires but does not include N8145017	N8145140
PTFE Sampling Probe Assembly with Carbon Fiber Support and Male Screw Fitting, 0.55 mm i.d. Requires but does not include N8145018*	N2024132*
PTFE sampling probe for FAST systems, carbon fiber support 0.8 mm i.d. (blue marker), 1/4-28 fitting, (1.2 m long). Requires but does not include B0172053	N0777434*
PTFE sampling probe for oneFAST system, carbon fiber support 1.0 mm i.d. (grey marker), 1/4-28 fitting, (100 cm long).	N2024131*

\* Note: Keep spare ferrules on hand, N8145237 Qty. 10



PTFE Sampling Probe Assembly with Carbon Fiber Support and Male Screw Fitting, 0.55 mm i.d. (N2024132)

### PerkinElmer HTS Sample Introduction Probes and Fittings

Description	Part No.
1.0 mm i.d. Corrosion Resistant, FEP Covered Carbon-Fiber Probe with Attached Capillary Tubing (2 blue bands)	N0811956
0.8 mm i.d. Corrosion Resistant, FEP Covered Carbon-Fiber Probe with Attached Capillary Tubing (red band)	N0811953
0.5 mm i.d. Corrosion Resistant, FEP Covered Carbon-Fiber Probe with Attached Capillary Tubing (blue band)	N0777547
Filtered Stainless Steel Autosampler Probe for Oils Applications	N0771529
Probe Fitting Kit - Includes one barb fitting (tan), one 1/16" flangeless fitting (blue) and two 1/16" CTFE ferrules (clear). To be purchased with autosampler probes N0811956, N0811953 and N0777547	N0830022

## Autosampler Tubes

### Round Bottom Tubes



Graduated 15 mL Autosampler Tube without Caps (N0777099)

Description	Type	Size (mm)	Pack	Qty	Part No.
8 mL Tubes without Caps	Standard	13 x 100	Bulk	1200	N0777156
8 mL Tubes without Caps	Standard	13 x 100	250/Bag	1000	N0777159
8 mL Tubes without Caps	Standard	13 x 100	Bulk	1000	B0508901
14 mL Tubes without Caps	Standard	17 x 100	Bulk	1000	N0777940
14 mL Tubes with Dual-Position Polyethylene Caps Included	Standard	17 x 100	Bulk	1000	N0777941
14 mL Tubes without Caps	Standard	17 x 100	250/Bag	1000	N0777168
15 mL Tubes without Plugs	Standard	17 x 100	Bulk	1200	N0777167
Plugs for 15 mL N0777167 Tubes, (Blue) Plugs Only	Standard		Bulk	1200	N0777599
15 mL Graduated Tubes without Plugs	Standard	17 x 100	Bulk	2000	N0777099
16 mL Tubes without Caps	Standard	17 x 100	250/Bag	1000	N9301205

### Conical Bottom Tubes with Graduation

Description	Type	Size (mm)	Pack	Qty	Part No.
15 mL Tubes with Red Flat Caps	Standard	17 x 120	50/bag	500	B0193233
15 mL Sterile Tubes with Teal Flat Caps Attached	Sterile	17 x 118	50/Bag	500	N0774095
15 mL Metal-Free Sterile Tubes with White Flat Caps	Sterile, Metal-Free	17 x 118	50/Bag	500	N0776117
15 mL Metal-Free Sterile Tubes Racked with White Flat Caps	Sterile, Metal-Free	17 x 118	25/Rack	500	N0776118
15 mL Tubes with Teal Flat Caps Attached	SuperClear	17 x 118	50/Bag	500	N0777701
15 mL Tubes with Teal Flat Caps Attached	SuperClear	17 x 118	Racked 50/Bag	500	N0777702
15 mL Tubes with Teal Flat Caps Attached	SuperClear	17 x 118	Bulk	500	N0777703
15 mL Tubes with Teal Plug Caps	SuperClear	17 x 118	Bulk	500	N0777706
50 mL Tubes with Red Flat Caps Attached, Freestanding	Standard	28 x 115	25/Bag	500	B0193234
50 mL Tubes with Teal Flat Caps Attached, Freestanding	PerformR	30 x 116	Bulk	500	N0777698
50 mL Tubes with Teal Plug Caps Included, Freestanding	PerformR	30 x 116	Bulk	500	N0777700
15 mL Metal-Free Sterile Tubes with White Flat Caps	Sterile, Metal-Free	17 x 118	50/Bag	500	N0776117
50 mL Metal-Free Sterile Tubes with White Flat Caps	Sterile, Metal-Free	29 x 115	50/Bag	500	N0776115
50 mL Metal-Free Sterile Tubes Racked with White Flat Caps	Sterile, Metal-Free	30 x 116	25/Rack	500	N0776116
15 mL Metal-Free Sterile Tubes Racked with White Flat Caps	Sterile, Metal-Free	17 x 118	25/Rack	500	N0776118
50 mL Tubes with Teal Flat Caps Included	SuperClear	29 x 115	Bulk	500	N0777693

### PFA and Non-PFA Bottles

Description	Material	Qty	Part No.
250 mL, 61 mm o.d. Polyethylene Wide Mouth Bottle	Polyethylene	12	N0777818
125 mL, 50 mm o.d. Polyethylene Bottle with Screw Cap	Polyethylene	12	N0777820
20 mL, 28 mm o.d. PFA Bottle and Cap	PFA	5	N0777406

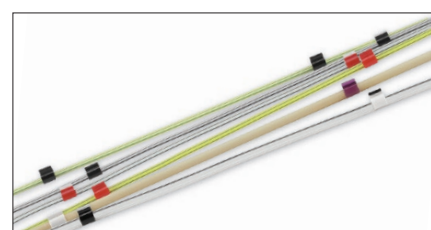
## Pump Tubing

### Peristaltic Pump Tubing for Standard Pump

Peripump Tubing Kit (Starter Pack of Tubing, includes 4 of each commonly used sizes)				Part No. N8150060
I.D. (mm)	Stop Colors	Quantity	Type	
0.19	Orange/Red	4	PVC Pure	
0.44	Green/Yellow	4	PVC Pure	
0.51	Yellow/Orange	4	PVC Pure	
1.30	Gray/Gray	4	Santoprene	

### Peristaltic Pump Tubing for PerkinElmer HTS System

I.D. (mm)	Stop Colors	PVC Pure (Pkg. 12)
0.19	Orange/Red	N8152401
0.44	Green/Yellow	N8152404
0.51	Yellow/Orange	N8152405
1.30	Santoprene Gray/Gray	N8152415



Pump Tubing

### Standard Pump Tubing

I.D. (mm)	Stop Colors	PVC Pure (Pkg. 12)	Solvent Flex (Pkg. 12)	Santoprene (Pkg. 12)
0.19	Orange Red	N8152401	N8152408	
0.25	Orange Blue	N8152402		
0.38	Orange Green	N8152403	N8152409	
0.44	Green Yellow	N8152404	N8152410	
0.51	Orange Yellow	N8152405		
0.64	Orange White	N8152406	N8152411	
0.76	Black Black	N8152407		
1.14	Red Red		N8152413	
1.30	Gray Gray			N8152415

## Consumables and Supplies Reference Guide

### MP<sup>2</sup> Two-Stop Peristaltic Pump Tubing

i.d. (mm)	Stop Colors	Slope (L/min per RPM)	Non-Flared			Flared	
			PVC 2-Stop	Santoprene 2-Stop	Solva 2-Stop	PVC 2-Stop	Solva 2-Stop
0.19	Orange Red	1.3				N8145195	
0.27	Orange Blue	2.7				N8145196	
0.38	Orange Green	4.7	N8145148	N8145171		N8145197	N8145211
0.44	Green Yellow	7.6	N8145149			N8145198	
0.51	Orange Yellow	9.5	N8145150			N8145199	
0.64	Orange White	14				N8145201	
0.76	Black Black	19	N8145153			N8145202	N8145216
0.89	Orange Orange	24	N8145154				
1.02	White White	31	N8145156				
1.09	White Red	33				N8145206	
1.14	Red Red	35	N8145158		N8145181	N8145207	
1.3	Grey Grey	47	N8145160	N8145173	N8145183		
1.42	Yellow Yellow	50	N8145161				
1.52	Yellow Blue	51	N8145162	N8145174			
3.17	Black White	70	N8145170	N8145175			

## Additional Tubing

Description	Part No.
Sample Uptake Tubing (per cm)	WE024375
Drain Tubing (per foot)	02506495
Torch Box Drain Assembly	N8140506

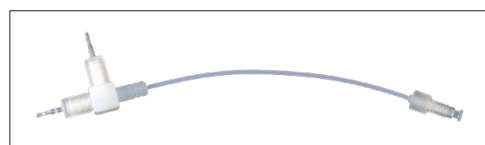
## Internal Standard Components

### Internal Standard Spares for Concentric Nebulizers

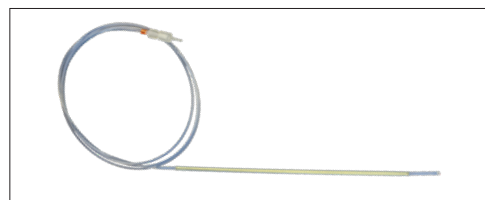
Description	Part No.
PFA Tee	N8152386
Nebulizer Fitting	N8152387
Internal Standard Probe	N8152452

### Internal Standard Spares for PFA Nebulizer

Description	Part No.
PFA Tee	N8152423
Internal Standard Probe	N8152452



Internal Standard Addition Tee (N8152386)



Internal Standard Probe (N8152452)

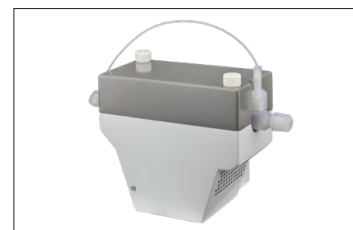
## Temperature Controlled Spray Chambers

The PC<sup>3x</sup> is a compact Peltier heated or cooled inlet system that controls the spray chamber temperature and improves long-term stability. In cooling mode, water or solvent is removed from the aerosol, reducing water or solvent loading on the plasma. In heating mode, sensitivity is increased and volatility-related sample transport effects can be reduced or eliminated.

PCH is a Peltier cooler/heater for the high-performance PFA spray chamber. It is a temperature-controlled inlet system that offers the advantages of low memory effects, rapid washout, and high sample transport efficiency for HF-containing samples. In heating mode, sensitivity can be improved for low-flow applications. Temperature control is available from -10 °C to +80 °C.



PC<sup>3x</sup> Sample Introduction (N8152382)






PCH Peltier Cooler/Heater Thermally Stabilized Inlet System (N8150445)

### PC<sup>3</sup> Peltier Coolers

Description	Part No.
PC <sup>3x</sup> Sample Introduction Kit (Spray Chamber not included)	N8152382
PCH Peltier Cooler/Heater Thermally Stabilized Inlet System with PFA Spray Chamber and PFA-ST nebulizer	N8150446
PCH Peltier Cooler/Heater Thermally Stabilized Inlet System with PFA Spray Chamber (Nebulizer not included)	N8150445
Cyclonic Peltier Cooler for NexION 1100	N8171067

## Sample Introduction Kits

NexION SMARTintro sample introduction solutions are tested as a complete unit to ensure intra- and inter-laboratory performance. Each system utilizes specific combinations of sample introduction components and the kits are color-coded according to the application of the user.






SMARTintro Sample Introduction Kit	Description	Part No.
	<b>PFA-Platinum HF-Resistant Sample Introduction Kit (Platinum) – Kit Includes:</b> PFA Spray Chamber with PFA End Cap (N8150450), Cyclonic Spray Chamber Adapter (N8152551), Demountable SiIQ Torch (N8152448), Platinum 2 mm Injector (N8152429), Plasma Light Shield (N8152377), Cassette Only Platinum (N8152434), Gas Line Addition (N8152449), Spray Chamber Drain Line 1/4-28 Fitting (N8152456), PFA-200 PFA MicroFlow Nebulizer (N8152420), Torch Injector Nut Removal Tool (N8152443), Torch Injector Nut (N8152451)	<b>N8150033</b>
	<b>Organics Sample Introduction Kit (Orange) – Kit Includes:</b> Quartz Nebulizer Type C (N8152372), Cyclonic Quartz Spray Chamber (N8152383), Quartz 1.5 mm Organics Torch One-Piece (N8152616), Nebulizer Sample Line (N8145503), Plasma Light Shield (N8152377), Internal Standard Nebulizer Tee (N8152386), Spray Chamber Light Shield (N8152425), Gas Line Addition (N8152426), Cassette Only Orange (N8152435), Torch Injector Nut Removal Tool (N8152443), Torch Injector Nut (N8152451), Internal Standard Probe Qty. 2 (N8152452), PeriPump Union Male Fitting (N8152453), Spray Chamber Drain Line 1/4-28 Fitting (N8152456)	<b>N8150031</b>
	<b>Geological Samples Introduction Kit (Gold) – Kit Includes:</b> PeriPump Tubing Starter Kit (N8150060), PFA Spray Chamber with PFA End Cap (N8150450), Plasma Light Shield (N8152377), PFA Nebulizer (N8152378), Quartz Torch (N8152379), Internal Standard Addition Tee for PFA-ST Nebulizer (N8152423), Torch Injector Nut Removal Tool (N8152443), Sapphire 2 mm Injector (N8152447), Gas Line Addition (N8152449), Torch Injector Nut (N8152451), Internal Standard Probe Qty. 2 (N8152452), Sample Introduction Line for PFA-ST Nebulizer (N8152455), Spray Chamber Drain Line 1/4-28 Fitting (N8152456), PC3 Spray Chamber Light Shield (N8152475), Cassette Only Gold (N8152532)	<b>N8150040</b>

## Consumables and Supplies Reference Guide

SMARTintro Sample Introduction Kit	Description	Part No.
	<p><b>Sample Introduction Kit for NexION 1000 ICP-MS (Purple) – Kit Includes:</b></p> <p>Nebulizer Sample Line (N8145503), PeriPump Tubing Starter Kit (N8150060), Glass Nebulizer Type C (N8152373), Matrix Gas Line 9 in. (N8152374), Plasma Light Shield (N8152377), Glass Cyclonic C3 High Sensitivity Spray Chamber with Matrix Port (N8152389), Spray Chamber Light Shield (N8152425), Torch Injector Nut Removal Tool (N8152443), Torch Injector Nut (N8152451), Internal Standard Probe Qty. 2 (N8152452), PeriPump Union Male Fitting (N8152453), Spray Chamber Drain Line 1/4-28 Fitting (N8152456), One Piece 2.0 mm Quartz Torch (N8152472), Cassette Only Purple (N8152533)</p>	<b>N8150038</b>
	<p><b>Sample Introduction Kit for NexION 1100/2200 (Dark Green) – Kit Includes:</b></p> <p>Glass Nebulizer Type C (N8152373), Glass Cyclonic Spray Chamber with Matrix Gas Port (N8152389), One Piece 2.0 mm Quartz Torch (N8152472), Nebulizer Sample Line (N8145503), PeriPump Tubing Starter Kit (N8150060), Plasma Light Shield (N8152377), Matrix Gas Line (N8152374), Spray Chamber Light Shield (N8152425), Cassette Only Dark Green (N8171005), Torch Injector Nut Removal Tool (N8152443), Internal Standard Probe Qty. 2 (N8152452), PeriPump Union Male Fitting (N8152453), Spray Chamber Drain Line 1/4-28 Fitting (N8152456), Torch Injector Nut (N8152451)</p>	<b>N8170020</b>
	<p><b>Sample Introduction Kit for NexION 2000B ICP-MS (Blue) – Kit Includes:</b></p> <p>Glass Nebulizer Type C (N8152373), Glass Cyclonic Spray Chamber (N8152389), One Piece 2.0 mm Quartz Torch (N8152472), Nebulizer Sample Line (N8145503), Matrix Gas Line 9 in. (N8152374), Plasma Light Shield (N8152377), Internal Standard Nebulizer Tee (N8152386), Spray Chamber Light Shield (N8152425), Cassette Only Blue (N8152430), Torch Injector Nut Removal Tool (N8152443), Internal Standard Probe Qty. 2 (N8152452), PeriPump Union Male Fitting (N8152453), Spray Chamber Drain Line 1/4-28 Fitting (N8152456), Torch Injector Nut (N8152451)</p>	<b>N8150034</b>
	<p><b>Sample Introduction Kit for NexION 2000C ICP-MS (Green) – Kit Includes:</b></p> <p>Glass Nebulizer Type C (N8152373), Glass Cyclonic C3 High Sensitivity Spray Chamber with Matrix Port (N8152389), Quartz Torch (N8152473), Nebulizer Sample Line (N8145503), PeriPump Tubing Starter Kit (N8150060), Matrix Gas Line 9 in. (N8152374), Plasma Light Shield (N8152377), PC3 Peltier Cooler (N8152382), Internal Standard Nebulizer Tee (N8152386), Spray Chamber Light Shield (N8152425), Cassette Only Green (N8152431), Torch Injector Nut Removal Tool (N8152443), Internal Standard Probe Qty. 2 (N8152452), PeriPump Union Male Fitting (N8152453), Spray Chamber Drain Line 1/4-28 Fitting (N8152456), PC3 Spray Chamber Light Shield (N8152475), Torch Injector Nut (N8152451)</p>	<b>N8150035</b>
	<p><b>Sample Introduction Kit for NexION 2000P ICP-MS (Black) – Kit Includes:</b></p> <p>PFA Nebulizer (N8152378), Quartz Torch (N8152379), Glass Cyclonic C3 High Sensitivity Spray Chamber with Matrix Port (N8152389), Cyclonic Peltier Cooler (N8152382), CTFE Liquid Fitting Female (N8145018), FAST 1 Valve (N8152359), MP2 Micro Peristaltic Pump with Single FAST Valve (N8152364), Matrix Gas Line 9 in. (N8152374), 2 mm Demountable Quartz Injector (N8152376), Plasma Light Shield (N8152377), FAST Pump Tubing Starter Kit (N8152380), Spray Chamber Light Shield (N8152425), Cassette Only Black (N8152433), Torch Injector Nut Removal Tool (N8152443), PFA-ST Nebulizer Sample Introduction Line (N8152455), Spray Chamber Drain Line 1/4-28 Fitting (N8152456), Spray Chamber Light Shield (N8152475), Nuts/Ferrules 1/4-28 for HF Valve Qty.10 (N0777283), Sample Loop for Ports 1 and 4, 0.5 mL (N0777289), HF Ferrule/Nut Set, 1/4-28 Qty.10 (N0777297), 0.8 mm i.d. Sample Loop PTFE, 300 µL (N0777430), 2 mL Sample Loop (N0777479), 4 mL Sample Loop (N0777480), 1 mL Loop 1 mm i.d. (N0777664), 1.5 mL Loop 1 mm i.d. (N0777665), 0.5 mL Loop 1 mm i.d. (N0777666), L 3 mL Loop 1 mm i.d. (N0777667), Holder with FAST Bottles (N0777777), RS485 Serial Cable – Power/Communication (N2024191), CTFE Female Barb Fitting (N8145018), Waste Line for SC-FAST (N8145536), FAST 1 Valve with P7 (N8152359), Torch Injector Nut (N8152451)</p>	<b>N8150036</b>



## Consumables and Supplies Reference Guide

SMARTintro Sample Introduction Kit	Description	Part No.
	<p><b>Sample Introduction Kit for NexION 2000S ICP-MS (White) – Kit Includes:</b></p> <p>PFA Nebulizer (N8152378), Quartz SiIQ Spray Chamber (N8152424), Quartz Fixed Torch (N8152428), Cyclonic Peltier Cooler (N8152382), CTFE Liquid Fitting Female (N8145018), PeriPump Tubing Starter Kit (N8150060), Plasma Light Shield (N8152377), Internal Standard Addition Tee for PFA ST Nebulizer (N8152423), Cassette Only White (N8152432), Torch Injector Nut Removal Tool (N8152443), Internal Standard Probe Qty. 2 (N8152452), PFA-ST Nebulizer Sample Introduction Line (N8152455), Spray Chamber Drain Line 1/4-28 Fitting (N8152456), Spray Chamber Light Shield (N8152475)</p>	<p><b>N8150037</b></p>
	<p><b>Sample Introduction Kit for NexION 5000B ICP-MS (Green/White Stripe) - Kit Includes:</b></p> <p>Glass Low Volume Type C Nebulizer (N8152373), Sample Line for Low Volume Type C Nebulizer (N8145503), Kit-Peristaltic Pump Starter Tubing (N8150060), Gas Line-Matrix (N8152374), Shield-Plasma Light (N8152377), High Sensitivity Glass Cyclonic Spray Chamber with Matrix Gas Port C3 (N8152389), Light Shield-Spray Chamber (N8152425), Tool-Torch Injector Nut Removal (N8152443), Nut-Torch Injector (N8152451), 0.5 mm i.d. Internal Standard Probe (N8152452), Peripump Union with Male Barb CTFE Fitting (N8152453), Drain Line-1/4-28 fitting 0.8 mm i.d. PTFE (N8152456), One-Piece Quartz Torch with 2.0 mm i.d. Injector (N8152472), Cassette-Green with White Stripe without Torch (N8161044)</p>	<p><b>N8160023</b></p>
	<p><b>Sample Introduction Kit for NexION 5000S ICP-MS (White/Green Stripe) – Kit Includes:</b></p> <p>CTFE Barbed Fittings, Small Female Qty.1 (N0777829), PeriPump Tubing Starter Kit (N8150060), Plasma Light Shield (N8152377), PFA Nebulizer (N8152378), Spray Chamber Light Shield (N8152425), Quartz Fixed Torch (N8152428), Torch Injector Nut Removal Tool (N8152443), Torch Injector Nut (N8152451), PFA-ST Nebulizer Sample Introduction Line (N8152455), Spray Chamber Drain Line 1/4-28 Fitting (N8152456), 0.25 mm i.d. Self-aspirating Probe 80 cm (N8152512), 0.30 mm i.d. Self-aspirating Probe 80 cm (N8152513), SiIQ Spray Chamber with Matrix Gas Port (N8152539), White Torch Cassette with Green Stripe (N8161009)</p>	<p><b>N8160020</b></p>
	<p><b>Sample Introduction Kit for NexION 5000G ICP-MS (Light Blue/Green Stripe) - Kit Includes:</b></p> <p>PFA Nebulizer w/Integrated Gas Line (N8152378), Union FEM Pump Fluoropolymer Peristaltic (N0777829), Fitting-Male Barb CTFE (N8145017), Peristaltic Pump Tubing Starter Kit (N8150060), Shield-Plasma Light (N8152377), Light Shield-Spray Chamber (N8152425), Torch-Fixed 2.0 mm Ultra-Pure Quartz (N8152428), Tool-Torch Injector Nut Removal (N8152443), Nut-Torch Injector (N8152451), Line-Sample Introduction PFA-ST Neb (N8152455), Drain Line-1/4-28 Fitting 0.8mm i.d. PTFE (N8152456), Self Aspirating Probe-with ST Neb Fit 0.25 mm i.d. (N8152512), Self-Aspirating Probe with ST Neb Fit 0.30 mm i.d. (N8152513), Spray Chamber SiIQ with Gas Port (N8152539), Cassette-Light Blue with Green Stripe without Torch (N8991016)</p>	<p><b>N8990040</b></p>
	<p><b>Sample Introduction Kit for NexION 1000G, 2000G, and 2200G ICP-MS (Light Blue) – Kit Includes:</b></p> <p>Nebulizer Sample Line (N8145503), Peripump Starter Tubing Kit (N8150060), Glass Low Volume Type C Nebulizer (N8152373), Matrix Gas Line (N8152374), Plasma Light Shield (N8152377), Glass Cyclonic Spray Chamber with AMS (N8152389), UV Light Shield (N8152425), Torch Injector Nut Removal Tool (N8152443), Torch Injector Nut (N8152451), Internal Standards Probe (N8152452), Peripump Union Male (N8152453), Threaded Spray Chamber Drain Line (N8152456), Fixed 2.0 mm Injector Quartz Torch Blue (N8152472), Light Blue Cassette (N8991010)</p>	<p><b>N8990039</b></p>

## SMARTintro Sample Introduction Modules and Cassettes

SMARTintro sample introduction modules and cassettes include cassettes, torch, PTFE torch light shield and torch injector removal tool.

### Sample Introduction Modules

Color		Part No.
	Blue Colored Cassette, One-Piece Quartz Torch, 2.0 mm Quartz Injector	N8152437
	Black Colored Cassette, Demountable Quartz Torch, 2.0 mm Quartz Injector	N8152440
	Purple Colored Cassette, One-piece Quartz Torch, 2.0 mm Quartz Injector	N8152536
	Gold Colored Cassette, One-Piece Quartz Torch, 2.0 mm Quartz Injector	N8152535
	White/Green Stripe Colored Cassette, One-Piece SilQ Torch, 2.0 mm Quartz Injector	N8161013
	Dark Green Colored Cassette, One-Piece Quartz Torch, 2.0 mm Quartz Injector	N8171004
	Green/White Stripe Colored Cassette, One-Piece Quartz Torch, 2.0 mm Quartz Injector	N8161016

### Sample Introduction Cassettes Only

Color		Part No.
	Blue (Included with NexION 2000B ICP-MS)	N8152430
	Green (Included with NexION 2000C ICP-MS)	N8152431
	Black (Included with NexION 2000P ICP-MS)	N8152433
	Platinum (High Purity/HF Resistant)	N8152434
	White/Green Stripe (Organic Solvents)	N8161009
	Dark Green (Included with NexION 1100/2200 ICP-MS)	N8171005

*All sample introduction modules and sample introduction cassettes are compatible with NexION 1000/2000/2200/5000 instruments*

## Reference Standards

### Multi-Element Instrument Calibration Standards

Matrix	Contents	Vol.	Part No.
<b>NexION 1000/2000 Solution Kit</b>			<b>N8140504</b>
Solution Kit	NexION Wash Solution	250 mL	N8145050
	NexION Setup Solution	500 mL	N8145051
	NexION Standard/Reaction Mode Detection Limit Blank Solution	100 mL	N8145055
	NexION Standard/Reaction Mode Detection Limit Solution	100 mL	N8145056
	NexION Dual Detector Solution	100 mL	N8145059
	NexION Collision Setup Solution	250 mL	N8145052
	NexION ICP-MS Cell Stability Solution	500 mL	N8145054
	NexION Collision Mode Detection Limit Blank Solution	100 mL	N8145057
	NexION Collision Mode Detection Limit Solution	100 mL	N8145058
<b>NexION 1100/2200 Solution Kit</b>			<b>N8170112</b>
Solution Kit	NexION 2200 STD/DRC Mode Detection Limit Standard Solution	100 mL	N8171046
	NexION Rinse Solution	250 mL	N8145050
	NexION Setup Solution	500 mL	N8145051
	NexION KED Setup Solution	250 mL	N8145052
	NexION Cell Stability Solution	500 mL	N8145054
	NexION Detection Limit STD/DRC Blank Solution	100 mL	N8145055
	NexION Detection Limit KED Blank Solution	100 mL	N8145057
	NexION Detection Limit KED Standard Solution	100 mL	N8145058
	NexION Dual Detector Solution	125 mL	N8145059
<b>NexION 1100/2200 STD/DRC Mode Detection Limit Standard Solution</b>			
0.5% HNO <sub>3</sub>	1 µg/mL: Be, Bi, Ca, Fe, In	100 mL	N8171046
<b>NexION 5000 Solution Kit</b>			<b>N8160029</b>
Solution Kit	NexION Wash Solution	250 mL	N8145050
	NexION 5000 Setup Solution	500 mL	N8162301
	NexION 5000 Stability Solution	500 mL	N8162302
	NexION 5000 DRC NH <sub>3</sub> and Standard Detection Limit	125 mL	N8162303
	NexION 5000 DRC O <sub>2</sub> Detection Limit Solution	125 mL	N8162304
	Water Blank ASTM® Type I Water, 18 megohm	250 mL x 2	N9303813
<b>NexION 5000 Setup Solution</b>			
1% HNO <sub>3</sub>	200 ng/L Be, Ce, Fe, In, Li, Mg, Pb, U	500 mL	N8162301
<b>NexION 5000 Stability Solution</b>			
1% HNO <sub>3</sub>	200 ng/L Cd, Cr, Fe, Li, Mg, Pb 1 µg/L Cu	500 mL	N8162302
<b>NexION 5000 DRC NH<sub>3</sub> and Standard Detection Limit</b>			
1% HNO <sub>3</sub>	200 ng/L Al, Be, Ca, Cr, Cu, Fe, In, K, Mn, Ni, U, Zn	125 mL	N8162303
<b>NexION 5000 DRC O<sub>2</sub> Detection Limit Solution</b>			
H <sub>2</sub> O	200 ng/L Al, Be, Ca, Cr, Cu, Fe, In, K, Mn, Ni, U, Zn	125 mL	N8162304
<b>NexION Setup Solution</b>			
1% HNO <sub>3</sub>	1 µg/L: Be, Ce, Fe, In, Li, Mg, Pb, U	500 mL	N8145051

Matrix	Contents	Vol.	Part No.
<b>NexION Collision Setup Solution</b>			
1% HCl	10 µg/L: Co	250 mL	N8145052
	1 µg/L: Ce	100 mL	N8145056
<b>NexION Non-Cell Stability Solution</b>			
1% HNO <sub>3</sub>	1 µg/L: Cd, Cu, Mg, Pb	500 mL	N8145053
<b>NexION Cell Stability Solution</b>			
1% HNO <sub>3</sub>	1 µg/L: Cd, Cr, Fe, In, Mg, Pb	500 mL	N8145054
	10 µg/L: Co, Cu, Se (without Se N8151033)		
<b>NexION Standard/Reaction Mode Detection Limit Blank Solution</b>			
0.5% HNO <sub>3</sub>		100 mL	N8145055
<b>NexION Standard/Reaction Mode Detection Limit Solution</b>			
0.5% HNO <sub>3</sub>	1 µg/L: Be, Ca, Co, Fe, In, U	100 mL	N8145056
<b>NexION Collision Mode Detection Limit Blank Solution</b>			
1% HCl		100 mL	N8145057
<b>NexION Collision Mode Detection Limit Solution</b>			
1% HCl	10 µg/L: V, As, Se	100 mL	N8145058
<b>NexION 1000/1100/2000/2200 Dual Detector Solution</b>			
2% HNO <sub>3</sub>	200 µg/L: Al, Ba, Ce, Co, Cu, In, Li, Mg, Mn, Ni, Pb, Tb, U, Zn	100 mL	N8145059
<b>NexION 5000 Dual Detector Solution</b>			
2% HNO <sub>3</sub>	20 µg/L of Al, Ba, Ce, Co, In, Li, Mg, Mn, Ni, Pb, Tb, U, Zn	100 mL	N8162305
<b>NexION AFT Single-Element Solution</b>			
2% HNO <sub>3</sub>	2 µg/L: Fe	100 mL	N8145060
<b>NexION AFT Multi-Element Solution</b>			
2% HNO <sub>3</sub>	2 µg/L: Ag, Al, As, Ba, Be, Bi, Ca, Cd, Co, Cr, Cs, Cu, Fe, Ga, In, K, Li, Mg, Mn, Na, Ni, Pb, Rb, Se, Sr, Ti, U, V, Zn	100 mL	N8145061
<b>NexION AFT Single-Element Solution (H/He)</b>			
2% HNO <sub>3</sub>	10 µg/L Se	125 mL	N8151037
<b>NexION AFT Multi-Element Solution (H/He)</b>			
2% HNO <sub>3</sub>	10 µg/L Ag, Al, AS, Ba, Be, Bi, Ca, Cd, Co, Cr, Cs, Cu, Fe, Ga, In, K, Li, Mg, Mn, Na, Ni, Pb, Rb, Se, Sr, Ti, U, V, Zn	125 mL	N8151038
<b>NexION Wash Solution</b>			
1% HNO <sub>3</sub>		250 mL	N8145050
<b>Internal Standard Mix</b>			
5-10% HNO <sub>3</sub>	10 µg/mL: 6Li, Sc, Ge, Y, In, Tb, Bi	125 mL	N9303832
<b>Multi-Element Internal Standard</b>			
2% HNO <sub>3</sub>	10 µg/mL: Bi, Ho, In, 6Li, Sc, Tb, Y	125 mL	N9303834
<b>Quality Control Standard – 21 Elements</b>			
5% HNO <sub>3</sub> /tr Tartaric Acid/tr HF	100 µg/mL: As, Be, Ca, Cd, Co, Cr, Cu, Fe, Li, Mg, Mn, Mo, Ni, Pb, Sb, Se, Sr, Ti, Tl, V, Zn	125 mL	N9300281
<b>Instrument Calibration Standard 2</b>			
5% HNO <sub>3</sub> /tr Tartaric Acid/tr HF	100 mg/L: Ag, Al, As, Ba, Be, Ca, Cd, Co, Cr, Cu, Fe, K, Mg, Mn, Mo, Na, Ni, Pb, Sb, Se, Sn, Sr, Ti, Tl, V, Zn	125 mL	N9301721
<b>Cold Plasma Solution</b>			
0.1% HNO <sub>3</sub>	1 µg/L Li, Co, Rb, In, Cs, Pb	250 mL	N8151032

## Reference Standards

### Single-Element Reference Standards, 1,000 µg/mL – Pure Plus Grade 125 mL

Element	Symbol	Matrix	Part No.
Aluminum	Al	2% HNO <sub>3</sub>	N9303726
Antimony	Sb	H <sub>2</sub> O/0.6% Tart. Acid/Tr. HNO <sub>3</sub>	N9303750
Arsenic	As	2% HNO <sub>3</sub>	N8151035
Barium	Ba	2% HNO <sub>3</sub>	N9303726
Beryllium	Be	2% HNO <sub>3</sub>	N8151035
Bismuth	Bi	2% HNO <sub>3</sub>	N9303726
Cadmium	Cd	2% HNO <sub>3</sub>	N8151035
Calcium	Ca	2% HNO <sub>3</sub>	N9303726
Chromium	Cr	2% HNO <sub>3</sub>	N8151035
Cobalt	Co	2% HNO <sub>3</sub>	N9303726
Copper	Cu	2% HNO <sub>3</sub>	N8151035
Iron	Fe	2% HNO <sub>3</sub>	N9303726
Lead	Pb	2% HNO <sub>3</sub>	N8151035
Magnesium	Mg	2% HNO <sub>3</sub>	N9303726
Manganese	Mn	2% HNO <sub>3</sub>	N8151035
Mercury	Hg	10% HNO <sub>3</sub>	N9303726
Molybdenum	Mo	H <sub>2</sub> O	N8151035
Nickel	Ni	2% HNO <sub>3</sub>	N9303726
Potassium	K	2% HNO <sub>3</sub>	N8151035
Selenium	Se	2% HNO <sub>3</sub>	N9303726
Silver	Ag	2% HNO <sub>3</sub>	N8151035
Sodium	Na	2% HNO <sub>3</sub>	N9303726
Tin	Sn	1% HNO <sub>3</sub> /1% HF	N8151035
Thallium	Tl	2% HNO <sub>3</sub>	N9303726
Thorium	Th	2% HNO <sub>3</sub>	N8151035
Titanium	Ti	H <sub>2</sub> O/0.24% F	N9303726
Uranium	U	2% HNO <sub>3</sub>	N8151035
Vanadium	V	2% HNO <sub>3</sub>	N9303726
Zinc	Zn	2% HNO <sub>3</sub>	N8151035

### Single-Element Pure Plus ICP-MS Grade Standards – 1 µg/mL

Element	Symbol	Matrix	Pure Plus Grade 125 mL Part No.
Aluminum	Al	2% HCl	N9304200
Aluminum	Al	2% HNO <sub>3</sub>	N9304201
Antimony	Sb	H <sub>2</sub> O/Tr. HNO <sub>3</sub> /Tr. Tart	N9304203
Arsenic	As	2% HCl	N9304204
Arsenic	As	2% HNO <sub>3</sub>	N9304205
Barium	Ba	2% HNO <sub>3</sub>	N9304207
Beryllium	Be	2% HNO <sub>3</sub>	N9304208
Bismuth	Bi	2% HNO <sub>3</sub>	N9304209
Boron	B	H <sub>2</sub> O	N9304210
Cadmium	Cd	2% HNO <sub>3</sub>	N9304211
Cadmium	Cd	2% HNO <sub>3</sub>	N9304212
Calcium	Ca	5% HCl	N9304213
Calcium	Ca	2% HNO <sub>3</sub>	N9304214
Cerium	Ce	2% HNO <sub>3</sub>	N9304216
Cesium	Cs	2% HNO <sub>3</sub>	N9304217
Chromium	Cr	2% HCl	N9304218
Chromium	Cr	5% HNO <sub>3</sub>	N9304219
Cobalt	Co	2% HCl	N9304221
Cobalt	Co	2% HNO <sub>3</sub>	N9304222
Copper	Cu	2% HCl	N9304223
Copper	Cu	2% HNO <sub>3</sub>	N9304224
Gadolinium	Gd	2% HNO <sub>3</sub>	N9304228
Gallium	Ga	2% HNO <sub>3</sub>	N9304229
Germanium	Ge	H <sub>2</sub> O	N9304230
Gold	Au	1% HNO <sub>3</sub> /3% HCL	N9304231
Iridium	Ir	2% HCl	N9304235
Iron	Fe	2% HCl	N9304236
Iron	Fe	2% HNO <sub>3</sub>	N9304237
Lanthanum	La	2% HNO <sub>3</sub>	N9304238
Lead	Pb	2% HNO <sub>3</sub>	N9304239
Lithium	Li	2% HNO <sub>3</sub>	N9304241
Magnesium	Mg	2% HNO <sub>3</sub>	N9304244
Manganese	Mn	2% HNO <sub>3</sub>	N9304245

## Reference Standards

Single-Element Pure Plus ICP-MS Grade Standards –  
1 µg/mL *continued...*

Element	Symbol	Matrix	Pure Plus Grade 125 mL Part No.
Potassium	K	2% HNO <sub>3</sub>	N9304248
Molybdenum	Mo	H <sub>2</sub> O	N9304249
Niobium	Nb	H <sub>2</sub> O/tr HF	N9304250
Nickel	Ni	2% HNO <sub>3</sub>	N9304252
Palladium	Pd	2% HCl	N9304253
Phosphorus	P	H <sub>2</sub> O	N9304254
Platinum	Pt	2% HCl	N9304255
Rubidium	Rb	2% HNO <sub>3</sub>	N9304257
Rhenium	Re	H <sub>2</sub> O	N9304258
Ruthenium	Ru	2% HCl	N9304260
Scandium	Sc	2% HNO <sub>3</sub>	N9304262
Selenium	Se	2% HNO <sub>3</sub>	N9304263
Silicon	Si	H <sub>2</sub> O/tr. F-	N9304264
Silver	Ag	2% HNO <sub>3</sub>	N9304265
Sodium	Na	2% HCl	N9304266
Strontium	Sr	2% HNO <sub>3</sub>	N9304269
Sulfur	S	H <sub>2</sub> O	N9304270
Tantalum	Ta	H <sub>2</sub> O/tr HF	N9304271
Tellurium	Te	5% HCl	N9304272
Thallium	Tl	2% HNO <sub>3</sub>	N9304275
Thorium	Th	2% HNO <sub>3</sub>	N9304276
Thulium	Tm	2% HNO <sub>3</sub>	N9304277
Tin	Sn	2% HNO <sub>3</sub> /tr HF	N9304279
Titanium	Ti	H <sub>2</sub> O/tr HF	N9304280
Tungsten	W	2% HNO <sub>3</sub> /tr HF	N9304281
Uranium	U	2% HNO <sub>3</sub>	N9304283
Vanadium	V	2% HNO <sub>3</sub>	N9304285
Ytterbium	Yb	2% HNO <sub>3</sub>	N9304286
Yttrium	Y	2% HNO <sub>3</sub>	N9304287
Zinc	Zn	2% HNO <sub>3</sub>	N9304289
Zirconium	Zr	2% HNO <sub>3</sub>	N9304291

Single-Element Pure Plus ICP-MS Grade Standards –  
10 µg/mL

Element	Symbol	Matrix	Pure Plus Grade 125 mL Part No.
Bismuth	Bi	2% HNO <sub>3</sub>	N9303731
Germanium	Ge	H <sub>2</sub> O tr HF	N9303739
Indium	In	2% HNO <sub>3</sub>	N9303741
Rhodium	Rh	2% HCl	N9303749
Scandium	Sc	2% HNO <sub>3</sub>	N9303751
Terbium	Tb	2% HNO <sub>3</sub>	N9303753
Yttrium	Y	2% HNO <sub>3</sub>	N9303757

## Environmental Analysis Standards

We offer a variety of environmental analysis standards for all your environmental application needs.

Matrix	Contents	Vol.	Part No.
<b>Environmental Standard Kit for NexION Standard Mode</b> Includes 1 bottle of each:			<b>N9307118</b>
5% HNO <sub>3</sub>	1,000 µg/mL: Ca, K, Mg, Na	125 mL	N9307805
5% HNO <sub>3</sub>	1,000 µg/mL: Al, Fe	125 mL	N9307806
5% HNO <sub>3</sub> / trace Tartaric Acid/ trace HF	100 µg/mL: Ag, Al, As, Ba, Be, Ca, Cd, Co, Cr, Cu, Fe, K, Mg, Mn, Mo, Na, Ni, Pb, Sb, Se, Sn, Sr, Ti, Tl, V, Zn	125 mL	N9301721
2% HNO <sub>3</sub>	100 µg/mL: B, Th, U	125 mL	N9307807
5% HNO <sub>3</sub>	10 µg/mL: Hg	125 mL	N9300253
2% HNO <sub>3</sub>	50 µg/mL: Sc, Ge 10 µg/mL: Rh, In, Tb	125 mL	N9308591
<b>Environmental Standards for NexION KED/DRC Mode</b>			
5% HNO <sub>3</sub>	1,000 µg/mL: Ca, K, Mg, Na	125 mL	N9307805
5% HNO <sub>3</sub>	1,000 µg/mL: Al, Fe	125 mL	N9307806
5% HNO <sub>3</sub> / trace Tartaric Acid/ trace HF	100 µg/mL: Ag, Al, As, Ba, Be, Ca, Cd, Co, Cr, Cu, Fe, K, Mg, Mn, Mo, Na, Ni, Pb, Sb, Se, Sn, Sr, Ti, Tl, V, Zn	125 mL	N9301721
2% HNO <sub>3</sub>	100 µg/mL: B, Th, U	125 mL	N9307807
5% HNO <sub>3</sub>	10 µg/mL: Hg	125 mL	N9300253
2% HNO <sub>3</sub>	100 µg/mL: Sc 50 µg/mL: Ge 1 µg/mL: Rh, In, Tb	125 mL	N9308592
<b>Additional Internal Standard Mixes</b>			
5% HNO <sub>3</sub> / trace HCL	50 µg/mL: Sc 20 µg/mL: Ge 10 µg/mL: In, Ir, Li6, Rh, Tb, Y	125 mL	N9307808
5% HNO <sub>3</sub>	10 µg/mL: Bi, Ge, In, Li6, Sc, Tb, Y	125 mL	N9303832
5% HNO <sub>3</sub>	10 µg/mL: Bi, Ge, In, Li6, Sc, Tb, Y	500 mL	N9303833
2% HNO <sub>3</sub>	10 µg/mL: Bi, Ho, In, Li6, Sc, Tb, Y	125 mL	N9303834
2% HNO <sub>3</sub>	100 µg/L: Li6	125 mL	N9303955
<b>NexION 5000 Setup Solution</b>			
1% HNO <sub>3</sub>	200 ng/L Be, Ce, Fe, In, Li, Mg, Pb, U	500 mL	N8162301
<b>NexION 5000 Stability Solution</b>			
1% HNO <sub>3</sub>	200 ng/L Cd, Cr, Fe, Li, Mg, Pb 1 µg/L Cu	500 mL	N8162302
<b>NexION 5000 DRC NH<sub>3</sub> and Standard Detection Limit</b>			
1% HNO <sub>3</sub>	200 ng/L Al, Be, Ca, Cr, Cu, Fe, In, K, Mn, Ni, U, Zn	125 mL	N8162303

Matrix	Contents	Vol.	Part No.
<b>Contract Lab Program Modification Standards</b>			
2% HNO <sub>3</sub> / 5% HCl	10 µg/mL: Ba, Be, Ce, Co, In, Li, Mg, Pb, Rh, Tl, U, Y	125 mL	N9303843
2% HNO <sub>3</sub> / trace Tartaric Acid	20 µg/mL: Ag, Al, As, Ba, Be, Cd, Co, Cr, Cu, Mn, Mo, Ni, Pb, Sb, Se, Th, Tl, U, V, Zn	125 mL	N9303816
5% HNO <sub>3</sub> / trace Tartaric Acid/trace HF	100 µg/mL: Ag, Al, As, Ba, Be, Ca, Cd, Co, Cr, Cu, Fe, K, Mg, Mn, Mo, Na, Ni, Pb, Sb, Se, Sn, Sr, Ti, Tl, V, Zn	125 mL	N9301721
2% HNO <sub>3</sub> / trace Tartaric Acid/trace HF	10 µg/mL: Ag, Al, As, Ba, Be, Cd, Co, Cr, Cu, Mn, Ni, Pb, Sb, Se, Tl, V, Zn	125 mL	N9303821
5% HNO <sub>3</sub>	10 µg/mL: Hg	125 mL	N9300253
2% HNO <sub>3</sub>	200 µg/mL: Ca, Fe, K, Mg, Na	125 mL	N9303822
2% HNO <sub>3</sub> / trace HF	10 µg/mL: Mo, Sn, Sr, Ti	125 mL	N9303824
5% HNO <sub>3</sub> / trace HF	21,215 µg/mL: Cl 3,000 µg/mL: Ca 2,500 µg/mL: Fe, Na 2,000 µg/mL: C 1,000 µg/mL: Al, K, Mg, P, S 20 µg/mL: Mo, Ti	125 mL	N9303827
5% HNO <sub>3</sub> / trace Tartaric Acid/ trace HF	500 µg/mL: Fe 250 µg/mL: Ba, Zn 100 µg/mL: Co, Cr, Cu, Mn, Ni, Sb, V 50 µg/mL: As, Pb 25 µg/mL: Ag, Be, Cd, Se, Tl	125 mL	N9303839
5% HNO <sub>3</sub> / trace Tartaric Acid/ trace HF	250 µg/mL: Ba, Cr, Cu, Zn 150 µg/mL: V 125 µg/mL: Ni 100 µg/mL: Co, Pb, Sb 50 µg/mL: As, Cd 5 µg/mL: Ag, Be, Se, Tl	125 mL	N9303840
5% HNO <sub>3</sub>	1,000 µg/mL: Al, Ca, Fe, K, Mg, Na 20 µg/mL: Ag, As, Ba, Be, Cd, Co, Cr, Cu, Mn, Ni, Pb, Se, Tl, V, Zn	125 mL	N9303835



## Pharmaceutical Analysis Standards

### USP Standards

As a part of our United States Pharmacopeia (USP) compliant offering, we provide calibration and check standards for the analysis of metals in drugs, pharmaceutical substances and raw materials to verify Oral Daily Dose PDE, Parenteral Component Limit and Parenteral Daily Dose PDE.

Our trace metal USP standards can be used to calibrate arsenic, cadmium, lead and mercury, the four toxic metals as part of the USP regulations. When you couple our accurate and reliable USP standards with your ICP-MS, you will remain compliant with the new and changing regulations.

### ICH Standards

In conjunction with our United States Pharmacopeia (USP) calibration standards, we offer six calibration standards for the analysis of metals in pharmaceutical materials and products following the International Conference on Harmonization (ICH) guidelines.

The ICH standards verify elemental impurities, precious metals impurities and parenteral elemental impurities. Couple our accurate and reliable ICH standards with your ICP-MS and you will remain compliant with the new and changing regulations.

### USP and ICH Standards



Matrix	Content	Volume	Part No.
ICH Elemental Impurities Standards Kit – Oral PDEs Kit Includes:			N9304361
5% HNO <sub>3</sub> / 2% HCl	ICH Class 1 Elements + TI – Oral PDE: 30 µg/mL Hg 15 µg/mL As 8 µg/mL TI 5 µg/mL Cd, Pb	125 mL	N9304362
2% HNO <sub>3</sub>	ICH Class 2A Elements – Oral and Parenteral PDEs: 200 µg/mL Ni 100 µg/mL V 50 µg/mL Co	125 mL	N9304363
2% HCl	100 µg/mL Au	125 mL	N9303728
15% HCl	ICH Class 2B Precious Metal Elements (no Os) – All PDEs: 100 µg/mL Ir, Pd, Pt, Rh, Ru	125 mL	N9304364
2% HNO <sub>3</sub>	ICH Class 2B Se and Ag – Oral PDE: 150 µg/mL Ag, Se	125 mL	N9304365
5% HNO <sub>3</sub> /trace Tartaric Acid/ trace HF	ICH Class 3 Elements – Oral PDEs: 1100 µg/mL Cr 600 µg/mL Sn 300 µg/mL Cu, Mo 140 µg/mL Ba 120 µg/mL Sb 55 µg/mL Li	125 mL	N9304366
2% HNO <sub>3</sub>	Internal Standards – All PDEs: 50 µg/mL Ge 20 µg/mL Ga 1 µg/mL In, Tb	125 mL	N9304367

Matrix	Content	Volume	Part No.
ICH Elemental Impurities Standards – Parenteral PDEs			
5% HNO <sub>3</sub> / 2% HCl	ICH Class 1 Elements + TI – Parenteral PDE: 15 µg/mL As 8 µg/mL TI 5 µg/mL Pb 3 µg/mL Hg 2 µg/mL Cd	125 mL	N9304369
2% HNO <sub>3</sub>	ICH Class 2A Elements – Oral and Parenteral PDEs: 200 µg/mL Ni 100 µg/mL V 50 µg/mL Co	125 mL	N9304363
2% HCl	100 µg/mL Au	125 mL	N9303728
15% HCl	ICH Class 2B Precious Metal Elements (no Os) – All PDEs: 100 µg/mL Ir, Pd, Pt, Rh, Ru	125 mL	N9304364
2% HNO <sub>3</sub>	ICH Class 2B Se and Ag – Parenteral PDE: 80 µg/mL Se 10 µg/mL Ag	125 mL	N9304372
2% HNO <sub>3</sub> /trace Tartaric Acid/ trace HF	ICH Class 3 Elements – Parenteral PDEs: 150 µg/mL Mo 110 µg/mL Cr 60 µg/mL Sn 30 µg/mL Cu 70 µg/mL Ba 25 µg/mL Li 9 µg/mL Sb	125 mL	N9304373
2% HNO <sub>3</sub>	Internal Standards – All PDEs: 50 µg/mL Ge 20 µg/mL Ga 1 µg/mL In, Tb	125 mL	N9304367



## Consumables and Supplies Reference Guide

### USP and ICH Standards *continued...*

Matrix	Content	Volume	Part No.
ICH Elemental Impurities Standards – Inhalational PDEs			
2% HNO <sub>3</sub>	Internal Standards – All PDEs: 50 µg/mL Ge 20 µg/mL Ga 1 µg/mL In, Tb	125 mL	N9304367
2% HCl	100 µg/mL Au	125 mL	N9303728
15% HCl	ICH Class 2B Precious Metal Elements (no Os) – All PDEs: 100 µg/mL Ir, Pd, Pt, Rh, Ru	125 mL	N9304364

Matrix	Content	Volume	Part No.
ICH Class 2B Precious Metal Elements – All PDEs			
15% HCl	ICH Class 2B Precious Metal Elements – All PDEs: 100 µg/mL Ir, Os, Pd, Pt, Rh, Ru	125 mL	N9304382
Osmium			
15% HCl	100 µg/mL Os	125 mL	N9304383

## Gold Nanoparticle Standards

Fully characterized gold nanospheres designed for Single Particle ICP-MS. These particle standards have a nearly perfect spherical shape, narrow size (CV < 5%) and particle number distributions (CV < 5%).

Description	Diameter (nm)	Volume (mL)	Part No.
Super Uniform Gold Nanospheres Particle Standard 1e7 particles/mL, PEG Carboxyl 0.8 kDa	30	25 mL	N8151034
Super Uniform Gold Nanospheres Particle Standard 1e7 particles/mL, PEG Carboxyl 0.8 kDa	50	25 mL	N8151035
Super Uniform Gold Nanospheres Particle Standard 1e7 particles/mL, PEG Carboxyl 0.8 kDa	100	25 mL	N8151036



Super Uniform Gold Nanospheres

## Other Replacement Parts

Description	Part No.	Description	Part No.
Metal Spill Tray for Vacuum Pump	N8145022	Helium Cell Gas Filter Assembly with Ultra-Torr Fittings	N8146004
Vacuum Fluid Pump – Fomblin® 1 L	N8145003	Helium Cell Gas Filter Assembly with VCR Fittings	N8150302
Instrument Filter (Left) for NexION 1000 and 2000 ICP-MS	N8152167	Replacement Filters for Helium Cell Gas Filtration System	N8146005
Instrument Filter (Right) for NexION 1000, 2000 and 2200 ICP-MS	N8152136	Roller head replacement part for Peristaltic Pump Assembly	N8152770
NexION Sample Spill Tray for NexION 1000 and 2000 ICP-MS	N8150361	Green Coolant for GreenCT Cooling System	N8171229
NexION Sample Spill Tray for NexION 2200 and 5000 ICP-MS	N8160361	ICP-MS Chiller Coolant Mix	WE016558
		Air Filter with Frame for Whispercool Chiller, 13 in. x 14 in.	N0777095
		Passive Air Filter for Durachill Chiller	N0772058

## NexION 1100/2200/5000 ICP-MS Consumable Kits

Typical one year's supply of consumables. Consists of the following parts:

Configurations			1100	2200	5000S
Description	Qty.	Part No.	N8170104	N8170105	N8160027
Fixed 2.0 mm Injector Quartz Torch (Blue/Purple Mark)	2	N8152472	✓	✓	
Fixed 2.0 mm Injector Ultra-Pure Quartz Torch (White Mark)	2	N8152428			✓
Glass Cyclonic High Sensitivity Spray Chamber with Matrix Gas Port	1	N8152389	✓	✓	
SilQ Cyclonic Spray Chamber with Matrix Gas Port	1	N8152539			✓
Threaded Spray Chamber Drain Line	2	N8152456	✓	✓	✓
Nickel Skimmer Cone	2	W1026356	✓		
Nickel Skimmer Cone	2	N8161117			
Nickel Skimmer Cone	2	N8171142		✓	
Platinum Skimmer Cone for NexION 5000	2	N8161041			✓
Nickel Sampler Cone	2	W1033612	✓	✓	
Platinum Sampler Cone	2	N8161140			✓
Hyper-Skimmer Cone for NexION 1000/1100/2000	1	W1033995	✓		
Hyper-Skimmer Cone for NexION 2200/5000	1	N8160120		✓	✓
Sampler Cone Gasket	3	W1040148	✓	✓	✓
Hyper-Skimmer Cone O-Ring (0.614 i.d. 0.070WD)	2	09902123	✓		
Hyper-Skimmer O-Ring (0.926 i.d. 0.070WD VITON 60 DUR)	3	09211389		✓	✓
Gray/Gray Pump Tubing	2	N8152415	✓	✓	✓
Flared PVC 0.51 mm Orange/Yellow Pump Tubing	4	N8152405	✓	✓	✓
Drain Tubing (feet)	6	02506495	✓	✓	✓
Sample Uptake Tubing	12	WE024375	✓	✓	✓
Sample Introduction Line for PFA-ST Nebulizer	1	N8152455			✓
Sample Introduction Line for Concentric Nebulizer	1	N8145503	✓	✓	
Female Barb CTFE Fitting	4	N8152388	✓	✓	✓
Male Barb CTFE Fitting	4	N8152453	✓	✓	✓
Instrument Filter (Right)	2	N8152136	✓	✓	✓
Instrument Filter (Rear)	2	N8152465	✓	✓	✓
Torch Injector Nut	2	N8152451	✓	✓	✓
Self-Aspirating Probe, Purple Marker, 0.25 mm i.d.	1	N8152512			✓
Self-Aspirating Probe, Purple Marker, 0.3 mm i.d.	1	N8152513			✓
Internal Standards Probe, 0.5 mm i.d.	2	N8152452			

## NexION 1000/2000 ICP-MS Consumable Kits

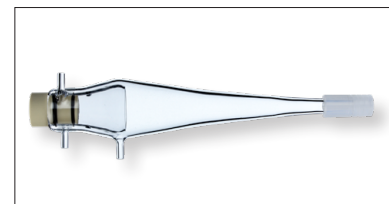
Typical one year's supply of consumables. Consists of the following parts:

Configurations			1000	2000B	2000C	2000P	2000S
Description	Qty.	Part No.	N8150058	N8150052	N8150055	N8150056	N8150057
Fixed 2.0 mm Injector Quartz Torch (Blue/Purple Mark)	2	N8152472	✓	✓			
Fixed 2.5 mm Injector Quartz Torch (Green Mark)	2	N8152473			✓		
Demountable Quartz Torch PVDF	2	N8152379				✓	
2.0 mm Demountable Quartz Injector	2	N8152376				✓	
Fixed 2.0 mm Injector UHP Quartz Torch	2	N8152428					✓
Glass Cyclonic High Sensitivity Spray Chamber with Matrix Gas Port	1	N8152389	✓	✓	✓	✓	
SilQ Cyclonic High Sensitivity Spray Chamber	1	N8152424					✓
SilQ Cyclonic Spray Chamber with Matrix Gas Port	1	N8152539					✓
Threaded Spray Chamber Drain Line	2	N8152456	✓	✓	✓	✓	✓
Nickel Skimmer Cone	2	W1026356	✓	✓	✓	✓	
Nickel Sampler Cone	2	W1033612	✓	✓	✓	✓	
Platinum Skimmer Cone	2	W1026907					✓
Platinum Sampler Cone	2	N8161140					✓
Hyper-Skimmer Cone for NexION 1000/1100/2000	1	W1033995	✓	✓	✓	✓	✓
Sampler Cone Gasket	3	W1040148	✓	✓	✓	✓	✓
Hyper-Skimmer Cone O-Ring (0.614 i.d. 0.070WD)	2	09902123	✓	✓	✓	✓	✓
Gray/Gray Pump Tubing	2	N8152415	✓	✓	✓		✓
Gray/Gray Pump Tubing MP2	2	N8145173				✓	
Flared PVC 0.51 mm Orange/Yellow Pump Tubing	4	N8152405	✓	✓	✓		✓
Flared PVC 0.51 mm Orange/Yellow Pump Tubing MP2	4	N8145150				✓	
Drain Tubing (feet)	6	02506495	✓	✓	✓	✓	✓
Sample Uptake Tubing	12	WE024375	✓	✓	✓	✓	✓
Sample Introduction Line for PFA-ST Nebulizer	1	N8152455				✓	✓
Sample Introduction Line for Concentric Nebulizer	1	N8145503	✓	✓	✓		
Female Barb CTFE Fitting	4	N8152388	✓	✓	✓	✓	✓
Male Barb CTFE Fitting	4	N8152453	✓	✓	✓	✓	✓
Instrument Filter (Left)	2	N8152167	✓	✓	✓	✓	✓
Instrument Filter (Right)	2	N8152136	✓	✓	✓	✓	✓
Instrument Filter (Rear)	2	N8152465	✓	✓	✓	✓	✓
Torch Injector Nut	2	N8152451	✓	✓	✓	✓	✓

## Single Cell Sample Introduction Kit

The Single Cell Application Module within PerkinElmer's Syngistix™ for ICP-MS software is the ideal tool for laboratories analyzing metal content in cells, either from uptake of metals or metallic nanoparticles in solution or measuring the intrinsic metal content of the cell. This unique software module allows the differentiation and quantification between the dissolved and cellular fractions of the same analyte. In a single analysis, cell metal or particulate content, cell-containing metal concentration, mass per cell, and mass distribution can all be determined without the need for subsequent data processing. Coupled with the NexION ICP-MS series, the Syngistix Single Cell Application Module is the world's first single cell ICP-MS dedicated software, delivering speed, flexibility, automation, and ease-of-use.

Description	Qty.	Part No.
Single Cell Kit for NexION 1000/2000/2200/5000		N8150032
<b>Kit Includes:</b>		
Asperon™ Spray Chamber	1	N8152493
Spray Chamber Drain Line	1	N8152494
Spray Chamber Addition Gas Tee 1/8 Fitting	1	N8152495
Nebulizer HEN	1	N8152496
Sample Line for HEN Nebulizer	1	N8152489
UV Light Shield	1	N8152497
Plasma Light Shield	1	N8152377
Fixed 2.0 mm Injector UHP Quartz Torch	1	N8152428
Torch Cassette White	1	N8152432
Torch Injector Nut	1	N8152451
Syngistix Single Cell Software	1	N8150321
CTFE Male Barb Fitting	2	N8145017
CTFE Female Barb Fitting	2	N8145018
Orange/Red 0.19 mm i.d. PVC Pkg.12	1	N8152401
MP2 PVC Flared 0.19 mm Orange/Red Pkg.12	1	N8145195
Description	Part No.	
Single Cell Automation Kit for NexION 1000/2000/2200/5000	N8150039	
<b>Kit Includes:</b>		
Includes Single Cell Micro DX Autosampler. Does not include Single Cell Kit (N8150032)		



Asperon™ Spray Chamber



Single Cell Automation Kit for NexION 1000/2000/2200/5000 (N8150039)

[SHOP ONLINE](#)

[www.perkinelmer.com/supplies](http://www.perkinelmer.com/supplies)

PerkinElmer U.S. LLC  
710 Bridgeport Ave.  
Shelton, CT 06484-4794 USA  
(+1) 855-726-9377  
[www.perkinelmer.com](http://www.perkinelmer.com)



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright ©2024, PerkinElmer U.S. LLC. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer U.S. LLC. All other trademarks are the property of their respective owners.

84114 (45246)