

Inorganic Aqueous Standards



Whatever your application, you can depend on PerkinElmer Pure and Pure Plus Standards to yield reliable, accurate results.

Table of Contents

	Page
Single-Element Standards.....	2
Mixed Calibration Standards	3
Initial Calibration Verification Standards.....	3
Matrix Blanks	3
Quality Control Standards	3
Instrument Calibration Standards for CLP	3
GFAAS Mixed Standard.....	4
AA Test Mix.....	4
Contract Required Detection Limits (CRDL)	4
Matrix Modifiers for Graphite Furnace AA	4
Reagents.....	4
Instrument Check Standards	4
Interference Check Standards.....	4-5
Universal Data Acquisition Standards Kit.....	5
Water Pollutant Standards.....	5
Instrument Setup Solutions	5
Spike Sample Analysis.....	5-6
Internal Standard Solutions	6
Isotope Standard	6
Wash Standards.....	6
Multi-Element Standards.....	6-7
Instrument Calibration Standards	7-8
SmartTune Standards	8
Environmental Method Sets	8-10
Toxicity Characteristic Leachate Procedure (TCLP)	10
Performance Verification Standards	11
USP Standards Kit.....	11
Part Number Index.....	12

PerkinElmer offers a complete selection of atomic spectroscopy aqueous standards. Each solution is supplied with a comprehensive Certificate of Analysis that documents the quality and reliability.

Pure Grade Standards for AA and ICP-OES

- Analyzed by ICP-OES
- Analyzed by Classical Wet Assay
- 32 trace impurities analyzed by ICP-MS of the final solution and reported on the certificate
- Impurities reported at ppm level
- All Standards are prepared and certified under ISO Guide 34 and ISO 17025 – certified by A2LA



Pure Plus Grade Standards for ICP-MS

- Analyzed by ICP-OES
- Analyzed by Classical Wet Assay
- 67 trace impurities analyzed by ICP-MS of the final solution and reported on the certificate
- Impurities reported at ppb level
- All Standards are prepared and certified under ISO Guide 34 and ISO 17025 – certified by A2LA

PerkinElmer Pure Inorganic Aqueous Standards are compatible with other manufacturers' instruments.

Single-Element Standards – 1,000 mg/L

Element	Symbol	Matrix	Pure Grade 125 mL Part No.	Pure Grade 500 mL Part No.	Pure Plus Grade 125 mL Part No.
Aluminum	Al	2% HNO ₃	N9300184	N9300100	N9303726
Antimony	Sb	2% HNO ₃	N9300207	N9300101	N9303750
Arsenic	As	2% HNO ₃	N9300180	N9300102	N9303727
Barium	Ba	2% HNO ₃	N9300181	N9300103	N9303729
Beryllium	Be	2% HNO ₃	N9300172	N9300104	N9303730
Bismuth	Bi	10% HNO ₃	N9303761	N9300105	N9303731 ¹
Boron	B	H ₂ O	N9303760	N9300106	
Cadmium	Cd	2% HNO ₃	N9300176	N9300107	N9303734
Calcium	Ca	2% HNO ₃	N9303763	N9300108	N9303733
Carbon	C	H ₂ O	N9303762	N9300109	
Cerium	Ce	2% HNO ₃	N9303765	N9300110	
Cesium	Cs	2% HNO ₃	N9303767	N9300111	
Chromium	Cr	2% HNO ₃	N9300173	N9300112	N9303736
Cobalt	Co	2% HNO ₃	N9303766	N9300113	N9303735
Copper	Cu	2% HNO ₃	N9300183	N9300114	N9303737
Dysprosium	Dy	2% HNO ₃	N9303768	N9300115	
Erbium	Er	2% HNO ₃	N9303769	N9300116	
Europium	Eu	2% HNO ₃	N9303770	N9300117	
Gadolinium	Gd	2% HNO ₃	N9303773	N9300118	
Gallium	Ga	2% HNO ₃	N9303772	N9300119	
Germanium	Ge	H ₂ O/0.16% F-	N9303774	N9300120	N9303739 ¹
Gold	Au	10% HCl	N9303759	N9300121	N9303728 ²
Hafnium	Hf	2% HCl	N9303775	N9300122	
Holmium	Ho	2% HNO ₃	N9303776	N9300123	
Indium	In	2% HNO ₃	N9303777	N9300124	N9303741 ¹
Iridium	Ir	10% HCl	N9303778	N9300125	
Iron	Fe	2% HNO ₃	N9303771	N9300126	N9303738
Lanthanum	La	2% HNO ₃	N9303780	N9300127	
Lead	Pb	2% HNO ₃	N9300175	N9300128	N9303748
Lithium	Li	2% HNO ₃	N9303781	N9300129	
Lutetium	Lu	2% HNO ₃	N9303782	N9300130	
Magnesium	Mg	2% HNO ₃	N9300179	N9300131	N9303743
Manganese	Mn	2% HNO ₃	N9303783	N9300132	N9303744
Mercury	Hg	10% HNO ₃	N9300174	N9300133	N9303740 ¹
Molybdenum	Mo	H ₂ O	N9303784	N9300134	N9303745
Neodymium	Nd	2% HNO ₃	N9303787	N9300135	
Nickel	Ni	2% HNO ₃	N9300177	N9300136	N9303747
Niobium	Nb	H ₂ O/0.4% HF	N9303786	N9300137	
Palladium	Pd	10% HCl	N9303789	N9300138	
Phosphorus	P	H ₂ O	N9303788	N9300139	
Platinum	Pt	10% HCl	N9303791	N9300140	
Potassium	K	2% HNO ₃	N9303779	N9300141	N9303742
Praseodymium	Pr	2% HNO ₃	N9303790	N9300142	
Rhenium	Re	H ₂ O	N9303793	N9300143	
Rhodium	Rh	10% HCl	N9303794	N9300144	N9303749 ¹
Rubidium	Rb	2% HNO ₃	N9303792	N9300145	
Ruthenium	Ru	10% HCl	N9303795	N9300146	
Samarium	Sm	2% HNO ₃	N9303800	N9300147	
Scandium	Sc	2% HNO ₃	N9303798	N9300148	N9303751 ¹
Selenium	Se	2% HNO ₃	N9300182	N9300149	N9303752
Silicon	Si	H ₂ O	N9303799	N9300150	
Silver	Ag	2% HNO ₃	N9300171	N9300151	N9303725
Sodium	Na	2% HNO ₃	N9303785	N9300152	N9303746
Strontium	Sr	2% HNO ₃	N9303802	N9300153	
Sulfur	S	H ₂ O	N9303796	N9300154	
Tantalum	Ta	H ₂ O/0.8% HF	N9303803	N9300155	
Tellurium	Te	5% HNO ₃	N9303805	N9300156	
Terbium	Tb	2% HNO ₃	N9303804	N9300157	N9303753 ¹
Tin	Sn	20% HCl	N9303801	N9300161	N9303838
Thallium	Tl	2% HNO ₃	N9300170	N9300158	N9303755
Thorium	Th	2% HNO ₃	N9303842		
Thulium	Tm	2% HNO ₃	N9303807	N9300160	
Titanium	Ti	H ₂ O/tr 0.24% F-	N9303806	N9300162	N9303754
Tungsten	W	H ₂ O	N9303809	N9300163	
Uranium	U	2% HNO ₃		N9303844	
Vanadium	V	2% HNO ₃	N9303808	N9300165	N9303756
Ytterbium	Yb	2% HNO ₃	N9303811	N9300166	
Yttrium	Y	2% HNO ₃	N9303810	N9300167	N9303757 ¹
Zinc	Zn	2% HNO ₃	N9300178	N9300168	N9303758
Zirconium	Zr	2% HNO ₃	N9303812	N9300169	

¹ 10 mg/L

² 100 mg/L

Mixed Calibration Standards

Matrix	Contents	Vol.	Part No.
Mixed Calibration Standard			
2% HNO ₃	50 mg/L: As, K 10 mg/L: La, Li, Mn, Ni, Sr, Zn 1 mg/L: Ba, Mg	500 mL	N0691579
Mixed Calibration Standard 1			
2% HNO ₃	500 mg/L: Pb 200 mg/L: Se 150 mg/L: Cd, Zn 100 mg/L: Mn 50 mg/L: Be	125 mL	N9300200
Mixed Calibration Standard 2			
5% HNO ₃	10,000 mg/L: Fe 100 mg/L: Ba, Co, Cu, V	125 mL	N9300201
Mixed Calibration Standard 3			
2% HNO ₃ /tr HF	500 mg/L: As 100 mg/L: Mo, Si	125 mL	N9300202
Mixed Calibration Standard 4			
5% HNO ₃	1,000 mg/L: Ca 400 mg/L: K 200 mg/L: Al, Na 20 mg/L: Cr, Ni	125 mL	N9300203
Mixed Calibration Standard 5			
5% HNO ₃ /tr Tartaric Acid/tr HF	1,000 mg/L: Mg 200 mg/L: Sb, Tl 100 mg/L: B 50 mg/L: Ag	125 mL	N9300204

Initial Calibration Verification Standards

Matrix	Contents	Vol.	Part No.
5% HNO ₃	500 mg/L: Ca, K, Mg, Na 200 mg/L: Al, Ba 100 mg/L: Fe 60 mg/L: Sb 50 mg/L: Co, V 40 mg/L: Ni 25 mg/L: Cu 20 mg/L: Zn 15 mg/L: Mn 10 mg/L: Ag, As, Cr, Tl 5 mg/L: Cd, Se 3 mg/L: Pb	500 mL	N9300224
5% HNO ₃	500 mg/L: Ca, K, Mg, Na 200 mg/L: Al, Ba 100 mg/L: Fe 60 mg/L: Sb 50 mg/L: Co, V 40 mg/L: Ni 25 mg/L: Cu 20 mg/L: Zn 15 mg/L: Mn 10 mg/L: Ag, As, Cr, Tl 5 mg/L: Cd, Se 3 mg/L: Pb	125 mL	N9303953

Matrix Blanks

Matrix	Contents	Vol.	Part No.
Water Blank			
ASTM® Type I Water	18 megohm	125 mL	N9300237
ASTM® Type I Water	18 megohm	250 mL	N9303813
Hydrochloric Acid Blank			
2% HCl in ASTM® Type I Water		125 mL	N9300238
2% Nitric Acid Blank			
2% HNO ₃ in ASTM® Type I Water		125 mL	N0773120
1% Nitric Acid Blank			
1% HNO ₃ in ASTM® Type I Water		125 mL	N9303732

Quality Control Standards

Matrix	Contents	Vol.	Part No.
Quality Control Standard, 21 Elements Pure (Pure XVI)			
5% HNO ₃ /tr Tartaric Acid/tr HF	100 mg/L: As, Be, Ca, Cd, Co, Cr, Cu, Fe, Li, Mg, Mn, Mo, Ni, Pb, Sb, Se, Sr, Ti, Tl, V, Zn	125 mL	N9300281
Quality Control Standard, 21 Elements Pure Plus			
5% HNO ₃ /tr Tartaric Acid/tr HF	100 mg/L: As, Be, Ca, Cd, Co, Cr, Cu, Fe, Li, Mg, Mn, Mo, Ni, Pb, Sb, Se, Sr, Ti, Tl, V, Zn	125 mL	N9303837
Quality Control Standard, 7A Elements			
5% HNO ₃ /tr HF	1000 mg/L: K 500 mg/L: Si 100 mg/L: Al, B, Ba, Na 50 mg/L: Ag	125 mL	N9300280

Instrument Calibration Standards for CLP

Matrix	Contents	Vol.	Part No.
Instrument Calibration Standard 1			
5% HNO ₃	5,000 mg/L: Ca, K, Mg, Na	125 mL	N9300218
Instrument Calibration Standard 2			
5% HNO ₃	400 mg/L: Ni 200 mg/L: Zn 150 mg/L: Mn 100 mg/L: Ag, Cr	125 mL	N9300219
Instrument Calibration Standard 3			
5% HNO ₃	2,000 mg/L: Al, Ba 1,000 mg/L: Fe 500 mg/L: Co, V 250 mg/L: Cu 50 mg/L: Be	125 mL	N9300220
Instrument Calibration Standard 4			
5% HNO ₃	100 mg/L: As, Tl 50 mg/L: Cd, Se 30 mg/L: Pb	125 mL	N9300221

GFAAS Mixed Standard

Matrix	Contents	Vol.	Part No.
5% HNO ₃ w/ trace of HF	100 mg/L: Al, As, Pb, Sb, Se, Tl 50 mg/L: Ba, Co, Cu, Ni 20 mg/L: Cr, Fe, Mn 10 mg/L: Ag 5 mg/L: Be, Cd	125 mL	N9300244

AA Test Mix

Matrix	Contents	Vol.	Part No.
2% HCl	50 mg/L: Ca, Cr, Cu, Fe, Ni 20 mg/L: K 10 mg/L: Na, Zn	125 mL	02900540

Contract Required Detection Limits (CRDL)

Matrix	Contents	Vol.	Part No.
5% HNO ₃ /tr Tartaric Acid/tr HF	120 mg/L: Sb 100 mg/L: Co, V 80 mg/L: Ni 50 mg/L: Cu 40 mg/L: Zn 30 mg/L: Mn 20 mg/L: Ag, As, Cr, Tl 10 mg/L: Be, Cd, Se 6 mg/L: Pb	125 mL	N9300225

Matrix Modifiers for Graphite Furnace AA

Matrix	Contents	Vol.	Part No.
Mg(NO ₃) ₂	1% Mg (NO ₃) ₂ (as nitrate)	100 mL	B0190634
Pd	1% Pd (as nitrate)	50 mL	B0190635
NH ₄ H ₂ PO ₄	10% NH ₄ H ₂ PO ₄	100 mL	N9303445

Reagents

Matrix	Contents	Vol.	Part No.
Triton® X-100 Wetting Agent		100 mL	N9300260
Antifoaming Silicone Emulsion		500 mL	B0507226
Glycerol		1 L	B3141064

Instrument Check Standards

Matrix	Contents	Vol.	Part No.
Instrument Check Standard 1			
2% HNO ₃ /tr Tartaric Acid/tr HF	10 mg/L: Ag, Al, As, Ba, Be, Cd, Co, Cr, Cu, Mn, Ni, Pb, Sb, Se, Tl, V, Zn	125 mL	N9303821
Instrument Check Standard 3			
2% HNO ₃	200 mg/L: Ca, Fe, K, Mg, Na	125 mL	N9303822
Instrument Check Standard 4			
2% HNO ₃	10 mg/L: Mo, Th, U	125 mL	N9303823
Instrument Check Standard 5			
2% HNO ₃ /tr HF	10 mg/L: Mo, Sn, Sr, Ti	125 mL	N9303824

Interference Check Standards

Matrix	Contents	Vol.	Part No.
Interference Check Standard 1			
H ₂ O/tr HNO ₃ /0.6% Tartaric Acid	1,000 mg/L: Sb	125 mL	N9300207
Interference Check Standard 5			
5% HNO ₃	6,000 mg/L: Ca 5,000 mg/L: Fe 3,000 mg/L: Mg 1,200 mg/L: Al 1,000 mg/L: Na	125 mL	N9300208
Interference Check Standard 18			
5% HNO ₃	20,000 mg/L: K 1,000 mg/L: As, Pb, Tl 500 mg/L: Se 300 mg/L: Ag, Ba, Cd, Co, Cr, Cu, Ni, V, Zn 200 mg/L: Mn 100 mg/L: Be, Hg*	125 mL	N9300205

*Supplied in separate bottle.

Interferents A

5% HNO ₃	5,000 mg/L: Al, Ca, Mg 2,000 mg/L: Fe	500 mL	N9300226
---------------------	--	--------	-----------------

Alternate Interferents A

5% HNO ₃	1,000 mg/L: Cr, Cu, Mn, Ni, Ti, V	500 mL	N9300228
---------------------	--------------------------------------	--------	-----------------

Analytes B

5% HNO ₃ /tr Tartaric Acid/tr HF	100 mg/L: Cd, Ni, Zn 60 mg/L: Sb 50 mg/L: Ba, Be, Co, Cr, Cu, Mn, V 20 mg/L: Ag 10 mg/L: As, Tl 5 mg/L: Pb, Se	125 mL	N9300227
---	---	--------	-----------------

Alternate Analytes B

5% HNO ₃ /tr Tartaric Acid/tr HF	100 mg/L: Al, As, B, Mo, Na, Sb, Se, Tl 10 mg/L: Ca, Fe, Mg, Si	125 mL	N9300229
---	--	--------	-----------------

Interference Check Standards

H ₂ O/tr HNO ₃ /0.6% Tartaric Acid	1000 mg/L: Sb	500 mL	N9303797
--	---------------	--------	-----------------

Interference Check Standards, *continued*

Matrix	Contents	Vol.	Part No.
Interference Check Solution 1 (for SW-846 & ILM 05.2)			
5% HNO ₃ /tr HF	10,000 mg/L: Cl 2000 mg/L: C 1000 mg/L: Al, Ca, Fe, K, Mg, Na, P, S 20 mg/L: Mo, Ti	125 mL	N9303828
Interference Check Solution 2 (for SW-846)			
2% HNO ₃	10 mg/L: Ag, As, Cd, Co, Cr, Cu, Mn, Ni, Zn	125 mL	N9303830
Analytes C (for ILM 05.2)			
2% HNO ₃ /tr Tartaric Acid/tr HF	2 mg/L: Ag, As, Ba, Be, Cd, Co, Cr, Cu, Hg*, Mn, Ni, Pb, Sb, Se, Tl, V, Zn	125 mL	N9303831

*Supplied in separate bottle.

Universal Data Acquisition Standards Kit

Matrix	Contents	Vol.	Part No.
Solution Kit (Includes items listed below)			N9306225
5% HNO ₃	10 mg/L: Ce, Dy, Er, Eu, Gd, Ho, La, Lu, Nd, Pr, Sm, Sc, Tb, Th, Tm, Y, Yb	125 mL	N9300232
5% HNO ₃	10 mg/L: Al, Ag, 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	N9300233
10% HCl/ 1% HNO ₃	10 mg/L: Au, Hf, Ir, Pd, Pt, Rh, Ru, Sb, Sn, Te	125 mL	N9300234
5% HNO ₃	10 mg/L: Hg	125 mL	N9300253
H ₂ O/tr HF/tr HNO ₃	10 mg/L: B, Ge, Mo, Nb, P, Re, S, Si, Ta, Ti, W, Zr	125 mL	N9300235

Water Pollutant Standards

Matrix	Contents	Vol.	Part No.
Primary Drinking Water Metals			
2% HNO ₃	100 mg/L: Ba 10 mg/L: Ag, As, Cr, Hg*, Pb 5 mg/L: Cd, Se	125 mL	N9300216
Secondary Drinking Water Metals			
2% HNO ₃	500 mg/L: Zn 100 mg/L: Cu 30 mg/L: Fe 5 mg/L: Mn	125 mL	N9300217
Trace Metals I			
5% HNO ₃	500 mg/L: Al 250 mg/L: V 100 mg/L: As, Be, Co, Cr, Cu, Fe, Mn, Ni, Pb, Zn 25 mg/L: Cd, Se 10 mg/L: Hg*	125 mL	N9300211

*Supplied in separate bottle.

Water Pollutant Standards, *continued*

Matrix	Contents	Vol.	Part No.
Trace Metals II			
2% HNO ₃	20 mg/L: Sb, Tl 10 mg/L: Ag	125 mL	N9300212
Trace Metals III			
2% HNO ₃	500 mg/L: Ba, Ca, Mo, Na 100 mg/L: K, Mg	125 mL	N9300213
Alternate Metals I			
2% HNO ₃	20 mg/L: Al, Fe, V 10 mg/L: Co, Cu, Mn, Ni, Zn 5 mg/L: Be, Sb, Tl	125 mL	N9300214
Alternate Metals II			
2% HNO ₃	500 mg/L: Ca, Na 100 mg/L: K, Mg	125 mL	N9300215
2% HNO ₃	500 mg/L: Ca, Na 100 mg/L: K, Mg	500 mL	N9303952

Instrument Setup Solutions

Matrix	Contents	Vol.	Part No.
VIS Wavecal Solution			
2% HNO ₃	50 mg/L: K 10 mg/L: La, Li, Mn, Na, Sr 1 mg/L: Ba, Ca	250 mL	N9302946
UV Wavecal Solution			
5% HCl	100 mg/L: K, P, S 20 mg/L: As, La, Li, Mn, Mo, Na, Ni, Sc	250 mL	N0681470
5% HCl	100 mg/L: K, P, S 20 mg/L: As, La, Li, Mn, Mo, Na, Ni, Sc 1 mg/L: Ca	500 mL	N0582152
Low UV Standard			
2% HNO ₃	10 mg/L: Al, P, S	250 mL	N0691580
Calcium Stray Light Standard			
H ₂ O	10,000 mg/L: Ca	125 mL	N0691581

Spike Sample Analysis

Matrix	Contents	Vol.	Part No.
Spike Sample Standard I			
5% HNO ₃ /tr Tartaric Acid/tr HF	200 mg/L: Al, As, Ba, Se, Tl 100 mg/L: Fe 50 mg/L: Co, Mn, Ni, Pb, Sb, V, Zn 25 mg/L: Cu 20 mg/L: Cr 5 mg/L: Ag, Be, Cd	125 mL	N9300230

Spike Sample Analysis, *continued*

Matrix	Contents	Vol.	Part No.
Spike Sample Standard 1 (water)			
5% HNO ₃ /tr Tartaric Acid/tr HF	500 mg/L: Fe 250 mg/L: Ba, Zn 100 mg/L: Co, Cr, Cu, Mn, Ni, Sb, V 50 mg/L: As, Pb 25 mg/L: Ag, Be, Cd, Se, Tl	125 mL	N9303839
Spike Sample Standard 2 (soil)			
5% HNO ₃ /tr Tartaric Acid/tr HF	250 mg/L: Ba, Cr, Cu, Zn 150 mg/L: V 125 mg/L: Ni 100 mg/L: Co, Pb, Sb 50 mg/L: As, Cd 25 mg/L: Ag, Be, Se, Tl	125 mL	N9303840
Spike Sample Standard 3 (for ILM 05.2)			
5% HNO ₃ /tr Tartaric Acid/tr HF	200 mg/L: Al, Ba 50 mg/L: Co, Mn, Ni, V, Zn 25 mg/L: Cu 20 mg/L: Cr 10 mg/L: Sb 5 mg/L: Ag, Be, Cd, Tl 4 mg/L: As 2 mg/L: Pb 1 mg/L: Se	125 mL	N9303841

Internal Standard Solutions

Matrix	Contents	Vol.	Part No.
Multi-Element Internal Standard			
2% HNO ₃	10 mg/L: Bi, Ho, In, ⁶ Li, Sc, Tb, Y	125 mL	N9303834
Internal Standard Mix			
5-10% HNO ₃	10 mg/L: Bi, Ge, In, ⁶ Li, Sc, Tb, Y	125 mL	N9303832
Internal Standard Mix			
5-10% HNO ₃	10 mg/L: Bi, Ge, In, ⁶ Li, Sc, Tb, Y	500 mL	N9303833

Isotope Standard

Matrix	Contents	Vol.	Part No.
Lithium 6 Standard			
2% HNO ₃	100 mg/L: ⁶ Li	125 mL	N9303955

Wash Standards

Matrix	Vol.	Part No.
ELAN DRC Wash Solution		
0.5% HNO ₃	250 mL	N8125033
ELAN 9000/6X00/DRC-e Wash Solution		
1% HNO ₃	1000 mL	N8122038
NexION Wash Solution		
1% HNO ₃	250 mL	N8145050
Water Blank		
ASTM® Type I Water, 18 megohm	125 mL	N9303814

Multi-Element Standards

Matrix	Contents	Vol.	Part No.
Multi-Element Solution			
5% HNO ₃	1000 mg/L: Al, Ca, Fe, Mg	500 mL	N9307113
Multi-Element Solution			
5% HNO ₃	1000 mg/L: K, Na, P	500 mL	N9307114
Multi-Element Solution			
2-5% HNO ₃	1000 mg/L: Mo, Sb, Sn, W, Zr	500 mL	N9307115
Multi-Element Solution			
5% HNO ₃	1000 mg/L: As, Ba, Be, Cd, Cr, Co, Cu, La, Pb, Li, Mn, Ni, Sc, Sr, V, Y, Zn	500 mL	N9307116
Multi-Element Solution 1			
2% HNO ₃	10 mg/L: Be, Bi, Ce, Co, In, Mg, Ni, Pb, U	125 mL	N9300231
Multi-Element Solution 2			
5% HNO ₃	10 mg/L: Ce, Dy, Er, Eu, Gd, Ho, La, Lu, Nd, Pr, Sc, Sm, Tb, Th, Tm, Y, Yb	125 mL	N9300232
Multi-Element Solution 3			
5% HNO ₃	10 mg/L: Ag, Al, As, Ba, Be, Bi, Ca, Cd, Co, Cr, Cs, Cu, Fe, Ga, Hg*, In, K, Li, Mg, Mn, Na, Ni, Pb, Rb, Se, Sr, Tl, U, V, Zn	125 mL	N9300233
<i>*Supplied in separate bottle.</i>			
Multi-Element Solution 3 without Mercury			N9301720
Multi-Element Solution 4			
10% HCl/1% HNO ₃	10 mg/L: Au, Hf, Ir, Pd, Pt, Rh, Ru, Sb, Sn, Te	125 mL	N9300234
Multi-Element Solution 5			
H ₂ O/tr HF/tr HNO ₃	10 mg/L: B, Ge, Mo, Nb, P, Re, S, Si, Ta, Ti, W, Zr	125 mL	N9300235
Set of Multi-Element Solutions			
Includes: (1 bottle of each)			N9300236
	Multi-Element Solution 2	125 mL	N9300232
	Multi-Element Solution 3	125 mL	N9300233
	Multi-Element Solution 4	125 mL	N9300234
	Multi-Element Solution 5	125 mL	N9300235
	Water Blank	125 mL	N9300237
	Hydrochloric Acid Blank	125 mL	N9300238
	2% Nitric Acid Blank	125 mL	N9300239
PerkinElmer Pure I			
5% HNO ₃	400 mg/L: Tl 200 mg/L: Bi, In, Pb 150 mg/L: Ga 100 mg/L: Al 50 mg/L: Ag, Ni 25 mg/L: Cr 20 mg/L: Cd, Co, Cu, Zn 15 mg/L: B, Fe 5 mg/L: Ba, Mn 1 mg/L: Be, Sr	125 mL	N9303940

Multi-Element Standards, *continued*

Matrix	Contents	Vol.	Part No.
PerkinElmer Pure IV (Quality Control Standard 23)			
10% HNO ₃	1000 mg/L: Ag, Al, B, Ba, Bi, Ca, Cd, Co, Cr, Cu, Fe, Ga, In, K, Li, Mg, Mn, Na, Ni, Pb, Sr, Tl, Zn	125 mL	N9303941
PerkinElmer Pure VIII			
5% HNO ₃	100 mg/L: Al, B, Ba, Be, Bi, Ca, Cd, Co, Cr, Cu, Fe, Ga, K, Li, Mg, Mn, Na, Ni, Pb, Se, Sr, Te, Tl, Zn	125 mL	N9303942
PerkinElmer Pure IX			
5% HNO ₃	100 mg/L: As, Be, Pb, Cd,	125 mL	N9303943
PerkinElmer Pure X			
2% HNO ₃	35,000 ug/L: Ca 15,000 ug/L: Mg 8,000 ug/L: Na 3,000 ug/L: K 100 ug/L: B, Fe, Mo, Sr 50 ug/L: As, Ba, Ni, V, Zn 30 ug/L: Mn 25 ug/L: Co, Pb 20 ug/L: Be, Cd, Cr, Cu 10 ug/L: Bi, Se, Tl	125 mL	N9303944
8 mg/L Mercury in 5% HNO₃			
5% HNO ₃	8 mg/L: Hg	125 mL	N9303954
PerkinElmer Pure XI			
5% HNO ₃	2500 mg/L: Zn 900 mg/L: Cr, Pb 800 mg/L: Cu 200 mg/L: Ni 10 mg/L: Cd	125 mL	N9303945
5 mg/L Mercury in 5% HNO₃			
5% HNO ₃	5 mg/L: Hg in 5% HNO ₃	125 mL	N9303949
PerkinElmer Pure XIII			
5% HNO ₃	500 mg/L: Al 250 mg/L: V 100 mg/L: As, Be, Co, Cr, Cu, Fe, Mn, Ni, Pb, Zn 25 mg/L: Cd, Se	125 mL	N9303946
PerkinElmer Pure XVI (Quality Control Standard, 21 Elements)			
5% HNO ₃ /tr Tartaric Acid/tr HF	100 mg/L: As, Be, Ca, Cd, Co, Cr, Cu, Fe, Li, Mg, Mn, Mo, Ni, Pb, Sb, Se, Sr, Ti, Tl, V, Zn	125 mL	N9300281
PerkinElmer Pure XVII			
15% HCl/trace HF	100 mg/L: Hf, Ir, Sb, Sn, Ta, Tl, Zr	125 mL	N9303948

Instrument Calibration Standards

Matrix	Contents	Vol.	Part No.
Instrument Calibration Standard 1			
5% HNO ₃ /tr Tartaric Acid	20 mg/L: Ag, Al, As, Ba, Be, Cd, Co, Cr, Cu, Mn, Mo, Ni, Pb, Sb, Se, Th, Tl, U, V, Zn	125 mL	N9303816
Instrument Calibration Standard 2			
5% HNO ₃ /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
Instrument Calibration Standard 3			
5% HNO ₃	1,000 mg/L: Ca, Fe, K, Mg, Na	125 mL	N9303818
Initial Calibration Verification Standard 1			
5% HNO ₃ /tr Tartaric Acid	1,000 mg/L: Ca, Fe, K, Mg, Na, Sr 10 mg/L: Ag, Al, As, Ba, Be, Cd, Co, Cr, Cu, Mn, Mo, Ni, Pb, Sb, Se, Tl, V, Zn, Th, U	125 mL	N9303825
Initial Calibration Verification Standard 2			
2% HNO ₃ /tr HF	10 mg/L: Sn, Ti	125 mL	N9303826
ELAN 6100 DRC Setup/Stab/Masscal Solution			
0.5% HNO ₃	10 µg/L: Ba 1 µg/L: Al, Cd, Ce, Cr, Cu, In, Pb, Mg, Mn, Rh, Th	1,000 mL	N8125035
ELAN DRC/DRCplus/DRC II Solution Kit			
Solution Kit (Includes items listed below)			N8120541
2 x 1,000 mL: Setup/Stability/Masscal Solution			N8125035
2 x 250 mL: Wash Solution			N8125033
1 x 250 mL: Sensitivity/Detection Limit Solution			N8125034
1 x 125 mL: Methanol Blank Solution			N8125037
1 x 125 mL: Chromium in Methanol Solution			N8125038
ELAN 9000/6100 Solution Kit			
Solution Kit (Includes items listed below)			N8120522
2 x 1,000 mL: Setup/Stability/Masscal Solution			N8125030
1 x 125 mL: Dual-Detector Calibration Solution			N8125010
1 x 1000 mL: Wash Solution			N8122038
1 x 250 mL: Detection Limit Solution			N8125031
ELAN 9000/6100 Setup/Stability/Masscal Solution			
1% HNO ₃	10 µg/L: Ba, Cd, Ce, Cu, In, Pb, Mg, Rh, U	1,000 mL	N8125030
ELAN 9000/6X00 Dual-Detector Calibration Solution			
1% HNO ₃	200 µg/L: Cd, Cu, Pb, Mg, Rh	1,000 mL	N8125032
ELAN 6000/5000 Plasma Setup Solution			
2% HNO ₃	10 µg/L: Ba, Cd, Ce, Cu, Ge, Pb, Mg, Rh, Sc, Tb, Tl	1,000 mL	N8122014

Instrument Calibration Standards, *continued*

Matrix	Contents	Vol.	Part No.
ELAN 5000 Detection Limit Solution			
2% HNO ₃	10 µg/L: Be, Co, Ge, In, Tl, U	100 mL	N8122017

SmartTune Standards

Matrix	Contents	Vol.	Part No.
SmartTune Solution for Standard ELANs/DRC-e			
1% HNO ₃	10 µg/L: Ba, Be, Ce, Co, In, Pb, Mg, Rh, U	1,000 mL	N8125040
SmartTune Solution for DRC/DRCplus/DRC II			
0.5% HNO ₃	10 µg/L: Ba 1 µg/L: Be, Ce, Co, In, Fe, Pb, Mg, Th, U	1,000 mL	N8125041
Tuning Solution 1			
2% HNO ₃ /5% HCl	10 mg/L: Ba, Be, Ce, Co, In, Li, Mg, Pb, Rh, Tl, U, Y	125 mL	N9303843
NexION Setup Solution			
1% HNO ₃	1 µg/L: Be, Ce, Fe, In, Li, Mg, Pb, U	500 mL	N8145051

Environmental Method Sets

Description	Contents	Vol.	Part No.
Environmental Standard Kit for non-DRC/Standard ICP-MS Instruments			
1 bottle each			
	1,000 mg/L: Ca, K, Mg, Na	125 mL	N9307805
	1,000 mg/L: Al, Fe	125 mL	N9307806
	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
	100 mg/L: B, Th, U	125 mL	N9307807
	10 mg/L: Hg	125 mL	N9300253
	Internal Standard Mix	125 mL	N9307808
	50 mg/L: Sc		
	20 mg/L: Ge		
	10 mg/L: In, Ir, ⁶ Li, Rh, Tb, Y		
Environmental Standard Kit for DRC Instruments			
1 bottle each			
	1,000 mg/L: Ca, K, Mg, Na	125 mL	N9307805
	1,000 mg/L: Al, Fe	125 mL	N9307806
	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
	100 mg/L: B, Th, U	125 mL	N9307807
	10 mg/L: Hg	125 mL	N9300253
	Internal Standard Mix	125 mL	N9307738
	200 mg/L: Sc		
	20 mg/L: Ga		
	10 mg/L: In, Ir, Rh, Tm		

Environmental Method Sets, *continued*

Description	Contents	Vol.	Part No.
Contract Lab Program Modification Set			
1 bottle each			
2% HNO ₃ /5% HCl	10 mg/L: Ba, Be, Ce, Co, In, Li, Mg, Pb, Rh, Tl, U, Y	125 mL	N9303843
2% HNO ₃ /tr Tartaric Acid	20 mg/L: 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 ₃ /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
5% HNO ₃ /tr Tartaric Acid	500 mg/L: Ca, K, Mg, Na 20 mg/L: Al, Ba 10 mg/L: Fe 6 mg/L: Sb 5 mg/L: Co, V 4 mg/L: Ni 2.5 mg/L: Cu 2 mg/L: Zn 1.5 mg/L: Mn 1 mg/L: Ag, As, Cr, Tl 0.5 mg/L: Be, Cd, Se 0.3 mg/L: Pb	125 mL	N9303819
2% HNO ₃ /tr Tartaric Acid/tr HF	10 mg/L: Ag, Al, As, Ba, Be, Cd, Co, Cr, Cu, Mn, Ni, Pb, Sb, Se, Tl, V, Zn	125 mL	N9303821
5% HNO ₃	10 mg/L: Hg	125 mL	N9300253
2% HNO ₃	200 mg/L: Ca, Fe, K, Mg, Na	125 mL	N9303822
2% HNO ₃	10 mg/L: Mo, Th, U	125 mL	N9303823
2% HNO ₃ /tr HF	10 mg/L: Mo, Sn, Sr, Ti	125 mL	N9303824
5% HNO ₃ /tr HF	21215 mg/L: Cl 3000 mg/L: Ca 2500 mg/L: Fe, Na 2000 mg/L: C 1000 mg/L: Al, K, Mg, P, S 20 mg/L: Mo, Ti	125 mL	N9303827
2% HNO ₃	20 mg/L: Co, Cr, Cu, Mn, Ni, V 10 mg/L: As, Cd, Se, Zn 5 mg/L: Ag	125 mL	N9303829
5% HNO ₃ /tr Tartaric Acid/tr HF	500 mg/L: Fe 250 mg/L: Ba, Zn 100 mg/L: Co, Cr, Cu, Mn, Ni, Sb, V 50 mg/L: As, Pb 25 mg/L: Ag, Be, Cd, Se, Tl	125 mL	N9303839
5% HNO ₃ /tr Tartaric Acid/tr HF	250 mg/L: Ba, Cr, Cu, Zn 150 mg/L: V 125 mg/L: Ni 100 mg/L: Co, Pb, Sb 50 mg/L: As, Cd 25 mg/L: Ag, Be, Se, Tl	125 mL	N9303840

Environmental Method Sets, *continued*

Description	Contents	Vol.	Part No.
Contract Lab Program Modification Set, cont.			
5% HNO ₃	1000 mg/L: Al, Ca, Fe, K, Mg, Na 20 mg/L: Ag, As, Ba, Be, Cd, Co, Cr, Cu, Mn, Ni, Pb, Se, Tl, V, Zn	125 mL	N9303835
H ₂ O/tr HF	7200 mg/L: Cl 2000 mg/L: C 1000 mg/L: P, S 20 mg/L: Mo, Sb, Ti	125 mL	N9303836
Calibration Standards Set for Method 6010			N9307104
1 bottle each			
5% HNO ₃	5000 mg/L: Ca, K, Mg, Na	125 mL	N9300218
5% HNO ₃	400 mg/L: Ni 200 mg/L: Zn 150 mg/L: Mn 100 mg/L: Ag, Cr	125 mL	N9300219
5% HNO ₃	2000 mg/L: Al, Ba 1000 mg/L: Fe 500 mg/L: Co, V 250 mg/L: Cu 50 mg/L: Be	125 mL	N9300220
5% HNO ₃	100 mg/L: As, Tl 50 mg/L: Cd, Pb, Se	125 mL	N9300221
H ₂ O/tr HF	10 mg/L: B, Ge, Mo, Nb, P, Re, S, Si, Ta, Ti, W, Zr	125 mL	N9300235
2% HNO ₃	1000 mg/L: Li	500 mL	N9300129
2% HNO ₃	1000 mg/L: Sb	500 mL	N9300101
2% HNO ₃	1000 mg/L: Sr	500 mL	N9300153
20% HCl	1000 mg/L: Sn	500 mL	N9300161
2% HNO ₃	1000 mg/L: Cr	500 mL	N9300112
2% HNO ₃	1000 mg/L: Cu	500 mL	N9300114
2% HNO ₃	1000 mg/L: Mn	500 mL	N9300132
2% HNO ₃	1000 mg/L: Ni	500 mL	N9300136
H ₂ O/tr HF	1000 mg/L: Ti	500 mL	N9300162
2% HNO ₃	1000 mg/L: V	500 mL	N9300165
2% HNO ₃	1000 mg/L: Al	500 mL	N9300100
2% HNO ₃	1000 mg/L: Ca	500 mL	N9300108
2% HNO ₃	1000 mg/L: Fe	500 mL	N9300126
2% HNO ₃	1000 mg/L: Mg	500 mL	N9300131
Internal Standard Set for Method 6010 & 200.7			N9307105
1 bottle each			
2% HNO ₃	1000 mg/L: Y	500 mL	N9300167
2% HNO ₃	1000 mg/L: Sc	500 mL	N9300148

Environmental Method Sets, *continued*

Description	Contents	Vol.	Part No.
Interference Solutions for Method 6010			N9307106
1 bottle each			
2% HNO ₃	1000 mg/L: Al	125 mL	N9300184
2% HNO ₃	1000 mg/L: Ca	500 mL	N9300108
2% HNO ₃	1000 mg/L: Mg	125 mL	N9300179
2% HNO ₃	1000 mg/L: Fe	500 mL	N9300126
2% HNO ₃	1000 mg/L: K	500 mL	N9300141
2% HNO ₃	1000 mg/L: Na	500 mL	N9300152
2% HNO ₃	1000 mg/L: Cr	500 mL	N9300112
2% HNO ₃	1000 mg/L: Cu	500 mL	N9300114
2% HNO ₃	1000 mg/L: Mn	500 mL	N9300132
2% HNO ₃	1000 mg/L: Ni	500 mL	N9300136
H ₂ O/tr HF	1000 mg/L: Ti	500 mL	N9300162
2% HNO ₃	1000 mg/L: V	500 mL	N9300165
5% HNO ₃	5000 mg/L: Al, Ca, Mg 2000 mg/L: Fe	500 mL	N9300226
5% HNO ₃ / tr Tart-HF	100 mg/L: Cd, Ni, Zn 60 mg/L: Sb 50 mg/L: Ba, Be, Co, Cr, Cu, Mn, V 20 mg/L: Ag 10 mg/L: As, Tl 5 mg/L: Pb, Se	125 mL	N9300227
5% HNO ₃	1000 mg/L: Cr, Cu, Mn, Ni, Ti, V	500 mL	N9300228
5% HNO ₃ /tr Tartaric Acid/tr HF	100 mg/L: Al, As, B, Mo, Na, Sb, Se, Tl 10 mg/L: Ca, Fe, Mg, Si	125 mL	N9300229
Calibration Standards for Method 200.7			N9307107
5% HNO ₃	5000 mg/L: Ca, K, Mg, Na	125 mL	N9300218
5% HNO ₃	400 mg/L: Ni 200 mg/L: Zn 150 mg/L: Mn 100 mg/L: Ag, Cr	125 mL	N9300219
5% HNO ₃	2000 mg/L: Al, Ba 1000 mg/L: Fe 500 mg/L: Co, V 250 mg/L: Cu 50 mg/L: Be	125 mL	N9300220
5% HNO ₃	100 mg/L: As, Tl 50 mg/L: Cd, Pb, Se	125 mL	N9300221
H ₂ O/tr HF	10 mg/L: B, Ge, Mo, Nb, P, Re, S, Si, Ta, Ti, W, Zr	125 mL	N9300235
2% HNO ₃	1000 mg/L: Li	500 mL	N9300129
2% HNO ₃	1000 mg/L: Sb	500 mL	N9300101
2% HNO ₃	1000 mg/L: Sr	500 mL	N9300153
20% HCl	1000 mg/L: Sn	500 mL	N9300161

Environmental Method Sets, *continued*

Description	Contents	Vol.	Part No.
Interference Solutions for Method 200.7			N9307108
2% HNO ₃	1000 mg/L: Al	125 mL	N9300184
2% HNO ₃	1000 mg/L: Ca	500 mL	N9300108
2% HNO ₃	1000 mg/L: Mg	125 mL	N9300179
2% HNO ₃	1000 mg/L: Fe	500 mL	N9300126
2% HNO ₃	1000 mg/L: K	500 mL	N9300141
2% HNO ₃	1000 mg/L: Na	500 mL	N9300152
5% HNO ₃	5000 mg/L: Al, Ca, Mg 2000 mg/L: Fe	500 mL	N9300226
5% HNO ₃ /tr Tart-HF	100 mg/L: Cd, Ni, Zn 60 mg/L: Sb 50 mg/L: Ba, Be, Co, Cr, Cu, Mn, V 20 mg/L: Ag 10 mg/L: As, Tl 5 mg/L: Pb, Se	125 mL	N9300227
Environmental EPA Set 1			N9307110
2% HNO ₃	500 mg/L: Pb 200 mg/L: Se 150 mg/L: Cd, Zn 100 mg/L: Mn 50 mg/L: Be	125 mL	N9300200
5% HNO ₃	10,000 mg/L: Fe 100 mg/L: Ba, Co, Cu, V	125 mL	N9300201
2% HNO ₃ /tr HF	500 mg/L: As 100 mg/mL: Mo, Si	125 mL	N9300202
5% HNO ₃	1000 mg/L: Ca 400 mg/L: K 200 mg/L: Al, Na 20 mg/L: Cr, Ni	125 mL	N9300203
5% HNO ₃ / tr Tart-HF	1000 mg/L: Mg 200 mg/L: Sb, Tl 100 mg/L: B 50 mg/L: Ag	125 mL	N9300204
5% HNO ₃	20,000 mg/L: K 1000 mg/L: As, Pb, Tl 500 mg/L: Se 300 mg/L: Ag, Ba, Cd, Co, Cr, Cu, Ni, V, Zn 200 mg/L: Mn 100 mg/L: Be	125 mL	N9300205
5% HNO ₃	100 mg/L: Hg	125 mL	N9300223
5% HNO ₃	6000 mg/L: Ca 5000 mg/L: Fe 3000 mg/L: Mg 1200 mg/L: Al 1000 mg/L: Na	125 mL	N9300208
H ₂ O/tr HNO ₃ / 0.6% Tartaric Acid	1000 mg/L: Sb	125 mL	N9300207
5% HNO ₃	ASTM® Type I Water	500 mL	N9307809
5% HCl	ASTM® Type I Water	500 mL	N9307810

Environmental Method Sets, *continued*

Description	Contents	Vol.	Part No.
Environmental EPA Set 2			N9307109
2% HNO ₃	500 mg/L: Pb 200 mg/L: Se 150 mg/L: Cd, Zn 100 mg/L: Mn 50 mg/L: Be	125 mL	N9300200
5% HNO ₃	10,000 mg/L: Fe 100 mg/L: Ba, Co, Cu, V	125 mL	N9300201
2% HNO ₃ /tr HF	500 mg/L: As 100 mg/L: Mo, Si	125 mL	N9300202
5% HNO ₃	1000 mg/L: Ca 400 mg/L: K 200 mg/L: Al, Na 20 mg/L: Cr, Ni	125 mL	N9300203
5% HNO ₃ /tr Tart-HF	1000 mg/L Mg 200 mg/L: Sb, Tl 100 mg/L: B 50 mg/L: Ag	125 mL	N9300204
5% HNO ₃	20,000 mg/L: K 1000 mg/L: As, Pb, Tl 500 mg/L: Se 300 mg/L: Ag, Ba, Cd, Co, Cr, Cu, Ni, V, Zn 200 mg/L: Mn 100 mg/L: Be	125 mL	N9300205
5% HNO ₃	6000 mg/L: Ca 5000 mg/L: Fe 3000 mg/L: Mg 1200 mg/L: Al 1000 mg/L Na	125 mL	N9300208
H ₂ O/tr HNO ₃ / 0.6% Tartaric Acid	1000 mg/L: Sb	125 mL	N9300207
5% HNO ₃	ASTM® Type I Water	500 mL	N9307809
5% HCl	ASTM® Type I Water	500 mL	N9307810

Toxicity Characteristic Leachate Procedure (TCLP) Standard

Description	Contents	Vol.	Part No.
TCLP Standard 1			
2% HNO ₃	500 mg/L: Ba 25 mg/L: Ag, As, Cr, Pb 100 mg/L: Hg* 5 mg/L: Cd, Se	500 mL	N9300241

*Supplied in separate bottle.

Performance Verification Standards

Matrix	Contents	Vol.	Part No.
Methanol Blank Solution			
1% Semiconductor Grade		125 mL	N8125037
Chromium in Methanol Solution			
1% Semiconductor Grade Methanol		125 mL	N8125038
	10 mg/L: Cr		
Selenium Solution			
5% HNO ₃	10 mg/L: Se	250 mL	N8125039
NexION 300Q Solution Kit – Non-Cell			N8140503
Solution Kit (Includes items listed below)			
	1 x 250 mL NexION Wash Solution		N8145050
	1 X 500 mL NexION Setup Solution		N8145051
	1 X 500 mL NexION 300Q Non-cell Stability Solution		N8145053
	1 x 100 mL NexION Standard/DRC Mode Detection Limit Blank Solution		N8145055
	1 X 100 mL NexION Standard/DRC Mode Detection Limit Solution		N8145056
	1 x 100 mL NexION Dual Detector Solution		N8145059
NexION 300 X/D/S Solution Kit – Cell Instruments			N8140504
Solution Kit (Includes items listed below)			
	1 x 250 mL NexION Wash Solution		N8145050
	1 X 500 mL NexION Setup Solution		N8145051
	1 x 250 mL NexION KED Setup Solution		N8145052
	1 x 500 mL NexION 300X/D/S Cell Stability Solution		N8145054
	1 x 100 mL NexION Standard/DRC Mode Detection Limit Blank Solution		N8145055
	1 x 100 mL NexION Standard/DRC Mode Detection Limit Solution		N8145056
	1 x 100 mL NexION KED Mode Detection Limit Blank Solution		N8145057
	1 x 100 mL NexION KED Mode Detection Limit Solution		N8145058
	1 x 100 mL NexION Dual Detector Solution		N8145059
NexION Setup Solution			
1% HNO ₃	1 µg/L: Be, Ce, Fe, In, Li, Mg, Pb, U	500 mL	N8145051
NexION KED Setup Solution			
1% HCl	10 µg/L: Co 1 µg/L: Ce	250 mL	N8145052
NexION 300Q Non-cell Stability Solution			
1% HNO ₃	1 µg/L: Cd, Cu, Mg, Pb	500 mL	N8145053
NexION 300X/D/S Cell Stability Solution			
1% HNO ₃	1 µg/L: Cd, Cr, Fe, In, Mg, Pb 10 µg/L: Co, Cu, Se	500 mL	N8145054

Performance Verification Standards, *continued*

Matrix	Contents	Vol.	Part No.
NexION Standard/DRC Mode Detection Limit Blank Solution			
0.5% HNO ₃		100 mL	N8145055
NexION Standard/DRC Mode Detection Limit Solution			
0.5% HNO ₃	1 µg/L: Be, Ca, Co, Fe, In, U	100 mL	N8145056
NexION KED Mode Detection Limit Blank Solution			
1% HCl		100 mL	N8145057
NexION KED Mode Detection Limit Solution			
1% HCl	10 µg/L: V, As, Se	100 mL	N8145058
NexION Dual Detector Solution			
2% HNO ₃	50 µg/L: Al, Ba, Ce, Co, In, Li, Mg, Mn, Ni, Pb, Tb, U, Zn	100 mL	N8145059
NexION AFT Single-Element Solution			
2% HNO ₃	2 µg/L: Fe	100 mL	N8145060
NexION AFT Multi-Element Solution			
2% HNO ₃	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, Tl, U, V, Zn	100 mL	N8145061

USP Elemental Impurities Kit

Matrix	Contents	Vol.	Part No.
	Includes: 1 bottle each		N9303957
	Solution 1		
5% HNO ₃	2500 mg/kg: Cu, Mn 250 mg/kg: Cr, Mo, Ni, V 15 mg/kg: As, Hg 10 mg/kg: Pb 5 mg/kg: Cd	125 mL	
	Solution 2		
15% HCl	100 mg/kg: Ir, Os, Pd, Pt, Rh, Ru	125 mL	

Part Number Index

Part No.	Page	Part No.	Page	Part No.	Page	Part No.	Page	Part No.	Page
02900540	4	N9300121	2	N9300184	2, 9, 10	N9303740	2	N9303803	2
B0190634	4	N9300122	2	N9300200	3, 10	N9303741	2	N9303804	2
B0190635	4	N9300123	2	N9300201	3, 10	N9303742	2	N9303805	2
B0507226	4	N9300124	2	N9300202	3, 10	N9303743	2	N9303806	2
B3141064	4	N9300125	2	N9300203	3, 10	N9303744	2	N9303807	2
N0582152	5	N9300126	2, 9, 10	N9300204	3, 10	N9303745	2	N9303808	2
N0681470	5	N9300127	2	N9300205	4, 10	N9303746	2	N9303809	2
N0691579	3	N9300128	2	N9300207	2, 4, 10	N9303747	2	N9303810	2
N0691580	5	N9300129	2, 9	N9300208	4, 10	N9303748	2	N9303811	2
N0691581	5	N9300130	2	N9300211	5	N9303749	2	N9303812	2
N8120522	7	N9300131	2, 9	N9300212	5	N9303750	2	N9303813	3
N8120541	7	N9300132	2, 9	N9300213	5	N9303751	2	N9303814	6
N8122014	7	N9300133	2, 8	N9300214	5	N9303752	2	N9303816	7, 8
N8122017	8	N9300133	2	N9300215	5	N9303753	2	N9303818	7
N8122038	6, 7	N9300134	2	N9300216	5	N9303754	2	N9303819	8
N8125010	7	N9300135	2	N9300217	5	N9303755	2	N9303821	4, 8
N8125030	7	N9300136	2, 9	N9300218	3, 9	N9303756	2	N9303822	4, 8
N8125031	7	N9300137	2	N9300219	3, 9	N9303757	2	N9303823	4, 8
N8125032	7	N9300138	2	N9300220	3, 9	N9303758	2	N9303824	4, 8
N8125033	6, 7	N9300139	2	N9300221	3, 9	N9303759	2	N9303825	7
N8125034	7	N9300140	2	N9300223	10	N9303760	2	N9303826	7
N8125035	7	N9300141	2, 9, 10	N9300224	3	N9303761	2	N9303827	8
N8125037	7, 11	N9300142	2	N9300225	4	N9303762	2	N9303828	5
N8125038	7, 11	N9300143	2	N9300226	4, 9, 10	N9303763	2	N9303829	8
N8125039	11	N9300144	2	N9300227	4, 9, 10	N9303765	2	N9303830	5
N8125040	8	N9300145	2	N9300228	4, 9	N9303766	2	N9303831	5
N8125041	8	N9300146	2	N9300229	4, 9	N9303767	2	N9303832	6
N8140503	11	N9300147	2	N9300230	5	N9303768	2	N9303833	6
N8140504	11	N9300148	2, 9	N9300231	6	N9303769	2	N9303834	6
N8145050	6, 11	N9300149	2	N9300232	5, 6	N9303770	2	N9303835	9
N8145051	8, 11	N9300150	2	N9300233	5, 6	N9303771	2	N9303836	9
N8145052	11	N9300151	2	N9300234	5, 6	N9303772	2	N9303837	3
N8145053	11	N9300152	2, 9, 10	N9300235	5, 6, 9	N9303773	2	N9303838	2
N8145054	11	N9300153	2, 9	N9300236	6	N9303774	2	N9303839	6, 8
N8145055	11	N9300154	2	N9300237	3, 6	N9303775	2	N9303840	6, 8
N8145056	11	N9300155	2	N9300238	3, 6	N9303776	2	N9303841	6
N8145057	11	N9300156	2	N9300239	6	N9303777	2	N9303842	2
N8145058	11	N9300157	2	N9300241	10	N9303778	2	N9303843	8
N8145059	11	N9300158	2	N9300244	4	N9303779	2	N9303844	2
N8145060	11	N9300160	2	N9300253	5, 8	N9303780	2	N9303955	6
N8145061	11	N9300161	2, 9	N9300260	4	N9303781	2	N9303957	11
N9300100	2, 9	N9300162	2, 9	N9300280	3	N9303782	2	N9306225	5
N9300101	2, 9	N9300163	2	N9300281	3, 7	N9303783	2	N9307103	8
N9300102	2	N9300165	2, 9	N9301720	6	N9303784	2	N9307104	9
N9300103	2	N9300166	2	N9301721	7, 8	N9303785	2	N9307105	9
N9300104	2	N9300167	2, 9	N9302946	5	N9303786	2	N9307106	9
N9300105	2	N9300168	2	N9303445	4	N9303787	2	N9307107	9
N9300106	2	N9300169	2	N9303725	2	N9303788	2	N9307108	10
N9300107	2	N9300170	2	N9303726	2	N9303789	2	N9307109	10
N9300108	2, 9, 10	N9300171	2	N9303727	2	N9303790	2	N9307110	10
N9300109	2	N9300172	2	N9303728	2	N9303791	2	N9307111	8
N9300110	2	N9300173	2	N9303729	2	N9303792	2	N9307112	8
N9300111	2	N9300174	2	N9303730	2	N9303793	2	N9307113	6
N9300112	2, 9	N9300175	2	N9303731	2	N9303794	2	N9307114	6
N9300113	2	N9300176	2	N9303732	3	N9303795	2	N9307115	6
N9300114	2, 9	N9300177	2	N9303733	2	N9303796	2	N9307116	6
N9300115	2	N9300178	2	N9303734	2	N9303797	4	N9307738	8
N9300116	2	N9300179	2, 9, 10	N9303735	2	N9303798	2	N9307805	8
N9300117	2	N9300180	2	N9303736	2	N9303799	2	N9307806	8
N9300118	2	N9300181	2	N9303737	2	N9303800	2	N9307807	8
N9300119	2	N9300182	2	N9303738	2	N9303801	2	N9307808	8
N9300120	2	N9300183	2	N9303739	2	N9303802	2		

All of our Standards are guaranteed compatible with PerkinElmer and other manufacturers' instruments.

Order today at: www.perkinelmer.com/supplies

Every day, you can count on PerkinElmer to provide you with solutions that deliver reliable performance, control operating costs and maximize operational time. Our complete portfolio of consumables, parts, supplies, training and service helps you meet both routine and demanding measurement challenges. We invest heavily in testing and validating our products to ensure you receive guaranteed compatibility and performance – on-time, for every instrument in your laboratory.

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright © 2011, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

009515B_01