

HUMAN HEALTH

ENVIRONMENTAL HEALTH



FINALLY

REPRODUCIBILITY
COMES TO SPE



Supra-Clean® and Supra-Poly®
Solid Phase Extraction (SPE) Solutions


PerkinElmer®
For the Better

SUPERIOR REPRODUCIBILITY SUPERIOR RESULTS

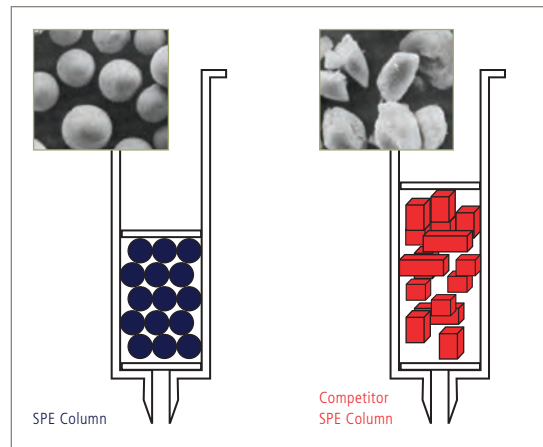


Imagine how much faster, easier and more efficient your sample preparation would be if your Solid Phase Extraction (SPE) solutions were precisely manufactured for guaranteed reproducibility. Not just column to column. Not just batch to batch. But lot to lot—no matter how many days, weeks or years between production times.

That's exactly the kind of confidence and convenience you can expect with the innovative line of Supra-Clean™ silica-based and Supra-Poly™ polymer-based SPE cartridges and columns from PerkinElmer.

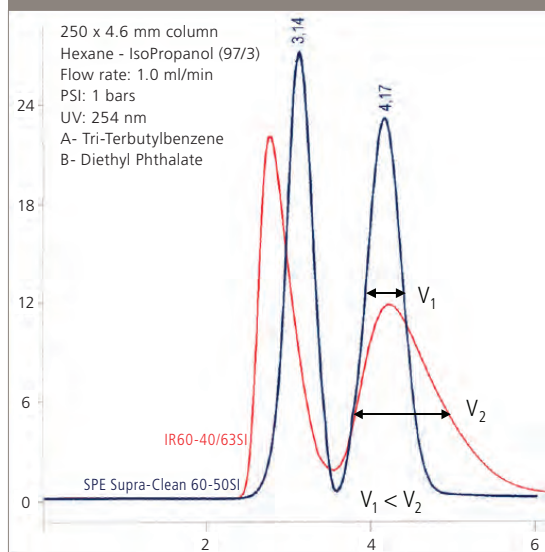
Our Precise-Bed Technology™ provides precision and manufacturing tolerances of $\pm 1\%$, ensuring guaranteed repeatability and optimized recovery reproducibility. The distribution and size of particles is controlled through stringent manufacturing processes, and each finished product is delivered with an individual quality certificate.

Experience the difference for yourself. Or—to be more exact—experience the uniformity.



Spherical media and consistent particle distribution enable smaller elution volumes and better, more reproducible extraction purification, concentration and recovery.

Performance Comparison of PerkinElmer Spherical SPE Media vs. Leading Irregular SPE Silica.



PerkinElmer's smaller, more homogeneous, spherical media deliver sharper, narrower peaks for faster, more accurate sample analysis.

Consistently. Excellent. Performance.

The unsurpassed reproducibility of SPE solutions is the result of a unique bed-packing technique and spherical media. Columns are evenly and consistently filled with particles sized for optimum distribution, delivering a level of performance unattainable with the irregularly shaped silica used in competitive solutions.

PerkinElmer's homogeneous filling and mono-dispersed spherical media not only provide unparalleled reproducibility, they also deliver superior performance. Sample diffusion through the media is better than in other systems, improving flow efficiency and avoiding channeling and clogging due to fines. This enables smaller sample sizes to be used without compromising recovery levels, dramatically reducing solvent volumes, analytical costs and processing times.

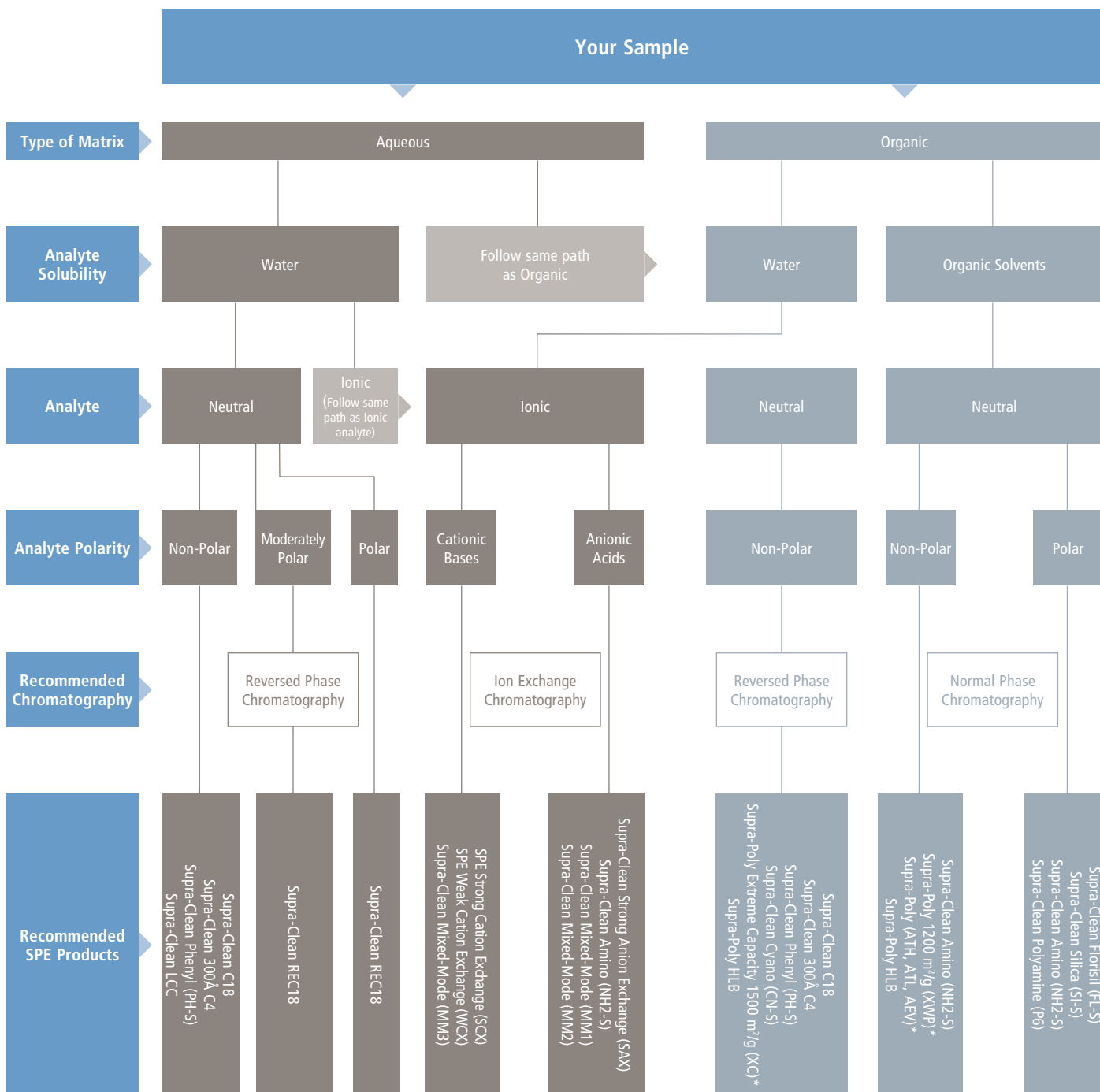
With PerkinElmer SPE solutions, you can work faster and more cost-effectively, all while experiencing the complete recovery, reproducibility and reliability you need for consistently excellent results.

Answering Your Needs. Empowering Your Lab.

Ideal for a broad array of analytes and matrices, SPE solutions are available in a variety of formats including Large Reservoir Capacity (LRC) columns, Polypropylene (PP) columns and cartridges, 96 well plate format, and glass columns. Each technology is offered with a wide selection of polymer and silica sorbents. Varied column capacities and bed weights allow you to perform scalable analyses depending upon your required detection limits.

Choose the Best SPE Solution For Your Method

To find the ideal SPE column for your particular application, simply follow the flow diagram below, selecting your path based on the options at each junction.



*AEV = Advanced Environmental
ATH = Hydrophobic/Hydrophilic
ATL = Lipophilic

HLB = Hydrophilic Lipophilic Balanced
XC = Extreme Capacity
XWP = Extra Wide Particle

A WIDE ARRAY OF COLUMNS FOR A BROAD RANGE OF APPLICATIONS

SUPRA-CLEAN SPHERICAL SILICA

Phase	Mechanism	Interaction mode	Compounds	Matrix
C18-S	Hydrophobic	Reversed Phase	Polar to Non-Polar compounds	Biological fluids, aqueous samples
High Recovery REC18	Hydrophobic	Reversed Phase	Non-polar and mid-polar compounds including 100% water solvents	Biological fluids, aqueous samples toxins in food
Phenyl (PH-S)	Hydrophobic	Reversed Phase	Non-polar to mid-polar aromatic compounds	Biological fluids
Silica (SI-S)	Hydrophilic	Normal Phase	Polar compounds	Non-polar organics, oils, lipids
Amino (NH2-S)	Hydrophilic	Normal Phase	Polar to Mid-Polar aromatic compounds	Biological fluids, aqueous samples, buffered organics
Strong Cation Exchange (SCX)	Ion Exchange	Ion Exchange	Basic compounds	Biological fluids, aqueous samples, buffered organics
Weak cation Exchange (WCX)	Ion Exchange	Ion Exchange	Strong basic compounds	Biological fluids, aqueous samples
Strong Anion Exchange (SAX)	Ion Exchange	Ion Exchange	Acidic compounds	Biological fluids, aqueous samples
Cyano (CN-S)	Hydrophilic	Normal Phase	Polar to Mid-Polar compounds	Non-polar organics, oils, lipids
Florisil (FL-S)	Hydrophilic	Normal Phase	Polar compounds	Ideal for polar compounds in non-polar matrix
Florisil Pesticide (FL-S)	Hydrophilic	Normal Phase	Polar compounds	Ideal for polar compounds in non-polar matrix
Polyamine (P6)	Hydrophilic	Reversed Phase	Carboxylic acids, phenolics and nitroaromatics	Aqueous and mid-polar matrices
300 A (C4)	Hydrophobic	Reversed Phase	Non-polar to mid-polar compounds	Biological Samples
LCC	Hydrophobic	Reversed Phase	Non-polar to mid-polar compounds	Biological fluids, aqueous samples
Mixed mode (MM1)	Ion Exchange/ Hydrophobic	Reversed Phase / SCX	Basic compounds	Biological samples
Mixed mode (MM2)	Ion Exchange/ Hydrophobic	Reversed Phase / WCX	Very basic compounds	Biological samples
Mixed mode (MM3)	Ion Exchange/ Hydrophobic	Reversed Phase / SAX	Acidic compounds	Biological samples

SUPRA-POLY SPHERICAL POLYMER

Extreme Capacity (XC)	Hydrophobic	Reversed Phase	Polar and non-polar	Aqueous or organic
Extreme Capacity Wide Pore (XWP)	Hydrophobic	Reversed Phase	Polar and non-polar	Biological and viscous samples
Hydrophilic (ATH)	Hydrophilic	Reversed Phase	Mid to non-polar compounds	Aqueous or organic
Lipophilic (ATL)	Lipophilic	Reversed Phase	Mid to non-polar compounds	Crude samples
Environmental (AEV)	Hydrophilic/ Hydrophobic	Reversed Phase	Mid to non-polar compounds	Aqueous or Organic
HLB	Hydrophilic/lipophilic balanced	Reversed Phase	Mid to non-polar compounds	Aqueous or organic

Typical Applications	pH Range	End-capping	Pore Size (Å)	Surface Area (m ² /g)	Particle Size (µm)	Comments
Drugs and drug metabolites in biological matrices, trace organic material in water, toxins in food	2-8	Yes	60	500	50	18% Carbon Load (CL)
Drugs and drug metabolites in biological matrices, trace organic material in water,	2-8	Yes	NA	NA	50	High capacity and better recovery especially for high aqueous conditions. 15% CL
Benzodiazepines in biological matrices, extraction of aromatic compounds	2-8	No	60	500	50	9% CL
Aldehydes, amines, pesticides, herbicides, carotenoids, fat soluble vitamins, aflatoxins, fatty acids, and phospholipids	2-8	No	60	500	50	Bare Silica
Basic compounds, polar amine compounds, carbohydrates	2-8	No	60	500	50	5% CL
Cations, antibiotics, drugs, amino acids, catecholamines, herbicides, nucleic acid bases, nucleosides, and surfactants	2-8	No	60	450	60	Strong Acid - Sulfonic acid; Exchange capacity 0.70 meq/g
Cations, amines, antibiotics, drugs, amino acids, catecholamines, nucleic acid bases, nucleosides, and surfactants	2-8	No	60	450	60	Weak Acid - Carboxylic acid; Exchange capacity 0.22 meq/g
Acidic food pigments, organic acids, phenol compounds, nucleic acids, nucleotides, surfactants	2-8	No	60	450	60	Strong Base - quaternary amine; Exchange capacity 0.30 meq/g
Polar compounds in hexane and oil	2-8	Yes	60	500	50	8% CL; Mid-range polarity between silica and C18
Pesticides, Polychlorinated Biphenyls (PCB)	2-8	No	NA	NA	200	Standard grade. Alternative to silica for viscous matrices due to large particle size. Granular shape.
Pesticides	2-8	No	NA	NA	200	High purity pesticide grade. Alternative to silica for viscous matrices. Granular shape.
Aromatic and natural products; Flavones, Chalkones, Anthraquinones	2-8	No	NA	NA	100	Nylon 6
Hydrophobic peptides and polypeptides	2-8	No	300	-	-	Large pore size for isolation of large biomolecules
Non-polar compounds in aqueous solution	2-8	Yes	60	500	50	10% CL; Lower carbon load than C18-S and REC18
Drugs and drug metabolites	2-8	No	60	450	60	Exchange capacity 0.09 meq/g
Drugs and drug metabolites	2-8	No	60	450	60	Exchange capacity 0.10 meq/g
Drugs and drug metabolites	2-8	No	60	450	60	Exchange capacity 0.14 meq/g
Drugs and drug metabolites biological fluids	0-14	No	NA	1500	70	High capacity polystyrene-divinylbenzene (PSDVB)
Drugs and drug metabolites biological fluids	0-14	No	Wide Pore	1200	90	High capacity PSDVB for large biomolecules and viscous matrices
Mid-polar and non-polar compounds in aqueous and organic solvents	1-13	No	70	800	75	Mixed hydrophilic/hydrophobic interactions
Lipids	0-14	No	70	800	100	PSDVB; Alternative to high flow silica for mid-polar to non-polar compounds (<3000D) in crude samples
Aqueous environmental compounds that are not retained on C18	1-12	No	70	800	75	Advanced environmental; Polystyrene-co-2-hydroxyethyl methacrylate (PSHEMA)
Mid-polar and non-polar compounds in aqueous and organic solvents	0-14	No	80	850	30 & 60	Hydrophilic-lipophilic-balanced reversed-phase sorbent for acids, bases and neutrals

PerkinElmer offers a comprehensive selection of SPE products that have been precisely manufactured for guaranteed reproducibility. Not just column to column. Not just batch to batch. But lot to lot — no matter how many days, weeks or years between production times. Supra-Clean and Supra-Poly SPE products are available in a range of media weights, volumes and formats, including easy to order selection kits and application packs.

SUPRA-CLEAN COLUMNS AND 96 WELL PLATES

- Pure spherical silica
- Pore size 60 - 120 Å
- 20 chemistries with pH range 2-8

Supra-Clean C18

Columns

Media Weight	Volume	Part No.	Quantity
50 mg	1 ml	N9306476	50
100 mg	1 ml	N9306478	100
100 mg	3 ml	N9306523	50
200 mg	3 ml	N9306462	50
500 mg	3 ml	N9306438	50
500 mg	3 ml*	N9306642	50
200 mg	6 ml	N9306634	30
500 mg	6 ml	N9306448	30
500 mg	6 ml*	N9306640	30
1 g	6 ml	N9306422	30
2 g	6 ml	N9306430	30
2 g	15 ml	N9306479	20
2 g	25 ml	N9306475	20

* Not end-capped

96 Well Plates

Media Weight	Volume	Part No.	Quantity
25 mg	2 ml	N9306566	1
50 mg	2 ml	N9306567	1
100 mg	2 ml	N9306568	1

Supra-Clean C18 Cartridges

Media Weight	Volume	Part No.	Quantity
390 mg	--	N9306587	50
910 mg	--	N9306588	50
1690 mg	--	N9306589	50

Supra-Clean REC18

Columns

Media Weight	Volume	Part No.	Quantity
50 mg	1 ml	N9306519	50
100 mg	1 ml	N9306520	100
100 mg	3 ml	N9306455	50
200 mg	3 ml	N9306521	50
500 mg	3 ml	N9306522	50
200 mg	6 ml	N9306633	30
500 mg	6 ml	N9306457	30
1 g	6 ml	N9306491	30

96 Well Plates

Media Weight	Volume	Part No.	Quantity
25 mg	2 ml	N9306563	1
50 mg	2 ml	N9306564	1
100 mg	2 ml	N9306565	1

Supra-Clean Strong Anion Exchange (SAX)

Columns

Media Weight	Volume	Part No.	Quantity
50 mg	1 ml	N9306553	50
100 mg	1 ml	N9306471	100
100 mg	3 ml	N9306554	50
200 mg	3 ml	N9306482	50
500 mg	3 ml	N9306555	50
500 mg	6 ml	N9306556	30

96 Well Plates

Media Weight	Volume	Part No.	Quantity
25 mg	2 ml	N9306581	1
50 mg	2 ml	N9306582	1
100 mg	2 ml	N9306583	1

Supra-Clean 300Å C4

Columns

Media Weight	Volume	Part No.	Quantity
50 mg	1 ml	N9306590	50
100 mg	1 ml	N9306591	100
100 mg	3 ml	N9306592	50
200 mg	3 ml	N9306593	50

Supra-Clean Strong Cation Exchange (SCX)

Columns

Media Weight	Volume	Part No.	Quantity
50 mg	1 ml	N9306536	50
100 mg	1 ml	N9306432	100
100 mg	3 ml	N9306537	50
200 mg	3 ml	N9306538	50
500 mg	3 ml	N9306539	50
500 mg	6 ml	N9306540	30

96 Well Plates

Media Weight	Volume	Part No.	Quantity
25 mg	2 ml	N9306575	1
50 mg	2 ml	N9306576	1
100 mg	2 ml	N9306577	1

Supra-Clean Weak Cation Exchange (WCX)

Columns

Media Weight	Volume	Part No.	Quantity
50 mg	1 ml	N9306544	50
100 mg	1 ml	N9306545	100
100 mg	3 ml	N9306546	50
200 mg	3 ml	N9306547	50
500 mg	3 ml	N9306420	50
500 mg	6 ml	N9306407	30

Supra-Clean Mixed-Mode (MM1)

Columns

Media Weight	Volume	Part No.	Quantity
50 mg	1 ml	N9306541	50
100 mg	1 ml	N9306542	100
100 mg	3 ml	N9306419	50
200 mg	3 ml	N9306543	50
500 mg	3 ml	N9306481	50
500 mg	6 ml	N9306416	30
200 mg	15 ml	N9306713	20

96 Well Plates

Media Weight	Volume	Part No.	Quantity
25 mg	2 ml	N9306578	1
50 mg	2 ml	N9306579	1
100 mg	2 ml	N9306580	1

Supra-Clean Mixed-Mode (MM2)

Columns

Media Weight	Volume	Part No.	Quantity
50 mg	1 ml	N9306548	50
100 mg	1 ml	N9306549	100
100 mg	3 ml	N9306550	50
200 mg	3 ml	N9306551	50
500 mg	3 ml	N9306411	50
500 mg	6 ml	N9306552	30

Supra-Clean Mixed-Mode (MM3)

Columns

Media Weight	Volume	Part No.	Quantity
500 mg	6 ml	N9306649	30

Supra-Clean Florisil* (FL-S)

Columns

Media Weight	Volume	Part No.	Quantity
200 mg	3 ml	N9306511	50
500 mg	3 ml	N9306512	50
500 mg	6 ml	N9306494	30
1 g	6 ml	N9306413	30
2 g	6 ml	N9306513	20
2 g	15 ml	N9306514	20
2 g	25 ml	N9306515	20

* Granular shape.

Supra-Clean Florisil* (FL-S) Pesticide Grade

Columns

Media Weight	Volume	Part No.	Quantity
200 mg	3 ml	N9306516	50
500 mg	3 ml	N9306400	50
500 mg	6 ml	N9306517	30
1 g	6 ml	N9306436	30
2 g	6 ml	N9306470	30
2 g	15 ml	N9306443	20
2 g	25 ml	N9306447	20

* Granular shape.

Supra-Clean Silica (SI-S)

Columns

Media Weight	Volume	Part No.	Quantity
100 mg	3 ml	N9306532	50
200 mg	3 ml	N9306444	50
500 mg	3 ml	N9306402	50
500 mg	6 ml	N9306466	30
1 g	6 ml	N9306404	30
2 g	6 ml	N9306533	20
2 g	15 ml	N9306534	20
2 g	25 ml	N9306535	20

Cartridges

Media Weight	Volume	Part No.	Quantity
300 mg	--	N9306584	50
700 mg	--	N9306585	50
1300 mg	--	N9306586	50

Supra-Clean Cyano (CN-S)

Columns

Media Weight	Volume	Part No.	Quantity
500 mg	3 ml	N9306645	50
500 mg	6 ml	N9306644	30

Supra-Clean Amino (NH2-S)

Columns

Media Weight	Volume	Part No.	Quantity
50 mg	1 ml	N9306528	50
100 mg	1 ml	N9306410	100
100 mg	3 ml	N9306529	50
500 mg	3 ml	N9306414	50
200 mg	6 ml	N9306530	50
500 mg	6 ml	N9306531	30

96 Well Plates

Media Weight	Volume	Part No.	Quantity
25 mg	2 ml	N9306572	1
50 mg	2 ml	N9306573	1
100 mg	2 ml	N9306574	1

Supra-Clean Polyamine (P6)

Columns

Media Weight	Volume	Part No.	Quantity
500 mg	3 ml	N9306518	50
500 mg	6 ml	N9306434	30

Supra-Clean Phenyl (PH-S)

Columns

Media Weight	Volume	Part No.	Quantity
50 mg	1 ml	N9306401	50
100 mg	1 ml	N9306524	100
100 mg	3 ml	N9306525	50
200 mg	3 ml	N9306490	50
500 mg	3 ml	N9306421	50
500 mg	6 ml	N9306526	30
1 g	6 ml	N9306527	30

96 Well Plates

Media Weight	Volume	Part No.	Quantity
25 mg	2 ml	N9306569	1
50 mg	2 ml	N9306570	1
100 mg	2 ml	N9306571	1

Supra-Clean LCC

Columns

Media Weight	Volume	Part No.	Quantity
500 mg	3 ml	N9306643	50
500 mg	6 ml	N9306641	30

SUPRA-POLY HLB COLUMNS AND 96 WELL PLATES

- Contains macro-porous polymers with ultra-pure, pharmaceutical grade spherical particles
- Shorter analysis times, greater load capacity and reduced solvent usage
- Ideal for high throughput assays

Supra-Poly HLB

30 µm Columns

Media Weight	Volume	Part No.	Quantity
30 mg	1 ml	N9306650	50
50 mg	1 ml	N9306655	50
60 mg	1 ml	N9306656	50
100 mg	1 ml	N9306657	50
30 mg	3 ml	N9306651	50
60 mg	3 ml	N9306658	50
100 mg	3 ml	N9306659	50
200 mg	3 ml	N9306660	50
500 mg	3 ml	N9306661	30
100 mg	6 ml*	N9306672	30
150 mg	6 ml	N9306662	30
200 mg	6 ml	N9306663	30
200 mg	6 ml*	N9306673	30
500 mg	6 ml*	N9306674	30
500 mg	6 ml	N9306664	30
500 mg	15 ml	N9306665	20
1 g	15 ml	N9306666	20
30 mg	15 ml**	N9306668	50
60 mg	15 ml**	N9306669	50
100 mg	15 ml**	N9306670	50
200 mg	15 ml**	N9306671	50
1 g	25 ml	N9306667	20

60 µm Columns

Media Weight	Volume	Part No.	Quantity
30 mg	1 ml	N9306652	50
50 mg	1 ml	N9306675	50
60 mg	1 ml	N9306676	50
100 mg	1 ml	N9306677	50
30 mg	3 ml	N9306653	50
60 mg	3 ml	N9306678	50
100 mg	3 ml	N9306679	50
200 mg	3 ml	N9306680	50
500 mg	3 ml	N9306681	30
100 mg	6 ml*	N9306692	30
150 mg	6 ml	N9306682	30
200 mg	6 ml	N9306683	30

Media Weight	Volume	Part No.	Quantity
200 mg	6 ml*	N9306693	30
500 mg	6 ml	N9306684	30
500 mg	15 ml	N9306685	20
500 mg	6 ml*	N9306694	30
1 g	15 ml	N9306686	20
30 mg	15 ml**	N9306688	50
60 mg	15 ml**	N9306689	50
100 mg	15 ml**	N9306690	50
200 mg	15 ml**	N9306691	50
1 g	25 ml	N9306687	20

* Glass columns ** LRC columns

30 µm 96 Well Plates

Media Weight	Volume	Part No.	Quantity
30 mg	2 ml	N9306698	1
50 mg	2 ml	N9306699	1
60 mg	2 ml	N9306700	1

60 µm 96 Well Plates

Media Weight	Volume	Part No.	Quantity
30 mg	2 ml	N9306695	1
50 mg	2 ml	N9306696	1
60 mg	2 ml	N9306697	1

SUPRA-POLY COLUMNS AND 96 WELL PLATES

- Contains macro-porous polymers with ultra-pure, pharmaceutical grade spherical particles
- Shorter analysis times, greater load capacity and reduced solvent usage
- Ideal for high throughput assays

Supra-Poly Extreme Capacity 1500 m²/g (XC)

Columns

Media Weight	Volume	Part No.	Quantity
30 mg	1 ml	N9306441	50
50 mg	1 ml	N9306500	50
60 mg	1 ml	N9306501	50
100 mg	1 ml	N9306403	50
60 mg	3 ml	N9306502	50
100 mg	3 ml	N9306440	50
200 mg	3 ml	N9306428	50
200 mg	6 ml	N9306635	30
500 mg	6 ml	N9306405	30
1 g	15 ml	N9306503	20

96 Well Plates

Media Weight	Volume	Part No.	Quantity
30 mg	2 ml	N9306557	1
50 mg	2 ml	N9306558	1
60 mg	2 ml	N9306559	1

Supra-Poly Extra Wide Particle 1200 m²/g (XWP)

Columns

Media Weight	Volume	Part No.	Quantity
30 mg	1 ml	N9306504	50
50 mg	1 ml	N9306427	50
60 mg	1 ml	N9306505	50
100 mg	1 ml	N9306506	50
60 mg	3 ml	N9306507	50
100 mg	3 ml	N9306508	50
200 mg	3 ml	N9306509	50
500 mg	6 ml	N9306418	30
1 g	15 ml	N9306510	20

96 Well Plates

Media Weight	Volume	Part No.	Quantity
30 mg	2 ml	N9306560	1
50 mg	2 ml	N9306561	1
60 mg	2 ml	N9306562	1

Supra-Poly Environmental (AEV)

Columns

Media Weight	Volume	Part No.	Quantity
100 mg	3 ml	N9306648	50

Supra-Poly Hydrophilic (ATH)

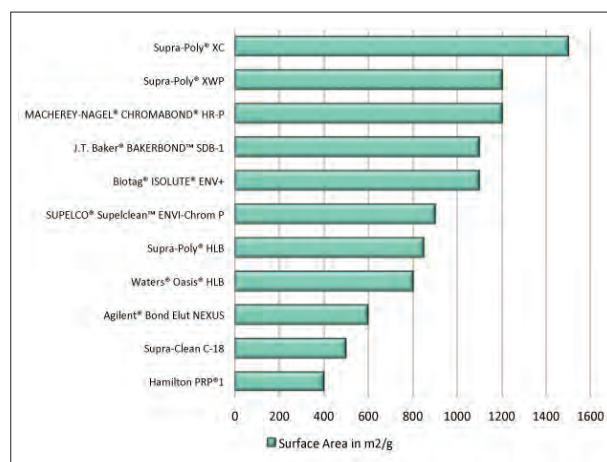
Columns

Media Weight	Volume	Part No.	Quantity
100 mg	3 ml	N9306646	50
200 mg	3 ml	N9306638	50
200 mg	6 ml	N9306636	30

Supra-Poly Lipophilic (ATL)

Columns

Media Weight	Volume	Part No.	Quantity
100 mg	3 ml	N9306647	50
200 mg	3 ml	N9306639	50
200 mg	6 ml	N9306637	30



Compared to other industry leading SPE materials, our Extreme Capacity (XC) and Extra Wide Particles (XWP) lead market in surface area. This allows for higher capacities at lower bed weights.

SPE SELECTION KITS: Enables quick column selection for development of reproducible and repeatable SPE methods.

Description	Media Weight	Volume	Part No.	Quantity
Pre-Concentration of Hydrophobic Compounds from Aqueous Matrix	200 mg	6 ml	N9306594	50
	200 mg	3 ml	N9306595	50
Extraction of Hydrophobic Compounds from Aqueous Matrix	500 mg	6 ml	N9306596	50
	500 mg	3 ml	N9306597	50
Pre-Concentration of Hydrophilic Compounds	500 mg	6 ml	N9306598	30
	500 mg	3 ml	N9306599	30
Removal of Polar Compounds from Aqueous and Organic Matrix	500 mg	6 ml	N9306600	30
	500 mg	3 ml	N9306601	30
Extraction of Acidic Basic and Neutral Compounds from Aqueous or Organic Matrix	100 mg	3 ml	N9306602	50
Extraction of Carboxylic Acids and Strong Bases from Aqueous Matrix	500 mg	6 ml	N9306603	40
Extraction of Weak Bases from Aqueous Matrix	500 mg	6 ml	N9306604	30

SPE APPLICATION PACKS: Ideal for extraction of known entities from a variety of matrices, our packs are expertly tailored to meet your application needs and are designed to support major EPA methods and applications.

Description	Media Weight	Volume	Part No.	Quantity
Extraction of Basic Drugs from Biological Fluids	200 mg	3 ml	N9306605	50
Extraction of Bisphenol A from Aqueous Matrix	1 g	6 ml	N9306613	30
Extraction of Oil and Grease from Aqueous Matrix-EPA 1664	500 mg	3 ml	N9306612	50
	1 g	6 ml	N9306611	30
Extraction of PAH from Soil and Oil	1.5 g	6 ml	N9306609	30
Extraction of PAH from Soil and Oil (Glass Straight Column)	1.5 g	6 ml	N9306610	30
Extraction of PAH from Water Containing Humic Acids	1.5 g	6 ml	N9306608	30
Extraction of PAH from Water or Soil	4 g	6 ml	N9306606	30
Extraction of PAH from Water or Soil (Glass Straight Column)	4 g	6 ml	N9306607	30
Extraction of PCB from Oil	1 g	6 ml	N9306617	30
	1 g	3 ml	N9306616	50
Extraction of Pesticides and Herbicides from Aqueous Matrix	500 mg	3 ml	N9306614	50
Extraction of Steroids from Biological Fluids	500 mg	6 ml	N9306615	30
Extraction of SVOC from Water-EPA 525	1 g	6 ml	N9306618	30

SPE VACUUM PUMPS, MANIFOLDS AND ACCESSORIES: SPE Vacuum Manifolds accommodate either 12 or 24 cartridges; 1, 3, 6, 15, and 25 mL columns can be used. Manifolds are equipped with a vacuum port to connect a standard laboratory vacuum pump. Vacuum pulls the sample through the stationary phase, metered by the stopcocks, to control the speed of the extraction process and the sample flow. Waste and wash solvents are discarded and analytes are collected in sample tubes below the manifold completing the extraction.

Description	12 Position Part No.	Quantity	24 Position Part No.	Quantity	
Vacuum Manifold	N9306619	1 each	N9306626	1 each	
Replacement Chamber (Glass)	N9306620	1 each	N9306627	1 each	
Cover Gasket - 12 Position	N9306621	1 each	N9306628	1 each	
Stopcocks - 12 Position	N9306624	12 pack	N9306631	24 pack	
Needles - Polypropylene	N9306622	12 pack	N9306629	24 pack	
Needles - Stainless Steel	N9306623	12 pack	N9306630	24 pack	
Drying Attachment	N9306625	1 each	N9306632	1 each	
Description	20 L/min 115V	60 L/min 115V	17 L/min 230V	58 L/min 230V	Quantity
Vacuum Pumps	N9308035	N9308063	N9308331	N9308332	1 each

THE PERFECT PARTNERSHIP FROM START TO FINISH



At PerkinElmer, we understand that sample preparation is one of the most critical steps in the analytical process. Often accounting for 60% or more of your timetable, it has a fundamental impact on a wide range of operational parameters. Any errors within this process undermine the quality of your data at all subsequent stages of your analysis.

Solid Phase Extraction helps avoid potential errors in sample preparation, reducing re-runs and dramatically increasing productivity. As one of the most cost-effective and flexible tools within the laboratory environment, SPE also provides efficient sample concentration and purification prior to many of today's most popular analytical techniques, including HPLC, LC/MS, GC and GC/MS.

SPE columns are part of PerkinElmer's complete offering of analytical solutions—an industry-leading portfolio that encompasses instruments, accessories, consumables, supplies, training and service. It's a breadth of capabilities that enables us to offer the ease and convenience of having a single supplier for all your needs at every stage of your workflow. So you can benefit from greater responsiveness, superior reliability, and dramatic cost savings.

Couple SPE With Leading Instrumentation For Ultimate Laboratory Efficiency

Clarus® 680/580/480 GC

Clarus SQ 8 GC/MS

Flexar™ SQ 300 MS

Flexar LC and UHPLC

The right accessories, consumables, methods and application support are as integral to the success of your laboratory as your instrumentation. That's why we invest heavily in testing and validating our complete portfolio of solutions to ensure you receive accurate, repeatable results on time, every time.

Our OneSource Laboratory Services division even takes it a step further. With more than 1,500 trained and certified field service engineers and service personnel around the world, OneSource offers the most comprehensive portfolio of professional laboratory services in the industry, including complete care programs for virtually every technology and manufacturer.

So turn to PerkinElmer. For the experience. For the confidence. For the better.

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 ©2011-2012, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

009795A_01

Printed in USA