



Optimal Results.
Compliments of Agilent.

Thank you for trying Agilent SampliQ SPE products

Start getting more from your sample preparation, start using SampliQ SPE products today. Get more information about method development and guidelines for selecting the right SampliQ product for your analysis – visit www.agilent.com/chem/SampliQ

These high-quality and high-purity sorbents for your SPE method are available in a wide selection of phases for use in any analyses when consistent results are crucial. SampliQ SPE products improve and simplify separations, increase column lifetime and improve detection limits. To ensure product integrity, all SampliQ products are packaged with triple lamination and include a certificate of performance.



Agilent’s SampliQ SPE is the first step in getting optimal results. Count on our full line of high-quality columns and supplies to optimize your full analysis.

- Agilent J&W columns provide many options to meet your specific needs. For fast GC analyses, we recommend Agilent J&W High-Efficiency columns. For your most sensitive trace analysis, there is Agilent J&W Ultra Inert columns.
- For HPLC applications, Agilent Poroshell 120 and Agilent ZORBAX Eclipse Plus columns provide the most innovative technology for reproducible results.

Don’t wait for optimal results —
order your SampliQ SPE products today!

Contact your local Agilent sales office or Agilent Authorized Distributor.

For technical or product questions, contact Agilent Technical Support:
Call **1-800-227-9770 (U.S. and Canada)**. For outside U.S./Canada, please go to www.agilent.com/chem/contactus

SampliQ from Agilent

HIGH PERFORMANCE SOLID PHASE EXTRACTION PRODUCTS



The SampliQ line has polymers, silica and non-silica options to meet your specific SPE needs. Use the charts below to find the SampliQ product best suited for your sample preparation.

POLYMERS

	OPT	DVB	SCX	SAX	PS-DVB
1ml 30mg, 100/pk	5982-3013	5982-3113	5982-3213	5982-3313	
3ml 60mg, 50/pk	5982-3036	5982-3136	5982-3236	5982-3336	
6ml 150mg, 30/pk	5982-3067	5982-3167	5982-3267	5982-3367	
6ml 500mg, 30/pk					5982-3165
6ml 1000mg, 30/pk					5982-3160

SILICA

Reversed Phase (Non-Polar)					
	C18	C18EC	C8	C2	Phenyl
1ml 100mg, 100/pk	5982-1111	5982-1311	5982-1011	5982-1411	5982-1511
3ml 200mg, 50/pk	5982-1132	5982-1332	5982-1032	5982-1432	5982-1532
3ml 500mg, 50/pk	5982-1135	5982-1335	5982-1035	5982-1435	5982-1535
6ml 500mg, 30/pk	5982-1165	5982-1365	5982-1065	5982-1465	
6ml 1000mg, 30/pk	5982-1160	5982-1360		5982-1460	
Bulk 25g Bottle	5982-1182	5982-1382	5982-1082		5982-1582

Normal Phase (Polar)				
	Silica	Diol	Cyano	Amino
1ml 100mg, 100/pk	5982-2211		5982-1711	5982-1811
3ml 200mg, 50/pk	5982-2232			5982-1832
3ml 500mg, 50/pk	5982-2235	5982-1635		5982-1835
6ml 500mg, 30/pk	5982-2265		5982-1765	5982-1865
6ml 1000mg, 30/pk	5982-2260		5982-1760	5982-1860
Bulk 25g Bottle	5982-2282			5982-1882

Mixed Mode		
	C8/SCX	EVDX
1ml 100mg, 100/pk	5982-1911	
3ml 200mg, 50/pk	5982-1932	5982-2332
3ml 500mg, 50/pk	5982-1935	
6ml 500mg, 30/pk	5982-1965	5982-2364*
6ml 1000mg, 30/pk	5982-1960	
Bulk 25g Bottle		

Ion Exchange	
Si-SAX	Si-SCX
5982-2011	5982-2111
5982-2032	5982-2132
5982-2035	5982-2135
5982-2065	5982-2165
5982-2060	5982-2060
5982-2082	

*6ml/400mg

ADDITIONAL SORBENTS

	Alumina Acid	Alumina Basic	Alumina Neutral	Florisil	Carbon
1ml 100mg, 100/pk					5982-4411
3ml 200mg, 50/pk				5982-4332	
3ml 200mg, 50/pk	5982-4035	5982-4135	5982-4235	5982-4335	5982-4435*
6ml 500mg, 30/pk				5982-4365	5982-4465
6ml 1000mg, 30/pk	5982-4060	5982-4160	5982-4260	5982-4360	
Bulk 25g Bottle	5982-4082	5982-4182	5982-4282	5982-4382	5982-4482

*3ml/250mg

Agilent takes pride in manufacturing SPE products in the U.S. with the same exacting quality that is used to manufacture its industry-leading instruments and columns.





呈现最佳结果
来自安捷伦的问候

感谢您试用安捷伦的 SampliQ SPE 产品

要获得更好的样品制备结果，现在就开始使用 SampliQ SPE 产品吧。如需更多有关方法开发的知识和根据您的具体分析正确选择 SampliQ SPE 产品的方法，请访问 www.agilent.com/chem/SampliQ

这些用于 SPE 方法的高质量和高纯度的吸附剂选择范围广，可用于对结果的一致性要求非常严格的任何分析应用。SampliQ SPE 产品可以改进和简化分离过程，延长柱寿命，提高检测限。为了保证产品的完整性，所有 SampliQ SPE 产品使用三层叠加包装，并附有质量保证书。



安捷伦的 SampliQ SPE 是获得最佳结果的第一步。凭借我们全方位的高质量的色谱柱和消耗品可使您的整个分析过程得到优化

- Agilent J&W 色谱柱有许多选择，可满足您的特殊需要。我们建议您使用 Agilent J&W 高效色谱柱用于快速气相色谱分析。如果您要进行高灵敏痕量分析，可以使用 Agilent J&W 超高惰性色谱柱。
- 对于高效液相色谱应用，可以使用采用最新工艺制造的 Agilent Poroshell 120 和 ZORBAX Eclipse Plus 色谱柱，以获得重现性优异的结果。

即刻获得最佳结果——
现在购买三个 SampliQ SPE 产品，
免费获赠一个！
有效期限 12/31/2008

请与您当地的安捷伦分公司或安捷伦授权的代理商联系。
仅限在中国提供优惠活动。

如有技术和产品方面的问题，请拨打安捷伦技术支持电话 800-820-3278
或请访问：www.agilent.com/chem/cncontactus

SampliQ 安捷伦出品

高品质的固相萃取产品



SampliQ 产品系列有聚合物、硅胶和非硅胶固定相供选择以满足特殊 SPE 的需求，使用下列图表您可以找到最适合您样品制备需要的 SampliQ 产品。

聚合物

	OPT	DVB	SCX	SAX	PS-DVB
1ml 30mg, 100/包	5982-3013	5982-3113	5982-3213	5982-3313	
3ml 60mg, 50/包	5982-3036	5982-3136	5982-3236	5982-3336	
6ml 150mg, 30/包	5982-3067	5982-3167	5982-3267	5982-3367	
6ml 500mg, 30/包					5982-3165
6ml 1000mg, 30/包					5982-3160

硅胶

反相（非极性）					
	C18	C18EC	C8	C2	Phenyl
1ml 100mg, 100/包	5982-1111	5982-1311	5982-1011	5982-1411	5982-1511
3ml 200mg, 50/包	5982-1132	5982-1332	5982-1032	5982-1432	5982-1532
3ml 500mg, 50/包	5982-1135	5982-1335	5982-1035	5982-1435	5982-1535
6ml 500mg, 30/包	5982-1165	5982-1365	5982-1065	5982-1465	
6ml 1000mg, 30/包	5982-1160	5982-1360		5982-1460	
Bulk 25g Bottle	5982-1182	5982-1382	5982-1082		5982-1582

正相（极性）				
	Silica	Diol	Cyano	Amino
1ml 100mg, 100/包	5982-2211		5982-1711	5982-1811
3ml 200mg, 50/包	5982-2232			5982-1832
3ml 500mg, 50/包	5982-2235	5982-1635		5982-1835
6ml 500mg, 30/包	5982-2265		5982-1765	5982-1865
6ml 1000mg, 30/包	5982-2260		5982-1760	5982-1860
Bulk 25g Bottle	5982-2282			5982-1882

混合型		
	C8/SCX	EVDX
1ml 100mg, 100/包	5982-1911	
3ml 200mg, 50/包	5982-1932	5982-2332
3ml 500mg, 50/包	5982-1935	
6ml 500mg, 30/包	5982-1965	5982-2364*
6ml 1000mg, 30/包	5982-1960	
Bulk 25g Bottle		

离子交换	
Si-SAX	Si-SCX
5982-2011	5982-2111
5982-2032	5982-2132
5982-2035	5982-2135
5982-2065	5982-2165
5982-2060	5982-2060
5982-2082	

*6ml/400mg

其他吸附剂

	三氧化二铝 酸性	三氧化二铝 碱性	三氧化二铝 中性	Florisil	碳
1ml 100mg, 100/包					5982-4411
3ml 200mg, 50/包				5982-4332	
3ml 200mg, 50/包	5982-4035	5982-4135	5982-4235	5982-4335	5982-4435*
6ml 500mg, 30/包				5982-4365	5982-4465
6ml 1000mg, 30/包	5982-4060	5982-4160	5982-4260	5982-4360	
Bulk 25g Bottle	5982-4082	5982-4182	5982-4282	5982-4382	5982-4482

*3ml/250mg

安捷伦为在美国制造的 SPE 产品具有与领导行业潮流的安捷伦仪器和色谱柱一样的严格意义上的高质量而引以自豪。

