

TAKE CONTROL OF YOUR SAMPLES

Agilent Bond Elut Plexa Polymeric SPE

Complex biological samples, such as plasma, are highly variable. But with Agilent Bond Elut Plexa SPE, you can reduce variability and simplify your analysis.

Bond Elut Plexa SPE eliminates common matrix interferences with a unique combination of hydroxylated exterior, hydrophobic interior, and advanced polymerization. What's more, the lot-to-lot reproducibility of Bond Elut Plexa ensures optimal and consistent flow rates, recovery, and cleanliness, which improve performance and simplify workflows.

Choosing the right sorbent for your analytes

Analyte Type	Bond Elut Plexa Phase	Mechanism of Separation	Loading Solvent	Eluting Solvent
Acidic, basic, neutral Log P >1.5 pKa 3-6 (acidic load) pKa 6-10 (basic load)	Plexa	Polymeric Reversed-Phase (hydrophobic-hydrophilic)	Water or buffer	Intermediate P' e.g. MeOH, ACN
Ionic (ionizable), basic Log P > 0.8 pKa 6-10	Plexa PCX	Polymeric Cation Exchange (Strong)	Water or buffer (pH =pKa-2)	• Buffer (pH=pKa+2) • pH value where analyte is neutral • Buffer with high ionic strength
Ionic (ionizable), acidic Log P > 1.0 pKa < 5	Plexa PAX	Polymeric Anion Exchange (Strong)	Water or buffer (pH = pKa+2)	• Buffer (pH=pKa-2) • pH value where analyte is neutral • Buffer with high ionic strength

Confidently clean samples and elute analytes


Bond Elut Plexa Polymeric SPE Protocol

	ACIDS	NEUTRALS		BASES
Analyte	LogP > 1.0 pKa < 5	LogP > 1.5 pKa 3-6 pKa 6-10		LogP > 0.8 pKa 6-10
	PLEXA PAX	PLEXA <small>Acid Load Method</small>	PLEXA <small>Base Load Method</small>	PLEXA PCX
Sample Pre-treatment	2% NH ₄ OH	1% HCO ₂ H	2% NH ₄ OH	2% H ₃ PO ₄
Sorbent Condition	100% MeOH	100% MeOH		100% MeOH
Equilibration	100% H ₂ O	100% H ₂ O		100% H ₂ O
Load	Apply pre-treated sample			
Wash	100% H ₂ O	5% MeOH in H ₂ O		2% HCO ₂ H in H ₂ O
Elution 1/Wash 2	100% MeOH Neutrals	100% MeOH Neutrals		1:1 MeOH/ACN Acids, Neutrals
Elution 2	5% HCO ₂ H in MeOH Acids			5% NH ₃ in 1:1 MeOH/ACN Bases
Analysis	Prepare extracts for instrumental analysis			


Complete your pharmaceutical research workflow with these high-performing Agilent LC/MS systems



Agilent 6470 Triple Quadrupole LC/MS
Perform confident quantitation and streamline your analytical workflow with improved sensitivity, precision, and scan speed.



Agilent 6230 LC/MS TOF
Detect both targets and unknowns with greater accuracy, and use enhanced MassHunter Workstation software to mine and analyze data.



Agilent 6545 LC/MS Q-TOF
Obtain better answers, faster with resolving power and sensitivity on the order of five times higher than previous generations.

Bond Elut Plexa SPE Ordering Information

Bond Elut Plexa		
Description	Unit	Part No.
Straight Barrel Cartridges		
30 mg, 1 mL	100/pk	12109301
30 mg, 3 mL	50/pk	12109303
60 mg, 1 mL	100/pk	12109601
60 mg, 3 mL	50/pk	12109603
200 mg, 3 mL	50/pk	12109610
200 mg, 6 mL	30/pk	12109206
500 mg, 6 mL	30/pk	12258506
Bond Elut Jr		
200 mg	50/pk	12169610B
Mega Bond Elut Plexa		
500 mg, 12 mL	20/pk	327832
96-Well Plates		
10 mg, 1 mL	1/pk	A4969010
30 mg, 1 mL	1/pk	A4969030
10 mg, 2 mL	1/pk	A3969010
30 mg, 2 mL	1/pk	A3969030

Bond Elut Plexa PCX		
Description	Unit	Part No.
Straight Barrel Cartridges		
30 mg, 1 mL	100/pk	12108301
60 mg, 1 mL	100/pk	12108601
30 mg, 3 mL	50/pk	12108303
60 mg, 3 mL	50/pk	12108603
200 mg, 6 mL	30/pk	12108206
500 mg, 6 mL	30/pk	12258506
96-Well Plates		
10 mg, 1 mL	1/pk	A4968010
30 mg, 1 mL	1/pk	A4968030
10 mg, 2 mL	1/pk	A3968010
30 mg, 2 mL	1/pk	A3968030

Bond Elut Plexa PAX		
Description	Unit	Part No.
Straight Barrel Cartridges		
30 mg, 1 mL	100/pk	12107301
60 mg, 1 mL	100/pk	12107601
30 mg, 3 mL	50/pk	12107303
60 mg, 3 mL	50/pk	12107603
200 mg, 6 mL	30/pk	12107206
500 mg, 6 mL	30/pk	12257506
96-Well Plates		
10 mg, 1 mL	1/pk	A4967010
30 mg, 1 mL	1/pk	A4967030
10 mg, 2 mL	1/pk	A3967010
30 mg, 2 mL	1/pk	A3967030

The importance of SPE on instrumentation

Use SPE to remove common matrix background from your samples. This will minimize interfering peaks, reduce matrix effects to maximize sensitivity, reduce maintenance, and increase instrument lifetime by preventing contamination.

For Research Use Only. Not for use in diagnostic procedures. Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc. 2016
Printed in the U.S.A. June 22, 2016
5991-7020EN



Agilent Technologies