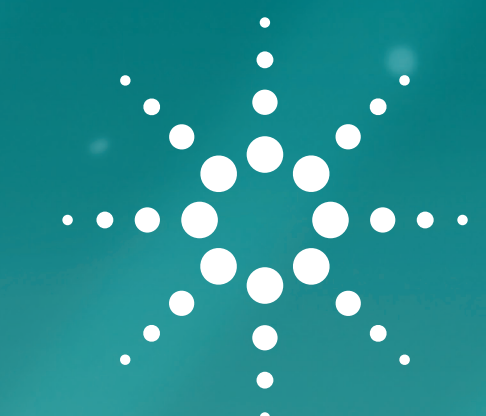


WORRY-FREE SAMPLE PREPARATION FOR FOOD AND ENVIRONMENTAL ANALYSIS

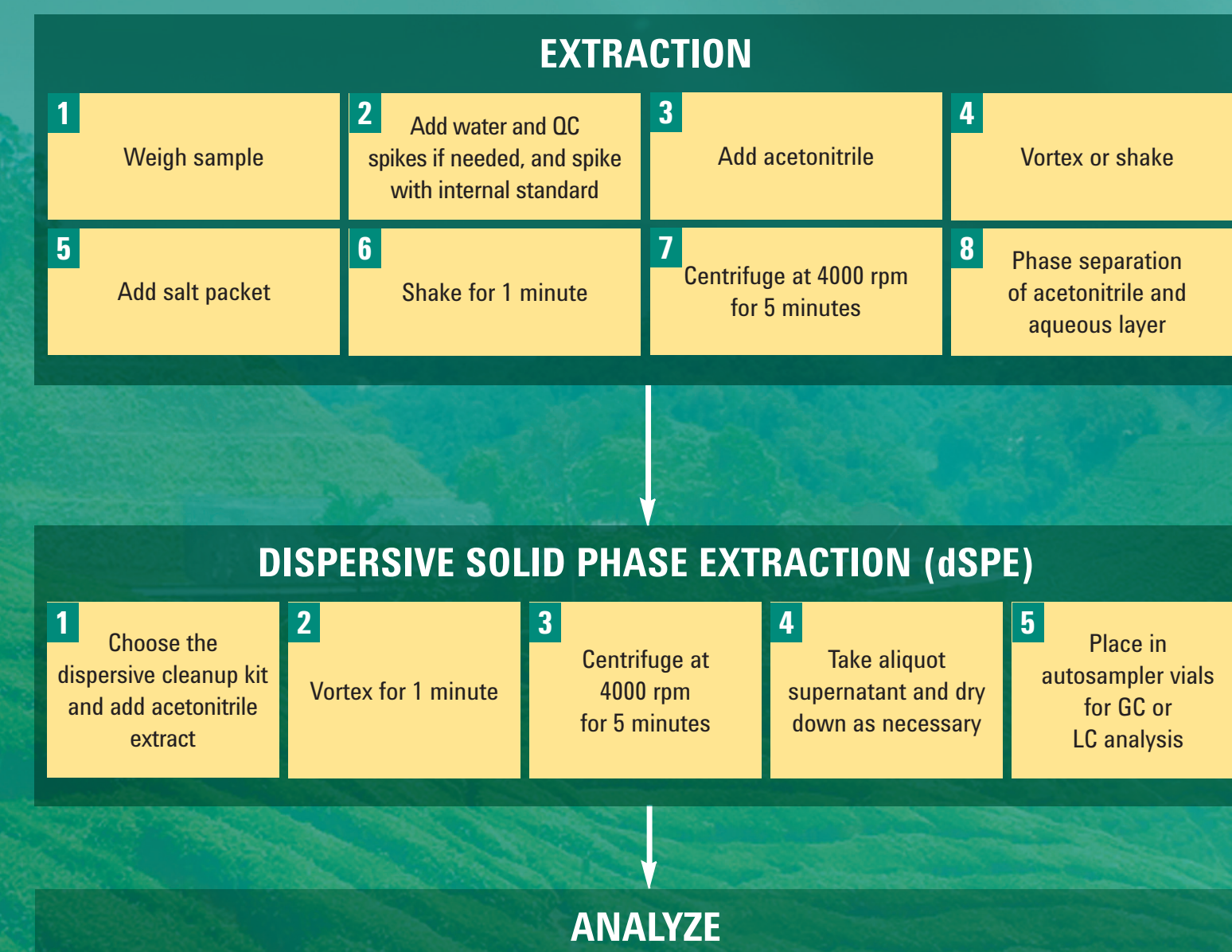
Agilent Bond Elut sample preparation ensures reproducible results right from the start



Save time and labor QuEChERS Kits

These easy-to-use kits remove interfering matrix compounds, allowing you to analyze samples for a broad range of pesticides, veterinary drugs, environmental contaminants, and other compounds of interest.

General Protocol for QuEChERS



QuEChERS application notes

- Food**
- 5990-4977EN Agilent Food Safety Applications Notebook: Volume 2 – Bond Elut QuEChERS
 - 5991-3638EN Sample Prep for Trace Analysis of Adulterants in Erectile Dysfunction Dietary Supplements
 - 5991-5553EN QuEChERS Combined with GC/MS/MS for Analysis of Over 200 Pesticides in Cereals
- Environmental**
- 5990-5452EN Analysis of Polycyclic Aromatic Hydrocarbons in Soil with Agilent Bond Elut HPLC-FLD

For more application notes please visit: www.agilent.com

Optimize productivity from sample prep to analysis and support

Consistent sample preparation

From SPE... to QuEChERS... to filtration... Agilent sample preparation products let you quickly produce aliquots that are representative, reproducible, and homogenous.

High-performing Agilent columns

Agilent JMW Ultra Inert GC columns deliver consistent column inertness and exceptionally low column bleed for lower LODs and more accurate data.

For the most efficient LC separations and reproducible results, choose InfinityLab Poroshell 120 LC columns.

Agilent QQQ and Q-TOF instruments

Multi-target screening: Agilent 6400 Series QQQ LC/MS and 7000C QQQ GC/MS

Non-targeted screening: Agilent 6500 Series Accurate-Mass Q-TOF and 7200 GC/Q-TOF

MassHunter Software

MassHunter enables you to produce high-quality MS data for targets and unknowns. The software helps your lab transform data into action.

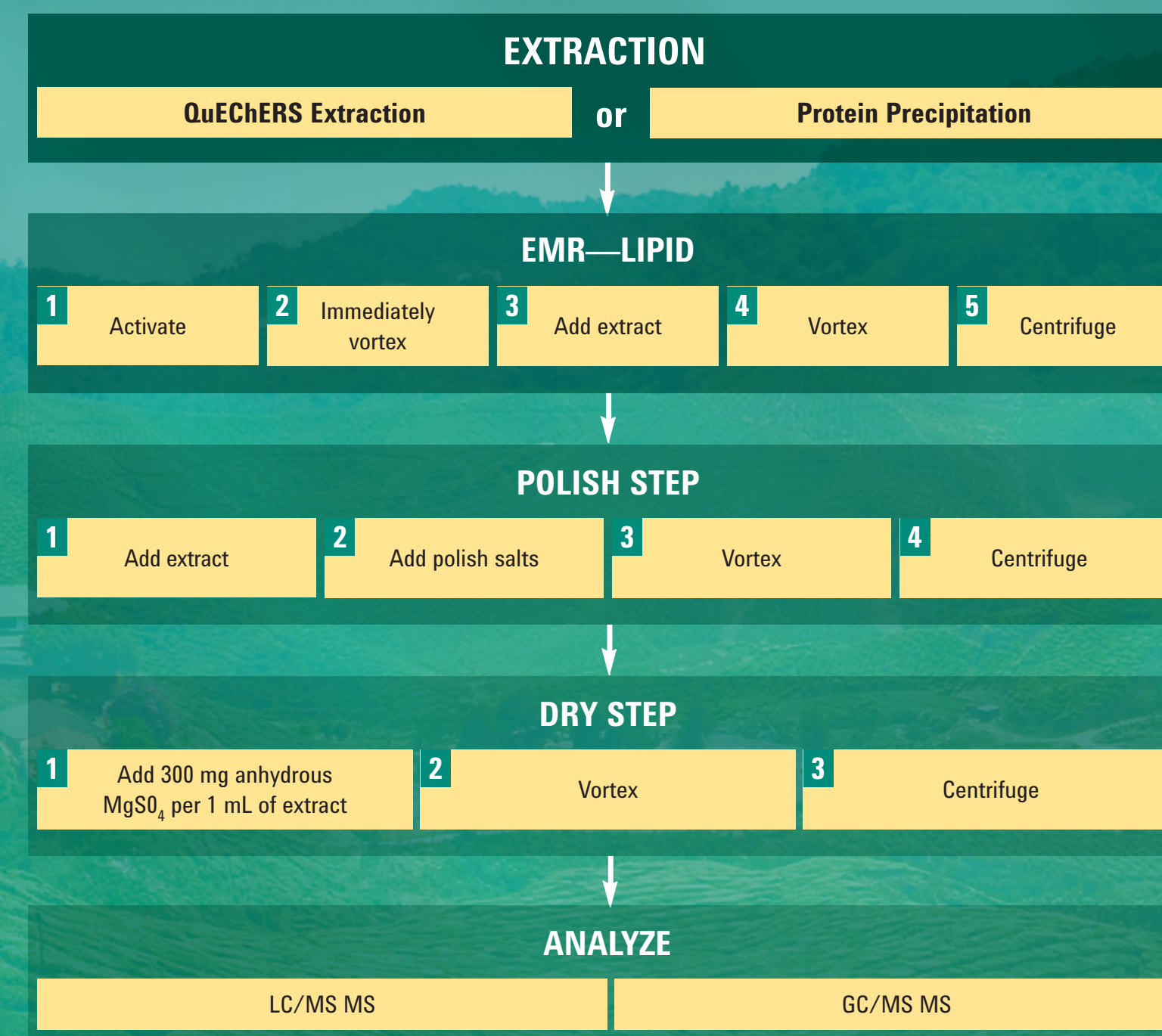
Agilent Services

Agilent CrossLab offers a comprehensive suite of enterprise services, learning and consultative solutions, as well as instrument services to maximize uptime, simplify administration, and protect the investments in your laboratory.

Lose the lipids, keep your analytes Enhanced Matrix Removal—Lipid

The innovative sorbent in Enhanced Matrix Removal—Lipid (EMR—Lipid) works with your current QuEChERS methodology to selectively remove lipids in complex matrices, without losing analytes of interest.

General Protocol for EMR—Lipid



EMR—Lipid application notes

- Food**
- 5991-6818EN Aflatoxin Analysis in Infant Formula with Enhanced Matrix Removal—Lipid by LC/MS/MS
 - 5991-6097EN Multiresidue Analysis of Pesticides in Avocado by GC/MS/MS with Bond Elut Enhanced Matrix Removal—Lipid
 - 5991-6096EN Multiresidue Analysis of Veterinary Drugs in Bovine Liver by LC/MS/MS with Bond Elut Enhanced Matrix Removal—Lipid
- Environmental**
- 5991-6088EN PAH Analysis in Salmon with Bond Elut Enhanced Matrix Removal—Lipid

Confidently clean samples and elute analytes Bond Elut SPE

The Bond Elut SPE family provides the largest choice of sorbent formats in the market today and includes polymeric SPE phases: Plexa (non-polar polymeric), Plexa PCX (cation exchanger mixed mode), and Plexa PAX (anion exchanger mixed mode).

General Protocol for Trouble-free SPE

Analyte	ACIDS	NEUTRALS	BASES
	LogP > 1.0 pKa < 5	LogP > 1.5 pKa 3-6 pKa 6-10	LogP > 0.8 pKa 6-10
	PLEXA PAX	PLEXA Acid Load Method PLEXA Base Load Method	PLEXA PCX
Sample Pre-treatment	2% NH ₄ OH	1% HCO ₂ H	2% NH ₄ OH
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		

SPE application notes

- Food**
- 5990-4917EN Agilent Food Safety Applications Notebook: Volume 1 – SPE, SPME, SLE, and Captiva Filtration
 - 5991-5414EN Analyze Quinolone Residues in Milk with an Agilent Poroshell 120 4 μm Column
 - 5991-5421EN Analyze Sulfonamide Residues in Pork with Agilent Poroshell 120 EC-C18, 4 μm Columns
- Environmental**
- 5991-5304EN Solid Phase Extraction of Organophosphorus Pesticides in Water with Agilent Bond Elut PPL
 - 5991-5490EN Monitoring of Bisphenol A and its Analogues in Environmental Matrices Using the Agilent 6550 Q-TOF LC/MS
 - 5991-5594EN EPA – Method 540: Selected Organic Contaminants Using Agilent Plexa Cartridges

Bond Elut QuEChERS, EMR—Lipid, and SPE Part Numbers

Method	Buffered	Contents	Part No.	Part No.	Part No.
			With Tubes	50/pk	Packets Only 200/pk
ADAC 2007.01	Yes	6 g MgSO ₄ , 1.5 g NaAcetate	5982-5755	5982-6755	5982-7755
Original (10 g samples)	No	4 g MgSO ₄ , 1 g NaCl	5982-5550	5982-6550	5982-7550
Original (15 g samples)	No	6 g MgSO ₄ , 1.5 g NaCl	5982-5555	5982-6555	5982-7555
EN 15682	Yes	4 g MgSO ₄ , 1 g NaCl, 1 g NaCitrate, 0.5 g disodium citrate sesquihydrate	5982-5650	5982-6650	5982-7650
Acrylamides	No	4 g MgSO ₄ , 0.5 g NaCl	5982-5850		
Veterinary Drugs	No	4 g Na ₂ SO ₄ , 1 g NaCl	5982-0032		

Kit	Size	Unit	AOAC 2007.01 Method	European Method EN 15662
			Kit Contents Part No.	Kit Contents Part No.
General fruits and vegetables	15 mL	50/pk	400 mg PSA, 1200 mg MgSO ₄ , 5982-5058, 5982-5058CH	150 mg PSA, 900 mg MgSO ₄ , 5982-5056, 5982-5056CH
Fruits and vegetables with fats and waxes	15 mL	50/pk	400 mg PSA, 400 mg C18EC, 1200 mg MgSO ₄ , 5982-5158, 5982-5158CH	150 mg PSA, 150 mg C18EC, 900 mg MgSO ₄ , 5982-5156, 5982-5156CH
Pigmented fruits and vegetables	15 mL	50/pk	400 mg PSA, 400 mg GCB, 1200 mg MgSO ₄ , 5982-5258, 5982-5258CH	150 mg PSA, 15 mg GCB, 885 mg MgSO ₄ , 5982-5256, 5982-5256CH
Highly pigmented fruits and vegetables	15 mL	50/pk		150 mg PSA, 45 mg GCB, 855 mg MgSO ₄ , 5982-5356, 5982-5356CH
Fruits and vegetables with pigments and fats	15 mL	50/pk	400 mg PSA, 400 mg GCB, 1200 mg MgSO ₄ , 400 mg C18EC, 5982-5456, 5982-5456CH	
Other food methods	15 mL	50/pk	150 mg C18, 900 mg MgSO ₄ , 5982-4956, 5982-4956CH	
All food types (universal)	15 mL	50/pk	400 mg PSA, 400 mg C18, 45 mg GCB, 1200 mg MgSO ₄ , 5982-0029, 5982-0029CH	

Part numbers ending in CH indicate tubes containing ceramic homogenizers. Agilent also offers many kits in 2 mL formats.

EMR—Lipid		
Description	Unit	Part No.
Bond Elut EMR—Lipid dSPE	50/pk	5982-1010
Bond Elut EMR—Lipid Final Polish Tube	50/pk	5982-0101
Bond Elut EMR—Lipid MgSO ₄ Polish Pouch	50/pk	5982-0102

Bond Elut Plexa			Bond Elut Plexa PCX			Bond Elut Plexa PAX		
Description	Unit	Part No.	Description	Unit	Part No.	Description	Unit	Part No.
Straight Barrel Cartridges			Straight Barrel Cartridges			Straight Barrel Cartridges		
30 mg, 1 mL	100/pk	12109301	30 mg, 1 mL	100/pk	12109301	30 mg, 1 mL	100/pk	12107301
30 mg, 3 mL	50/pk	12109303	30 mg, 1 mL	100/pk	12109601	60 mg, 1 mL	100/pk	12107901
60 mg, 1 mL	100/pk	12109601	30 mg, 3 mL	50/pk	12109303	30 mg, 3 mL	50/pk	12107903
60 mg, 3 mL	50/pk	12109603	60 mg, 3 mL	50/pk	12109603	60 mg, 3 mL	50/pk	12107903
200 mg, 3 mL	50/pk	12109610	200 mg, 6 mL	30/pk	12108206	200 mg, 6 mL	30/pk	12107206
200 mg, 6 mL	30/pk	12109206	500 mg, 6 mL	30/pk	12258506	500 mg, 6 mL	30/pk	12257506
500 mg, 6 mL	30/pk	12259506						
Bond Elut Jr			Bond Elut Jr			Bond Elut Jr		
200 mg	50/pk	121696108						
Mega Bond Elut Plexa			Mega Bond Elut Plexa			Mega Bond Elut Plexa		
500 mg, 12 mL	20/pk	327832						

Learn more about Agilent Sample Prep Solutions www.agilent.com/chem/sampleprep

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc. 2018
Printed in the U.S.A. May 23, 2018
5991-5996EN