

MAKE QuEChERS AS EASY AS 1-2-3 WITH AGILENT BOND ELUT QuEChERS KITS

Step 1 EXTRACT

Choose your salts

AOAC Method 2007.01:

6 g MgSO₄,
1.5 g Na acetate
(5982-5755)

or

EN Method 15662:

4 g MgSO₄, 1 g NaCl,
1 g Na citrate, 0.5 g disodium
citrate sesquihydrate
(5982-5650)

or

Original QuEChERS Extraction:

10 g sample (5982-5550):
4 g MgSO₄, 1 g NaCl
15 g sample (5982-5555):
6 g MgSO₄, 1.5 NaCl



Step 2 CLEAN

Choose a dispersive
SPE kit* specific to
your matrix

AOAC

5982-5058
400 mg PSA, 1,200 mg MgSO₄

5982-5258
400 mg PSA, 400 mg GCB,
1,200 mg MgSO₄

5982-5158
400 mg PSA, 400 mg C18EC,
1,200 mg MgSO₄

5982-5456
400 mg PSA, 400 mg GCB, 400
mg C18EC, 1,200 mg MgSO₄

5982-0029
400 mg PSA, 400 mg C18,
45 mg GCB, 1,200 mg MgSO₄

5982-4950
50 mg PSA, 150 mg C18EC,
900 mg Na₂SO₄ (2 mL)

5982-4956
150 mg C18, 900 mg MgSO₄

General Fruits and Vegetables

Removes polar organic acids,
some sugars and lipids

Pigmented Fruits and Vegetables

Removes polar organic acids,
some sugars and lipids, and
carotenoids and chlorophyll;
not for use with planar
pesticides

Highly Pigmented Fruits and Vegetables

Removes polar organic acids,
some sugars and lipids, plus
high levels of carotenoids and
chlorophyll; not for use with
planar pesticides

Fruits and Vegetables with Fats and Waxes

Removes polar organic acids,
some sugars, more lipids and
sterols

Fruits and Vegetables with Pigments and Fats

Removes polar organic acids,
some sugars and lipids, plus
carotenoids and chlorophyll;
not for use with planar
pesticides

High-Fat Matrices

Removes lipids and other
fatty interferences, including
straight-chain aliphatic
compounds; also effective in
removing some pigments

General Food

Removes all matrix-interfering
materials including polar
organic acids, lipids, sugars,
proteins, carotenoids, and
chlorophyll

Animal Origin Food

Removes matrix interferences
such as polar organic salts,
sugars, lipids and proteins

Other Foods

Removes biological matrix
interferences, including
hydrophobic substances (fats,
lipids) and proteins

EN

5982-5056
150 mg PSA,
900 mg MgSO₄

5982-5256
150 mg PSA,
15 mg GCB,
885 mg MgSO₄

5982-5356
150 mg PSA,
45 mg GCB,
855 mg MgSO₄

5982-5156
150 mg PSA,
150 mg C18EC,
900 mg MgSO₄

Enhanced Matrix Removal—Lipid

Bond Elut QuEChERS dSPE Enhanced Matrix Removal—Lipid (5982-1010) in combination with Bond Elut Final Polish for Enhanced Matrix Removal—Lipid (5982-0101 or 5982-0102) is specifically designed to remove lipids from high-fat samples. The sorbent removes interfering lipids from samples while leaving analytes behind.

Step 3 ANALYZE

with Poroshell LC Columns or Ultra Inert GC Columns
using the 6400 Series Triple Quadrupole LC/MS
or 7000C Triple Quadrupole GC/MS System for
targeted screening or the 6500 Series Accurate-Mass
Quadrupole Time-of-Flight (Q-TOF) or 7200 GC/Q-TOF
for non-targeted screening.



For more information on Agilent Bond Elut QuEChERS products visit www.agilent.com/chem/quechers

Agilent
UNIVERSITY

DID YOU KNOW?

We now offer a complete curriculum of flexible learning solutions from Agilent University.

Agilent University caters to all your individual and enterprise training needs by delivering a complete curriculum of flexible learning solutions. At Agilent University, you can find a wide range of learning options to meet the needs of your lab, helping optimize productivity on your new or existing equipment.

To learn more go to www.agilent.com/crosslab/university

*Shown here are kits in a 15 mL format unless otherwise noted. Agilent also offers many kits in 2 mL formats.

Information, descriptions, and specifications in this publication are subject to change without notice.
© Agilent Technologies, Inc. 2015 Printed in the U.S.A. November 20, 2015 5990-5324EN