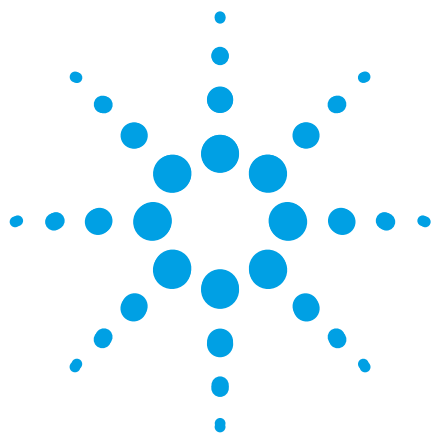




Screening PAHs in Soil Using RTL Database with Agilent Bond Elut QuEChERS Extraction Kits and Agilent 5975 LTM GC/MS

Application Note

This application note is temporarily unavailable.

www.agilent.com/chem

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

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

© Agilent Technologies, Inc., 2012
Published in the USA,
August 30, 2012
5990-6324EN



Agilent Technologies