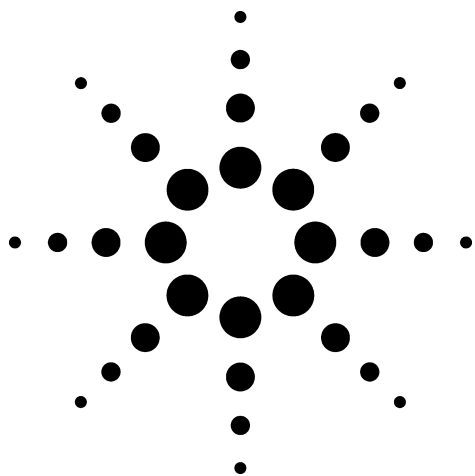
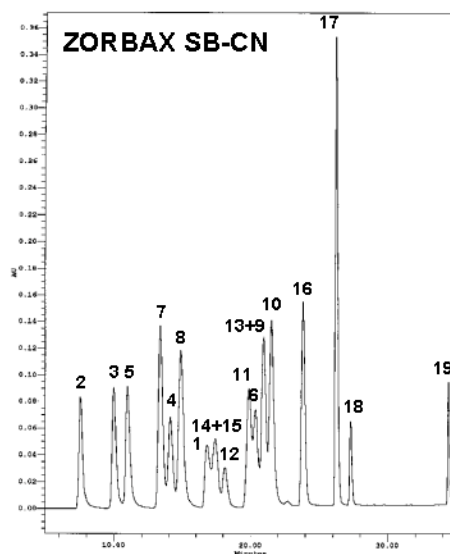
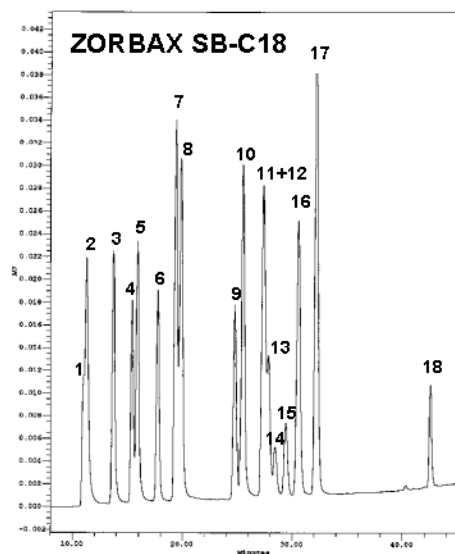




Qualitative and Quantitative Analysis of Explosives and Related Compounds Using Polar and Nonpolar HPLC Columns

Application
Environmental
Robert Ricker



Courtesy of Th. Renner, Hess. Landesanstalt f. Umwelt, Dez. III/2, P.O. Box 3209, 65022 Wiesbaden, Germany

Conditions:
ZORBAX® SB-C18, SB-CN (2.1 x 150 mm) (Agilent P/N: 883700.905, 883700.922)
Mobile Phase: A: ACN + 5% H₂O + 5 mM CF₃COONH₄
 B: H₂O + 5% ACN + 5 mM CF₃COONH₄
 pH 2.7 (CF₃COOH)
Inject: 10 µL of 19 nitromethanes in ACN:H₂O (20:80), 5mM CF₃COONH₄;
 0.23mL/min, 18°C, Detect. UV(210, 240, 360 nm)



Agilent Technologies

Gradient Shapes

ZORBAX[®] SB-C18 (2.1x150 mm) (P/N:883700.922)
Sample: 0.5ng / µL

Time [min]	Flow [mL/min]	% A [%]	% B [%]
0	0.23	20	80
2	0.23	20	80
10	0.23	30	70
20	0.23	35	65
25	0.23	40	60
35	0.23	70	30
40	0.23	70	30
42	0.23	20	80
45	0.23	20	80

ZORBAX[®] SB-CN (2.1x150 mm) (P/N:883700.905)
Sample: 5ng / µL

Time [min]	Flow [mL/min]	% A [%]	% B [%]
0	0.23	20	80
1	0.23	20	80
15	0.23	30	70
30	0.23	80	20
35	0.23	80	20
37	0.23	20	80
40	0.23	20	80

Robert Ricker is an application chemist
based at Agilent Technologies, Wilmington,
Delaware.

For more information on our products and
services, visit our website at:
www.agilent.com/chem

Table 1 - List of Analytes

	Compound	Rt [min] on Zorbax SB-C18	Rt [min] on Zorbax SB-CN	Quantitation on Zorbax SB
1	picric acid	11.02	16.82	CN (360 nm)
2	4-amino-2-nitrotoluene	11.31	7.57	C18 (240 nm)
3	2-amino-6-nitrotoluene	13.72	10.00	C18 (240 nm)
4	RDX	15.42	14.13	C18 (210 nm)
5	2-amino-4-nitrotoluene	15.92	11.02	C18 (240 nm)
6	HMX	17.76	20.37	C18 (210 nm)
7	1,3-dinitrobenzene	19.41	13.40	CN (240 nm)
8	1,3,5-trinitrobenzene	19.87	14.90	CN (240 nm)
9	2-amino-4,6-dinitrotoluene	24.79	20.98	C18 (240 nm)
10	2,4-dinitrotoluene	25.52	21.55	C18 (240 nm)
11	4-amino-2,6-dinitrotoluene	27.37	19.90	CN (240 nm)
12	2-nitrotoluene	27.41	18.13	CN (240 nm)
13	2,6-dinitrotoluene	27.84	20.87	C18 (240 nm)
14	4-nitrotoluene	28.47	17.39	C18 (240 nm)
15	3-nitrotoluene	29.46	17.43	C18 (240 nm)
16	2,4,6-trinitrotoluene	30.59	23.85	CN (240 nm)
17	tetryl	32.22	26.30	CN (240 nm)
18	diphenylamine	42.59	27.32	CN (210 nm)
19	hexyl	45.78	34.45	CN (360 nm)

Copyright© 2002 Agilent Technologies, Inc.
All Rights Reserved. Reproduction,
adaptation or translation without prior
written permission is prohibited, except as
allowed under the copyright laws.

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.

Printed in the USA

April 25, 2002

5988-6345EN



Agilent Technologies