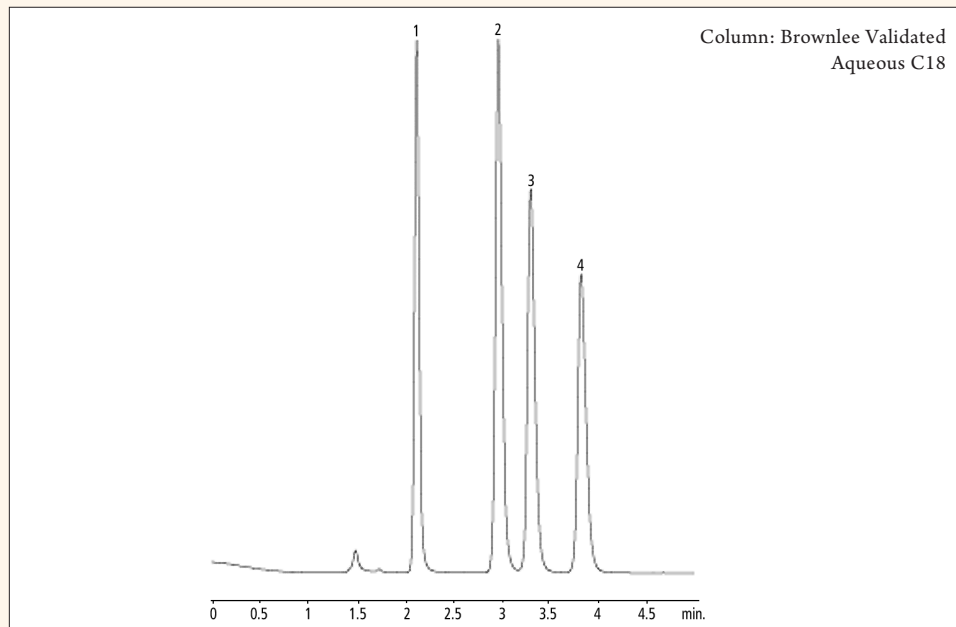


CARBOXYLIC ACIDS

HPLC

Biochemical: Carboxylic Acids on Brownlee Validated Aqueous C18



Peak List

Peak List	Concentration (mg/mL)
1. glycolic acid	2.0
2. malonic acid	1.8
3. lactic acid	3.3
4. acetic acid	2.1

Concentration (mg/mL)

Conditions

Mobile phase	50 mM potassium phosphate, pH 2.5
Flow	1.0 mL/min
Temperature	ambient
Detector	UV @ 210 nm

Column

Column name	Brownlee Validated Aqueous C18
Part number	N9303546
Dimensions	150 x 4.6 mm
Particle size	5 μ m
Pore size	100Å

Sample

Injection volume	10 μ L
Solvent	mobile phase

Notes

Recommended Guard Cartridge	N9303691
Recommended Guard Holder	N9303713

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright © 2006-2012 PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

007568A_07