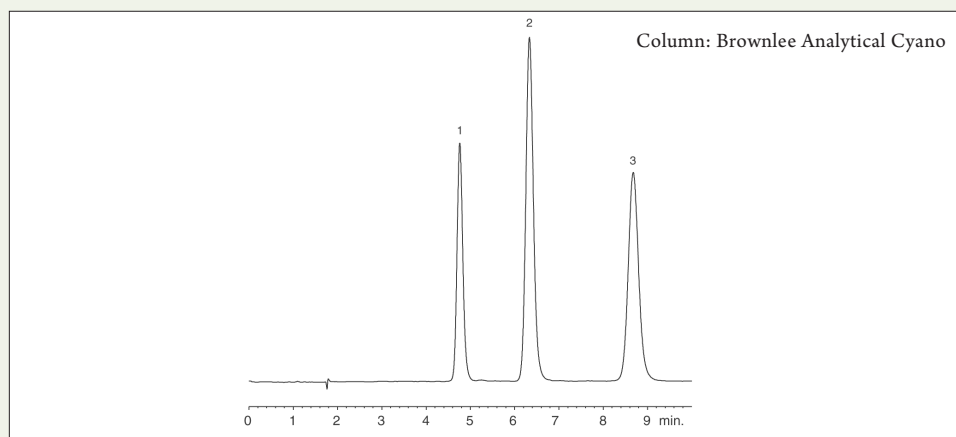




CORTICOSTEROIDS/ BRONCHODILATOR HPLC

Pharmaceutical: Corticosteroids on Brownlee Analytical Cyano (Reversed Phase)



Peak List

	Concentration (mg/mL)
1. cortisone	0.1
2. corticosterone	0.5
3. cortisone acetate	0.7

Concentration (mg/mL)

Conditions

Mobile phase:	water:methanol (65:35, v/v)
Flow:	1.0 mL/min
Temperature:	30 °C
Detector:	UV @ 254 nm
Pressure:	133 bar

Column

Column name:	Brownlee Analytical Cyano
Part number:	N9303522
Dimensions:	150 x 4.6 mm
Particle size:	5 µm
Pore size:	110Å

Sample

Injection volume:	5.0 µL
Solvent:	methanol

Notes

Recommended Guard Cartridge	N9303681
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


PerkinElmer[®]
For the Better

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

Copyright ©2006-2014 PerkinElmer, Inc. All rights reserved. PerkinElmer[®] is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

007573A_41