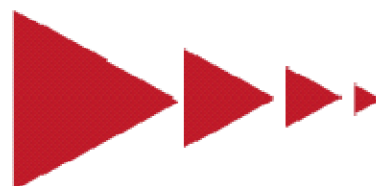


CHIRAL

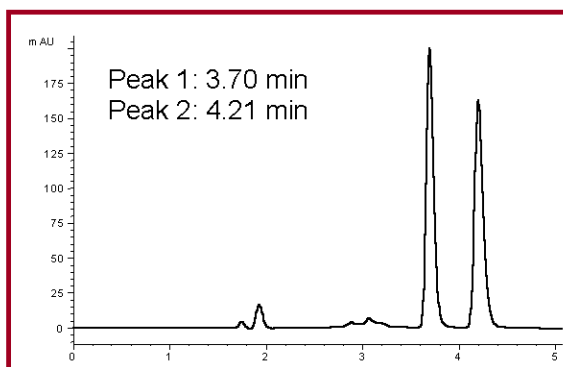
APPLICATION NOTE



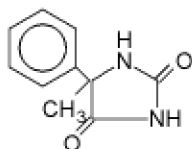
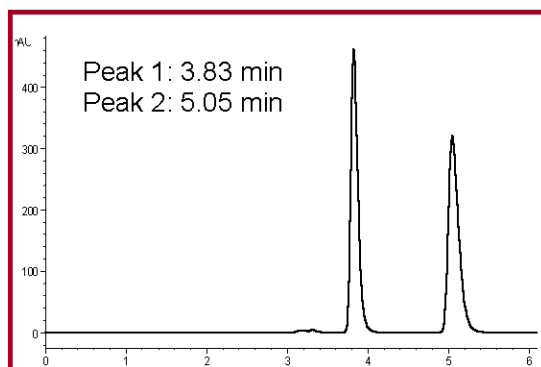
PERFORMANCE TESTING OF CHIROBIOTIC™ COLUMNS

To ensure the selectivity and performance of all CHIROBIOTIC™ LC columns, periodically test your columns. This can now be accomplished with a single compound for all the CHIROBIOTIC™ phases in a very simple mobile phase of 100% MeOH.

CHIROBIOTIC V



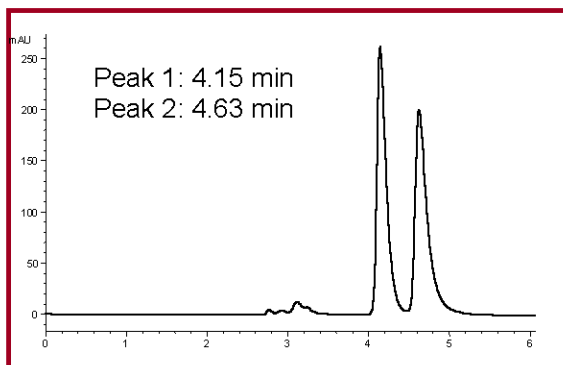
CHIROBIOTIC T



CONDITIONS FOR ALL COLUMNS:

Sample: 5-Methyl-5-Phenylhydantoin (Aldrich 18,082-3)
Column size: 250 x 4.6mm
Mobile phase: 100% MeOH
Flow rate: 1 mL/min
UV: 220nm

CHIROBIOTIC R



CHIROBIOTIC TAG

