

Application Report 140

Separation of Aromatic Compounds Using Ascentis™ RP-Amide

This application demonstrates the suitability of Ascentis RP-Amide in determining aromatic selectivity of amylbenzene, butylbenzene, o-terphenyl and triphenylene by HPLC.

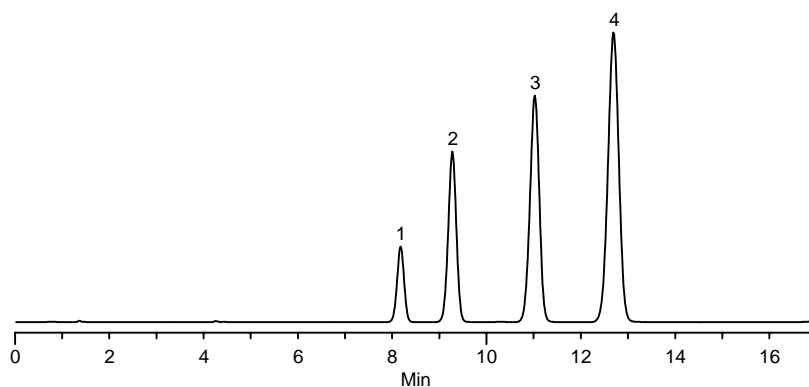
Key Words

Ascentis RP-Amide, 565324-U, aromatics, selectivity, amylbenzene 538-68-1, 11,317-4, butylbenzene 104-51-8, B9,020-3, o-terphenyl, 84-15-1, T2800, triphenylene, 217-59-4, T8,260-0

Author: Jacinth A.M. McKenzie

Acquisition System: Waters

Notebook Reference: 1550-27



G002607

Conditions

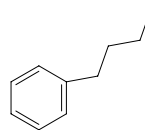
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)
mobile phase: 30:70, water:acetonitrile
temp.: 35 °C
flow rate: 1.0 mL/min.
det.: UV, 254 nm
injection: 2 µL
sample: as indicated in acetonitrile

Peak IDs

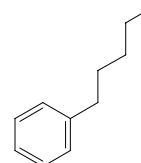
1. Butylbenzene (2.5 mL/mL)
2. o-terphenyl (150 mg/mL)
3. Amylbenzene (2.5 mL/mL)
4. Triphenylene (50 mg/mL) (2.5 mL/mL)

Structures

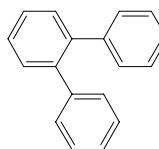
Butylbenzene - G002604



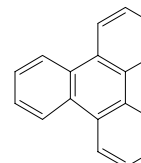
Amylbenzene - G002605



o-terphenyl - G002606



Triphenylene - G002608



Ascentis is a trademark of Sigma-Aldrich Co.