

Application Report 226

Phenolic and Benzyl Compounds Using Ascentis™ RP-Amide

This application demonstrates the suitability of Ascentis RP-Amide for the efficient separation of phenolic and benzyl compounds by HPLC.

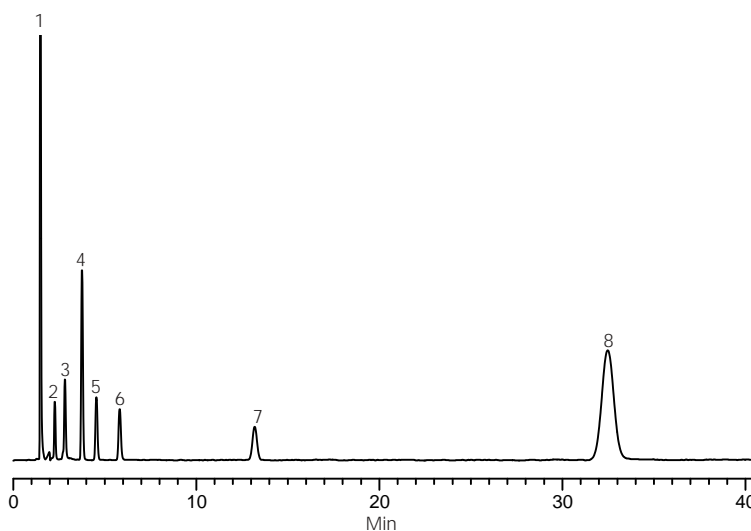
Key Words

Ascentis RP-Amide, 565324-U, uracil, 66-22-8, U0750, pyrogallol, 87-66-1, 254002, phloroglucinol, 6099-90-7, P38005, benzamide, 55-21-0, 135828, phenol, 108-95-2, 431516, resorcinol, 108-46-3, R406, catechol, 120-80-9, 135011, nitrobenzene, 98-95-3, 252379

Author: Hugh M. Cramer

Acquisition System: Shimadzu
Mongaup

Notebook Reference: 1551-81



G002899

Conditions

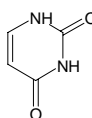
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)
mobile phase: 80:20, 10 mM ammonium acetate (pH 6.7 unadjusted):acetonitrile
flow rate: 1 mL/min.
temp.: 35 °C
det.: UV at 254 nm
injection: 10 µL
sample: as indicated in mobile phase

Peak IDs

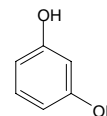
- | | |
|------------------------------|----------------------------|
| 1. Uracil (10 µg/mL) | 5. Resorcinol (50 µg/mL) |
| 2. Phloroglucinol (50 µg/mL) | 6. Catechol (50 µg/mL) |
| 3. Pyrogallol (50 µg/mL) | 7. Phenol (50 µg/mL) |
| 4. Benzamide (50 µg/mL) | 8. Nitrobenzene (50 µg/mL) |

Structures

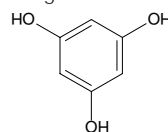
Uracil - G002739



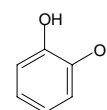
Resorcinol - G002594



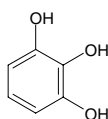
Phloroglucinol - G001274



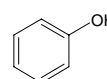
Catechol - G002590



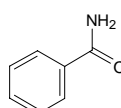
Pyrogallol - G002897



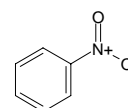
Phenol - G002896



Benzamide - G002894



Nitrobenzene - G002895



Ascentis is a trademark of Sigma-Aldrich Co.