

Application Report 143

Separation of Hydrophobic Bases Using Ascentis™ RP-Amide

This application demonstrates the suitability of Ascentis RP-Amide for the efficient separation of peptides with varying hydrophobicities.

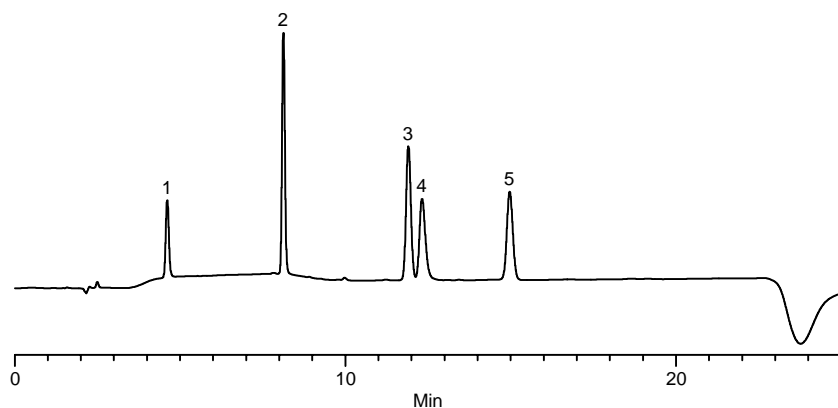
Key Words

Ascentis RP-Amide, 565324-U, peptides, angiotensin II, methionine enkephalin, leucine enkephalin, H2016

Author: Jacinth A.M. McKenzie

Acquisition System: Hitachi LC

Notebook Reference: 1550-22



G002613

Conditions

column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)
mobile phase A: 0.1 % TFA in water
mobile phase B: 0.1 % TFA in acetonitrile
flow rate: 1.0 mL/min.
temp.: 30 °C
det.: UV, 220 nm
injection: 5 µL
sample: as indicated in mobile phase A
gradient:

Min	%B
5	9-21
15	21-27

Peak IDs

1. Gly-Phe (62.5 mg/mL)
2. Val-Tyr-Val (250 mg/mL)
3. Methionine Enkephalin (250 mg/mL)
4. Angiotensin II (250 mg/mL)
5. Leucine Enkephalin (250 mg/mL)

Structures

Methionine Enkephalin - G002614

YGGFM

Leucine Enkephalin - G002615

YGGFL

Angiotensin II - G002616

RVYIHPF

Ascentis is a trademark of Sigma-Aldrich Co.