

Application Report 189

Analysis of Hydrophobic Peptide Probes Using Ascentis™ RP-Amide

Excellent resolution of these hydrophobic peptide probes is demonstrated on Ascentis RP-Amide, as is the excellent peak shape even in the absence of an ion-pairing reagent.

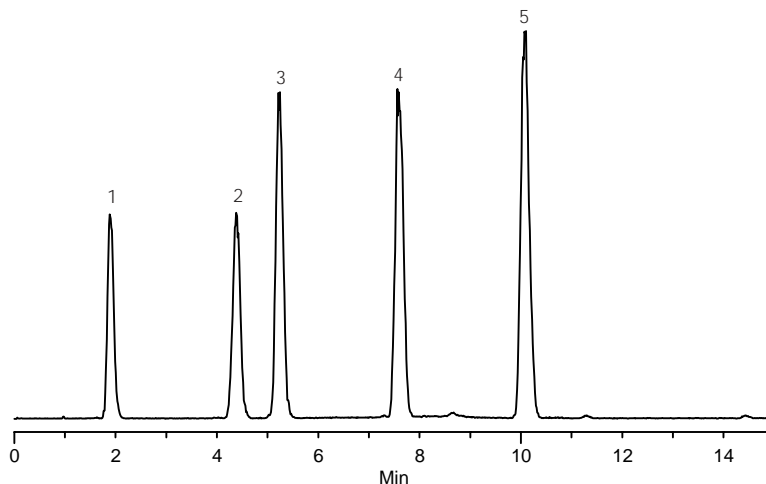
Key Words

hydrophobic peptides; peptides; Ascentis RP-Amide, 565304-U

Author: Hillel Brandes

Acquisition System: Micro-Tech Scientific UltraPlus II / Thermo LCQ Advantage

Notebook Reference: 1524



G002682

Conditions

column: Ascentis RP-Amide, 10 cm x 2.1 mm I.D., 5 µm particles (565304-U)
mobile phase A: 50:50, 50 mM formic acid titrated with ammonium hydroxide (pH 3.0):water
mobile phase B: 50:50, 50 mM formic acid titrated with ammonium hydroxide (pH 3.0):acetonitrile
flow rate: 0.2 mL/min.
temp.: ambient
det.: ESI (+)
injection: 4 µL
sample: as indicated in 10 mM formic acid
gradient:

Min	%A	%B
0	85	15
15	55	45

Peak IDs

1. Peptide 1 (20 mg/L)
2. Peptide 2 (20 mg/L)
3. Peptide 3 (20 mg/L)
4. Peptide 4 (20 mg/L)
5. Peptide 5 (20 mg/L)

Structures

Peptide 1 - G002683

RGAVGLGLGK-amide

Peptide 2 - G002684

ac-RGGGGLGLGK-amide

Peptide 3 - G002685

ac-RGAGGLGLGK-amide

Peptide 4 - G002686

ac-RGVGGLGLGK-amide

Peptide 5 - G002687

ac-RGVVGLGLGK-amide

Ascentis is a trademark of Sigma-Aldrich Co.