

Application Report 191

Analysis of Serum-derived Peptide Probes Using Ascentis™ RP-Amide

These naturally occurring peptide sequences vary in their number of positive charges at low pH from +2 to +4, and exhibit excellent peak shape on Ascentis RP-Amide even in the absence of ion-pairing reagents.

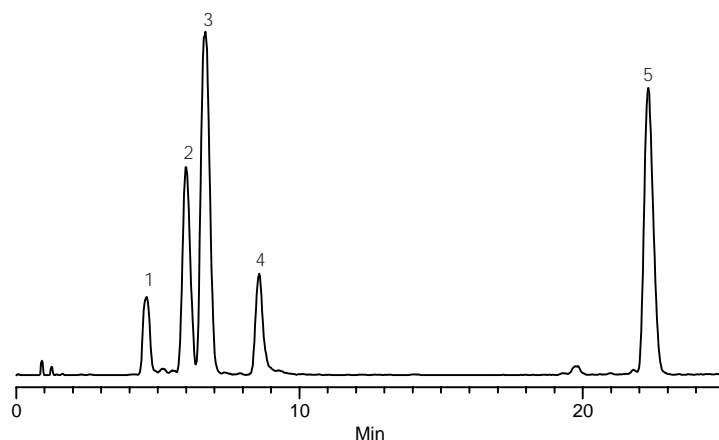
Key Words

serum peptides, cationic peptides, LC-MS, Ascentis RP-Amide, 565304-U

Author: Hillel Brandes

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

Notebook Reference: 1524



G002689

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	70	30
25	45	55

Peak IDs

1. Peptide 1, (M+H)⁺ = 1142.0, (M+2H)²⁺ = 572.6; (10 mg/L)
2. Peptide 2, (M+2H)²⁺ = 810.0; (10 mg/L)
3. Peptide 3, (M+2H)²⁺ = 649.5; (10 mg/L)
4. Peptide 4, (M+2H)²⁺ = 885.5; (10 mg/L)
5. Peptide 5, (M+2H)²⁺ = 977.8, (M+3H)³⁺ = 652.6; (10 mg/L)

Structures

Peptide 1 - G002690

HINIDQFVR

Peptide 2 - G002691

TEVSSNHVLIYLDK

Peptide 3 - G002692

DRVYIHPFHL

Peptide 4 - G002693

DNENVVNEYSSELEK

Peptide 5 - G002694

FTVDRPFLFLIYEHR

Ascentis is a trademark of Sigma-Aldrich Co.