

Application Report 120

LC-MS Analysis of Basic Peptide Probes Using Ascentis™ C18

Lack of severe tailing with these basic peptides, chromatographed in the absence of any strong anionic ion-pairing reagents, demonstrates the very low silanol activity of the silica surface of Ascentis C18.

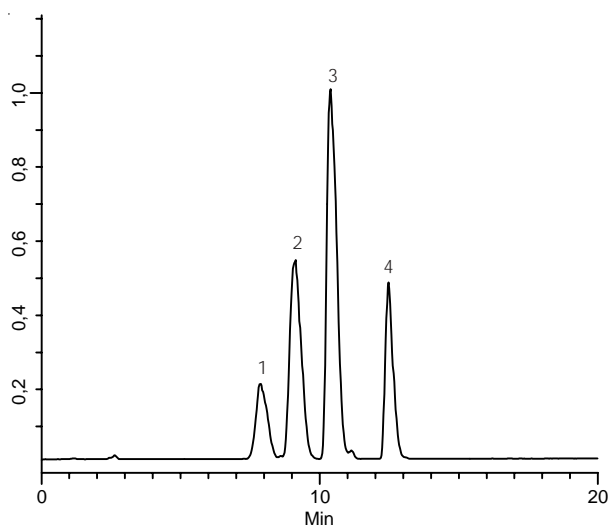
Key Words

basic peptides, cationic peptides, Ascentis C18; LC-MS; mass spectrometry, 581326-U

Author: Hillel Brandes

Acquisition System: Thermo Surveyor

Notebook Reference: 1498



G002444

Conditions

column: Ascentis C18, 10 cm x 2.1 mm I.D., 5 µm particles (581326-U)
mobile phase A: 20 mM formic acid in water
mobile phase B: 50:50, 20 mM formic acid in water:20 mM formic acid in acetonitrile
flow rate: 0.2 mL/min
temp.: 22 °C
det.: ESI (+), full scan
injection: 2 µL
gradient:

Min	%A	%B
0	80	20
20	40	60

Peak IDs

1. Peptide 1 (100 µg/mL)
2. Peptide 2 (100 µg/mL)
3. Peptide 3 (100 µg/mL)
4. Peptide 4 (100 µg/mL)

Structures

- Peptide 1
ac-GGGLGGAGGLK-amide
- Peptide 2
ac-KYGLGGAGGLK-amide
- Peptide 3
ac-GGAVKALKGLK-amide
- Peptide 4
ac-KYALKALKGLK-amide

Ascentis is a trademark of Sigma-Aldrich Co.