

# Application Report 190

## Analysis of Hydrophobic Peptide Probes Using Ascentis™ C18

Excellent resolution of these hydrophobic peptide probes is demonstrated on Ascentis C18, as is the excellent peak shape even in the absence of an ion-pairing reagent.

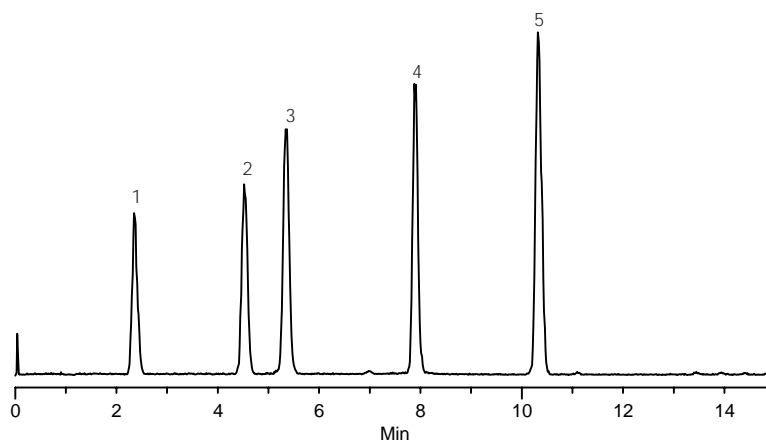
### Key Words

hydrophobic peptides; peptides; Ascentis C18, 581326-U

Author: Hillel Brandes

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

Notebook Reference: 1524



G002688

### Conditions

column: Ascentis C18, 10 cm x 2.1 mm I.D., 5 µm particles (581326-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	80	20
15	50	50

### 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.