

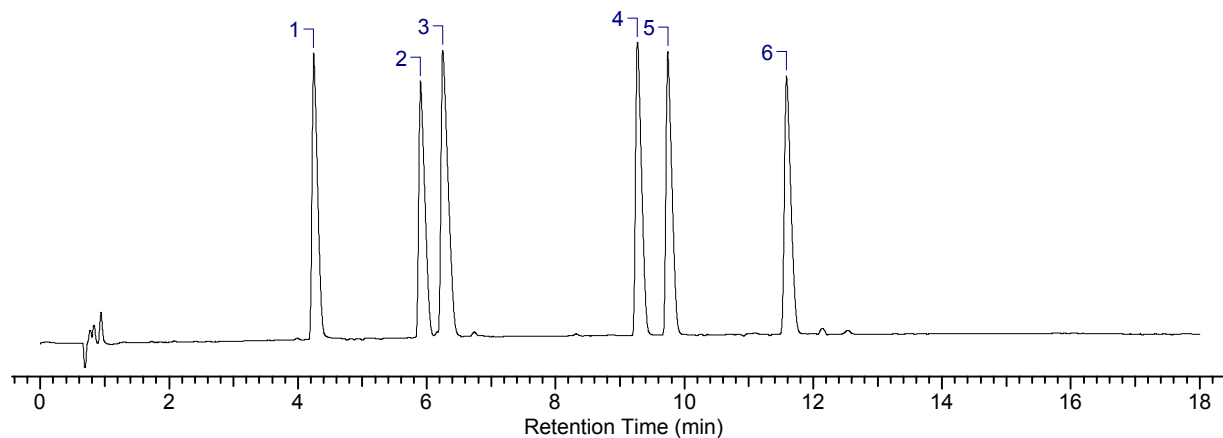
Application Report

Basic Peptides (Angiotensin analogs) on Ascentis Express Peptide ES-C18

Bonded Phase octadecyl
Column Name Ascentis Express Peptide ES-C18
Instrument Name Shimadzu LC-10ADvp

Key Words

Ascentis Express, angiotensin, basic peptide, 53313-U



Conditions

column: Ascentis Express Peptide ES-C18, 10 cm x 3 mm I.D.
2.7 μ m particles (53313-U)
mobile phase A: 0.1% (v/v) formic acid, pH 4 (with ammonium hydroxide)
mobile phase B: 50:50, [0.175% (v/v) formic acid, pH 4] : acetonitrile
flow rate: 0.6 mL/min.
initial pressure: 207 bar (3000 psi)
temp.: 35 °C
det.: UV 215 nm
injection: 2 μ L
sample: 0.5 g/mL ea. peptide
gradient:

min	% A	% B
0	80	20
15	50	50

Peak IDs

No.	Name	tR
1	IVYIHPR	4.2
2	KVYIHPI	5.9
3	RVYIHPI	6.2
4	ac-KVYIHPI	9.3
5	ac-RVYIHPI	9.7
6	ac-IVYIHPR	11.6