

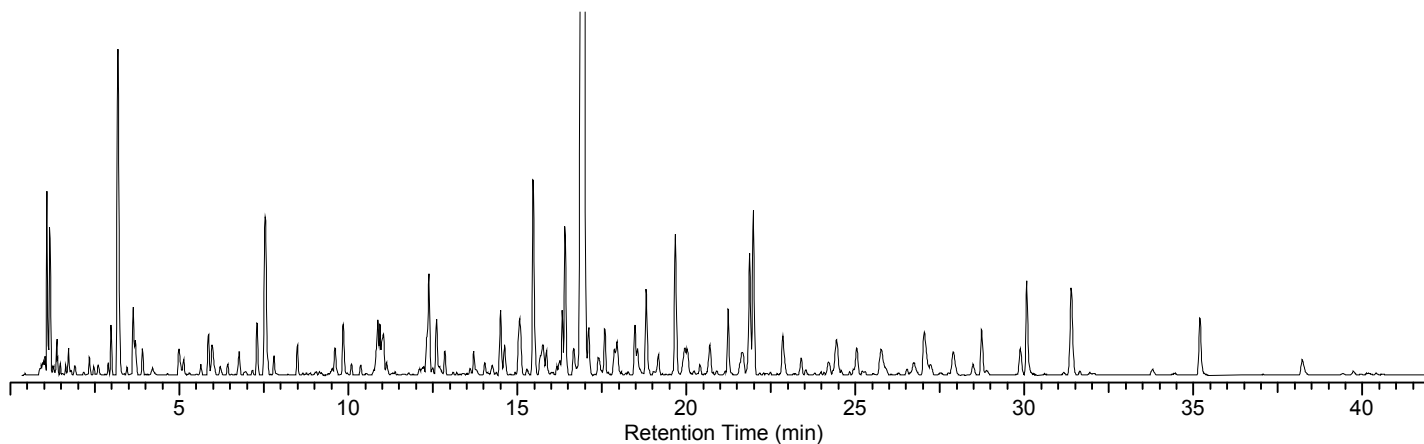
## Application Report

### Tryptic Digests of Protein Mixture on Ascentis Express Peptide ES-C18

**Bonded Phase** octadecyl  
**Column Name** Ascentis Express Peptide  
ES-C18  
**Instrument Name** Agilent 1200SL

#### Key Words

tryptic digest, proteolytic digest, trypsin, proteomics, peptide map, peptide mapping, 53307-U



#### Conditions

column: Ascentis Express Peptide ES-C18, 15 cm x 2.1 mm I.D.  
2.7  $\mu$ m particles (53307-U)  
mobile phase A: 0.1% formic acid in water  
mobile phase B: 25:75, (0.4% formic acid in water) : acetonitrile  
flow rate: 0.3 mL/min.  
temp.: 35 °C  
det.: ESI(+), TOF  
injection: 2  $\mu$ L  
sample: 10 pmol/L in 0.1% formic acid  
gradient: 

min	% A	% B
0	94	6
40	0	70