



## HPLC peptide standard mixture

# Dependence of peptide separations on sequence length and amino acid side chains

## Application Note

BioPharma

### Authors

Agilent Technologies, Inc.

### Introduction

The separation of a peptide standard mixture consisting of 5 peptides of varying length and composition was used to study this effect. The mix contained selected small peptides, e.g. a dipeptide, a tripeptide, 2 pentapeptides differing in one amino acid, and one 13-amino acid long peptide. This mixture was thus useful to study chromatographic differences arising due to a change in peptide length as well as monitor side-chain differences in peptides of the same length (e.g. methionine and leucine enkephalins). Distinct retention time differences are observed for the 5 peptides, a good base-line separation is achieved between methionine and leucine enkephalins. In contrast to other RP-HPLC columns, base-line separation between leucine enkephalin and angiotensin II is also obtained.



**Agilent Technologies**

## Conditions

Technique : Agilent Polaris C18-A, 4.6 x 150 mm, 5  $\mu$ m  
Part no. A2000150X046

Eluent : Solvent A: H<sub>2</sub>O + 0.1% TFA,  
Solvent B: CH<sub>3</sub>CN + 0.1% TFA

Gradient : 10% - 65% B in 10 min

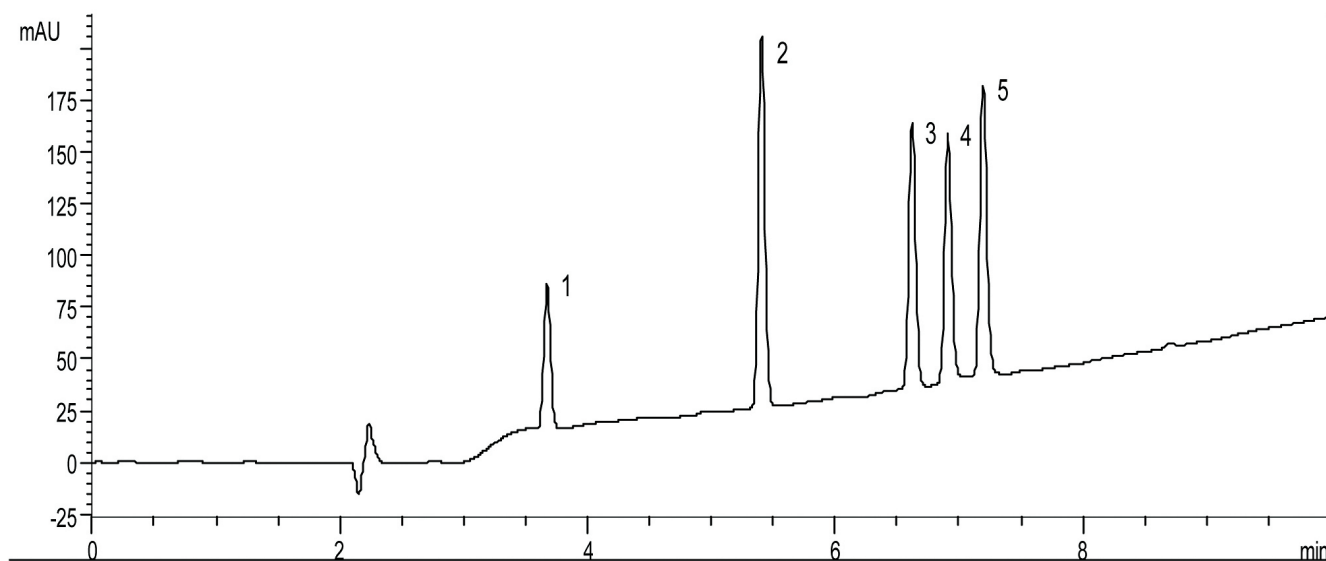
Flow : 1.0 mL/min

Temperature : ambient

Detection : 220 nm

## Peak identification

1. gly-tyr
2. vai-tyr-val
3. methionine enkephalin (tyr-gly-gly-phe-met)
4. leucine enkephalin (tyr-gly-gly-phe-leu)
5. angiotensin II (pglu-leu-tyr-glu-asn-lys-pro-arg-arg-pro-tyr-Ile-leu)



[www.agilent.com/chem](http://www.agilent.com/chem)

This information is subject to change without notice.

© Agilent Technologies, Inc. 2011

Printed in the USA

31 October, 2011

First published prior to 11 May, 2010

A02056



**Agilent Technologies**