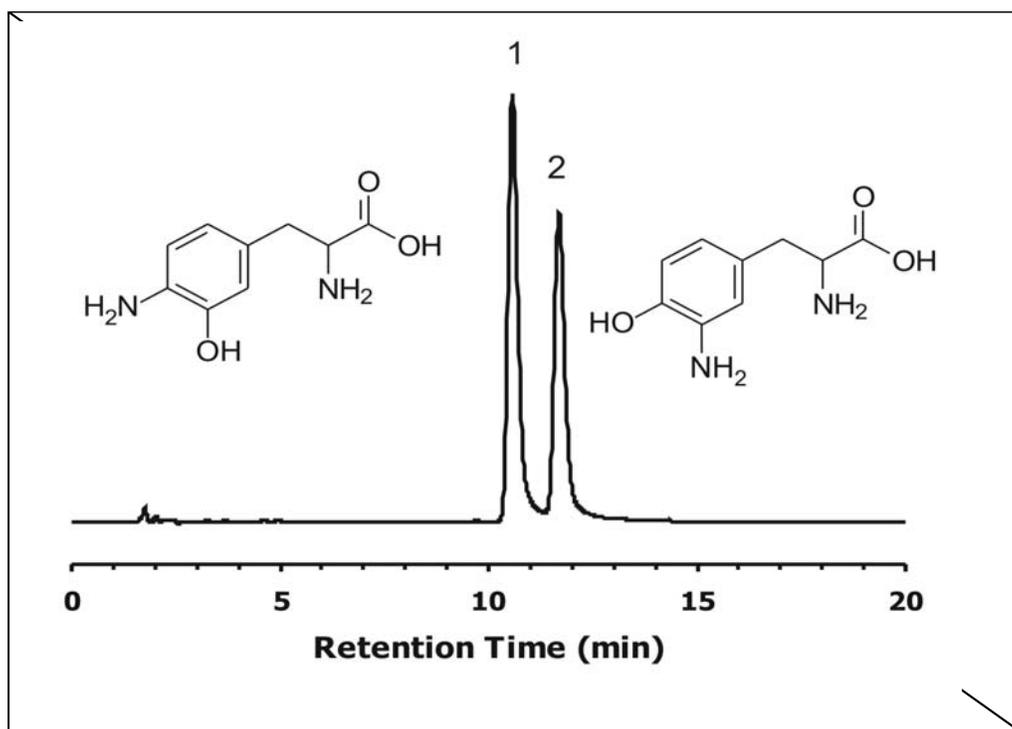


## ZIC<sup>®</sup>-HILIC Separation of 4-AHP and 3-AHP

### Chromatographic Conditions

Column: ZIC<sup>®</sup>-HILIC, PEEK 150 x 2.1 mm, 3.5  $\mu$ m, 100 Å (cat. no. 150442)  
 Injection: 5  $\mu$ L in mobile phase  
 Detection: UV @ 254 nm (UFS 1.0 V)  
 Flow Rate: 0.15 mL/min  
 Mobile Phase (v/v): 75%, Acetonitrile  
 25%, Ammonium Acetate 100 mM; pH 6.8 (total ionic strength 25 mM)



### Chromatographic Data

No.	Compound		Time (min)	Retention factor
	$t_0$	void volume	2.0	-
1	4-AHP	4-amino-3-hydroxyphenylalanine	10.6	4.3
2	3-AHP	3-amino-L-Tyrosine	11.7	4.8

A ZIC<sup>®</sup>-HILIC-ECD method for these analytes from melanin pigments can be found at: [Journal of Chromatography A, 1163 \(2007\) 70-79](#)