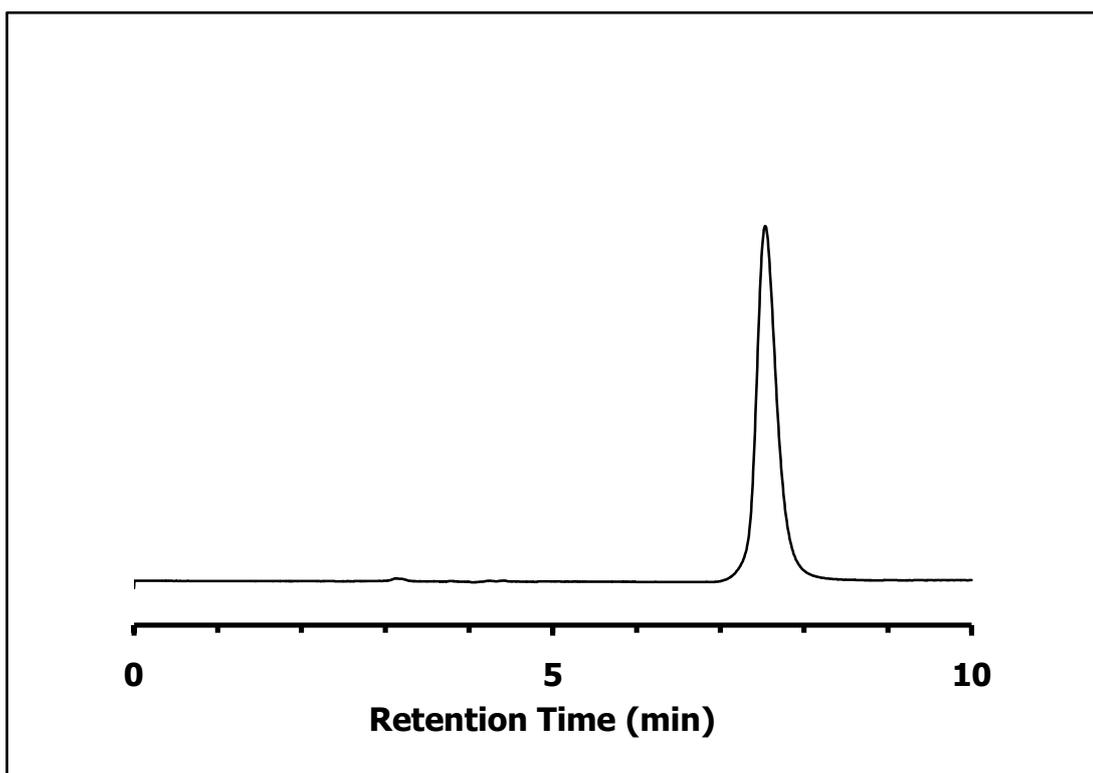


## ZIC®-HILIC Separation of Gly-His-Lys

### Chromatographic Conditions

Column: ZIC®-HILIC, PEEK 150 x 4.6 mm, 5 µm (cat. no. 150455)  
Injection: 2 µL of a 0.2 mg/ml standard solution in mobile phase  
Detection: UV @ 214 nm (UFS 1.0 V)  
Flow Rate: 0.5 mL/min  
Mobile Phase (v/v): 40%, Acetonitrile  
60%, Ammonium formate 100 mM, pH 6.2  
(overall ionic strength: 60 mM)



### Chromatographic Data

No.	Compound	Time (min)	Retention factor
	$t_0$ void volume	3,0	-
1	Gly-His-Lys	7,5	1,5