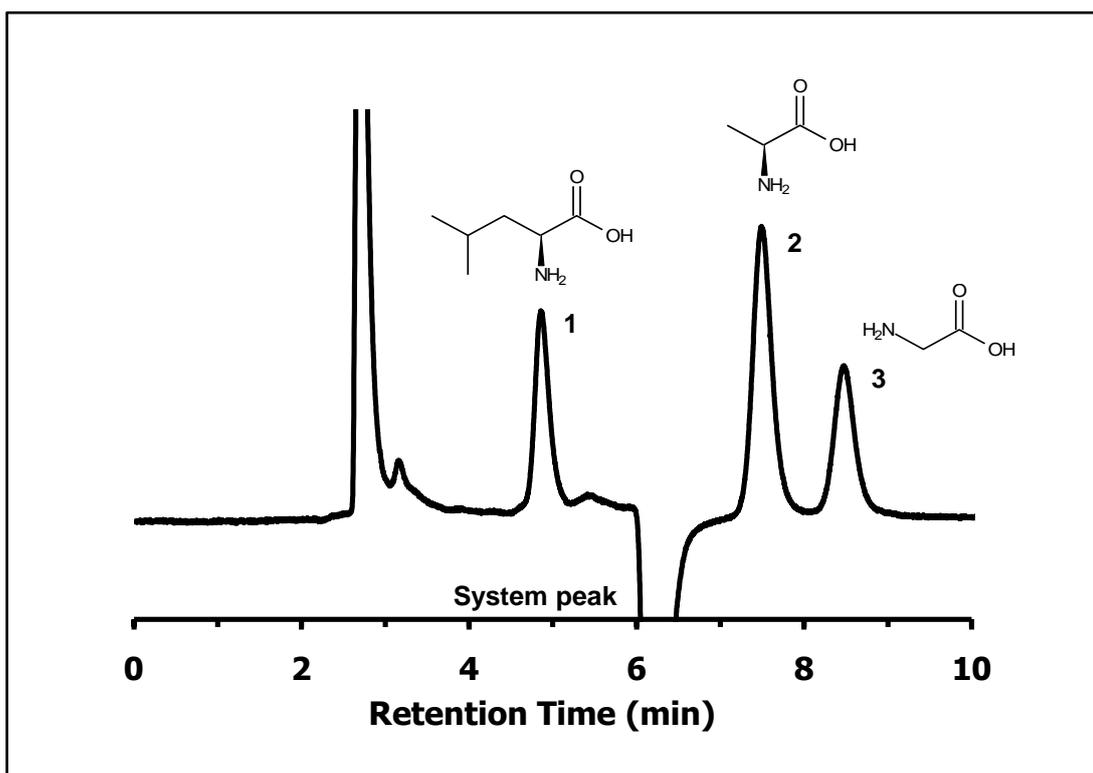


## ZIC®-pHILIC Separation of Aliphatic Amino Acids

### Chromatographic Conditions

Column: ZIC®-pHILIC, PEEK 150 x 4.6 mm, 5 µm, (cat. no. 150461)  
 Injection: 5 µL in mobile phase (with slightly lower buffer concentration)  
 Detection: UV @ 206 nm (UFS 1.0 V)  
 Flow Rate: 0.5 mL/min  
 Mobile Phase (v/v): 70%, Acetonitrile  
 30% Ammonium Formate, 20 mM, pH 6.2



### Chromatographic Data

No.	Compound	Time (min)	Retention factor
	void volume ( $t_0$ )	2,7	
1	Leucine	4,9	0,8
2	Alanin	7,5	1,8
3	Glycine	8,5	2,1