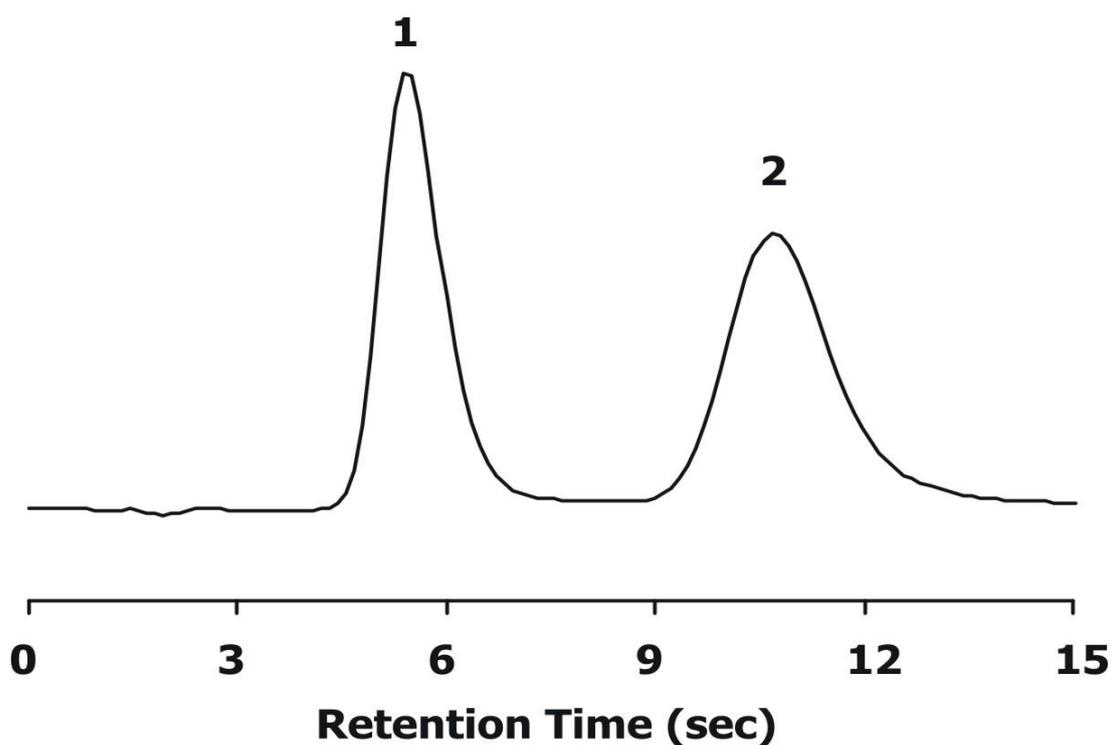


## ZIC®-HILIC Separation of AMP and ATP

### Chromatographic Conditions

Column: ZIC®-HILIC, PEEK 20 x 2.1 mm, 3.5 µm, 100 Å (cat. no. 150439)  
Injection: 750 nL in mobile phase  
Detection: UV @ 254 nm (UFS 1.0 V)  
Pressure Drop: 13.1 MPa (1886 psi)  
Flow Rate: 2.0 mL/min  
Mobile Phase (v/v): 70%, Acetonitrile  
30%, Ammonium Acetate 100 mM; pH 4.5 (total ionic strength 30 mM)



### Chromatographic Data

No.	Compound	Time (sec)	Retention factor	
1	t <sub>0</sub>	void volume	1,2	-
2	AMP	Adenosine monophosphate	5,4	3,5
3	ATP	Adenosine triphosphate	10,7	7,9