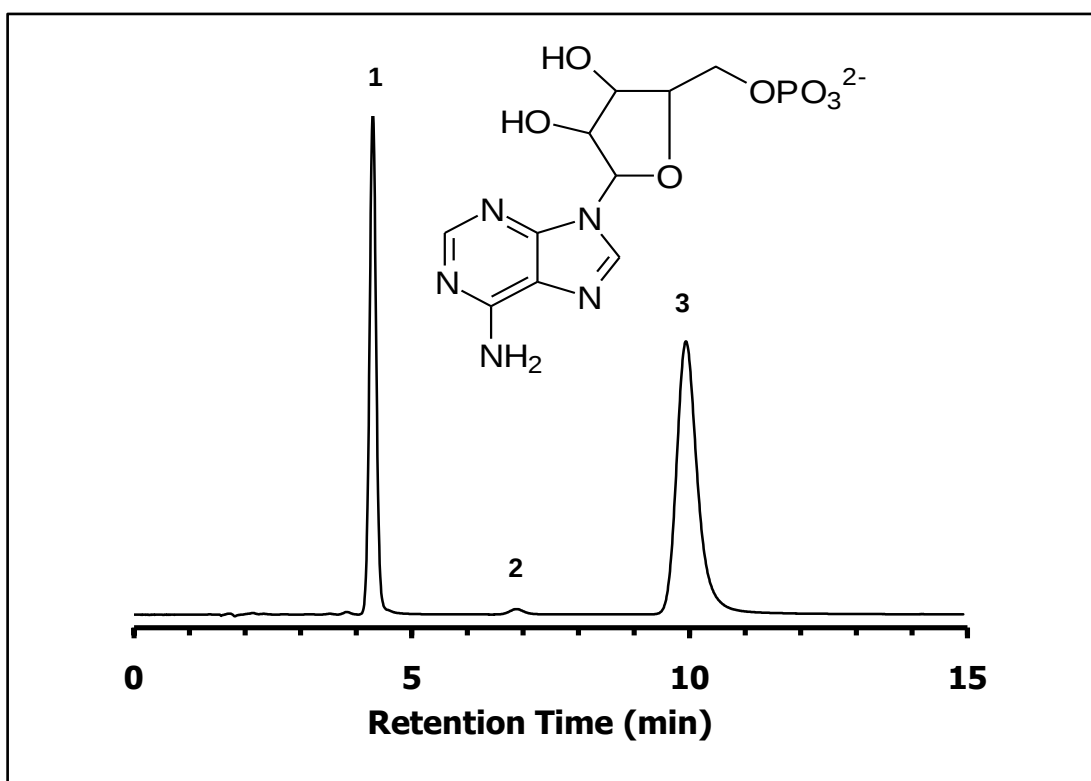


ZIC®-HILIC Separation of AMP, ADP and ATP

Chromatographic Conditions

Column: ZIC®-HILIC, PEEK 150 x 4.6 mm, 5 µm, 200 Å (cat. no. 150455)
Injection: 20 µL in mobile phase
Detection: UV @ 244 nm (UFS 1.0 V)
Pressure Drop: 6.3 MPa (913 psi)
Flow Rate: 1.0 mL/min
Mobile Phase (v/v): 70%, Acetonitrile
30%, Ammonium Acetate 100 mM; pH 4.0



Chromatographic Data

No.	Compound		Time (min)	Retention factor
	t_0	void volume	1,5	-
1	AMP	Adenosine monophosphate	4,3	1,9
2	ADP	Adenosine diphosphate	6,9	3,6
3	ATP	Adenosine triphosphate	9,9	5,6