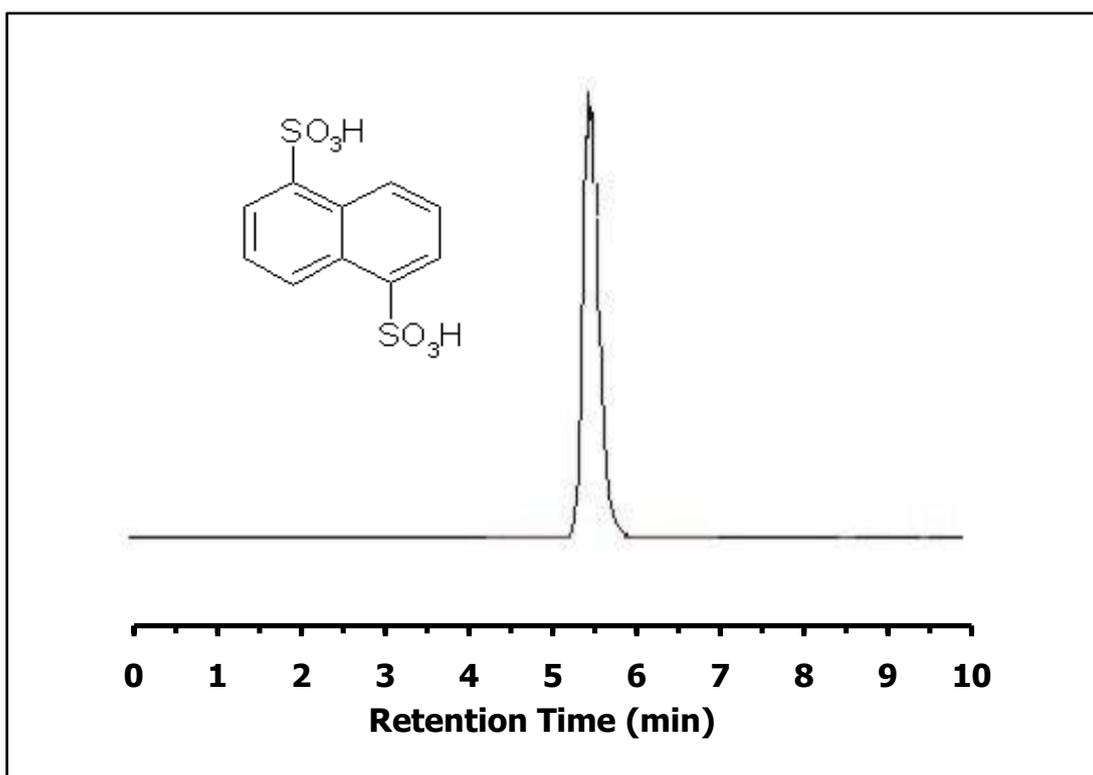


ZIC®-HILIC Separation of 1,5-naphthalenedisulphonic acid

Chromatographic Conditions

Column: ZIC®-HILIC, PEEK 150 x 4.6 mm, 5 µm, 200 Å (cat. no. 150455)
Injection: 10 µL in mobile phase
Detection: UV @ 288 nm (UFS 1.0 V)
Column Temperature: 40°C
Flow Rate: 0.5 mL/min
Mobile Phase (v/v): 80%, Acetonitrile
20%, Ammonium Acetate 10 mM, pH 6.8
(total ionic strength: 2 mM)



Chromatographic Data

No.	Compound	Time (min)	Retention factor
	t_0 void volume	3,0	-
1	1,5-naphthalenedisulphonic acid	5,5	0,6