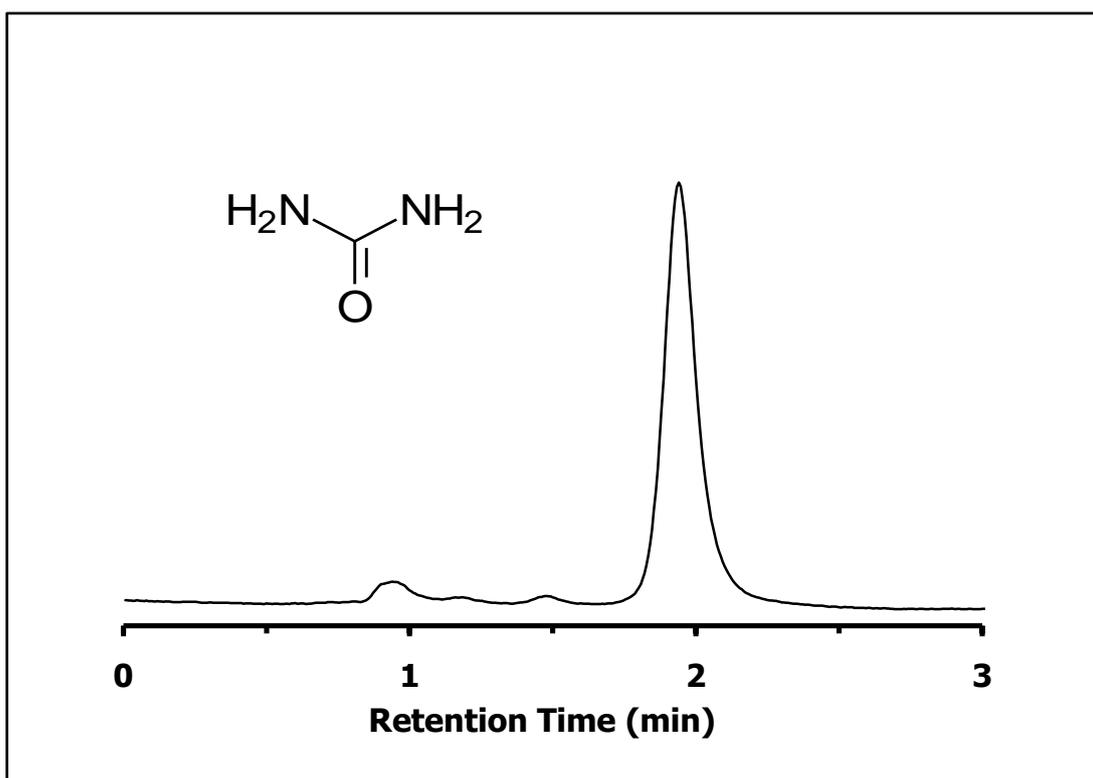


ZIC®-HILIC Separation of Urea

Chromatographic Conditions

Column: ZIC®-HILIC, PEEK 50 x 4.6 mm, 5 µm, 200 Å (cat. no. 150451)
Injection: 5 µL in mobile phase
Detection: UV @ 204 nm
Flow Rate: 0.5 mL/min
Mobile Phase (v/v): 90%, Acetonitrile
10%, Ammonium Acetate 5 mM, pH 6.8
(total ionic strength 0.5 mM)



Chromatographic Data

No.	Compound	Time (min)	Retention factor
	t_0 void volume	1,0	-
1	Urea (1.0 mg/mL)	1,9	0,9