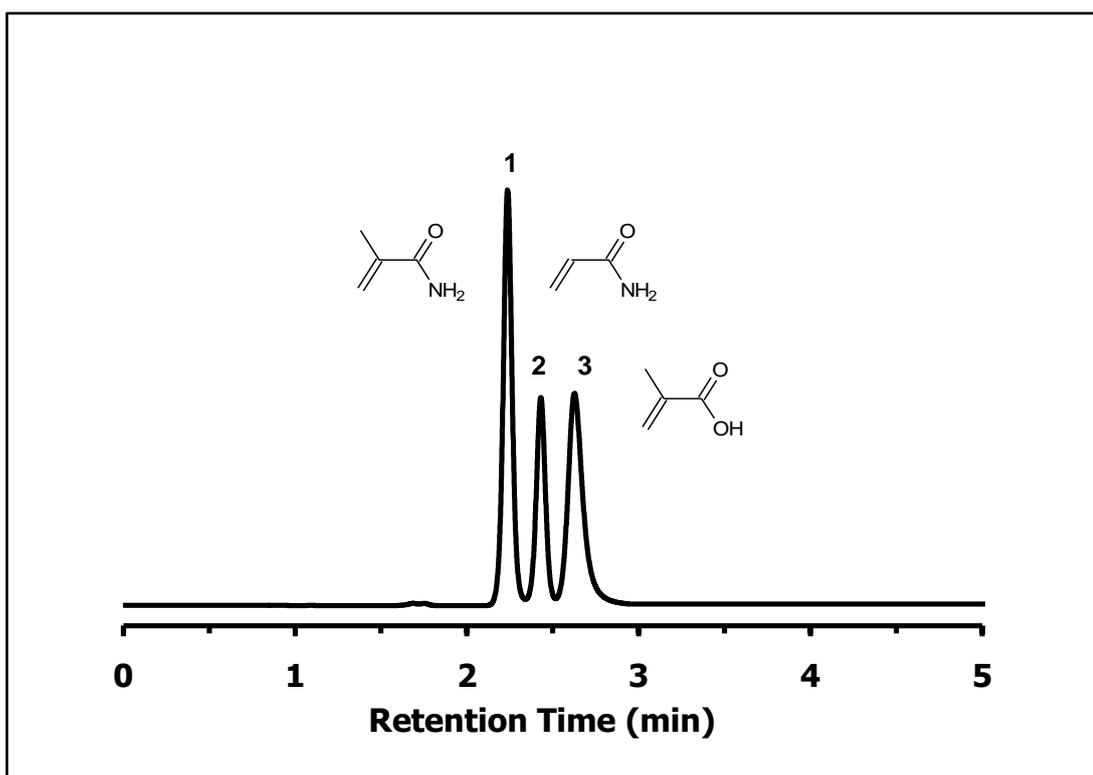


Separation of Methacrylamide, Acrylamide, Methacrylic Acid

Chromatographic Conditions

Column: ZIC®-HILIC, PEEK 150 x 4.6 mm, 5 µm, 200 Å (cat. no. 150455)
 Injection: 5.0 µL
 Detection: UV @ 210 nm (UFS 1.0 V)
 Pressure Drop: 8.0 MPa
 Flow Rate: 1.0 mL/min
 Mobile Phase (v/v): 95%, Acetonitrile
 5%, Acetic Acid 50 mM



Chromatographic Data

No.	Compound	Time (min)	Retention factor	Resolution
	void volume (t_0)	1,5	-	-
1	Methacrylamide	2,2	0,49	-
2	Acrylamide	2,4	0,62	2,00
3	Methacrylic acid	2,6	0,75	1,55