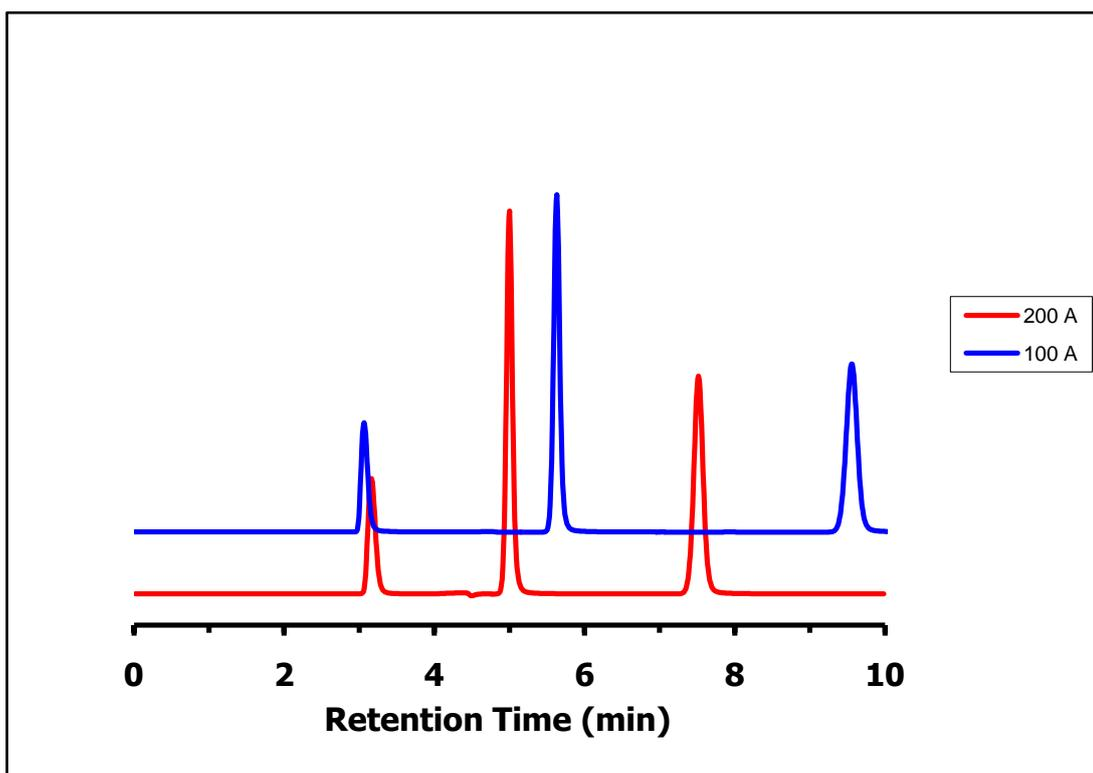


Retention Factor Vs Porosity (100 and 200 Å)

Chromatographic Conditions

Column: ZIC®-HILIC, PEEK 150 x 2.1 mm, 3.5 µm, (cat. no. 150448)
 Injection: 5 µL in mobile phase
 Detection: UV @ 254 nm (UFS 1.0 V)
 Flow Rate: 0.1 mL/min
 Mobile Phase (v/v): 80%, Acetonitrile
 20%, Ammonium Acetate 25 mM
 (total ionic strength: 5mM)



Chromatographic Data

	No.	Compound		Time (min)	Retention factor	Asymmetry
100 Å	1	Toluene	void volume (t_0)	3,1	-	-
	2	Uracil		5,6	0,88	1,10
	3	Cytosine		9,6	2,19	1,01
200 Å	1	Toluene	void volume (t_0)	3,2	-	-
	2	Uracil		5,0	0,67	1,10
	3	Cytosine		7,5	1,51	1,06