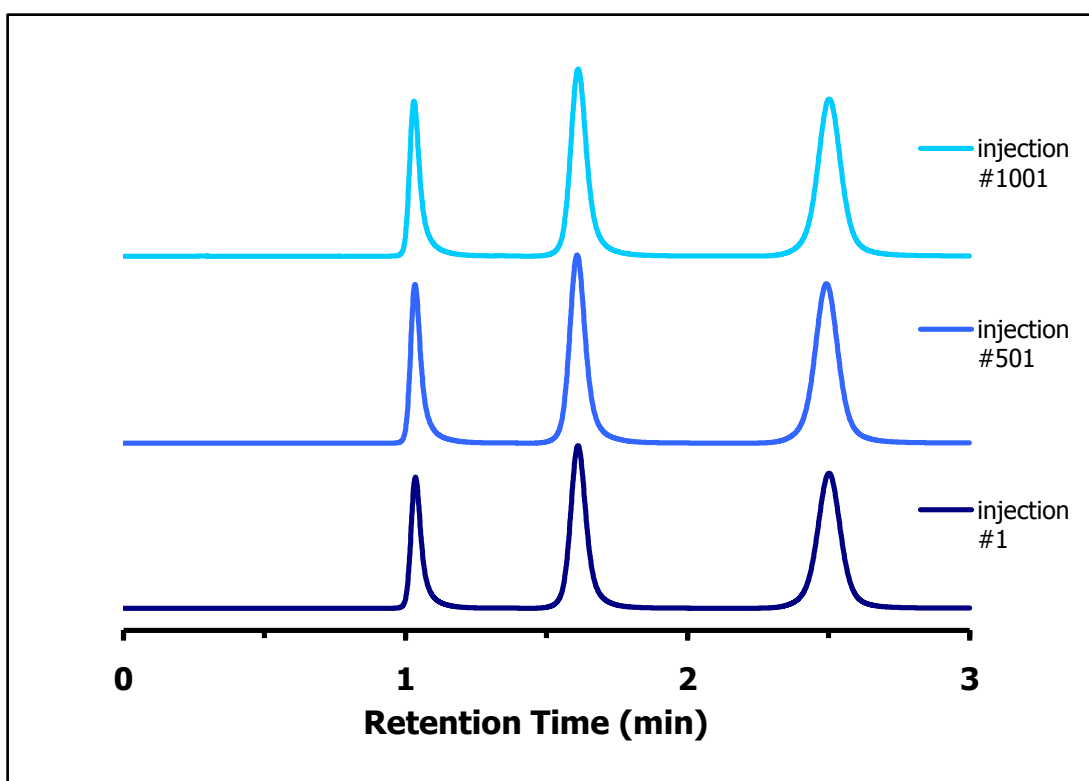


## Retention Time Stability on the ZIC®-HILIC Column

### 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 @ 254 nm (UFS 1.0 V)  
 Temperature: 24 °C  
 Pressure Drop: 1.1 MPa (158 psi)  
 Flow Rate: 0.5 mL/min  
 Mobile Phase (v/v): 80%, Acetonitrile  
 20%, Ammonium Acetate 5 mM



### Chromatographic Data

No.	Compound		Time (min)	Retention factor	Resolution
1	Toluene	void volume ( $t_0$ )	1.0	-	-
2	Uracil		1.6	0.61	6.6
3	Cytosine		2.5	1.50	6.7

The relative standard deviation for these data of injection #1, #501 and #1001 were 0.1-0.7%.