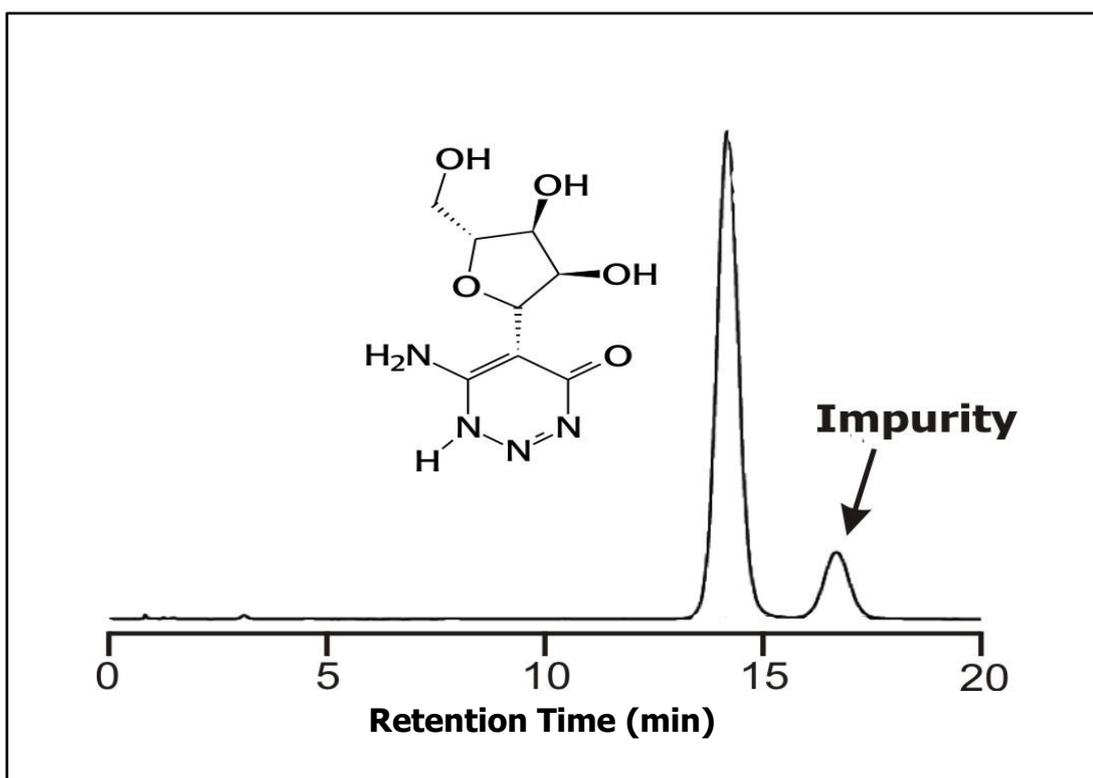


## ZIC®-HILIC Separation of Azacitidine

### Chromatographic Conditions

Column: ZIC®-HILIC, PEEK 150 x 4.6 mm, 5 µm, 200 Å (cat. no. 150455)  
 Injection: 10 µL of a 500 ppm solution diluted in mobile phase  
 Detection: UV @ 242 nm (UFS 1.0 V)  
 Column Temperature: Ambient  
 Flow Rate: 2.0 mL/min  
 Mobile Phase (v/v): 90%, Acetonitrile  
 10%, Ammonium Acetate 10 mM, pH 6.8  
 (total ionic strength: 1 mM)



### Chromatographic Data

No.	Compound	Time (min)	Retention factor
$t_0$	void volume	0,8	-
1	Azacitidine (LogP: -1.7)	14,2	25,2
2	impurity	16,7	29,6

Data provided by Gora Sharangi, Merck Specialities Pvt. Ltd, India.  
 For more details, contact [chromatography@merck.co.in](mailto:chromatography@merck.co.in)