

# Application Report 310

## Analysis of Antivirals Using Ascentis™ C8

This application demonstrates the suitability of the Ascentis C8 for the analysis of antivirals. Structures along with the optimized chromatogram obtained on the Ascentis C8 are presented.

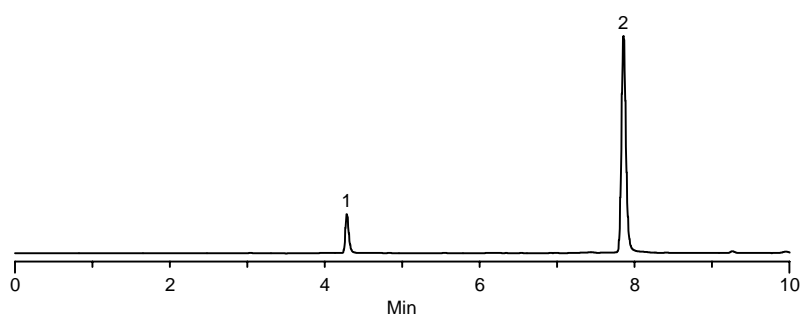
### Key Words

valacyclovir, 124832-27-5, imiquimod, 99011-02-6, Ascentis C8, 581402-U

Author: Hugh M. Cramer

Acquisition System: Shimadzu  
Barkaboom Mountain

Notebook Reference: 1558



G003201

### Conditions

column: Ascentis C8, 15 cm x 2.1 mm I.D., 3  $\mu$ m particles (581402-U)  
mobile phase A: 10 mM ammonium phosphate (pH 2.5 with phosphoric acid)  
mobile phase B: acetonitrile  
flow rate: 0.2 mL/min.  
temp.: 35 °C  
det.: UV at 254 nm  
injection: 2  $\mu$ L  
sample: as indicated in mobile phase  
gradient: 

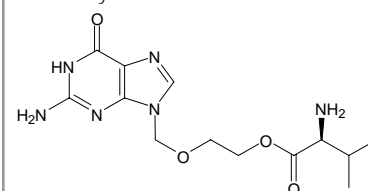
Min	%A	%B
0	95	5
10	20	80

### Peak IDs

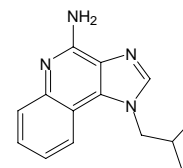
1. Valacyclovir (50  $\mu$ g/mL)
2. Imiquimod (50  $\mu$ g/mL)

### Structures

Valacyclovir - G003199



Imiquimod - G003200



Ascentis is a trademark of Sigma-Aldrich Co.