

# Application Report 311

## Analysis of Antiarrhythmics Using C8 Ascentis™ C8

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

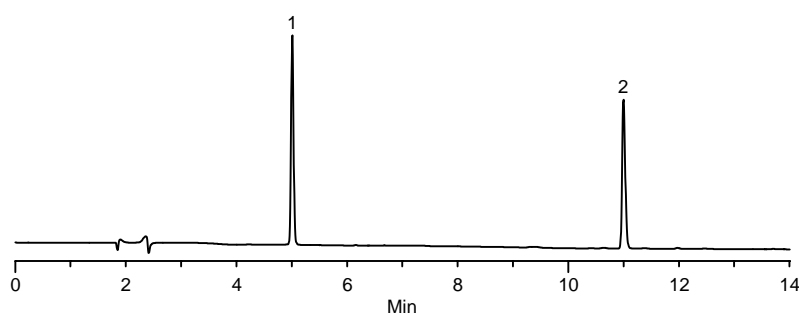
### Key Words

dofetilide, 115256-11-6, amiodarone, 1951-25-3, Ascentis C8, 581424-U, antiarrhythmics

Author: Hugh M. Cramer

Acquisition System: Hitachi Mt Fuji

Notebook Reference: 1558



G003204

### Conditions

column: Ascentis C8, 15 cm x 4.6 mm I.D., 5 µm particles (581424-U)  
mobile phase A: 25 mM ammonium phosphate (pH 3.0 with phosphoric acid)  
mobile phase B: acetonitrile  
flow rate: 1.0 mL/min.  
temp.: 35 °C  
det.: UV at 254 nm  
injection: 10 µL  
sample: as indicated in mobile phase

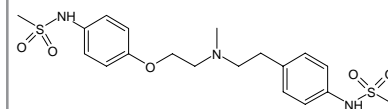
gradient:	Min	%A	%B
	0	85	15
	10	20	80
	15	20	80

### Peak IDs

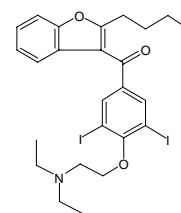
1. Dofetilide (50 µg/mL)
2. Amiodarone (50 µg/mL)

### Structures

Dofetilide - G003202



Amiodarone - G003203



Ascentis is a trademark of Sigma-Aldrich Co.