

# Application Report 101

## Analysis of Antipsychotics Using Ascentis™ C18

This application demonstrates the suitability of Ascentis C18 for the efficient separation of the antipsychotic drugs clozapine and thioridazine by HPLC.

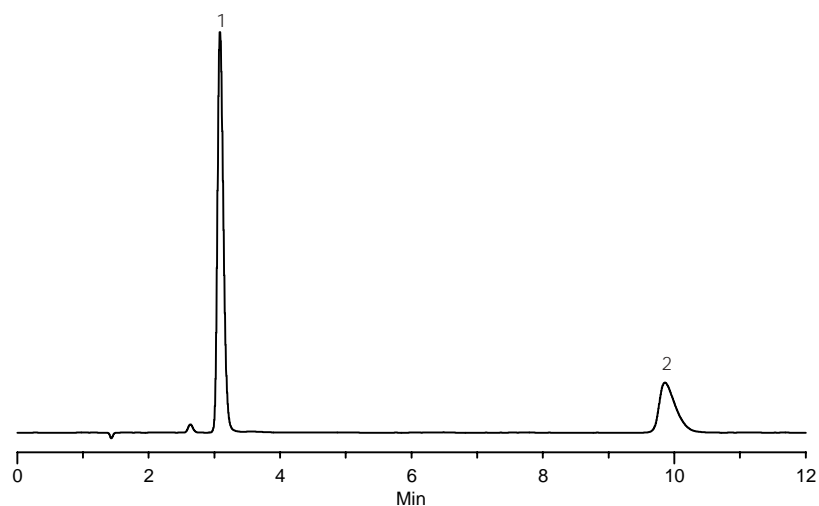
### Key Words

Ascentis C18, clozapine, 5786-21-0, C6305, thioridazine, 130-61-0, 581324-U

Author: Hugh Cramer

Acquisition System: Waters Alliance  
2690 ID 9371

Notebook Reference: 1486-86



G002400

### Conditions

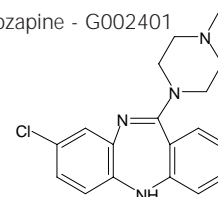
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)  
mobile phase: 40:60, 10 mM monobasic potassium phosphate (pH 3.0 with phosphoric acid):methanol  
flow rate: 1 mL/min.  
temp.: 35 °C  
det.: UV at 220 nm  
injection: 10 µL  
sample: 50 µg/mL each in mobile phase

### Peak IDs

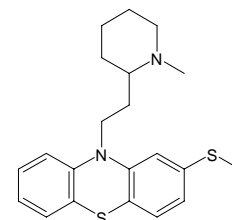
1. Clozapine
2. Thioridazine

### Structures

Clozapine - G002401



Thioridazine - G002402



Ascentis is a trademark of Sigma-Aldrich Co.