

Application Report 139

Analysis of Antipsychotics Using Ascentis™ RP-Amide

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

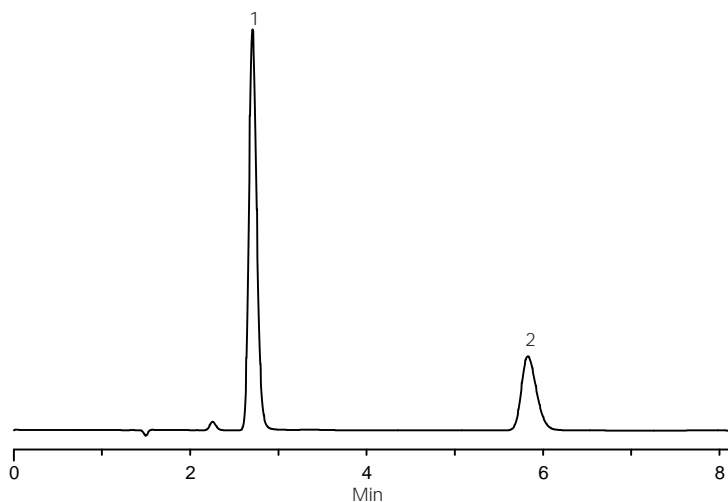
Key Words

Ascentis RP-Amide, clozapine, 5786-21-0, C6305, thioridazine, 130-61-0, 565324-U, antipsychotics

Author: Hugh Cramer

Acquisition System: Shimadzu
Mongaup

Notebook Reference: 1551



G002603

Conditions

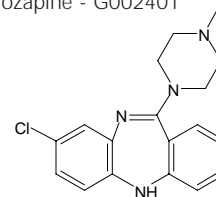
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)
mobile phase: 40:60, 10 mM monobasic potassium phosphate (pH 3.0 with phosphoric acid):methanol
flow rate: 1 mL/min.
temp.: 35 °C
det.: UV, 220 nm
injection: 10 µL
sample: 50 µg/mL each in mobile phase

Peak IDs

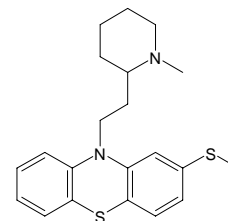
1. Clozapine
2. Thioridazine

Structures

Clozapine - G002401



Thioridazine - G002402



Ascentis is a trademark of Sigma-Aldrich Co.