

# Application Report 142

## Separation of Hydrophobic Bases Using Ascentis™ RP-Amide

This application demonstrates the suitability of Ascentis RP-Amide for the efficient separation of hydrophobic bases quinidine, diphenhydramine and fluoxetine by HPLC.

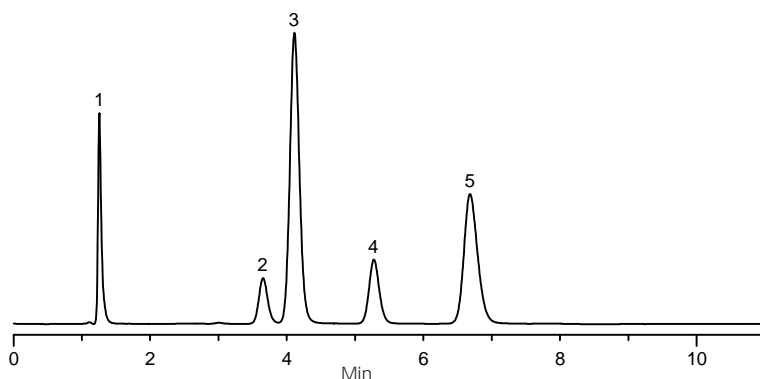
### Key Words

Ascentis RP-Amide, 565324-U, hydrophobic bases, N05158, quinidine HCl, 6151-40-2, 2-0750, diphenhydramine, 147-24-0, D3630, fluoxetine HCl, 59333-67-4, F132

Author: Jacinth A.M. McKenzie

Acquisition System: Hitachi LC

Notebook Reference: 1550-18



G002612

### Conditions

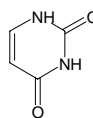
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)  
mobile phase: 35:65, 25 mM ammonium phosphate (pH 7):methanol  
temp. 35 °C  
flow rate: 1.2 mL/min.  
det.: UV, 230 nm  
injection: 10 µL  
sample: N05158

### Peak IDs

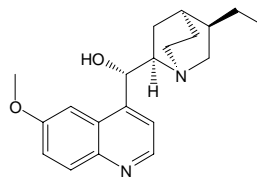
1. Uracil
2. Hydroquinidine
3. Quinidine
4. Diphenhydramine
5. Fluoxetine

### Structures

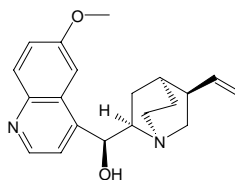
Uracil - G002739



Hydroquinidine - G002740

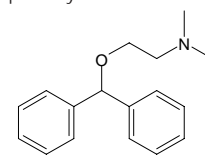


Quinidine - G002609

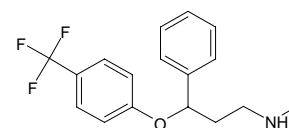


### Structures

Diphenhydramine - G002610



Fluoxetine - G002611



Ascentis is a trademark of Sigma-Aldrich Co.