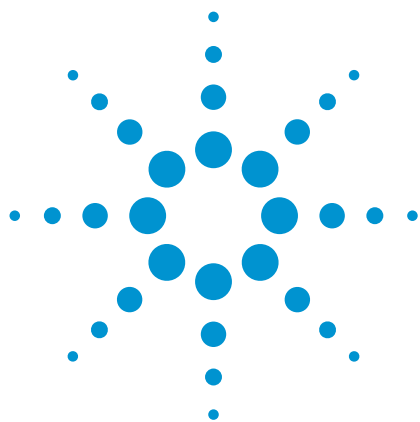




Tricyclic Antidepressants

Application Note

Pharmaceutical

Introduction

The selectivity of the Agilent Polaris C18-A column is demonstrated in the separation of eight tricyclic antidepressants (TCAs) from a KKG reference serum sample.



Agilent Technologies

TCAs according to the KKG¹ list (Psychofarmaca I and II with desmethylozapine).

Conditions

Sample	KKGT reference serum sample
Column	Agilent Polaris C18-A, 4.6 × 150 mm, 5 µm (p/n A2000150X046) (ChromSep cartridge system p/n A2000150C046)
Mobile phase	0.01 M KH ₂ PO ₄ , pH 3.0:acetonitrile, 64:36 (v/v)
Flow rate	1.3 mL/min
Injection volume	25 µL
Sample concentration	50-150 µg/L, clozapine and desmethylozapine > 200 µg/L (IS 250 µg/L)
Temperature	Ambient
Detector	UV, 254 nm

This is an example of the selectivity of tricyclic antidepressants on an Agilent Polaris C18-A column.

For More Information

These data represent typical results. For more information on our products and services, visit our Web site at www.agilent.com/chem.

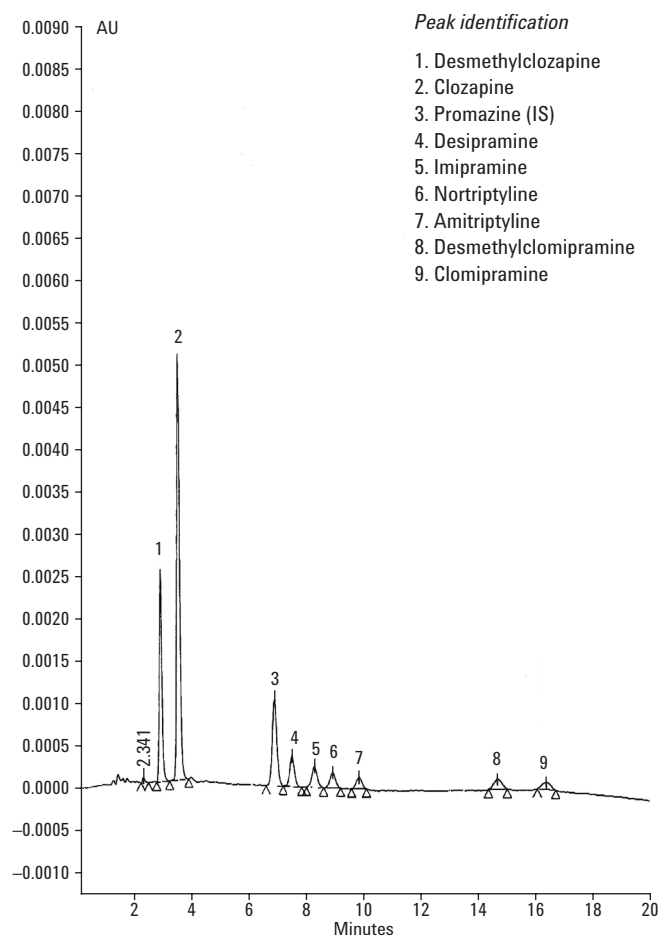


Figure 1. Tricyclic antidepressants.

www.agilent.com/chem

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice. This publication was originally published in 2009.

¹ Kwaliteitsbewaking Klinische Geneesmiddelenanalyse en Toxicologie (Association for Quality Assessment in Therapeutic Drug Monitoring and Clinical Toxicology).

