

Hydroxy-alprazolam in urine

Application Note

BioPharma

Authors

Agilent Technologies, Inc.

Introduction

The peak shape on a polymer SPE material was poor, as a coeluting substance was not removed during the washing step. This results in poor mass spectra, which were not sufficient for a library search. Using SPEC DAU and the low-bleed and high inertness of the Agilent VF-DA column, it was possible to determine the substance in Full Scan and to confirm this via library search.



Conditions

Technique : GC-capillary

Column : Agilent VF-DA, 0.20 mm x 12 m (df =optimized)

(Part no. CP8964)

Temperature : 70 °C, 1.2 min \rightarrow 200 °C, 20 °C/min, \rightarrow 270 °C,

7 °C/min, \rightarrow 320 °C, 20 °C/min

Carrier Gas : Helium, ca. 1.0 mL/min

Pressure Program : 58.7 kPa, 2.2 min, \rightarrow 97 kPa, 58 kPa/min \rightarrow

132 kPa, 3 kPa/min → 180 kPa, 12 kPa/min

Injector : Splitless

Detector : MS in SIM mode

Sample Size : 1 µL

Concentration Range : gram per μ L Solvent : methanol

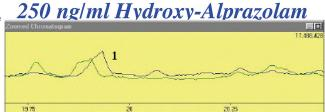
Courtesy : Joerg Szigan - Labor

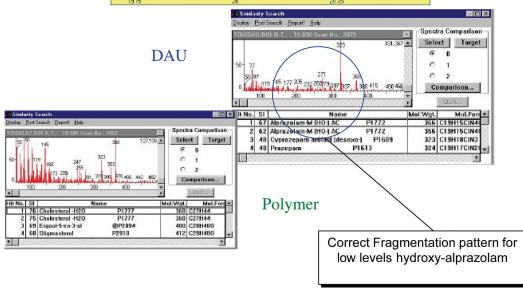
Dr. lembke, Dr. Lempfrid - Cologne

Peak identification

1. hydroxy-aiprazolam

DAU Polymer





www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2011

Printed in the USA
31 October, 2011

First published prior to 11 May, 2010

A01982

