

Application Report 04

Analysis of Amphetamines on Discovery C18

This application demonstrates the suitability of Discovery C18 for the analysis of amphetamines. Structures of Amphetamine, Methamphetamine, MDA and MDMA are shown. The optimized chromatogram obtained on Discovery C18 is presented in below.

Key Words

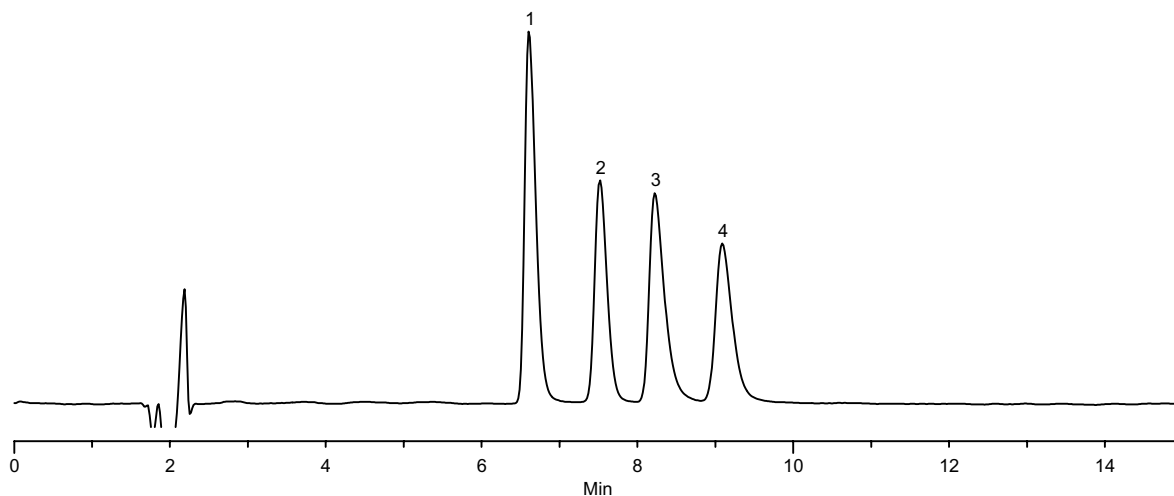
Amphetamine, Methamphetamine, MDA, 3,4-methylenedioxyamphetamine, MDMA, 3,4-methylenedioxymethamphetamine, Ecstasy, 504955

Author: Shawn R. Wyatt

Raw Data File Name:
Project "s_wyatt", Result ID #1623

Acquisition System:
Lab 35, Waters Alliance

Notebook Reference: 1445-22



G002031

Conditions

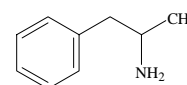
Column: Discovery C18, 15cm x 4.6mm ID, 5µm
Cat. No.: 504955
Mobile Phase: 10mM Ammonium Acetate in 90:10 Water:Acetonitrile, v/v
Flow Rate: 1.0mL/min
Temperature: 35°C
Detection: UV, 210nm
Injection Volume: 5µL
Sample: 50µg/mL each (Amphetamine, Methamphetamine, MDA and MDMA) in methanol

Peak IDs

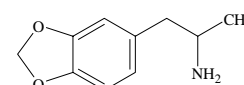
1. Amphetamine
2. MDA
3. Methamphetamine
4. MDMA

Structures

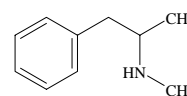
Amphetamine - G002037



MDA - G002038



Methamphetamine - G002035



MDMA - G002036

