

Application Report 31

Analysis of Methyl Paraben, Propyl Paraben and Potassium Sorbate on Discovery HS PEG

This application demonstrates the suitability of Discovery HS PEG for the analysis of methyl paraben, propyl paraben and potassium sorbate. Structures along with the optimized chromatogram obtained on Discovery HS PEG are presented below.

Key Words

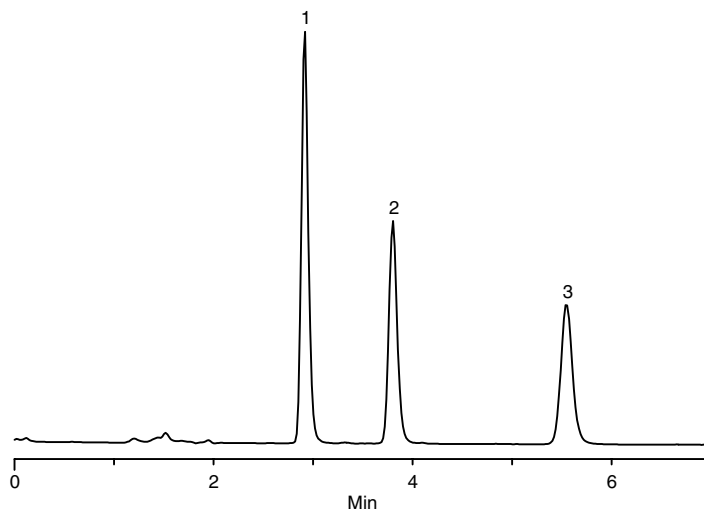
methyl paraben, p-hydroxybenzoic acid methyl ester, H6654, 99-76-3, propyl paraben, p-hydroxybenzoic acid propyl ester, n-propyl p-hydroxybenzoate, P5835, 94-13-3, potassium sorbate, S7420, 590-00-1, Discovery HS PEG, 567416-U

Author: Shawn R. Wyatt

Raw Data File Name: Project "s_wyatt", Result ID #4750

Acquisition System: Lab 35, Waters Alliance

Notebook Reference: 1445-39



G002083

Conditions

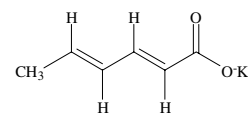
Column: Discovery HS PEG, 15cm x 4.6mm ID, 5µm
Cat. No.: 567416-U
Mobile Phase: (70:30) 10mM citric acid, pH 2.5:CH₃CN, v/v
Temperature: 35°C
Flow Rate: 1.0mL/min
Detection: UV, 254nm
Injection Volume: 5µL
Sample: 25µg/mL each (methyl paraben, propyl paraben and potassium sorbate) in 10mM citric acid, pH 2.5

Peak IDs

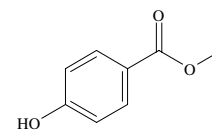
1. Potassium sorbate
2. Methyl paraben
3. Propyl paraben

Structures

Potassium sorbate - G002082



Methyl paraben - G000194



Propyl paraben - G000196

