

Application Report 05

Separation of Benzoic Acid, Benzyl Alcohol and Benzaldehyde on Discovery HS C18

This application demonstrates the efficient separation of the title benzyl analogs on Discovery HS C18.

Key Words

Benzoic Acid, Benzaldehyde, Benzyl Alcohol, Discovery HS C18, 569250-U, 48323, 442481

Author: David S. Bell

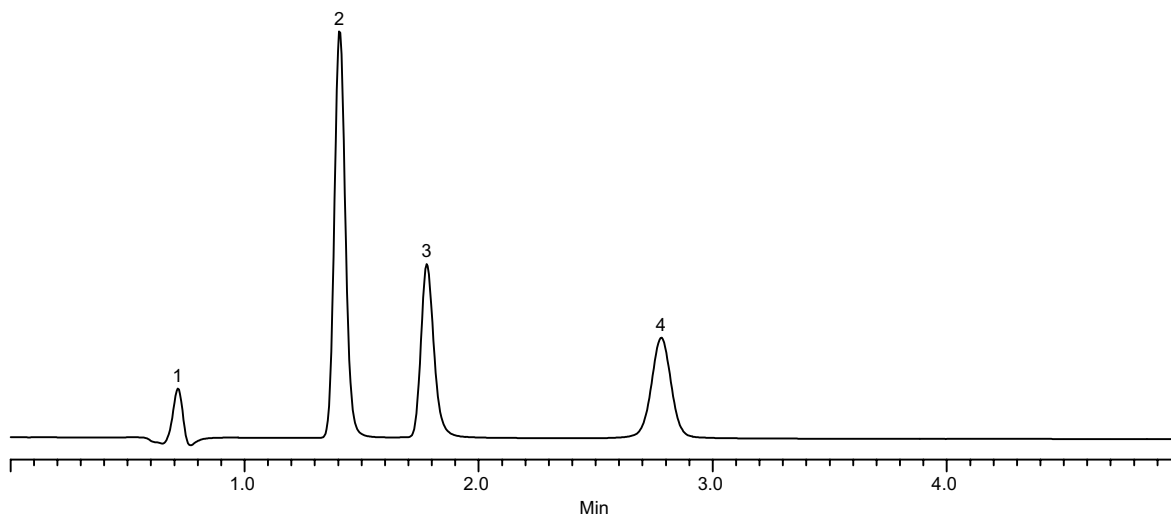
Raw Data File Name:

1462/13000013

Acquisition System:

SPE/HP1100/System2

Notebook Reference: 1462-13



G002032

Conditions

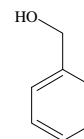
Column: Discovery HS C18, 5cm x 4.6mm ID, 3 μ m
Cat. No.: 569250-U
Mobile Phase: 10mM monobasic potassium phosphate,
pH to 3.0 w/ H₃PO₄:CH₃CN (70:30, v/v)
Flow Rate: 1mL/min
Temperature: 35°C
Detection: UV, 210nm
Injection Volume: 10 μ L
Sample: 10 μ g/mL benzoic acid, 15 μ g/mL each benzyl alcohol
and benzaldehyde

Peak IDs

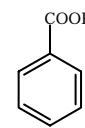
1. Void marker
2. Benzyl alcohol
3. Benzoic acid
4. Benzaldehyde

Structures

Benzyl alcohol - G002033



Benzoic acid - G001261



Benzaldehyde - G002034

