

# Application Report 29

## Analysis of Conjugated Flavonoids on Discovery HS C18

This application demonstrates the suitability of Discovery HS C18 for the analysis of disaccharide-conjugated flavonoids. Maysin, apimaysin and methoxymaysin structures along with the optimized chromatogram obtained on Discovery HS C18 are presented below.

### Key Words

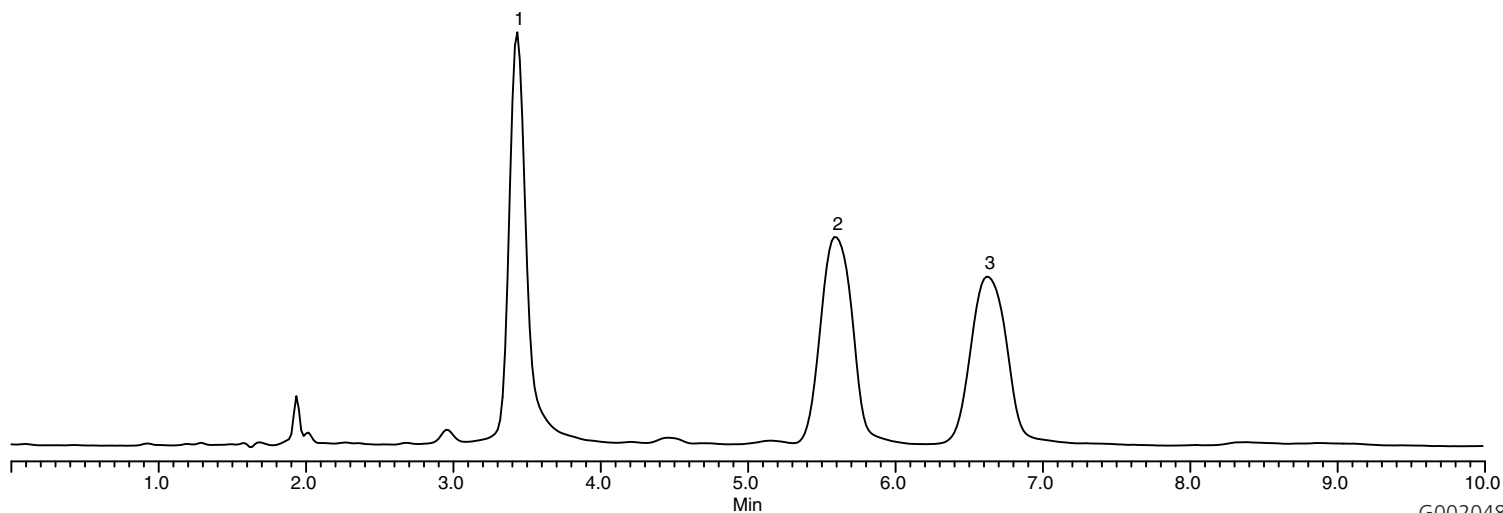
Maysin, Apimaysin, Methoxymaysin, Flavonoid, Flavone, Disaccharide, Aflatoxin, Maize, Corn, Cornsilk, Earworm, 569252-U, Discovery HS C18, Conjugated

**Author:** Shawn R. Wyatt

**Raw Data File Name:** Project "s\_wyatt", Result ID #4062

**Acquisition System:** Lab 35, Waters Alliance

**Notebook Reference:** 1445-29



G002048

### Conditions

Column: Discovery HS C18, 15cm x 4.6mm ID, 3µm  
Cat. No.: 569252-U  
Mobile Phase: (80:20) 10mM sodium citrate (pH 7.0 with 10mM citric acid):CH<sub>3</sub>CN, v/v  
Temperature: 60°C  
Flow Rate: 1.0mL/min  
Detection: UV, 340nm  
Injection Volume: 10µL  
Sample: 50µg/mL each (Maysin, Apimaysin, Methoxymaysin) in 10mM sodium citrate (pH 7.0 with 10mM citric acid)

### Peak IDs

1. Maysin
2. Apimaysin
3. Methoxymaysin

### Structures

