

# Application Report 127

## Analysis of Xanthines Using Ascentis™ C18

This application demonstrates the suitability of Ascentis C18 for the efficient separation of the xanthines 1-methylxanthine, 7-methylxanthine, 1,7-dimethylxanthine, and  $\beta$ -hydroxytheophylline by HPLC.

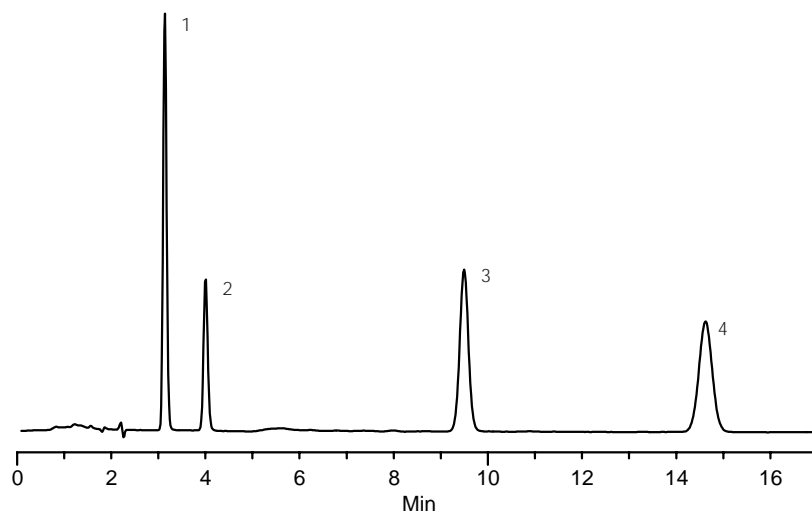
### Key Words

Ascentis C18, 1-methylxanthine, M3275, 6136-37-4, 7-methylxanthine, M9890, 552-62-5, 1,7-dimethylxanthine, D5385, 611-59-6,  $\beta$ -hydroxytheophylline, H9006, 519-37-9, 581324-U

Author: Hugh Cramer

Acquisition System: Waters Alliance  
2690 ID 9371

Notebook Reference: 1486-87



G002465

### Conditions

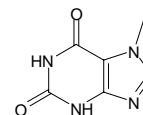
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5  $\mu$ m particles (581324-U)  
mobile phase: 94:6, 10 mM monobasic potassium phosphate (pH 3.0 with phosphoric acid):methanol  
flow rate: 1 mL/min.  
temp.: 35 °C  
det.: UV at 220 nm  
injection: 10  $\mu$ L  
sample: 50  $\mu$ g/mL each in mobile phase

### Peak IDs

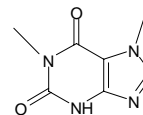
1. 7-methylxanthine
2. 1-methylxanthine
3. 1,7-dimethylxanthine
4.  $\beta$ -hydroxytheophylline

### Structures

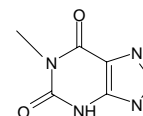
7-methylxanthine - G002466



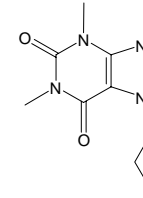
1-methylxanthine - G002467



1,7-dimethylxanthine - G002468



$\beta$ -hydroxytheophylline - G002469



Ascentis is a trademark of Sigma-Aldrich Co.