

# Application Report 206

## Analysis of Parabens Using Ascentis™ C18

This application demonstrates the suitability of the Ascentis C18 for the analysis of parabens.

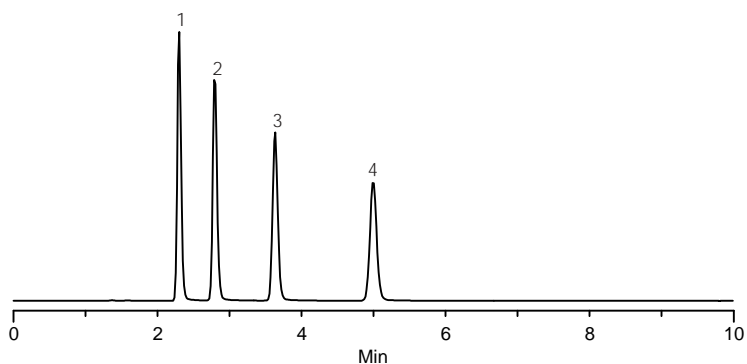
### Key Words

methyl paraben, 99-76-3, 47889, ethyl paraben, 120-47-8, 111988, propyl paraben, 94-13-3, P5835, butyl paraben, 94-26-8, H9503, Ascentis C18, 581324-U

**Author:** Hugh M. Cramer

**Acquisition System:** Cinnabar  
Waters 2690

**Notebook Reference:** 1551



G002849

### Conditions

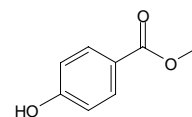
column: Ascentis C18, 15 cm x 4.6 mm I.D., 5 µm particles (581324-U)  
mobile phase: 45:55, water:acetonitrile  
flow rate: 1.0 mL/min.  
temp.: 35 °C  
det.: UV at 254 nm  
injection: 10 µL  
sample: as indicated in mobile phase

### Peak IDs

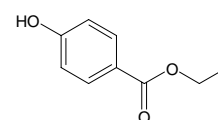
1. Methyl paraben (50 µg/mL)
2. Ethyl paraben (50 µg/mL)
3. Propyl paraben (50 µg/mL)
4. Butyl paraben (50 µg/mL)

### Structures

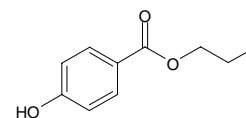
Methyl paraben - G000194



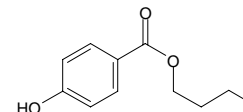
Ethyl paraben - G002626



Propyl paraben - G000196



Butyl paraben - G002848



Ascentis is a trademark of Sigma-Aldrich Co.