

# Application Report 08

## Furosemide and Indapamide by HPLC

This application demonstrates the suitability of Discovery C18 for the efficient separation of Lasix (furosemide) and indapamide by HPLC.

### Key Words

Discovery C18, Lasix, furosemide, indapamide, 504955

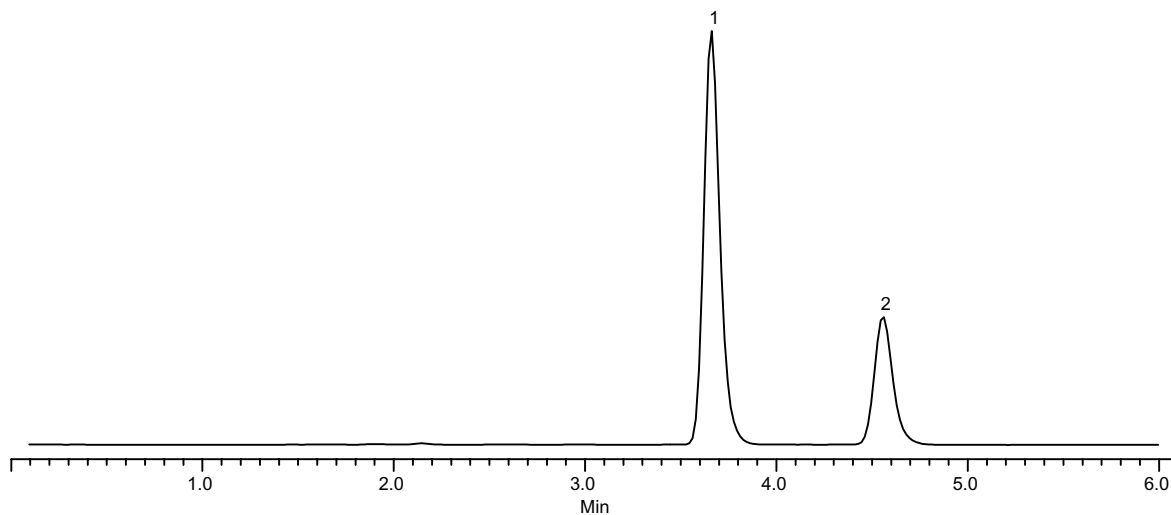
**Author:** Hugh Cramer

**Raw Data File Name:**

\\User\Public\Application  
Group\application reports\Application  
Report Furosemide and  
Indapamide.doc

**Acquisition System:** Waters Alliance  
2690 ID 9371

**Notebook Reference:** 1464-26



G002044

### Conditions

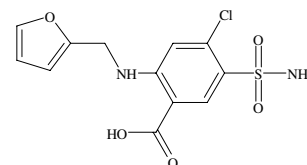
Column: Discovery C18, 15cm x 4.6mm, 5 $\mu$ m particles  
Cat. No.: 504955  
Mobile Phase: 40:60, 10mM monobasic potassium phosphate  
(pH 3.0 with H<sub>3</sub>PO<sub>4</sub>): CH<sub>3</sub>CN  
Flow Rate: 1mL/min  
Temp.: 35°C  
Det.: UV at 234nm  
Inj.: 10 $\mu$ L  
Sample: 10 $\mu$ g/mL each in mobile phase

### Peak IDs

1. Furosemide  
2. Indapamide

### Structures

Furosemide - G002045



Indapamide - G002046

