

Application Report 242

Analysis of Sulfonylurea Urea Pesticides Using Ascentis™ RP-Amide

This application demonstrates the suitability of the Ascentis RP-Amide for the analysis of sulfonylurea pesticides. Structures along with the optimized chromatogram obtained on the Ascentis RP-Amide are presented.

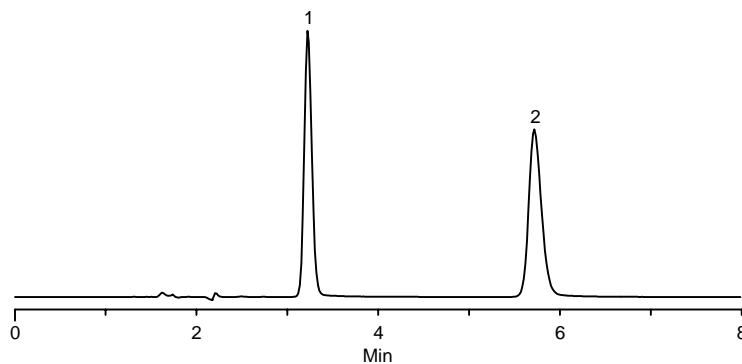
Key Words

nicosulfuron, 111991-09-4, chlorsulfuron, 64902-72-3, 34322, PS1065, Ascentis RP-Amide, 565324-U

Author: Hugh M. Cramer

Acquisition System: Cinnabar
Waters 2690

Notebook Reference: 1551



G002939

Conditions

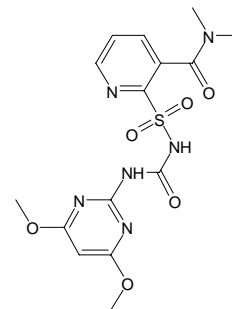
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)
mobile phase: 85:15, 10 mM potassium phosphate dibasic (pH 7.0 with phosphoric acid): acetonitrile
flow rate: 1.0 mL/min.
temp.: 35 °C
det.: UV at 254 nm
injection: 10 µL
sample: as indicated in 90:10 water:methanol

Peak IDs

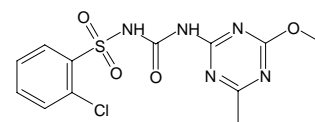
1. Nicosulfuron (50 µg/mL)
2. Chlorsulfuron (50 µg/mL)

Structures

Nicosulfuron - G002936



Chlorsulfuron - G002937



Ascentis is a trademark of Sigma-Aldrich Co.