

Application Report 169

Analysis of Bronchiodilator Terbutaline Using Ascentis™ RP-Amide

This application demonstrates the suitability of Ascentis RP-Amide for the efficient elution of terbutaline by HPLC.

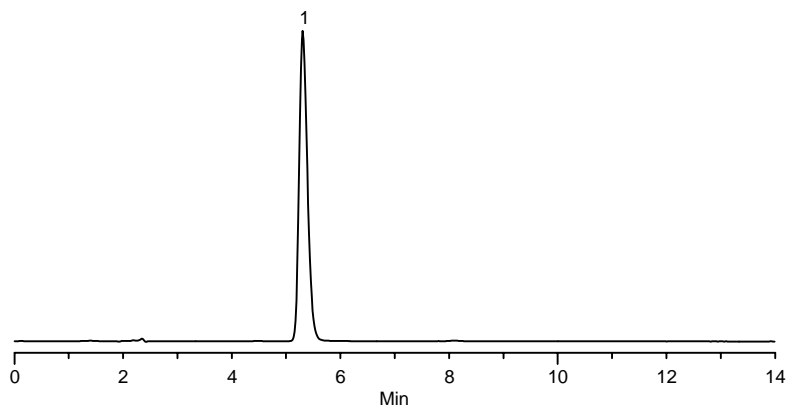
Key Words

Ascentis RP-Amide, terbutaline, 23031-32-5, T2528, 565324-U, bronchiodilator

Author: Hugh Cramer

Acquisition System: Waters Alliance
2690 Mt Hoyt

Notebook Reference: 1551



G002664

Conditions

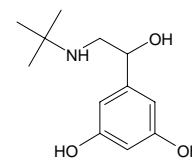
column: Ascentis RP-Amide, 15 cm x 4.6 mm I.D., 5 µm particles (565324-U)
mobile phase: 95:5, 10 mM monobasic potassium phosphate (pH 3.0 with phosphoric acid):acetonitrile
flow rate: 1 mL/min.
temp.: 35 °C
det.: UV, 220 nm
injection: 10 µL
sample: 50 µg/mL in mobile phase

Peak IDs

1. Terbutaline

Structures

Terbutaline - G002406



Ascentis is a trademark of Sigma-Aldrich Co.