

Application Report 128

Analysis of Calphostin C Using Ascentis™ C18

Calphostin C, naturally occurring perylenequinone antibiotic, is resolved from several contaminants, using Ascentis C18.

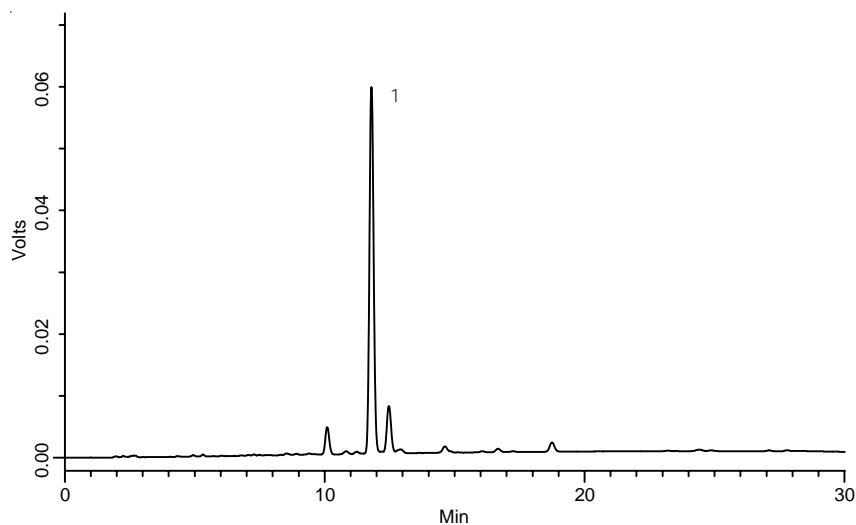
Key Words

Calphostin C, C6303, 121263-19-2, antibiotic, natural product; Ascentis C18, 581305-U

Author: Hillel Brandes

Acquisition System: Shimadzu LC-10AD System

Notebook Reference: #1396, p.72



G002470

Conditions

column: Ascentis C18, 25 cm x 2.1 mm I.D., 5 µm particles (581305-U)
mobile phase A: 50:50, 0.1% phosphoric acid in water:0.1% phosphoric acid in acetonitrile
mobile phase B: 0.1% phosphoric acid in acetonitrile
flow rate: 0.2 mL/min.
temp.: 24 °C
det.: UV at 254 nm
injection: 3 µL
sample: calphostin C 0.25 g/L in 40% methanol, 0.025% TFA

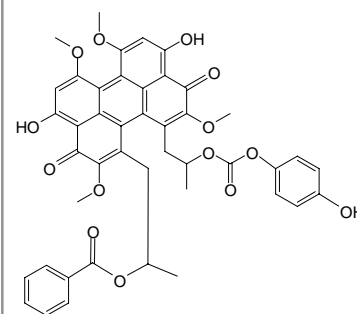
gradient:	Min	%A	%B
	0	70	30
	25	20	80
	30	20	80

Peak IDs

1. Calphostin C

Structures

Calphostin C - G002471



Ascentis is a trademark of Sigma-Aldrich Co.