

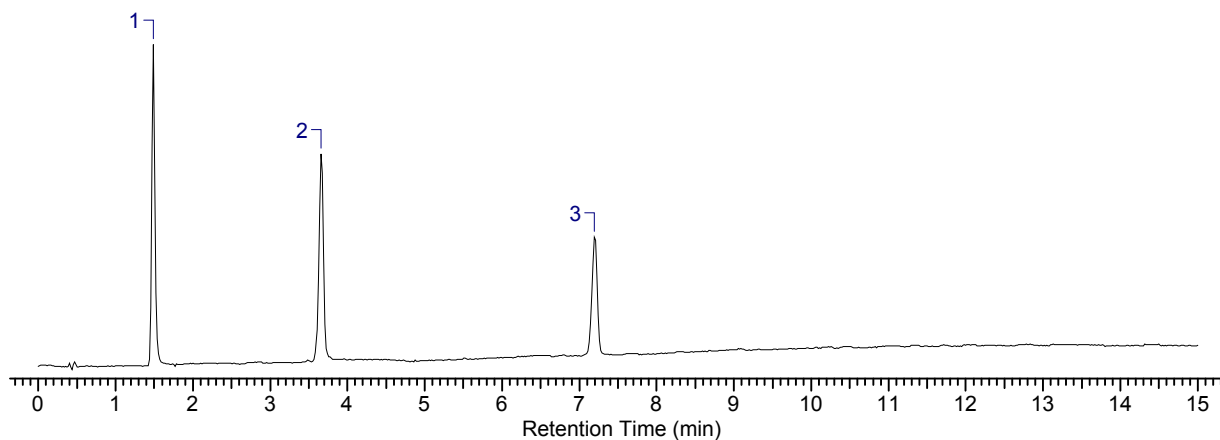
Application Report

Analysis of Bisphenol A and Related Compounds on Ascentis Express RP-Amide with UV Detection

Bonded Phase alkylamide
Column Name Ascentis Express RP-Amide
Instrument Name Waters 2690

Key Words

bisphenol A, dimethylbisphenol A, nonylphenol, 104-40-5, 80-05-7, 53922-U



Conditions

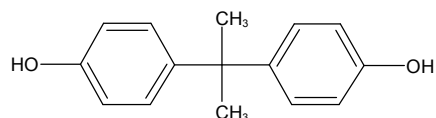
column: Ascentis Express RP-Amide, 5 cm x 4.6 mm I.D.
2.7 particles (53922-U)
mobile phase A: water
mobile phase B: acetonitrile
flow rate: 1.0 mL/min.
temp.: 35 °C
det.: UV 28 nm
injection: 10µL
sample: 50 mg/L ea. in 50:50, water: acetonitrile

gradient:	min	% A	% B
	0	50	50
	10	0	100

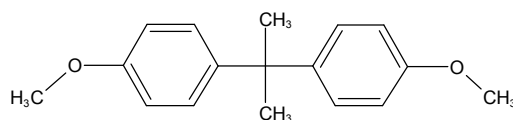
Peak IDs

No.	Name	tR
1	Bisphenol A	1.5
2	Dimethyl Bisphenol A	3.7
3	Nonylphenol	7.2

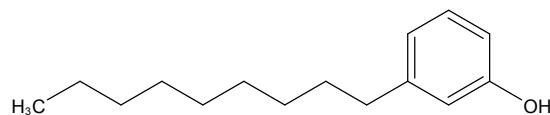
Structures



Bisphenol A



Dimethyl bisphenol A



Nonylphenol