

Liquid Chromatography

Epic Diphenyl Columns: Novel Aromatic Selectivity



The Epic™ Diphenyl bonded phase, with its unique selectivity and reduced hydrophobicity, exhibits increased speed and resolution for aromatic containing compounds. It utilizes dipole-dipole, hydrogen bonding, and π - π mechanisms for alternative selectivity of compounds containing double bonds or aromatic functional groups. In addition, the diphenyl arrangement of the phase can contribute to steric selectivity allowing for an additional chromatographic interaction. Epic Diphenyl may be used as a USP L11 column. Epic Diphenyl utilizes our proprietary high-density bonding technology enabling superior performance, durability and enhanced lot-to-lot reproducibility for each column.

As with all ES Industries phases, the Epic Diphenyl is available in both analytical and preparative media for maximum productivity and efficiency.

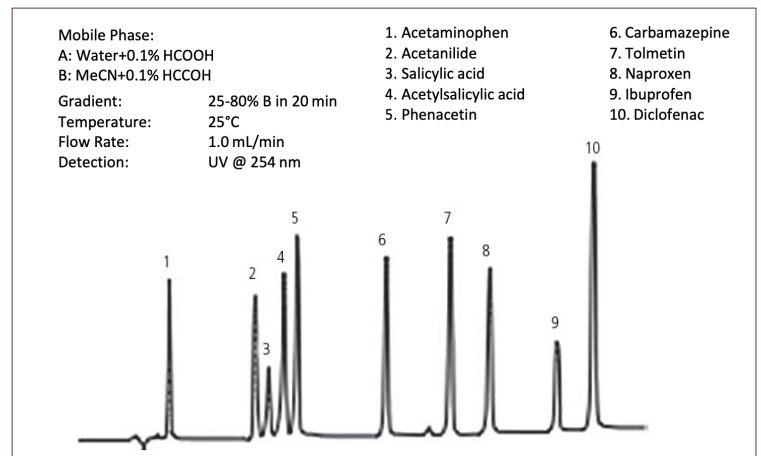


Figure 1. HPLC analysis of polar aromatic pharmaceutical compounds using Epic Diphenyl, 150 x 4.6 mm, 5 μ m.

Phase	Length (mm)	ID (mm)	Particle Size (µm)	Part No.
Epic Diphenyl	50	2.1	1.8	512A91-EDPH
Epic Diphenyl	50	2.1	3	112191-EDPH
Epic Diphenyl	50	2.1	5	112291-EDPH
Epic Diphenyl	100	2.1	1.8	522A91-EDPH
Epic Diphenyl	100	2.1	3	122191-EDPH
Epic Diphenyl	100	2.1	5	122291-EDPH
Epic Diphenyl	100	3.0	1.8	513A91-EDPH
Epic Diphenyl	100	4.6	3	125191-EDPH
Epic Diphenyl	150	2.1	3	132191-EDPH
Epic Diphenyl	150	2.1	5	132291-EDPH
Epic Diphenyl	150	4.6	3	135191-EDPH
Epic Diphenyl	150	4.6	5	135291-EDPH
Epic Diphenyl	250	4.6	5	155291-EDPH
Epic Diphenyl Prep	250	20	5	158291-EDPH
Epic Diphenyl Prep	250	30	5	15N291-EDPH
Epic Diphenyl Analytical Guard Cartridges (Pkg. 5)	10	3.0	5	500101-EDPH
Analytical Guard Cartridge Holder with integrated coupler	–	–	–	ES500100

Other column dimensions, particle sizes, and guard cartridges are available. Please enquire for more details at LCA.TechSupport@perkinelmer.com

PerkinElmer, Inc.
 940 Winter Street
 Waltham, MA 02451 USA
 P: (800) 762-4000 or
 (+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2021, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.