

Liquid Chromatography

Epic Polar HPLC Columns: Engineered for Polar Analytes



Epic™ Polar is a high density C18 packing specifically engineered for the retention of polar analytes such as small organic acids, water-soluble vitamins, catecholamines and other polar compounds. The retention characteristics of Epic Polar can be seen in the separation of 11 highly polar organic acids shown in Figure 1. This phase provides for superior hydrophobic performance even with 100% aqueous mobile phases. Under these highly aqueous conditions, our novel proprietary bonding chemistry allows the bonded chains to remain fully extended in the mobile phase, with no phase collapse.

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

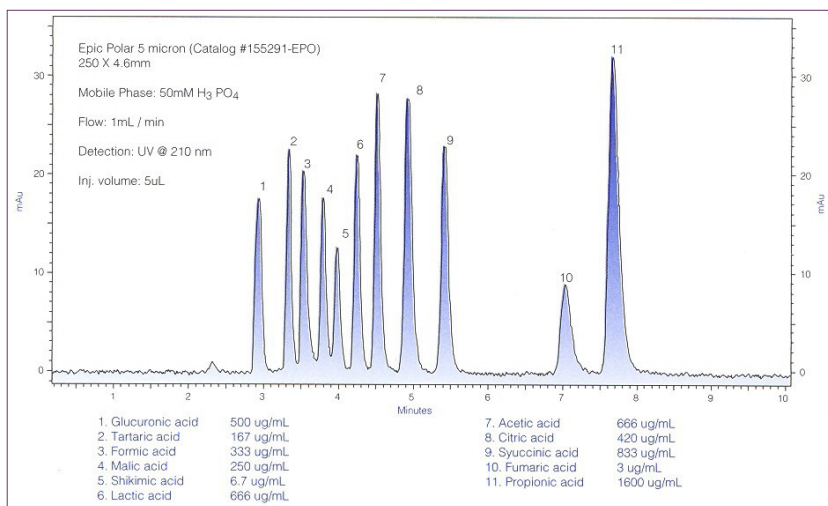


Figure 1. Low molecular weight polar organic acids on Epic Polar.

Phase	Length (mm)	ID (mm)	Particle Size (µm)	Part No.
Epic Polar	50	2.1	1.8	512A91-EPO
Epic Polar	50	2.1	3	112191-EPO
Epic Polar	50	2.1	5	112291-EPO
Epic Polar	50	4.6	3	115191-EPO
Epic Polar	50	4.6	5	115291-EPO
Epic Polar	100	2.1	1.8	522A91-EPO
Epic Polar	100	2.1	3	122191-EPO
Epic Polar	100	2.1	5	122291-EPO
Epic Polar	100	3.0	1.8	523A91-EPO
Epic Polar	100	3.0	3	123191-EPO
Epic Polar	100	4.6	3	125191-EPO
Epic Polar	150	2.1	1.8	532A91-EPO
Epic Polar	150	2.1	3	132191-EPO
Epic Polar	150	2.1	5	132291-EPO
Epic Polar	150	3.0	3	133191-EPO
Epic Polar	150	4.6	3	135191-EPO
Epic Polar	150	4.6	5	135291-EPO
Epic Polar	200	4.6	3	145191-EPO
Epic Polar	250	2.1	5	152291-EPO
Epic Polar	250	4.0	5	154291-EPO
Epic Polar	250	4.6	10	155391-EPO
Epic Polar	250	4.6	3	155191-EPO
Epic Polar	250	4.6	5	155291-EPO
Epic Polar	300	3.9	10	16e391-EPO
Epic Polar	300	4.0	5	164291-EPO
Epic Polar Prep	250	10	5	157291-EPO
Epic Polar Prep	250	20	10	158391-EPO
Epic Polar Prep	250	20	5	158291-EPO
Epic Polar Prep	250	30	5	15N291-EPO
Epic Polar Analytical Guard Cartridges (Pkg. 5)	10	3.0	5	500101-EPO
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.