

ProPac Evolution | 20 years of charge variant innovation

LC-MS/MS compatibility

ProPac Elite WCX LC Columns for maximum reproducibility

Resolution of biosimilars

Salt gradient

20 years of charge variant innovation brings ultimate versatility

pH gradients improve method transfer

Same selectivity with salt or pH

Part Number	Description	pH	Quantity
083273	Thermo Scientific™ CX-1 pH Gradient Buffer A	5.6	125 mL
083275	Thermo Scientific™ CX-1 pH Gradient Buffer B	10.2	125 mL
083274	CX-1 pH Gradient Buffer Kit; 125 mL Buffer A + 125 mL Buffer B	5.6 to 10.2	125 mL + 125 mL
085346	CX-1 pH Gradient Buffer A	5.6	250 mL
085348	CX-1 pH Gradient Buffer B	10.2	250 mL
085349	CX-1 pH Gradient Buffer Kit; 250 mL Buffer A + 250 mL Buffer B	5.6 to 10.2	250 mL + 250 mL
302779	CX-1 pH Gradient Buffer A	5.6	500 mL
302780	CX-1 pH Gradient Buffer B	10.2	500 mL
303274	CX-1 pH Gradient Buffer A	5.6	1 L
303275	CX-1 pH Gradient Buffer B	10.2	1 L
083381	CX-1 pH Gradient Starter Kit; 125 mL Buffer A + 125 mL Buffer B Thermo Scientific™ MAbPac SCX-10 Column (10 µm, 4 × 250 mm)	5.6 to 10.2	125 mL + 125 mL
083378	CX-1 pH Gradient High Throughput Kit; 125 mL Buffer A + 125 mL Buffer B MAbPac SCX-10 Column (5 µm, 4 × 50 mm)	5.6 to 10.2	125 mL + 125 mL

pH gradient buffers

Description	Particle Size (µm)	Part Number
Thermo Scientific™ ProPac™ Elite WCX columns		
ProPac Elite WCX, 2 × 50 mm	5	303028
ProPac Elite WCX, 2 × 150 mm	5	303027
ProPac Elite WCX, 2 × 250 mm	5	303026
ProPac Elite WCX, 4 × 50 mm	5	302973
ProPac Elite WCX, 4 × 150 mm	5	302972
ProPac Elite WCX, 4 × 250 mm	5	303025
ProPac Elite WCX, 4 × 150 mm, 3 Columns from 1 Lot	5	302976
ProPac Elite WCX, Analytical, 4 × 150 mm, 3 Columns from 3 Lots	5	302977
ProPac Elite WCX, Analytical, 4 × 250 mm, 3 Columns from 1 Lot	5	303061
ProPac Elite WCX, Analytical, 4 × 250 mm, 3 Columns from 3 Lots	5	303062

Versatility in column size formats

Speed with CX-1 and shorter column formats

Take charge of your variant analysis and see your separations in high resolution

For information please visit thermofisher.com/propac

For Research Use Only. Not for use in diagnostic procedures. © 2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. PO21984-EN 0919