

Consumables

Phase	Length (mm)	I.D. (mm)	µm	Part
Quasar C18	300	3.9	5	N9308800
Quasar C18	250	4.6	5	N9308801
Quasar C18	150	4.6	5	N9308802
Quasar C18	100	4.6	5	N9308803
Quasar C18	50	4.6	5	N9308804
Quasar C18	150	4.6	3	N9308805
Quasar C18	100	4.6	3	N9308806
Quasar C18	50	4.6	3	N9308807
Quasar C18	150	3.0	3	N9308808
Quasar C18	100	3.0	3	N9308809
Quasar C18	50	3.0	3	N9308810
Quasar C18	150	2.1	3	N9308811
Quasar C18	100	2.1	3	N9308812
Quasar C18	50	2.1	3	N9308813
Quasar C18	100	4.6	1.7	N9308814
Quasar C18	50	4.6	1.7	N9308815
Quasar C18	100	3.0	1.7	N9308816
Quasar C18	50	3.0	1.7	N9308817
Quasar C18	100	2.1	1.7	N9308818
Quasar C18	50	2.1	1.7	N9308819

Phase	Length (mm)	I.D. (mm)	µm	Part
Quasar C18 Guard Cartridge (3/pack)	10	3	5	N9308980
Quasar C18 Guard Cartridge (3/pack)	10	3	3	N9308981
Quasar Guard Cartridge Holder	-	-	-	N9306876
Quasar C18 Method Validation Kit, 250 x 4.6 mm (3 columns each from a different batch)	250	4.6	5	N9300940
Quasar C18 Method Validation Kit, 150 x 4.6 mm (3 columns each from a different batch)	150	4.6	5	N9300941

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 © 2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

014829_01

PKI

DRAFT