






# Clarus® 480/580/680 Capillary Split/Splitless Injector Inlet Supplies



## A INLET SEPTA (11 MM)

Product	Description	10-pk Part No.	50-pk Part No.
	BTO Orange Injection Port Septa (up to 400 °C)	N9306872	N9302972
	Green Injection Port Septa (up to 350 °C)	N9306218	N9306219
	Blue Injection Port Septa (up to 275 °C)	N9306873	N9306874

## B CAPILLARY INLET LINER O-RINGS

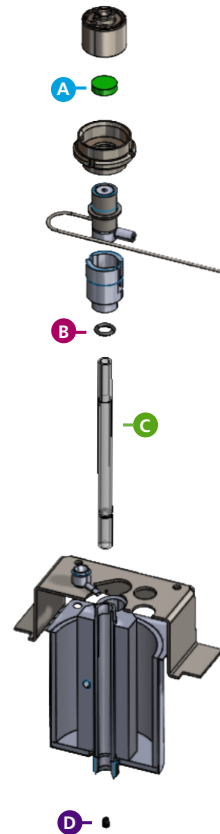
Product	Description	Quantity	Part No.
	O-ring, Viton® for Glass Liner (maximum injector temperature 300 °C)	10-pk	N9302783
	O-ring, Kalrez® for Glass Liner (maximum injector temperature 450 °C)	1-pk	N9302782

Visit our [Chromatography Consumables Catalog](#) for a complete product listing.




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

For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://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.



## C CAPILLARY SPLIT/SPLITLESS INLET LINERS\*

Product	Description (ID X OD X Length)	5-pk Part No.
	Capillary Split/Splitless Deactivated Glass Liners with Deactivated Wool (4 x 6.2 x 92.1 mm)	N9306233
	Capillary Split/Splitless Deactivated Glass Liners with Deactivated Wool and Tapered End (4 x 6.2 x 92.1 mm)	N9306235
	Capillary Split Deactivated Glass Liners with Deactivated Wool (4 x 6.2 x 92.1 mm)	N9306236

\*Additional liners are also available – see consumables catalogue for further details.

## D STANDARD FERRULES FOR CAPILLARY INJECTOR\*

Size	For use with Column ID (mm)	Ferrule ID (mm)	Quantity	Graphite Part No.	Graphite/Vespel Part No.
1/16 in.	≤ 0.25	0.4	10-pk	N9307245	09920104
1/16 in.	0.32	0.5	10-pk	09903700	09920105
1/16 in.	0.53	0.8	10-pk	09920141	09920107

\*2-hole ferrules are also available – see consumables catalogue for further details.

