

These new high quality and economical liners have been specially developed for the Clarus® gas chromatographs. They are color-coded to enable easy identification and selection for each injector type.

The liners are pre-packed with deactivated glass wool to ensure optimum sample flow path and inertness to deliver excellent performance even with critical applications.







- Engineered design for excellent vaporization
- Color coded with directional arrow per liner application
- Consistent reproducibility



INLET LINERS FOR CLARUS GC INSTRUMENTS

See reverse side for product details.

INLET LINERS FOR CLARUS GC INSTRUMENTS

Part No.	Color Code	Description	Applications	Qty.
N9306232	Yellow	PSS deactivated glass liners 2 mm i.d. x 4.0 mm o.d. x 86.2 mm with deactivated wool. Narrow bore and quartz wool increase volatilization and reproducibility. 	General purpose	5 per pack
N9306233	Green	Capillary splitless deactivated glass liners 4 mm i.d. x 6.2 mm o.d. x 92.1 mm with deactivated wool. 	General purpose Intermediate polarity	5 per pack
N9306234	Green	Capillary splitless deactivated glass liners 4 mm i.d. x 6.2 mm o.d. x 92.1 mm with deactivated wool. 	General purpose Intermediate polarity	1 per pack
N9306235	Red	Capillary splitless deactivated glass liners 4 mm i.d. x 6.2 mm o.d. x 92.1 mm with deactivated wool. 	Focus liner for split/ splitless operations	5 per pack
N9306236	Orange	Capillary splitless deactivated glass liners 4 mm i.d. x 6.2 mm o.d. x 92.1 mm with deactivated wool. Quartz wool is used to fully vaporize the sample. 	General purpose	5 per pack
N9306237	Blue	PSS Splitless deactivated glass liners 1.25 mm i.d. x 4.0 mm o.d. x 86.2 mm 	Low volume sample analysis Fast GC, gaseous samples (purge and trap)	5 per pack

For more details, call **1-800-762-4000** or visit www.perkinelmer.com/supplies



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

Copyright ©2010-2011, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

009478A_01