

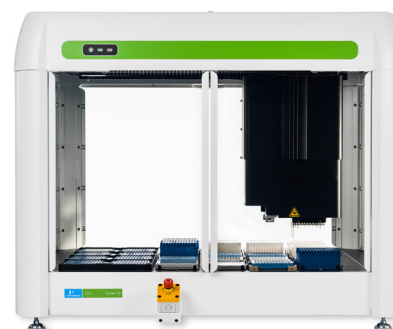


LESS HASSLE
GREATER
PRODUCTIVITY

Accessories for the Sciclone G3 and Zephyr G3 Automated Workstations

The Sciclone® G3 laboratory automation workstations provide labs with a broad range of highly accurate and precise liquid handling functions in a versatile, open-design format that adapts to almost any protocol. Each workstation is configured with your choice of liquid handling functionality and on-deck accessories, and are field reconfigurable to adapt as your research processes. Choose your deck configuration based on your application for high-throughput, consistent results.

The Zephyr® G3 workstation family includes NGS and SPE workstations. The NGS workstation is a benchtop liquid handler designed to automate preparation of libraries for next generation sequencing and to address the growing number of laboratories with requirements for the physical separation of pre-and post-PCR processes. The SPE workstation is a compact, robust and flexible liquid handler to automate the critical steps required for plate-based Solid Phase Extraction (SPE) purification.



Sciclone G3 Workstation



Zephyr G3 Workstation

DISPOSABLE TIPS

Automation pipette tips for 384-well microplates are available sterilized or non-sterilized and with either 10 trays or 35 trays per box. Our pipette tips are available in 96- and 384-well formats. These products are manufactured to precise specifications under stringent quality assurance standards.

All racks are designed for use with automation equipment, and the tips are designed to provide exceptional straightness. In addition, all tips are certified to be DNase and RNase free. Volume discounts are available for ordering 50 or more boxes. Please contact your local sales representative for details.

Part No.	Description	Quantity (per box)
125792	25 uL - 384 Rack Tips, Clear, Non-Sterilized	10
125791	25 uL - 384 well tips, Clear, Non-sterilized	35
125795	25 uL - 384 Rack Tips, Clear, Sterilized	10
125793	25 uL - 384 Rack Tips, Clear, Sterilized	35
105641	50 uL - 96 Rack Tips, Clear, Non-Sterilized	10
105643	50 uL - 96 Rack Tips, Clear, Non-Sterilized	30
105642	50 uL - 96 Rack Tips, Clear, Sterilized	10
105644	50 uL - 96 Rack Tips, Clear, Sterilized	30
105645	35 uL - 96 Rack Tips, Clear, Sterilized	10
109079	100 uL - 96 Rack Tips, Clear, Sterilized	10

Part No.	Description	Quantity (per box)
105647	100 uL - 96 Rack Tips, Clear, Non-Sterilized	30
109080	100 uL - 96 Rack Tips, Clear, Sterilized	10
105648	100 uL - 96 Rack Tips, Clear, Sterilized	30
111624	80 uL - 96 Rack Tips, Clear, Sterilized	10
109081	200 uL - 96 Rack Tips, Clear, Non-Sterilized	10
105649	200 uL - 96 Rack Tips, Clear, Non-Sterilized	30
109082	200 uL - 96 Rack Tips, Clear, Sterilized	10
105650	200 uL - 96 Rack Tips, Clear, Sterilized	30
111426	150 uL - 96 Rack Tips, Clear, Sterilized	10
111623	200 uL - 96 Rack Tips, Wide Bore, Clear	10

MICROPLATES FOR NEXT GENERATION SEQUENCING INSTRUMENTS

Microplates for next generation sequencing instruments from PerkinElmer are engineered to withstand the stresses of heat sealing, thermal cycling, and robotic handling. These plates are also engineered for stability, flatness and are warp-free which ensures

precise positioning for automation. Plates are available in 96- and 384-well formats. Below are recommended microplates for next generation sequencing instruments from PerkinElmer. To view all plates please visit www.perkinelmer.com/microplates

Part No.	Description	Quantity (per box)
6008700	Polypropylene deep-well reservoir, 12 column 21 mL	25
6008880	Polypropylene deep-well storage plate, 96 - 2 mL square well, v-bottom	25
6008870	Hard-Shell® PCR plate, 96-well	50
6008910	Hard-Shell® PCR plate, 384-well	50

Part No.	Description	Quantity (per box)
6008900	Polypropylene plate, 96 - 1 mL, round bottom	100
6008299	Polypropylene plate, 96 - 450 µL, v-bottom	200
6008890	Polypropylene low volume, 384 - 35 µL, round bottom	50

POPULAR ACCESSORIES

Part No.	Description	Quantity (per box)
6050174	TopSeal adhesive seal for PCR plates	100
6000030	Universal microplate lid	100

For more information visit www.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 © 2017, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.