

Polypropylene Autosampler Tubes

PerkinElmer translucent polypropylene tubes are designed to provide you with excellent chemical resistance. PerformR™ 15 mL and 50 mL conical tubes are designed for your routine everyday applications and storage needs. SuperClear™ tubes are made from a unique medical grade resin that is far more durable than commodity grade resins used in most other brands of tubes. PerformR and SuperClear tubes are available with either a flat cap with a two stage seal, or a plug style cap that includes a very deep sealing area.









| Round Bottom Tubes | | | | | | | | | |
|--------------------|--|----------|----------------|-----------|------|--|--|--|--|
| Part No. | Description | Туре | Size | Packaging | Qty. | | | | |
| B0193235 | 5 mL without caps | Standard | 12 mm x 75 mm | Bulk | 500 | | | | |
| N0777156 | 8 mL without caps | Standard | 13 mm x 100 mm | Bulk | 1200 | | | | |
| N0777159 | 8 mL without caps | Standard | 13 mm x 100 mm | 250/bag | 1000 | | | | |
| B0508901 | 8 mL without caps | Standard | 13 mm x 100 mm | Bulk | 1000 | | | | |
| N0777940 | 14 mL without caps | Standard | 17 mm x 100 mm | Bulk | 1000 | | | | |
| N0777941 | 14 mL with Dual Position Polyethylene Caps included | Standard | 17 mm x 100 mm | Bulk | 1000 | | | | |
| N0777942 | Dual Position Caps for 14 mL N0777940 Tubes (clear) Caps Only | Standard | | Bulk | 1000 | | | | |
| N0777168 | 14 mL without caps | Standard | 17 mm x 100 mm | 125/bag | 1000 | | | | |
| N0777167 | 15 mL without plugs | Standard | 17 mm x 100 mm | Bulk | 1200 | | | | |
| N0777599 | Plugs for 15 mL N0777167 Tubes (blue) Plugs Only | Standard | | Bulk | 1200 | | | | |
| N0777071 | 15 mL Graduated without plugs | Standard | 17 mm x 100 mm | Bulk | 2000 | | | | |
| N0777072 | Plugs for 15 mL N0777071 Tubes (white) Plugs Only | Standard | | Bulk | 1000 | | | | |
| N9301205 | 16 mL without caps | Standard | 17 mm x 100 mm | 125/bag | 1000 | | | | |
| | | | | | | | | | |



Figure 1. PerkinElmer S10 Autosampler













Figure 2. Flat and Plug style

| Conical | Bottom Tubes with Graduati | on | | | |
|----------|--|------------|----------------|---------------|------|
| Part No. | Description | Туре | Size | Packaging | Qty. |
| B0193233 | 15 mL with Red Flat Caps attached (Standard PerkinElmer S10 Tubes) | Standard | 17 mm x 120 mm | 50/bag | 500 |
| N0774095 | 15 mL Sterile Tubes, with Teal Flat Caps attached | PerformR | 17 mm x 118 mm | 50/bag | 500 |
| N0777701 | 15 mL with Teal Flat Caps attached | SuperClear | 17 mm x 118 mm | 50/bag | 500 |
| N0777702 | 15 mL with Teal Flat Caps attached | SuperClear | 17 mm x 118 mm | Racked 50/bag | 500 |
| N0777703 | 15 mL with Teal Flat Caps included | SuperClear | 17 mm x 118 mm | Bulk | 500 |
| N0777704 | 15 mL with Teal Plug Caps attached | SuperClear | 17 mm x 118 mm | 50/bag | 500 |
| N0777705 | 15 mL with Teal Plug Caps attached | SuperClear | 17 mm x 118 mm | Racked 50/bag | 500 |
| N0777706 | 15 mL with Teal Plug Caps | SuperClear | 17 mm x 118 mm | Bulk | 500 |
| B0193234 | 50 mL with Red Flat Caps attached (Standard PerkinElmer S10 Tubes), Freestanding | Standard | 28 mm x 115 mm | 25/bag | 500 |
| N0777697 | 50 mL with Teal Flat Caps attached, Freestanding | PerformR | 30 mm x 116 mm | 50/bag | 500 |
| N0777698 | 50 mL with Teal Flat Caps included, Freestanding | PerformR | 30 mm x 116 mm | Bulk | 500 |
| N0777699 | 50 mL with Teal Plug Caps attached, Freestanding | PerformR | 30 mm x 116 mm | 50/bag | 500 |
| N0777700 | 50 mL with Teal Plug Caps included, Freestanding | PerformR | 30 mm x 116 mm | Bulk | 500 |
| N0774096 | 50 mL Sterile Tubes with Teal Flat Caps attached | PerformR | 29 mm x 115 mm | 50/bag | 500 |
| N0777691 | 50 mL with Teal Flat Caps attached | SuperClear | 29 mm x 115 mm | 50/bag | 500 |
| N0777692 | 50 mL with Teal Flat Caps attached | SuperClear | 29 mm x 115 mm | Racked 50/bag | 500 |
| N0777693 | 50 mL with Teal Flat Caps included | SuperClear | 29 mm x 115 mm | Bulk | 500 |
| N0777694 | 50 mL with Teal Plug Caps attached | SuperClear | 29 mm x 115 mm | 50/bag | 500 |
| N0777695 | 50 mL with Teal Plug Caps attached | SuperClear | 29 mm x 115 mm | Racked 50/bag | 500 |
| N0777696 | 50 mL with Teal Plug Caps included | SuperClear | 29 mm x 115 mm | Bulk | 500 |

Every day you count on PerkinElmer to provide you with solutions that deliver reliable performance, control operating costs and maximize operational time. Our complete portfolio of consumables, parts, supplies, training and service helps you meet both routine and demanding measurement challenges. We invest heavily in testing and validating our products to ensure you receive guaranteed compatibility and performance — on-time, for every instrument in your laboratory.

For a complete listing of consumables, please visit www.perkinelmer.com/supplies

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

