

Samco Clicktainer Vials

Unique multi-click seal system to ensure sample protection and user safety

Thermo Scientific™ Samco™ Clicktainer™ Vials ordering information

| Description | Closure color | Label type | Product configuration | Quantity (pk x pks per case) | Catalog number |
|--|---------------|--------------------|-----------------------|------------------------------|----------------|
| 120mL/53mm sterile | | | | | |
| 120mL/53mm CLK SPC LBL OR CAP ST | Orange | Tabbed | Pre-assembled | 75x4 | 120ORG53-1000 |
| 120mL/53mm CLK O/R N/LBL OR CAP ST | Orange | None | Peel pouch | 1x100 | 120ORG53-4000 |
| 120mL/53mm CLK SPC LBL WT CAP ST | White | Tabbed | Pre-assembled | 75x4 | 120WHT53-1000 |
| 120mL/53mm CLK MSU, NO FUNNEL, ST | White | Tabbed | Mid-stream kit | 1x100 | 120WHT53-3000 |
| 120mL/53mm CLK O/R POUCH WT CAP ST | White | None | Peel pouch | 1x100 | 120WHT53-4000 |
| 120mL/53mm Non-sterile | | | | | |
| 120mL/53mm CLK SPC LBL OR CAP NS | Orange | Tabbed | Pre-assembled | 75x4 | 120ORG53-2000 |
| 120mL/53mm CLK SPC N/LBL WT CAP NS | White | None | Bagged separate | 300x1 | 120WHT53-2000 |
| 90mL/53mm sterile | | | | | |
| 90mL/53mm CLK SPC LBL OR CAP ST | Orange | Tabbed | Pre-assembled | 100x4 | 90ORG53-1000 |
| 90mL/53mm CLK SPC LBL OR CAP ST DISP | Orange | Tabbed | Dispenser box | 20x16 | 90ORG53-1001 |
| 90mL/53mm CLK SPC LBL WT CAP ST | White | Tabbed | Pre-assembled | 100x3 | 90WHT53-1000 |
| 90mL/53mm CLK SPC TSLBL WT CAP ST | White | Label w/temp strip | Pre-assembled | 100x3 | 90WHT53-1001 |
| 90mL/53mm CLK SPC TS WT CAP ST | White | Temp strip only | Pre-assembled | 100x4 | 90WHT53-1002 |
| 90mL/53mm CLK SPC LBL GN CAP ST | Green | Tabbed | Pre-assembled | 100x4 | 90GRN53-1000 |
| 90mL/53mm Non-sterile | | | | | |
| 90mL/53mm CLK SPC LBL WT CAP NS | White | Tabbed | Pre-assembled | 75x4 | 90WHT53-2000 |
| 90mL/53mm CLK SPC TSLBL WT CAP NS | White | Label w/temp strip | Pre-assembled | 300x1 | 90WHT53-2001 |
| 90mL/53mm CLK SPC N/LBL WT CAP NS | White | None | Bagged separate | 300x1 | 90WHT53-2002 |
| 90mL/53mm CLK SPC N/LBL WT CAP NS | White | None | Pre-assembled | 300x1 | 90WHT53-2003 |
| 90mL/53mm CLK SPC TS WT CAP NS | White | Temp strip only | Pre-assembled | 100x4 | 90WHT53-2004 |
| Non-sterile closures and containers | | | | | |
| 53mm CLK CAP ORANGE NS | Orange | None | Bulk closures only | 1500x1 | CAPORG53-1000 |
| 53mm CAP WHITE NS | White | None | Bulk closures only | 1500x1 | CAPWHT53-1000 |
| 53mm CLK CAP GREEN NS | Green | None | Bulk closures only | 1500x1 | CAPGRN53-1000 |
| 120mL/53mm CLK SPC N/LBL NO CAP NS | 120mL | None | Bulk container only | 300x1 | 12000-5300 |
| 90mL/53mm CLK SPC N/LBL NO CAP NS | 90mL | None | Bulk container only | 300x1 | 9000-5300 |

Note: Clicktainer Vial Mid-Stream Kit contains a container, 3 castile soap wipes, and an instruction sheet in a tear open bag

Find out more at thermofisher.com/clicktainer

ThermoFisher
SCIENTIFIC

Samco Clicktainer Non-Sterile Closures and Containers CAPORG53-1000, CAPWHT53-1000, CAPGRN53-1000, 12000-5300 and 9000-5300 are intended for General Lab Use Only. It's the customers responsibility to ensure that the performance of the product is suitable for customers' specific use or application. All other products listed are Class I Medical Devices. © 2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. BRLPECLICKTAVI 0919 Insert