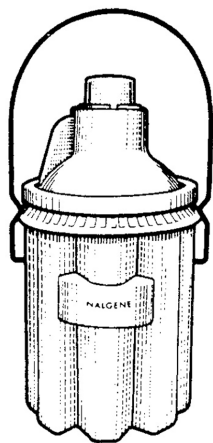


Thermo Scientific Nalgene Safety Bottle Carriers (low-density polyethylene with epoxy-coated steel handle, polycarbonate cap)

Cat. No. 6501-2500 – holds a 2.5 liter or 5 pint bottle

Cat. No. 6501-4000 – holds a 4 liter or 1 gallon bottle

These Nalgene® Safety Bottle Carriers cushion glass bottles to reduce the danger of injury from broken glass and spilled chemicals. The cover is easy to snap on and remove and the inside dimensions are large enough to accommodate plastic-coated glass bottles. A clear polycarbonate cap protects the bottle closure and lets you identify the color-coded closure on acid bottles. With the cap removed, you can pour from the bottle without removing it from the carrier. The handle locks the carrier cover in place for stability while carrying. Storage and transportation are made easier with the handle conveniently fitting over another carrier when stored side-by-side. This prevents accidental tipping.



To Remove Cover

1. Place the carrier on a firm, level surface and secure it firmly, minimizing movement. Make sure handle is in down position.
2. While holding handle firmly in one hand, grip cover under the beveled edge of the carrier with other hand. Gently pull up and out while sliding your fingers under and around the carrier lid (usually a minimum of 180°) until the cover comes loose.
3. Lift off the cover.

To Fasten Cover

1. Place the carrier on a firm, level surface and secure it firmly, minimizing movement.
2. Place the cover on the carrier.
3. Press down around the rim of the cover until it snaps down, over the beveled edge of the carrier.

NOTE: When using fingerloop bottles, make sure the slot in the cover is placed **over** the fingerloop.

To Lock Cover in Place

1. Lift the handle until it is in an upright position.

To Remove Plastic Cap to Pour

1. Twist cap firmly to the right and lift off.
2. To replace, line cap up with the notches in cover and turn firmly to the left until secure.

© 2011 Thermo Fisher Scientific Inc. All rights reserved. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.