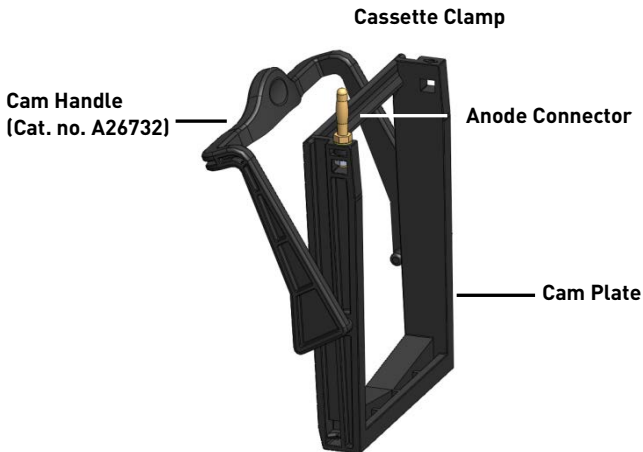


Instructions for replacing cam handles for the Bolt® Mini Gel Tank

Cat. Nos. A26732, B4477599

Pub. Part No. 100027558 Pub. No. MAN0010807 Rev. A

Upgrade your Bolt® Mini Gel Tank to accommodate the new 10 cm Bolt® mini gels (and existing NuPAGE® and Novex® Tris-Glycine mini gels) by replacing the current black cam handles on your cassette clamps with new gray cam handles.



Instructions

Two identical gray Cam Handles are provided. The replacement procedure is identical for Right and Left Cassette Clamps.

Note: Illustrations show the Right Cassette Clamp. The anode connector is in a different location on the Left Cassette Clamp.

View a video of these instructions at:

www.lifetechnologies.com/bolttanksupport

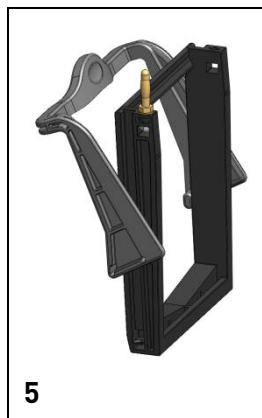
1. Carefully pull on one side of the black Cam Handle and separate the peg on the handle from the hole in the Cam Plate.
2. Slide the peg on the other side of the black Cam Handle out of the Cam Plate, and separate the two pieces.



3. Position the gray Cam Handle in the same orientation as the original black Cam Handle.

Note: each side of the gray Cam Handle has one round peg and one oval peg. Only the round peg fits into the Cam Plate.

4. Slide the peg on one side of the gray Cam Handle into the corresponding hole in the Cam Plate.
5. Carefully pull the other side of the gray Cam Handle, and slide the round peg into the hole in the Cam Plate.



Important licensing information

This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

DISCLAIMER: LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

© 2014 Thermo Fisher Scientific Inc. All rights reserved.
All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.
For support visit www.lifetechnologies.com/support or email techsupport@lifetech.com

www.lifetechnologies.com

11 July 2014

