

SureLock™ Tandem Midi Gel Tank

Cat. no. STM1001

Publication No. MAN0018517

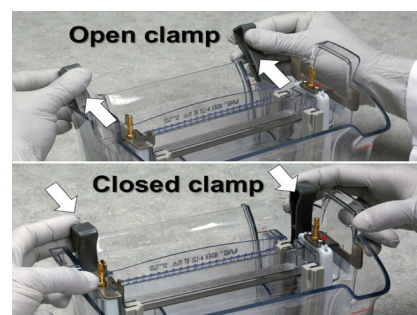
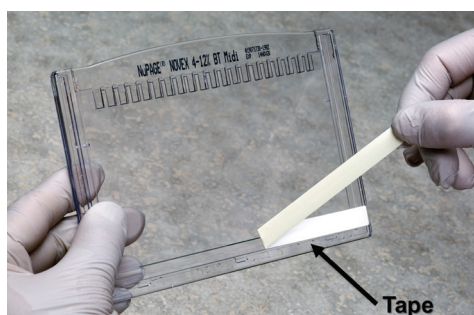
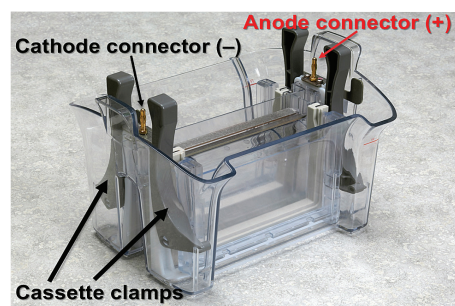
Rev. A.0

Instructions for using the SureLock™ Tandem Midi Gel Tank to perform electrophoresis are described below. For more detailed instructions, refer to the manual available from thermofisher.com.

Before Starting

- Ensure the power supply is adequate for the number of gels to be run.
- Select Invitrogen™ midi gels appropriate for your application.
- Prepare 1X running buffer appropriate for your Invitrogen™ midi gels. Each gel requires ~520 mL of buffer.

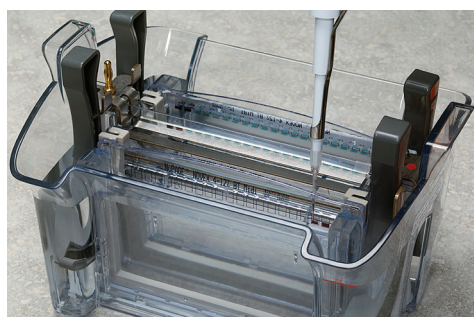
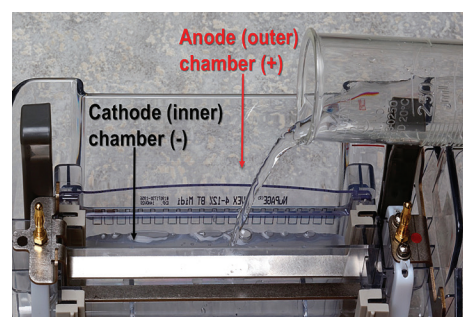
Procedure



1. Place the cassette clamp(s) into the chamber(s) with the flat side(s) facing the center of the tank.
The tank can run 2 gels at once with 1 gel per side.

2. Remove the comb and peel away the tape at the bottom of the gel cassette.
Rinse the wells 3 times with 1X running buffer.

3. Open the cassette clamps.
Place the gel cassette(s) inbetween the cassette clamp and the gray sealing gasket, with the cassette wells facing the tank center.
Lower the cassette so it rests on the bottom of the tank. Close the cassette clamps.



4. Fill the tank with running buffer. For each gel being run:
Add 1X buffer to the cathode (inner) chamber (-) such that the gel wells are submerged with buffer (~170 mL).
Add 1X buffer to the anode (outer) chamber (+) to the level of the red fill line printed on the outside of the tank (~350 mL).
WARNING! Do not overfill.

5. Load samples and protein standards.

6. Make sure the power supply is off.
WARNING! If only running one gel, ensure that the cassette clamp from the unused chamber is removed.
Place the lid on the tank and plug the electrode cords into the power supply.
Turn the power supply on to begin electrophoresis.

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 CORPORATION 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.

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.

Corporate entity: Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288

© 2019 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 thermofisher.com/support or email techsupport@lifetech.com

thermofisher.com

18 October 2019

ThermoFisher
S C I E N T I F I C