

SureLock™ Tandem Midi Blot Module

Cat. no. STM2001

Publication No. MAN0018516

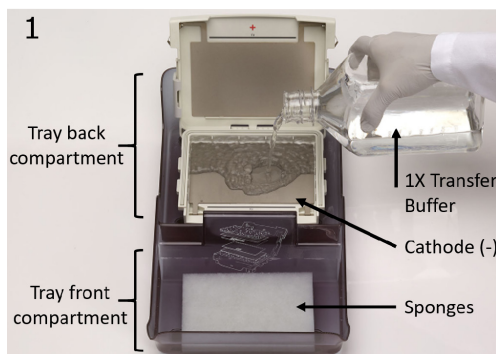
Rev. A.0

Instructions for using the SureLock™ Tandem Midi Blot Module to transfer proteins from a gel onto a membrane are described below. For more detailed instructions, refer to the SureLock™ Tandem Midi Blot Module User Guide available from thermofisher.com.

Before Starting

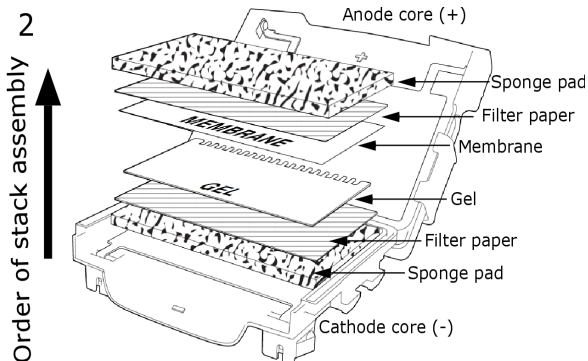
- Choose and prepare the transfer buffer compatible with the chemistry of the gel being transferred. Each transfer requires ~300 mL of 1X transfer buffer.
- Select and prepare a transfer membrane appropriate for your needs.
- Soak 2 sponge pads per transfer in 1X transfer buffer in the front compartment of the SureLock™ Tandem Tray (Cat. No. STM3001). Squeeze submerged pads or roll them using a Blotting Roller (Cat. No. LC2100) to ensure that all air bubbles are removed.
- Prepare the gel for transfer by trimming off the well fingers and foot from the gel.
- Equilibrate the gel:
 - If transferring a Novex™ Tris-Glycine Midi gel, soak the gel in 1X transfer buffer for 5 minutes prior to transfer.
 - If transferring a NuPAGE™ Bis-Tris or Tris-Acetate Midi gel, soaking is not recommended. Proceed directly to transfer.

Procedure



1. Place a Blot Module into the back compartment of the tray with the Cathode (-) side on the bottom.

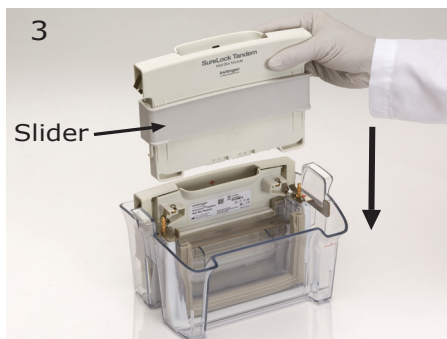
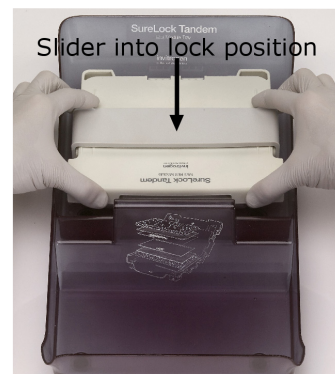
Add ~50 mL of 1X transfer buffer to the cathode shell prior to assembling the transfer stack.



2. Assemble the transfer stack in the Blot Module in the order shown.

- Soak filter paper briefly in 1X transfer buffer before use.
- Orient the gel with the wells toward the bottom of the Blot Module.
- Place the membrane on top of the gel.
- Use the Blotting Roller to remove any bubbles as each layer of the stack is assembled.

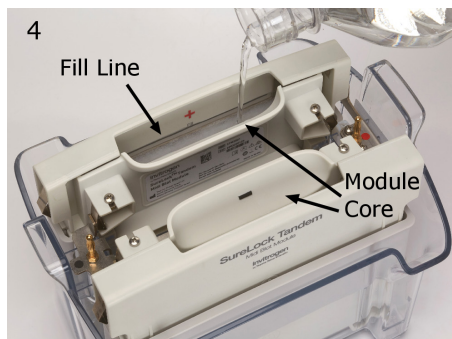
Close the Blot Module and shift the slider into the locked position.



3. Ensure that the cassette clamps used for gel electrophoresis are removed from the tank.

Note: The tank can run either 1 or 2 Blot Modules at a time.

Insert each Blot Module ensuring the slider is facing the outside of the tank and the Blot Module rests on the bottom of the tank.



4. Add 1X transfer buffer to the Fill Line shown on the inside of the Blot Module.

Note: Blot Modules do not need chilled buffer for transfer and there is no need to add water or transfer buffer in the space around the Blot Module.



5. Make sure the power supply is turned off.

Place the lid on the tank, plug the electrode cords into the power supply, then turn on the power supply.

Continued on reverse

For Research Use Only. Not for use in diagnostic procedures.

Continued from front page

Transfer Conditions

Recommended transfer conditions for Invitrogen™ Pre-Cast Midi Gels

Voltage	Time
25 V (constant)	30 minutes

CAUTION:

- Do not exceed 30 V.
- To prevent damaging the gel and the Blot Module, transfers should never be performed at 100 V, as may be required for other wet transfer devices.
- Refer to the SureLock™ Tandem Midi Blot Module User Guide for additional safety information.



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

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

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

© 2020 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

16 April 2020

ThermoFisher
SCIENTIFIC