

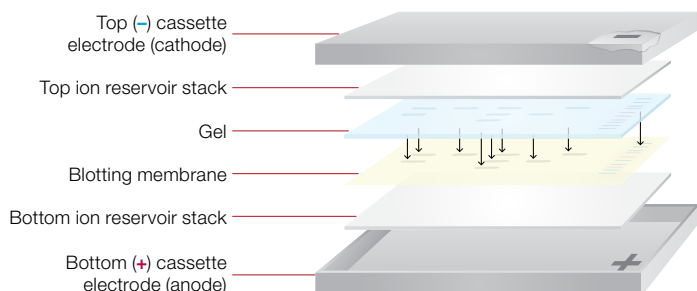
# BLOTTING

## Trans-Blot Turbo Transfer System

### Quick Start Guide

#### Instructions for Using the Trans-Blot Turbo System

**1. Assemble the Blotting Sandwich:** After gel electrophoresis, open the transfer pack and assemble the components on the cassette in the order shown below. For best results, use the roller to remove any air trapped between the layers.



If using a single mini or midi format sandwich, it should be placed in the center of the cassette bottom. If using two mini format gels, they should be placed on a midi stack with the foot of each gel toward the center.

**2. Insert Cassette into Instrument:** Place the lid on the cassette and lock by turning the knob clockwise, using the symbols on the lid as a guide. Slide the cassette into either bay. The upper bay is designated as bay A and the lower bay is designated as bay B. If running two cassettes simultaneously, refer to the table below for acceptable transfer pack combinations.

Combination	Acceptable*		Not Acceptable*	
	1	2	1	2
Upper bay A	1 mini gel	2 mini gels or 1 midi gel	1 mini gel	2 mini gels or 1 midi gel
Lower bay B	1 mini gel	2 mini gels or 1 midi gel	2 mini gels or 1 midi gel	1 mini gel

\* Conditions hold if trays are swapped.

**3. Start the Transfer:** With the cassette inserted into the instrument, press TURBO and select your gel type. Press A:RUN or B:RUN to begin the transfer. LIST will allow you to select from Bio-Rad's optimized protocols (see table below) or a user-defined protocol. NEW will allow you to create and run a new protocol.

Protocol Name	MW, kD	Time, min	2 Mini Gels or 1 Midi Gel	1 Mini Gel
STANDARD SD	Any	30	Up to 1.0 A; 25 V constant	
1.5 MM GEL	Any	10		
HIGH MW	>150	10	2.5 A constant; up to 25 V	1.3 A constant; up to 25 V
LOW MW	<30	5		
MIXED MW*	5–150	7		
1 Mini TGX*	5–150	3	N/A	2.5 A constant; up to 25 V

\* Also accessed via the TURBO navigation button.

**4. Downstream Processing:** At the end of the transfer, RUN COMPLETE will be displayed for the cassette slot.

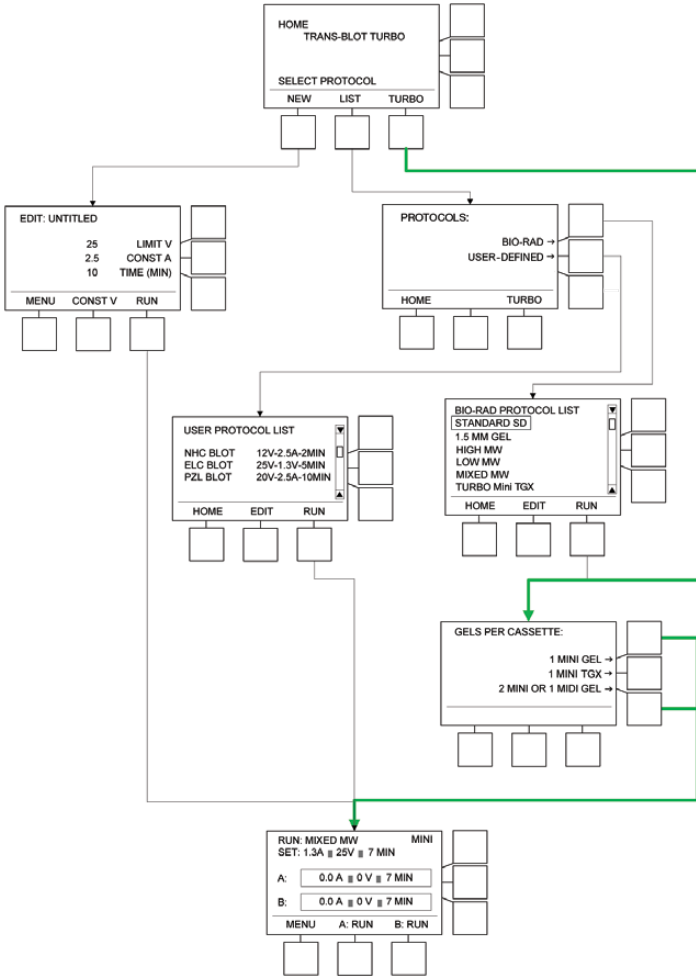
- Remove cassette from unit and unlock lid
- Disassemble blotting sandwich and extract membrane for downstream processing

**Note:** Clean and inspect the instrument and cassette after each use for damage. Contact Bio-Rad technical support if any damage is noted.

**BIO-RAD**

# Key Menu Screens

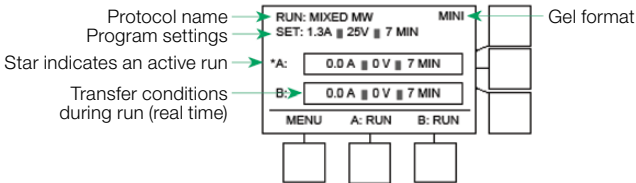
Refer to the Trans-Blot Turbo Blotting System Instruction Manual for a complete listing of menu screens.



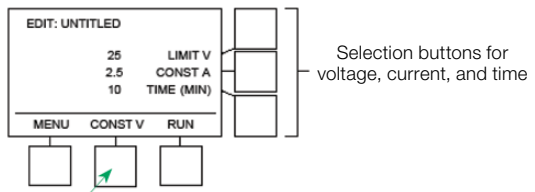
## Screen Features

### RUN Screen

MINI, MIDEI, MP TGX, or USER protocol



### EDIT Screen



Toggle between constant amps and constant volts

Bio-Rad and Trans-Blot are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.