

BLOTTING

Trans-Blot Turbo Transfer System

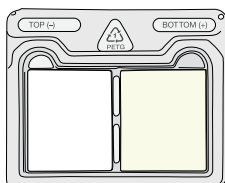
Transfer Pack

Quick Start Guide

Instructions for Using Trans-Blot Turbo Transfer Packs

Mini Transfer Packs

The anode stack and membrane are located in the right Bottom (+) well.



1. Place the membrane and bottom stack on the cassette base.*

- Use the blot roller to remove trapped air bubbles
- 1 mini or midi gel: place stack in the center of the cassette
- 2 mini gels on mini stacks: place stacks side by side in the center of the cassette
- 2 mini gels on midi stack: place midi stack in the center of the cassette

2. Place gel on top of membrane.

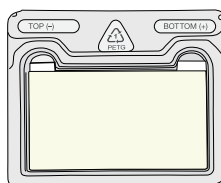
- Do not equilibrate the gel before transfer
- Remove any air bubbles with blot roller
- 2 mini gels on midi stack: place foot of gels toward the center

3. Place second wetted transfer stack on top of gel. This will serve as the top ion stack.

- Roll the assembled sandwich with blot roller to expel trapped air bubbles
- Please refrain from adding any extra transfer buffer to the cassette; saturated transfer stacks provide ample transfer buffer

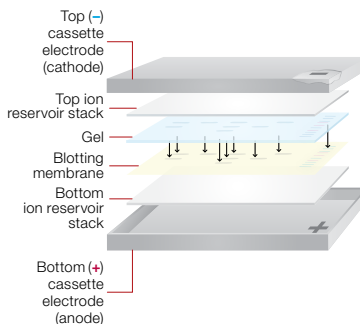
Midi Transfer Packs

The anode stack and membrane are located on top; use the right Bottom (+) finger tab to remove.



4. Close and lock cassette lid. Insert the cassette into the instrument and begin transfer.

* For detailed instructions, please refer to the Trans-Blot Turbo Blotting System Instruction Manual.



BIO-RAD



Please recycle transfer pack trays.

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.

Bio-Rad Preprogrammed Protocols

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	2.5 A constant; up to 25 V	1.3 A constant; up to 25 V
HIGH MW	>150	10		
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.

Notes

- Gels do not require equilibration and can be transferred immediately after electrophoresis
- To prevent the membrane from drying out, use transfer packs immediately after opening
- Transfer stacks will be warm after transfer. Avoid drying the membrane during stack disassembly
- Clean and inspect the instrument and cassette after each use for damage. Contact Bio-Rad technical support if any damage is noted

Call 1-800-4BIO-RAD (1-800-424-6723) or visit bio-rad.com for technical support.

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.