



BLOTTING

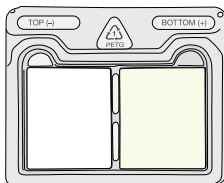
Trans-Blot® Turbo™ Transfer System

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: place midi stack in the center of the cassette

2. Place gel on top of membrane.

- Do not equilibrate the gel before transfer
- If needed, remove any air bubbles with blot roller
- 2 mini gels: place midi stack in the center of the cassette

3. Place foot of gel toward the center of the midi membrane.

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

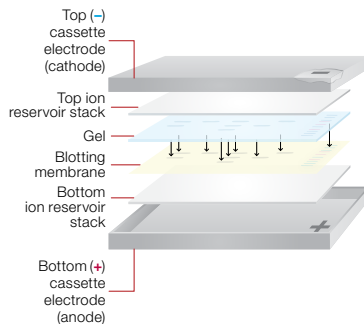
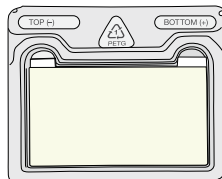
- Roll the assembled sandwich with blot roller to expel trapped air bubbles
- Please refrain from adding any extra transfer buffer to the cassette; saturated transfer stacks provide ample transfer buffer

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

* For detailed instructions, please refer to the Trans-Blot Turbo system instruction manual.

Midi Transfer Packs

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



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 for Efficient Transfer

- 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
- After transfer is complete, cassettes are immediately ready for another transfer; no cooling period is required

For technical support, call 1-800-4BIO-RAD (1-800-424-6723) or visit us at www.bio-rad.com.