



# Trans-Blot® Turbo™ Transfer System

Getting to the Finish Line Faster





The Trans-Blot® Turbo™ is a fast, efficient, and reproducible transfer system for transferring proteins from gels to membranes in as little as 3 minutes.





# Getting to the Finish Line Faster!



Bio-Rad introduces the Trans-Blot Turbo system — the next innovation in protein transfer. The Trans-Blot Turbo system reduces transfer protocols for gels to as little as 3 minutes, while maintaining high efficiency, throughput, and the flexibility to run turbo or traditional semi-dry protocols.



### Turbo Transfers with Trans-Blot Turbo Transfer Packs

- **3-minute protocol** — A single Mini-PROTEAN® TGX™ gel (for proteins with MW 5–150 kD) can be transferred in as little as 3 min
- **7-minute protocol** — Up to 4 mini or 2 midi gels with mixed-molecular weight proteins (MW 5–150 kD) can be efficiently transferred in 7 min
- **10-minute protocol** — Up to 4 mini or 2 midi gels with high-molecular weight proteins (MW 25–300+ kD) can be efficiently transferred in 10 min



3 min transfer



7 min transfer



10 min transfer

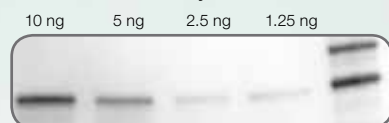
**Protein transferred using different protocols.** *E. coli* lysate (6 µg) was diluted two-fold. Samples were separated with Mini-PROTEAN TGX gels, transferred with the Trans-Blot Turbo system, stained with SYPRO Ruby and imaged on a VersaDoc™ 4000 MP system. Standards in lane 1 are Precision Plus Protein™ Unstained, with a top band of 250 kD.



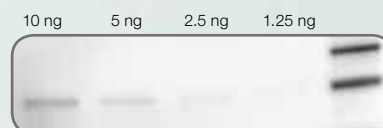
### Superior Transfer Efficiency

- Higher sensitivity and better transfer efficiency is seen with the Trans-Blot Turbo system in comparison to other blotting techniques. This data set demonstrates the successful transfer of the 1.25 ng protein band only with the Trans-Blot Turbo system.

#### A. Trans-Blot Turbo system



#### B. Tank blot



#### C. Semi-dry blot



#### D. iBlot system



**Superior transfer efficiency.** Serial dilutions of transferrin were separated on a Criterion™ TGX™ 4–20% gel and transferred using four different blotting techniques. **A**, Trans-Blot Turbo system (25 V for 7 min); **B**, Tank blotting (100 V for 30 min); **C**, Semi-dry blotting (25 V for 30 min); **D**, iBlot system (P3 protocol for 7 min).



**Throughput and Modularity**

- **High Throughput** — Up to 4 mini or 2 midi gels can be transferred simultaneously, doubling the throughput of our nearest competitor
- **Modular** — Assemble and run transfers independently with the two cassettes. A single unit and multiple cassettes can be purchased to satisfy a whole lab's blotting requirements



**System Flexibility**

- The Trans-Blot Turbo system accommodates both traditional semi-dry as well as rapid transfers



**Ready-to-use Prepacked Consumables**

- **Convenient, Green Transfer Packs** — Ready-to-use Trans-Blot Turbo transfer packs eliminate extra membrane, filter paper, and buffer preparation. Setup time is reduced to one minute from the opening of the gel cassette to the start of the transfer. Environmentally friendly packaging is recyclable and generates minimal waste; proprietary buffer contains no methanol



**Intuitive Interface**

- Provides customer confidence in protocol selection and execution. Select from optimized preloaded protocols or customize and save/recall up to 25 user-defined transfer protocols

Throughput Comparison based on Transfer Methodology

	Tank	Semi-dry	iBlot system	Trans-Blot Turbo system
# of Mini Blots	2	4	2	4
Transfer Time	30 min +	30 min +	7-10 min	3-10 min*

\* Transfer times are optimized for specific molecular weight ranges

System Flexibility

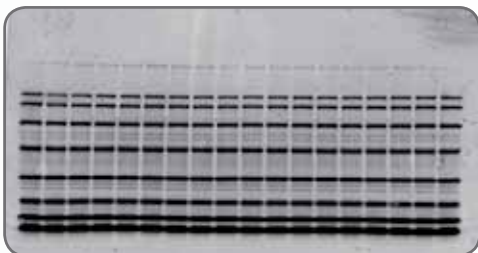
Current Method	Transfer Efficiency	Throughput	Speed
Tank transfer	●		
Semi-dry transfer		●	
Trans-Blot Turbo transfer	●	●	●



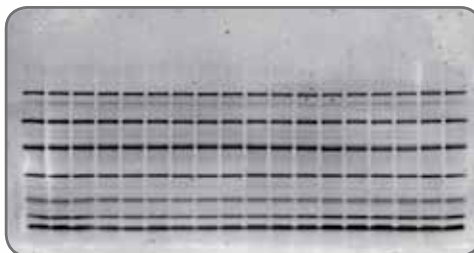
## Universal Rapid Transfer

The Trans-Blot Turbo system was developed to deliver the most uniform transfer for all proteins regardless of molecular weight, post-translational modifications, or protein pI

- **6x stronger signal intensity** — Signal intensities after the transfer were calculated to be 6x stronger with the Trans-Blot Turbo system compared to the iBlot system
- **50% decrease in CV** — CVs across a single blot were 50% lower with the Trans-Blot Turbo system than with the iBlot system



Trans-Blot Turbo system



iBlot system

### Intra-blot CV

Trans-Blot Turbo system

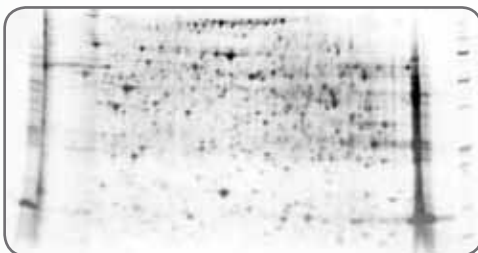
9%

iBlot system

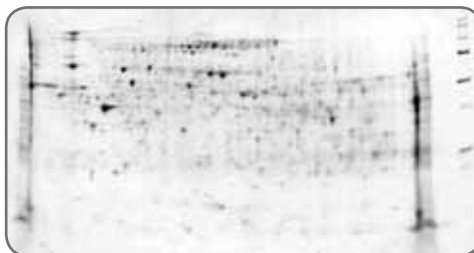
17%

**Reproducibility across blot.** Bio-Rad's SDS-PAGE Broad Range standards were separated on Criterion 4–20% gels, and transferred with the Trans-Blot Turbo and Invitrogen's iBlot systems, both using manufacturers' recommended 7 min protocol. The nitrocellulose membranes were subsequently stained with SYPRO Ruby and imaged on a VersaDoc™ 4000 MP system.

- **2x protein transfer** — Quantitation performed on equivalent 2-D gels transferred with the Trans-Blot Turbo and the iBlot systems demonstrated twice the number of proteins transferred and detected with the Trans-Blot Turbo system



Trans-Blot Turbo system



iBlot system

### 2-D Spot Quantitation

Trans-Blot Turbo system

1066

iBlot system

555

**Higher transfer efficiency using the Trans-Blot Turbo system.** Rat liver extract was focused on Bio-Rad's ReadyStrip™ 11 cm pH 5–8 IPG strips and separated on a Criterion™ TGX™ Any kD™ gel. Duplicate gels were transferred with the Trans-Blot Turbo and iBlot systems, both using manufacturers' recommended 7 min protocol. The nitrocellulose membranes were subsequently stained with SYPRO Ruby and imaged on a VersaDoc 4000 MP system.



## Ordering Information

Catalog #	Description
170-4155	<b>Trans-Blot Turbo Starter System</b>
170-4156	<b>Trans-Blot Turbo Transfer Pack, Mini, PVDF</b> , pkg of 10
170-4157	<b>Trans-Blot Turbo Transfer Pack, Midi, PVDF</b> , pkg of 10
170-4158	<b>Trans-Blot Turbo Transfer Pack, Mini, Nitrocellulose</b> , pkg of 10
170-4159	<b>Trans-Blot Turbo Transfer Pack, Midi, Nitrocellulose</b> , pkg of 10
170-4151	<b>Trans-Blot Turbo Cassette</b> , single
170-4152	<b>Trans-Blot Turbo Base</b> , no cassettes

Bio-Rad Laboratories, Inc. is licensed by Invitrogen Corporation to sell SYPRO products for research use only under U.S. Patent Number 5,616,502. iBlot is a trademark of Life Technologies. Precision Plus Protein standards are sold under license from Life Technologies Corporation, Carlsbad, CA for use only by the buyer of the product. The buyer is not authorized to sell or resell this product or its components.



**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699  
**Canada** 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00  
**France** 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300  
**Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666  
**New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04  
**Singapore** 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55  
**Taiwan** 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000