

MPX™ (MULTIPLEXER) BLOTTING SYSTEM

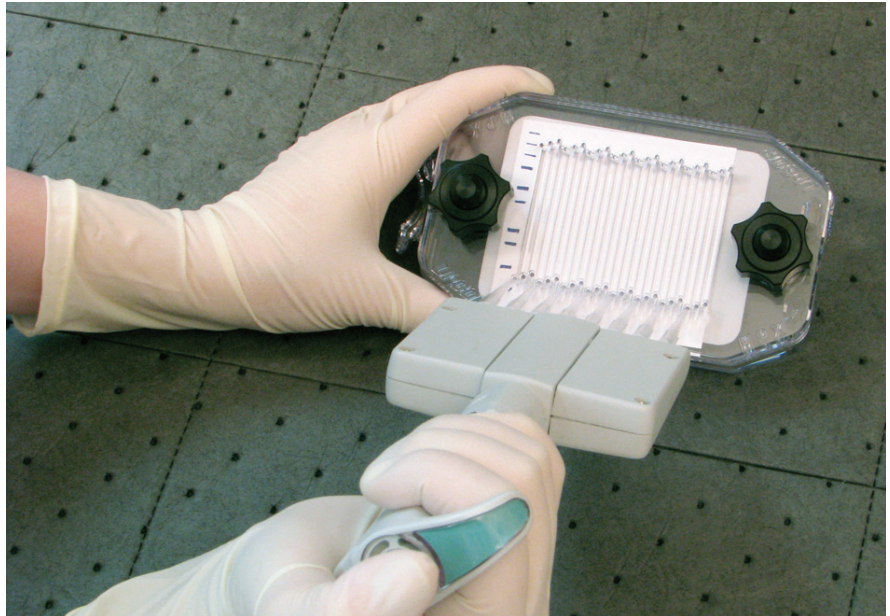
- **Screen multiple targets on the same blot.**
- **Accommodates most common Western blotting procedures.**
- **24 independent channels provide a maximum of 48 targets with infrared fluorescence and the Odyssey® Infrared Imaging System.**
- **Conserve precious antibodies and reagents with lower incubation volumes.**

Screen multiple targets all on the same blot. No cutting. No stripping. No cross-contamination. Adapt your existing Western blotting procedure for comprehensive antibody screening with the MPX Blotting System.

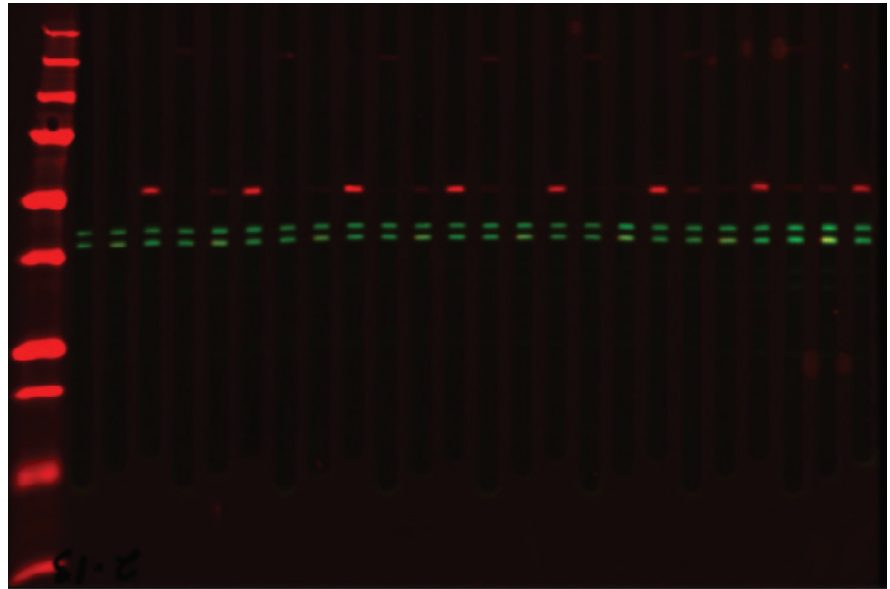
Ideal for any multiple-target Western blot application, the MPX Blotting System increases throughput, maximizes efficiency, and reduces relative overall cost. Quantitative analysis is possible with utilization of an Odyssey Infrared Imager and Odyssey optimized reagents.

Application Versatility

The MPX Blotting System can be incorporated into a variety of assays. Examples include primary and secondary antibody screening, monoclonal antibody screening, signal transduction pathway analysis, and time course studies. However, it is possible to adapt nearly all common Western blotting procedures that generate a PVDF or nitrocellulose blot of 7 x 8.5 cm to the MPX format.



MPX Blotting System



Representative MPX Western Blot with One Sample and Multiple Targets

Simple, Convenient Workflow

Channel ports are conveniently spaced, staggered, and beveled, creating a simple workflow for using both standard and multi-channel pipettes. The low-volume channel ports accommodate a

maximum volume of 160 μ L, thus conserving precious and costly antibody. The nesting dovetail joint design makes it easy to handle multiple-blot processing by allowing the continuous linkage of units.

Screen Up to 48 Targets on One Blot

Processing is as simple as clamping the blot into the MPX™ Blotting System to create up to 24 independent channels. The range of usable channels per sample is relative to comb size. Using LI-COR's selection of IRDye® secondary antibodies on an Odyssey® System provides a maximum of 48 targets on a single membrane – two per channel – and the option for quantitative analysis. Other detection methods, such as standard chemiluminescence and colorimetric methods can provide a maximum of 24 targets.

Compatibility

The MPX Blotting System is compatible with all common Western blotting procedures that generate a nitrocellulose or PVDF blot of 7 cm x 8.5 cm. It can be utilized with all standard chemiluminescence, colorimetric, and infrared protocols. The channel ports are spaced to accommodate most common multi-channel pipettes.

921-00000 MPX (Multiplex) Blotting System

Specifications

Overall Size: 15.5 cm x 8.5 cm x 3.5 cm (L x W x H)

Membrane Size: 7 cm x 8.5 cm

Channel Dimensions: 50 mm x 1.5 mm x 1.5 mm (L x W x H)

Number of Channels: 24

Total Volume per Channel Including Port: 160 µL

Related Products Optimized for use with MPX Blotting System:

921-00200	Single Marker/One Lane Comb
921-00202	Single Marker/Two Lane Comb
921-00204	Single Marker/Three Lane Comb
921-00206	Single Marker/Four Lane Comb
927-40000	Odyssey Blocking Buffer
Multiple P/Ns	IRDye Secondary Antibody Conjugates (IRDye 680RD/680LT/800CW)
926-31090	Odyssey Nitrocellulose Membrane (7 cm x 8.5 cm)
928-40000	Odyssey Protein Molecular Weight Marker
928-40006	IRDye (680/800) Protein Marker
928-40004	4X Protein Sample Loading Buffer
Multiple P/Ns	Running, Transfer, and Wash Buffers
921-11020	MPX Membrane Cushion

Western Blot Cost Comparison

Infrared Detection vs. Chemiluminescence

Reagents	IR Detection (2 Targets)	Chemiluminescence (1 Target)	Chemiluminescence (Strip and reprobe for second target) (2-Target Total)
Secondary Antibody (15 mL) Recommended Dilutions: (1:15,000 for IR#; 1:2,500 for Chemi)	\$0.68	\$0.33	\$0.66
Chemiluminescent Substrate (2 mL)	---	\$5.70 (2 mL)	\$11.40 (2 mL)
Film (2-4 pieces of film/blot)	---	\$7.68	\$15.36
Protein Markers Two-color Protein Marker for IR (2 µL) Standard Protein Marker for Chemi (10 µL)	\$1.16	\$4.68	No charge to reuse marker
Cost	\$1.84	\$18.39 (2 mL)	\$27.42 (2 mL)
Extra Cost Per Blot Compared to IR*		\$16.55	\$25.58

* IRDye 800CW and IRDye 680RD were used for IR calculations

* Based on GE Healthcare pricing, September 20, 2011. Assumes 10 x 10 cm blot.

Moving to digital imaging with Odyssey infrared imaging technology will save money for your lab.

Not only will you save on reagents, all Odyssey imaging systems eliminate the costs related to film and darkroom expenses.

LI-COR is an ISO 9001 registered company. ©2012 LI-COR, Inc. Specifications subject to change. LI-COR, Odyssey, MPX, and IRDye are trademarks or registered trademarks of LI-COR, Inc. in the United States and other countries. All other trademarks belong to their respective owners.



4647 Superior St. • P.O. Box 4000 • Lincoln, Nebraska 68504

LI-COR Biosciences North America: 800-645-4267 / 402-467-0700 • FAX: 402-467-0819 • Technical Support: 800-645-4260

LI-COR GmbH, Germany: Serving Europe, Africa, and the Middle East: +49 (0) 6172 17 17 71

LI-COR Ltd, UK: Serving UK, Ireland and Scandinavia: +44 (0) 1223 422104

In other countries, contact LI-COR Biosciences or a local LI-COR distributor: <http://www.licor.com/distributors>

www.licor.com/bio