

Technical Note

One-Blot Western Optimization Using the MPX™ Blotting System

Developed for:

**Aerius, and Odyssey®
Family of Imaging Systems**

*Please refer to your manual to confirm
that this protocol is appropriate for the
applications compatible with your
Odyssey Imager model.*



LI-COR®

Published May 2009. Revised May 2016.
The most recent version of this
Technical Note is posted at
<http://www.licor.com/bio/support>

Table of Contents

	Page
I. Introduction	2
II. Required Reagents	2
III. Gel Electrophoresis & Transfer	3
IV. Membrane Blocking	4
V. Alignment in MPX Blotter	5
VI. Primary & Secondary Antibody Application	5
VII. Imaging	8

I. Introduction

The independent channels of the LI-COR MPX (Multiplex) Blotter make it possible to optimize blocking buffer, primary antibody dilution, and secondary antibody dilution on a single Western blot. Western blotting procedures that generate a blot of 7.0 x 8.5 cm are easily adapted to the MPX format. The process fits into any laboratory's standard Western blot workflow. Both home-made and pre-cast gels can be used to generate blots.

Electrophoresis and transfer to nitrocellulose membrane are performed under standard conditions. Clamping the blot into the MPX Blotter creates up to 24 independent channels, allowing different conditions to be tested in each channel. The range of usable channels per sample is relative to comb size. For Western blot optimization, a single-well gel ("prep gel") is all that is needed. For this application, any detection method can be used, including near-infrared (NIR) fluorescence and chemiluminescence. This document presents general guidelines for use with the Odyssey® family of Infrared Imaging Systems.

II. Required Reagents

		LI-COR P/N
Sample Preparation	4X Protein Sample Loading Buffer	928-40004
Electrophoresis	Odyssey One-Color Protein Markers (Molecular Weight - 10 kDa to 250 kDa)	928-40000
	Chameleon® Duo Pre-stained Protein Ladder (8 kDa to 260 kDa)	928-60000
Blotting and Transfer	Tris Glycine	
	Odyssey Nitrocellulose (10 membranes; 7 x 8.5 cm) or	926-31090
	Odyssey Nitrocellulose (1 roll; 30 cm x 3 m)	926-31092

		LI-COR P/N
MPX Detection	Blocking Buffer Optimization Kit	927-40040
	• Odyssey® Blocking Buffer (TBS)	927-50000; 927-50100
	• Odyssey Blocking Buffer (PBS)	927-40000; 927-40100
	• Casein Blocking Buffer	927-40200; 927-40300
	• Blocking buffer of your choice (BSA, Milk, etc.)	
	IRDye® Labeled Secondary Antibodies	LI-COR
	NewBlot™ IR Stripping Buffer, 5X	928-40028
	• 10X PBS	
	• 10X TBS	
	Tween® 20	
	MPX Membrane Cushion	921-00120
Imaging	Odyssey Family Imager or Aeries Imager	

III. Gel Electrophoresis and Transfer

■ Gel Preparation

A wide variety of gel matrices are compatible with the MPX Blotter system.

If you are pouring your own gels, your gel casting system can be used with a single-well comb such as the LI-COR Single Marker/One Lane Comb (921-00200, 1 mm thickness).

Alternatively, pre-cast gels can be purchased and used. Table 1 lists several suggested types of pre-cast gels, and indicates the number of usable ports that each gel will provide for use with the MPX Blotter.

Table 1. Single-sample pre-cast gel options for use with the MPX Blotter.

Vendor	Well Designation	Sample #	MW Marker Well	Usable Ports
Invitrogen	2D	1	Yes	19
Bio-Rad	2D/Prep	1	Yes	21
C.B.S. Scientific	1 Well	1	No	23

■ Sample Preparation

When using a single-well gel, a larger volume of sample is required. Prepare your protein sample so that the sample volume and concentration is equivalent to running all the lanes on a standard 10-well gel. Example: 5 µg of lysate per lane = 50 µg in a total volume of 100 - 150 µL, including loading buffer.

The following procedure is suggested: Dilute the sample 1:4 in 4X Protein Sample Loading Buffer (LI-COR P/N 928-40004) with β -Mercaptoethanol. Heat the sample at 95 °C for 5 minutes.

■ Molecular Weight Marker

It is important to have a molecular weight marker that is visible to the eye because the marker is the primary tool used to align the blot in the MPX Blotter. Odyssey® One-Color Marker or Chameleon® Duo Protein Ladder is recommended.

■ Electrophoresis

IMPORTANT: The maximum length of the separating gel should not exceed 50 mm—the length of the channels on the MPX Blotter.

■ Transfer

- Always use clean forceps when handling membranes. Nitrocellulose membrane is recommended for this procedure.
- Once electrophoresis is complete, transfer proteins to Odyssey Nitrocellulose Membrane using standard transfer procedures.
- Mark the outside corners of the gel and sample wells with a pencil before separating the transferred gel from the membrane, as shown in Figure 1. The marks will help you correctly align the membrane when it is placed in the MPX Blotter.
- Allow the membrane to dry for a minimum of one hour before proceeding with detection. *IMPORTANT! Use pencil to mark the blot. Ink from most pens will fluoresce on the Odyssey Imager and cause increased membrane background.*

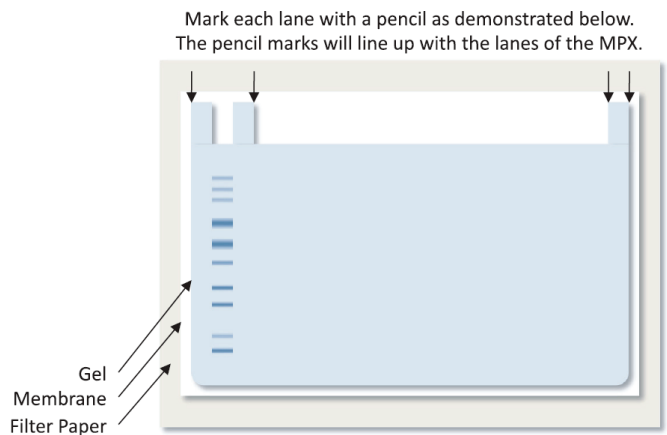


Figure 1. Mark the membrane with pencil, for later alignment into the MPX Blotter.

IV. Membrane Blocking

■ Membrane Preparation

- Cut the membrane into four individual blots, as shown in Figure 2.
- Each individual blot will be processed with a different blocking buffer, and that blocking buffer will also be used for dilution of antibodies. TBS or PBS buffer systems may be used for blocking. **During washing steps, rinse and wash each blot with an appropriate wash buffer that matches the buffer system used for blocking.**

Blot 1: Odyssey Blocking Buffer (TBS)

Blot 2: Odyssey Blocking Buffer (PBS)

Blot 3: Casein Blocking Buffer

Blot 4: Blocking buffer of your choice (milk, BSA, etc. in TBS or PBS)

- Pre-wet each membrane with TBS or PBS buffer as appropriate (see above).

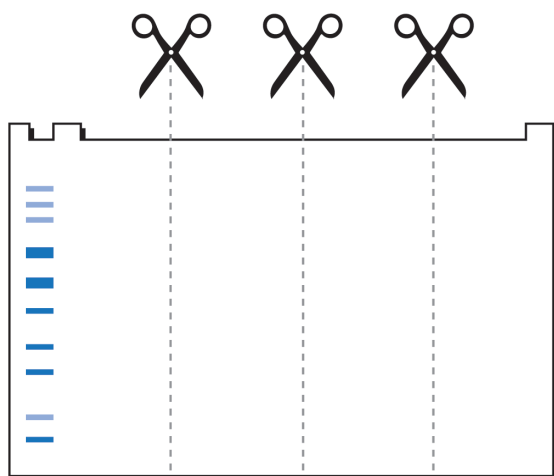


Figure 2. Cut the membrane into four individual blots for blocking buffer optimization.

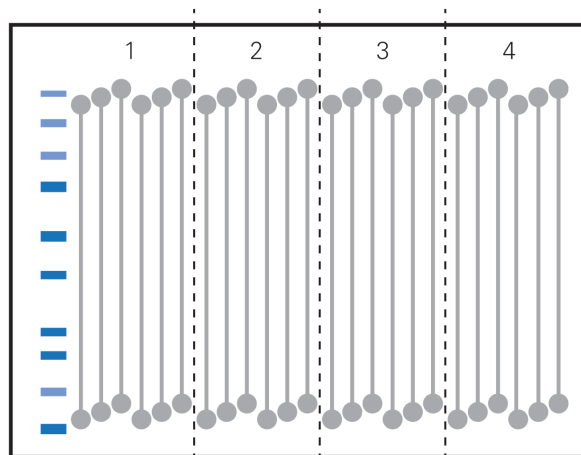


Figure 3. Place four individual blots into the MPX Blotter as shown.

■ Blocking

Place the membranes into 4 different incubation boxes. In each box, cover the entire membrane with blocking buffer (approximately 0.4 mL/cm²), using a different blocking buffer for each membrane. Block the membrane for 1 hour at room temperature with gentle shaking.

- Blot 1: Odyssey® Blocking Buffer (TBS)
- Blot 2: Odyssey Blocking Buffer (PBS)
- Blot 3: Casein Blocking Buffer
- Blot 4: Blocking buffer of your choice (milk, BSA, etc.)

V. Alignment in MPX Blotter

Detailed instructions for use of the MPX Blotter are found in the MPX Blotter Multiplex Western Blotting Accessory User Guide at <http://www.licor.com/mpxuserguide>

Place the four blocked membranes into the MPX Blotter so that there are at least 4 channels available for use on each membrane. See Figure 3.

VI. Primary & Secondary Antibody Application

■ Primary Antibody Preparation

Two dilutions of primary antibody should be made for each blocking buffer that is tested. Suggested starting dilutions are 1:500 and 1:1,000. You may wish to modify these dilutions, based on vendor recommendations.*

700 µL of each dilution will be needed. Dilute the primary antibody in the appropriate blocking buffer (see following table) with 0.2% Tween® 20.

**The correct working range for antibody dilution depends on the characteristics of your primary antibody. Start with the dilution recommended by the primary antibody vendor for Western blot applications.*

Blot	Blocker	Primary Antibody Dilutions	
1	Odyssey® Blocking Buffer (TBS)	1:500	1:1,000
2	Odyssey Blocking Buffer (PBS)	1:500	1:1,000
3	Casein Blocking Buffer	1:500	1:1,000
4	Blocking buffer of your choice (milk, BSA, etc.)	1:500	1:1,000

■ Primary Antibody Application

- Load the primary antibody/blocker dilutions into the MPX Blotter in the indicated locations for each blocking buffer you are testing.
 - Apply 2 replicates of each primary antibody dilution, as shown in Figure 4.
- Fill the unused channels with the appropriate corresponding blocking buffer.
- Incubate for 1 - 4 hours at room temperature.

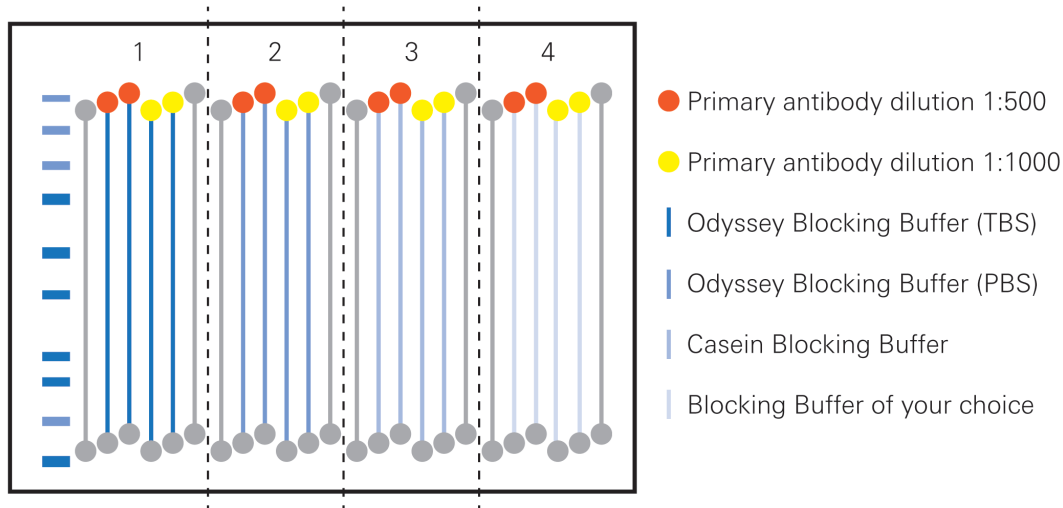


Figure 4. Placement of primary antibody dilutions in the channels of the MPX Blotter.

- Wash primary antibody from the channels thoroughly according to MPX Blotter manual instructions, using a buffer that matches the buffer system used for blocking. Wash buffers should contain 0.1% Tween® 20. **Do not remove blot from MPX blotting manifold during washing.**

Blot 1: TBS-T
 Blot 2: PBS-T
 Blot 3: PBS-T
 Blot 4: TBS-T or PBS-T, as appropriate

■ Secondary Antibody Preparation

Blot	Blocker	Secondary Antibody Dilutions	
1	Odyssey® Blocking Buffer (TBS)	1:5,000	1:10,000
2	Odyssey Blocking Buffer (PBS)	1:5,000	1:10,000
3	Casein Blocking Buffer	1:5,000	1:10,000
4	Blocking buffer of choice	1:5,000	1:10,000

Two dilutions of secondary antibody should be made for each blocking buffer that is tested. For IRDye® secondary antibodies, we recommend 1:5,000 and 1:10,000 as a starting point. Dilutions may be modified, based on vendor recommendations.

700 µL of each antibody will be needed. Dilute the secondary antibody in the appropriate blocking buffer with 0.2% Tween® 20.

■ Secondary Antibody Application

- Load the secondary antibody/blocker dilutions into the MPX Blotter in the indicated locations for each blocking buffer you are testing.
 - Add the secondary antibody dilutions to the channels previously stained with primary antibody, as shown in Figure 5.
- Fill the unused channels with the appropriate corresponding blocking buffer.
- Incubate 1 hour at room temperature. ***Protect from light during incubation.***

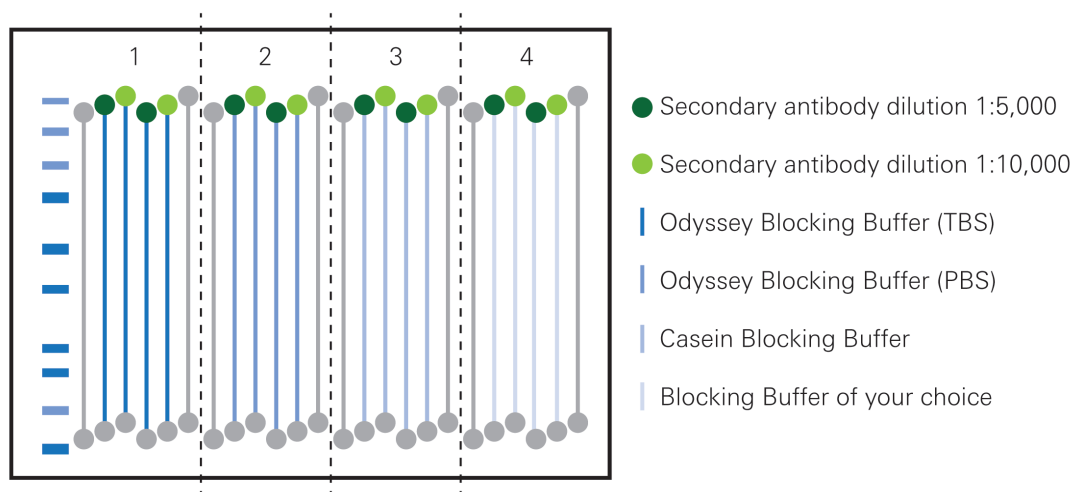


Figure 5. Placement of secondary antibody dilutions in the channels of the MPX Blotter.

VII. Imaging

Membranes can be imaged immediately.

- Image all four blots side-by-side, using standard Western blot imaging settings on any Odyssey® Family Imaging System.
- Visual inspection of images with Image Studio™ software or Odyssey application software can be used to determine which blocking buffer works best for the primary antibody you are testing.
 - View all blots together in a single image, with uniform image display settings, to compare membrane background levels and band intensity.
 - Individually adjust the image display settings for each blot to get the “best” image.
 - Evaluate non-specific banding in each blocking buffer condition.
- Look for blocking buffer conditions that provide strong signals for the expected band(s), low membrane background, and few non-specific background bands from the primary antibody.
 - Tradeoffs may be necessary. Blocking conditions that yield very strong bands might also have higher membrane background or non-specific banding.
- The “best” blocking conditions depend on the antigen-antibody pair you are using. Some primary antibodies are dramatically affected by blocking conditions. An inappropriate blocker can alter binding specificity, affecting the intensity of target bands and increasing non-specific banding. The pattern of non-specific bands may also be affected.
- Choose the blocking conditions that are most appropriate for the context and goals of your experiment.
- Quantitative analysis of specific bands on each blot will indicate if signal intensity (after background subtraction) is significantly different between blocker types.

© 2016 LI-COR, Inc. LI-COR, Odyssey, MPX, Image Studio, Chameleon, NewBlot, 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.

LI-COR is an ISO 9001 registered company.



4647 Superior Street • Lincoln, NE 68504
Toll free: 800-645-4267 • Fax: +1-402-467-0819
biosales@licor.com

LI-COR GmbH, Germany • Serving Europe, Africa, and the Middle East.
LI-COR Biosciences GmbH • Siemensstraße 25A • 61352 Bad Homburg • Germany
Phone: +49 (0) 6172 17 17 771 • Fax: +49 (0) 6172 17 17 799 • bio-eu@licor.com

LI-COR Ltd., United Kingdom • Serving Denmark, Ireland, Finland, Iceland, Norway, Sweden, and UK.
LI-COR Biosciences UK Ltd. • St. John's Innovation Centre • Cowley Road
Cambridge CB4 0WS • United Kingdom
Phone: +44 (0) 1223 422104 • Fax: +44 (0) 1223 422105 • bio-eu@licor.com