

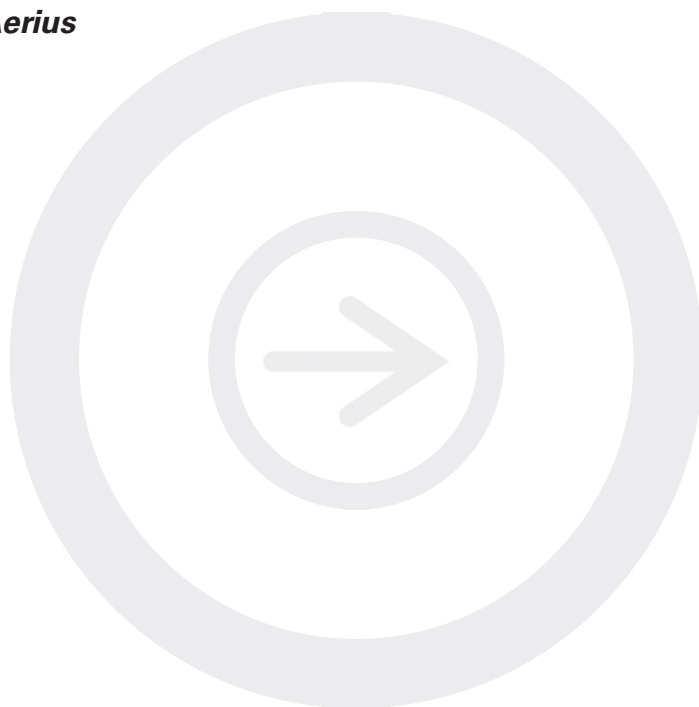
# Western Blot Analysis

---

Developed for:

**Aerius, and Odyssey® Family of Imagers**

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



Published 2003, revised January 2012. The most recent version of this document is posted at: <http://biosupport.licor.com>

## Table of Contents

		Page
I.	<b>Required Reagents</b> .....	.2
II.	<b>Western Detection Methods</b> .....	.2
III.	<b>Guidelines for Two-Color Detection</b> .....	.5
IV.	<b>Stripping the Membrane</b> .....	.6
V.	<b>Adapting Western Blotting Protocols for Odyssey® Detection</b> .....	.6
VI.	<b>General Tips</b> .....	.7
VII.	<b>Imaging of Coomassie-Stained Protein Gels</b> .....	.8
VIII.	<b>Troubleshooting Guide</b> .....	.9

### I. Required Reagents

- Blotted nitrocellulose (LI-COR, P/N 926-31090) or low-fluorescent PVDF membrane (LI-COR, P/N 926-31098)
- Odyssey Blocking Buffer (LI-COR, P/N 927-40000)
- Odyssey Pre-stained Molecular Weight Marker (LI-COR, P/N 928-40000)
- IRDye (680/800) Protein Markier (LI-COR, P/N 928-40006)
- Primary antibodies
- IRDye® secondary antibodies (LI-COR)
- Tween® 20
- PBS wash buffer (LI-COR, P/N 928-40018 or 928-40020)
- Ultrapure water
- Methanol for wetting of PVDF
- SDS (if desired)
- Other blocking buffers (if desired)
- NewBlot™ Stripping Buffer, if desired, for nitrocellulose (LI-COR, P/N 928-40030) or PVDF (LI-COR, P/N 928-40032) membranes

### II. Western Detection Methods

Nitrocellulose or PVDF membranes may be used for protein blotting. Pure cast nitrocellulose is generally preferable to supported nitrocellulose. Protein should be transferred from gel to membrane by standard procedures. Membranes should be handled only by their edges, with clean forceps.

After transfer, perform the following steps:

1. Wet the membrane in PBS for several minutes. If using a PVDF membrane that has been allowed to dry, pre-wet briefly in 100% methanol and rinse with ultrapure water before incubating in PBS.

**NOTES:**

- *Ink from most pens and markers will fluoresce on all Aeries or Odyssey Imagers. The ink may wash off and re-deposit elsewhere on the membrane, creating blotches and streaks. Pencil should be used to mark membranes. (The Odyssey pen doesn't fluoresce and can be used with nitrocellulose membranes, since the membrane will not be soaked in methanol, causing the ink to run.)*

- Block the membrane in Odyssey® Blocking Buffer for 1 hour. Be sure to use sufficient blocking buffer to cover the membrane (a minimum of 0.4 mL/cm<sup>2</sup> is suggested).

**NOTES:**

- *Membranes can be blocked overnight at 4°C if desired.*
- *Odyssey Blocking Buffer often yields higher and more consistent sensitivity and performance than other blockers. Nonfat dry milk or casein dissolved in PBS can also be used for blocking and antibody dilution, but be aware that milk may cause higher background on PVDF membranes. If using casein, a 0.1% solution in 0.2X PBS buffer is recommended (Hammersten-grade casein is not required).*
- *Milk-based reagents can interfere with detection when using anti-goat antibodies. They also deteriorate rapidly at 4°C, so diluted antibodies should not be kept and re-used.*
- *Blocking solutions containing BSA can be used, but in some cases they may cause high membrane background. **BSA-containing blockers are not generally recommended** and should be used only when the primary antibody requires BSA as blocker.*

- Dilute primary antibody in Odyssey Blocking Buffer. Optimum dilution depends on the antibody and should be determined empirically. A suggested starting range can usually be found in the product information from the vendor. To lower background, add 0.1 - 0.2% Tween® 20 to the diluted antibody before incubation. The optimum Tween 20 concentration will depend on the antibody.

**NOTES:**

- *Two-color detection requires careful selection of primary and secondary antibodies. For details, see III. Guidelines for Two Color Western Detection.*
- *The MPX™ Blotting System can be used to efficiently determine the optimum antibody concentration. For details, search for One Blot Western Optimization Using the MPX Blotting System at <http://biosupport.licor.com>.*

- Incubate blot in primary antibody for 60 minutes or longer at room temperature with gentle shaking. Optimum incubation times vary for different primary antibodies. Use enough antibody solution to completely cover the membrane.

- Wash membrane 4 times for 5 minutes each at room temperature in PBS + 0.1% Tween 20 with gentle shaking, using a generous amount of buffer.

- Dilute the fluorescently-labeled secondary antibody in Odyssey Blocking Buffer. **Recommended dilution can be found in the pack insert for the IRDye® conjugate.** Add the same amount of Tween 20 to the diluted secondary antibody as was added to the primary antibody.

**NOTES:**

- *Avoid prolonged exposure of the antibody vial to light.*
- *Be careful not to introduce contamination into the antibody vial.*
- *For best sensitivity and performance, use freshly diluted antibody solution.*
- *Adding 0.01% - 0.02% SDS to the diluted secondary antibody (in addition to Tween 20) will substantially reduce membrane background, particularly when using PVDF. However, DO NOT add SDS during blocking or to the diluted primary antibody. See **V. Adapting Western Blotting Protocols for Odyssey Detection** for more information about how and why to use SDS in the secondary antibody incubation.*
- *The MPX™ Blotting System can be used to efficiently determine the optimum antibody concentration. For details, search for One Blot Western Optimization Using the MPX Blotting System at <http://biosupport.licor.com>.*

7. Incubate blot in secondary antibody for 30-60 minutes at room temperature with gentle shaking. Protect from light during incubation.

**NOTES:**

- *Incubating more than 60 minutes may increase background.*
8. Wash membrane 4 times for 5 minutes each at room temperature in PBS + 0.1% Tween<sup>®</sup> 20 with gentle shaking. Protect from light.
  9. Rinse membrane with PBS to remove residual Tween 20. The membrane is now ready to scan.
  10. Image on Aerius, or Odyssey<sup>®</sup> Family of Imagers.
    - Scan in the appropriate channels.
    - Protect the membrane from light until it has been imaged.
    - Keep the membrane wet to strip and re-use it. Once a membrane has dried, stripping is ineffective.
    - Blots can be allowed to dry before scanning, if desired. Signal strength may be enhanced on a dry membrane. The membrane can also be re-wetted for imaging.
    - The fluorescent signal on the membrane will remain stable for several months, or longer, if protected from light. Membranes may be stored dry at 4°C.
    - If signal on membrane is too strong or too weak, re-image the membrane at a lower or higher scan intensity setting, respectively. Adjust image acquisition time for Odyssey Fc. AutoMode in Odyssey CLx may be used to improve the dynamic range of the image.

### ***Molecular Weight Marker***

If you loaded the Odyssey Prestained Molecular Weight Marker (LI-COR, P/N 928-40000), it will be visible in the 700 nm channel and also faintly visible in the 800 nm channel. If using IRDye (680/800) Protein Marker, it will be visible in both the 700 nm and 800 nm channels. Pre-stained blue molecular weight markers from other sources can also be used. Load 1/3 to 1/5 of the amount you would normally use for Western transfer. Too much marker can result in very strong marker bands that may interfere with visualization of sample lanes. If using multicolored markers, some bands may not be visualized.

### ***Optimization Tips***

1. ***Follow the protocol carefully.***
2. No single blocking reagent will be optimal for every antigen-antibody pair. Some primary antibodies may exhibit greatly reduced signal or different non-specific binding in different blocking solutions. If it is difficult to detect the target protein, changing the blocking solution may dramatically improve performance. If the primary antibody has worked well in the past using chemiluminescent detection, try that blocking solution for infrared detection.
3. Addition of detergent such as Tween<sup>®</sup> 20 can reduce membrane background and non-specific binding. Refer to **V. Adapting Western Blotting Protocols for Odyssey Detection** for details.
4. To avoid background speckles on blots, use ultrapure water for buffers and rinse plastic dishes well before and after use. Never perform Western incubations or washes in dishes that have been used for Coomassie staining.
5. Membranes should be handled only by their edges, with clean forceps.

6. After handling membranes that have been incubating in antibody solutions, clean forceps thoroughly with distilled water and/or ethanol. If forceps are not cleaned after being dipped in antibody solutions, they can cause spots or streaks of fluorescence on the membrane that are difficult to wash away.
7. Do not wrap the membrane in plastic when scanning.
8. If a Western blot will be stripped, do not allow the membrane to dry. Stripping is ineffective once a membrane has dried, or even partially dried.

### III. Guidelines for Two-Color Detection

Two different antigens can be detected simultaneously on the same blot using antibodies labeled with near-infrared dyes that are visualized in different fluorescence channels (700 and 800 nm). Two-color detection requires careful selection of primary and secondary antibodies.

The following guidelines will help design two-color experiments:

1. If the two primary antibodies are derived from different host species (for example, primary antibodies from mouse and chicken), IRDye® whole IgG secondary antibodies derived from the same host and labeled with different IRDye fluorophores must be used (for example, IRDye 800CW Donkey anti-Mouse and IRDye 680LT or IRDye 680RD Donkey anti-Chicken).
2. If the two primary antibodies are monoclonals (mouse) and are IgG<sub>1</sub>, IgG<sub>2a</sub>, or IgG<sub>2b</sub>, IRDye Subclass Specific secondary antibodies must be used. The same subclasses cannot be combined in a two-color Western blot (for example, two IgG<sub>1</sub> primary antibodies). For details, refer to *Western Blot and In-Cell Western™ Assay Detection Using IRDye Subclass Specific Antibodies*.
3. Before combining primary antibodies in a two-color experiment, always perform preliminary blots with each primary antibody alone to determine the expected banding pattern and possible background bands. Slight cross-reactivity may occur and can complicate interpretation of a blot, particularly if the antigen is very abundant. If cross-reactivity is a problem, load less protein or reduce the amount of antibody.
4. One secondary antibody must be labeled with a 700 nm channel dye, and the other with an 800 nm channel dye.
5. Always use highly cross-adsorbed secondary antibodies for two-color detection. Failure to use highly cross-adsorbed antibodies may result in cross-reactivity.
6. For best results, avoid using primary antibodies from mouse and rat together for a two-color experiment. It is not possible to completely adsorb away cross-reactivity because the species are so closely related. If using mouse and rat together, it is crucial to run single-color blots first with each individual antibody to be certain of expected band sizes.

When performing a two-color blot, use the standard Western blot protocol with the following modifications:

- Combine the two primary antibodies in the diluted antibody solution in step 3. Incubate simultaneously with the membrane (step 4).
- Combine the two dye-labeled secondary antibodies in the diluted antibody solution in step 6. Incubate simultaneously with the membrane (step 7).

#### **IV. Stripping the Membrane**

Typically, both PVDF and nitrocellulose membranes can be stripped up to three times. LI-COR® NewBlot™ Stripping Buffer is available under P/N 928-40030 for nitrocellulose or 928-40032 for PVDF. If a blot is to be stripped, DO NOT allow it to dry before, during, or after imaging (keep the blot as wet as possible). Complete usage instructions are given in the NewBlot Stripping Buffer pack insert that is shipped with the product. Before proceeding, read the instructions in the pack insert, including the frequently asked questions.

#### **V. Adapting Western Blotting Protocols for Detection with Odyssey® Systems**

When adapting Western blotting protocols for Odyssey detection or using a new primary antibody, it is important to determine the optimal antibody concentrations. Optimization will help achieve maximum sensitivity and consistency. Three parameters should be optimized: primary antibody concentration, dye-labeled secondary antibody concentration, and detergent concentration in the diluted antibodies.

##### ***Primary Antibody Concentration***

Primary antibodies vary widely in quality, affinity, and concentration. The correct working range for antibody dilution depends on the characteristics of the primary antibody and the amount of target antigen to detect. Suggested dilutions are 1:500, 1:1500, 1:5000, and 1:10,000 (start with the dilution factor normally used for chemiluminescent detection, or refer to the product information from the vendor). Use the MPX™ Blotting System to optimize the primary dilution to achieve maximum performance and conserve antibody (refer to *One Blot Western Optimization Using the MPX Blotting System* at <http://biosupport.licor.com>).

##### ***Secondary Antibody Concentration***

Optimal dilutions of dye-conjugated secondary antibodies should also be determined. Refer to the IRDye® conjugate pack insert for recommendations. The amount of secondary required varies depending on how much antigen is being detected – abundant proteins with strong signals require less secondary antibody. Use the MPX Blotting system to optimize (search for *One Blot Western Optimization Using the MPX Blotting System* at <http://biosupport.licor.com>).

##### ***Detergent Concentration***

Addition of detergents to diluted antibodies can significantly reduce background on the blot. Optimal detergent concentration will vary, depending on the antibodies, membrane type, and blocker used. Keep in mind that some primaries do not bind as tightly as others and may be washed away by too much detergent. Never expose the membrane to detergent until blocking is complete, as this may cause high membrane background.

**Tween® 20:**

- Add Tween 20 to both the primary antibody and secondary antibody solutions when the antibodies are diluted in blocking buffer. A final concentration of 0.1 - 0.2% is recommended for nitrocellulose membranes, and a final concentration of 0.1% is recommended for PVDF membranes (higher concentrations of Tween 20 may actually cause increased background on PVDF).
- Wash solutions should contain 0.1% Tween 20.

**SDS:**

- Adding 0.01 - 0.02% SDS to the diluted secondary antibody can dramatically reduce overall membrane background and also reduce or eliminate non-specific binding. It is critical to use only a very small amount, because SDS is an ionic detergent and can disrupt antigen-antibody interactions if too much is present at any time during the detection process.
- Addition of SDS is particularly helpful for reducing the higher overall background that is seen with PVDF membrane. When working with **IRDye® 680LT** conjugates on **PVDF membranes**, SDS (final concentration of 0.01 - 0.02%) and Tween 20 (final concentration of 0.1 - 0.2%) **must** be added during detection incubation step to avoid non-specific background staining.
- DO NOT add SDS during the blocking step or to the diluted primary antibody. Presence of SDS during binding of the primary antibody to its antigen may greatly reduce signal. Add SDS only to the diluted secondary antibody.
- Wash solutions should contain 0.1% Tween 20, but no SDS.
- Some antibody-antigen pairs may be more sensitive to the presence of SDS and may require even lower concentrations of this detergent (less than 0.01%) for best performance. Titrate the amount of SDS to find the best level for the antibodies used.
- If high background is seen when using BSA-containing blocking buffers, adding SDS to the secondary antibody may alleviate the problem.

## VI. General Tips

1. Milk-based blockers may contain IgG that can cross-react with anti-goat antibodies. This can significantly increase background and reduce antibody titer. Milk may also contain endogenous biotin or phospho-epitopes that can cause higher background.
2. Store the IRDye secondary antibody vial at 4°C in the dark. IRDye secondary antibodies may be aliquoted and frozen for long-term storage. Minimize exposure to light and take care not to introduce contamination into the vial. Dilute immediately prior to use. If particulates are seen in the antibody solution, centrifuge before use.
3. Protect membrane from light during secondary antibody incubations and washes.
4. Use the narrowest well size possible for the loading volume to concentrate the target protein.
5. The best transfer conditions, membrane, and blocking agent for experiments will vary, depending on the antigen and antibody. If there is high background or low signal level, a good first step is to try a different blocking solution.
6. Small amounts of purified protein may not transfer well. Adding non-specific proteins of similar molecular weight can have a “carrier” effect and substantially increase transfer efficiency.

7. For proteins <100 kDa, try blotting in standard Tris-glycine buffer with 20% methanol and no SDS. Addition of SDS to the transfer buffer can greatly reduce binding of transferred proteins to the membrane (for both PVDF and nitrocellulose).
8. Soak the gel in transfer buffer for 10-20 minutes before setting up the transfer. Soaking equilibrates the gel and removes buffer salts that will be carried over into the transfer tank.
9. To maximize retention of transferred proteins on the membrane, allow the membrane to air-dry completely after transfer (approximately 1-2 hours).
10. Do not over-block. Long blocking incubations, particularly with nonfat dry milk at 2% or higher, can cause loss of target protein from the membrane (*J. Immunol. Meth.* 122:129-135, 1989).
11. To enhance signal, try extended primary antibody incubation at room temperature or overnight incubation at 4°C. Avoid extended incubations in secondary antibody.

## VII. Imaging of Coomassie-Stained Protein

IRDye<sup>®</sup> Blue Protein Stain is a convenient, safe alternative for gel staining to provide confirmation of protein transfer to the membrane. Unlike traditional Coomassie Blue stains, which require methanol and acetic acid for staining and destaining, IRDye Blue Protein Stain is water-based and requires no hazardous solvents. This stain offers excellent detection sensitivity in the 700 nm channel of Aerius and Odyssey<sup>®</sup> imaging systems (< 5 ng of BSA can be detected). IRDye Blue Protein Stain is Coomassie-based and is provided as a ready-to-use 1X solution. Pre-washing and de-staining steps are performed in water.

1. Wash gels with ultrapure water for 15 minutes.
2. Submerge gel in IRDye Blue Stain for 1 hour.
3. Destain with ultrapure water for 30 minutes or overnight if needed.
4. Scan on an Aerius or Odyssey imaging system (Odyssey Classic, Odyssey CLx, Odyssey Sa, or Odyssey Fc) in the 700 nm channel only. *Please refer to the Tutorial Manual of each instrument for further information.*
  - If using Odyssey 3.0 software, select the Protein Gel scan preset.
  - If using Aerius and/or Odyssey Sa software, set the focus offset to 3.0 plus one-half the thickness of the gel.
  - In Image Studio, select Protein Gel, using the Custom setting, under the setup tab.

## VIII. Troubleshooting Guide

Problem	Possible Cause	Solution / Prevention	
High background, uniformly distributed.	BSA used for blocking.	Blocking solutions containing BSA may cause high membrane background. Try adding SDS to reduce background, or switch to a different blocker.	
	Not using optimal blocking reagent.	Compare different blocking buffers to find the most effective; try blocking longer.	
	Background on nitrocellulose.	Add Tween <sup>®</sup> 20 to the diluted antibodies to reduce background. Try adding SDS to diluted secondary antibody.	
	Background on PVDF.	Use low-fluorescent PVDF membrane. With IRDye <sup>®</sup> 680LT conjugates, always use SDS (0.01-0.02% final concentration) and Tween 20 (0.1-0.2% final) during the detection incubation step.	
	Antibody concentrations too high.	Optimize primary and secondary antibody dilutions using MPX <sup>™</sup> blotting system. For details, see One Blot Western Optimization Using the MPX Blotting System at <a href="http://biosupport.licor.com">http://biosupport.licor.com</a>	
	Insufficient washing.		Increase number of washes and buffer volume.
			Make sure that 0.1% Tween 20 is present in buffer and increase if needed. Note that excess Tween 20 (0.5-1%) may decrease signal.
	Cross-reactivity of antibody with contaminants in blocking		Use Odyssey Blocking Buffer instead of milk. Milk usually contains IgGs that cross-react with anti-goat secondary antibodies.
	Inadequate antibody volume used.		Increase antibody volume so entire membrane surface is sufficiently covered with liquid at all times (use heat-seal bags if volume is limiting). Do not allow any area of membrane to dry out.
Use agitation for all antibody incubations.			
Membrane contamination.		Always handle membranes carefully and with clean forceps. Do not allow membrane to dry. Use clean dishes, bags, or trays for incubations.	

<b>Problem</b>	<b>Possible Cause</b>	<b>Solution / Prevention</b>
Uneven blotchy or speckled background.	Blocking multiple membranes together in small volume.	If multiple membranes are being blocked in the same dish, ensure that blocker volume is adequate for all membranes to move freely and make contact with liquid.
	Membrane not fully wetted or allowed to partially dry.	Keep membrane completely wet at all times. This is particularly crucial if blot will be stripped and re-used.
		If using PVDF, remember to first pre-wet in 100% methanol.
	Contaminated forceps or dishes.	Always carefully clean forceps after they are dipped into an antibody solution, particularly dye-labeled secondary antibody. Dirty forceps can deposit dye on the membrane that will not wash away.
		Use clean dishes, bags or trays for incubations.
	Dirty scanning surface or silicone mat.	Clean scanning surface and mat carefully before each use. Dust, lint, and residue will cause speckles.
Incompatible marker or pen used to mark membrane.	Use only pencil or Odyssey® pen (nitrocellulose only) to mark membranes.	
Weak or no signal.	Not using optimal blocking reagent.	Primary antibody may perform substantially better with a different blocker.
	Insufficient antibody used.	Primary antibody may be of low affinity. Increase amount of antibody or try a different source.
		Extend primary antibody incubation time (try 4 - 8 hr at room temperature, or overnight at 4°C).
		Increase amount of primary or secondary antibody, optimizing for best performance.
		Try substituting a different dye-labeled secondary antibody.
Primary or secondary antibody may have lost reactivity due to age or storage conditions.		
Too much detergent present; signal being washed away.	Decrease Tween® 20 and/or SDS in diluted antibodies. Recommended SDS concentration is 0.01 - 0.02%, but some antibodies may require an even lower concentration.	

<b>Problem</b>	<b>Possible Cause</b>	<b>Solution / Prevention</b>
Weak or no signal. (Continued)	Insufficient antigen loaded.	Load more protein on the gel. Try using the narrowest possible well size to concentrate antigen.
	Protein did not transfer well.	Check transfer buffer choice and blotting procedure.
		Use pre-stained molecular weight marker to monitor transfer, and stain gel after transfer to make sure proteins are not retained in gel.
	Protein lost from membrane during detection.	Extended blocking times or high concentrations of detergent in diluted antibodies may cause loss of antigen from the blotted membrane.
	Proteins not retained on membrane during transfer.	Allow membrane to air dry completely (1 - 2 hr) after transfer. This helps make the binding irreversible.
		Addition of 20% methanol to transfer buffer may improve antigen binding. Note: Methanol decreases pore size of gel and can hamper transfer of large proteins.
SDS in transfer buffer may interfere with binding of transferred proteins, especially for low molecular weight proteins. Try reducing or eliminating SDS. Note: Presence of up to 0.05% SDS does improve transfer efficiency of some proteins.		
Non-specific or unexpected bands.	Antibody concentrations too high.	Reduce the amount of antibody used. Reduce antibody incubation times. Increase Tween® 20 in diluted antibodies.
		Add or increase SDS in diluted secondary antibodies.
	Not using optimal blocking reagent.	Choice of blocker may affect background bands. Try a different blocker.
	Cross-reactivity between antibodies in a two-color experiment.	Double-check the sources and specificities of the primary and secondary antibodies used (see III. Guidelines for Two-Color Detection).
Use only highly cross-adsorbed secondary antibodies.		

Cause	Possible Cause	Solution / Prevention
Non-specific or unexpected bands. <i>(Continued)</i>		There is always potential for cross-reactivity in two-color experiments. Use less secondary antibody to minimize.
		Always test the two colors on separate blots first so you know what bands to expect and where.
		Avoid using mouse and rat antibodies together, if possible. Because the species are so closely related, anti-mouse will react with rat IgG to some extent, and anti-rat with mouse IgG. Sheep and goat antibodies may exhibit the same behavior.
	Bleed through of signal from one channel into other channel.	Check the fluorescent dye used. Fluorophores such as Alexa Fluor® 750 may appear in both channels and are not recommended for use with the Odyssey® Imaging Systems.
		If signal in one channel is very strong (near or at saturation), it may generate a small amount of bleedthrough signal in the other channel. Minimize this by using a lower scan intensity setting in the problem channel. Try AutoMode on Odyssey CLx.
		Reduce signal by reducing the amount of protein loaded or antibody.

© 2012 LI-COR, Inc. LI-COR is an ISO 9001 registered company. LI-COR, Odyssey, IRDye, NewBlot, In-Cell Western, and MPX are trademarks or registered trademarks of LI-COR, Inc. in the United States and other countries. The Odyssey Imaging Systems and IRDye infrared dye-labeled biomolecules are covered by U.S. and foreign patents and patents pending. All other trademarks belong to their respective owners.



4647 Superior Street • P.O. Box 4000 • Lincoln, Nebraska 68504 USA  
 Technical Support: 800-645-4260 • North America: 800-645-4267  
 International: 402-467-0700 • Fax: 402-467-0819

LI-COR GmbH Germany, Serving Europe, Middle East, and Africa: +49 (0) 6172 17 17 771  
 LI-COR UK Ltd. UK, Serving UK, Ireland, and Scandinavia: +44 (0) 1223 422104  
 All other countries, contact LI-COR Biosciences or a local LI-COR distributor:  
<http://www.licor.com/distributors>

[www.licor.com/bio](http://www.licor.com/bio)

Doc# 988-12622  
 0112