

Consult the Odyssey Application Protocols Manual for complete guidelines and troubleshooting.

Sample Protocol

1. Wet membrane in 1X PBS for 5 minutes. (Pre-wet PVDF membrane in 100% methanol and rinse with ultrapure water before wetting.)
2. Block membrane with Odyssey Blocking Buffer for 1 hour at room temperature with gentle shaking.
3. Decant blocking buffer.
4. Add primary antibody using the vendor's recommended dilution for Western blot applications in Odyssey Blocking Buffer, plus 0.1 - 0.2% Tween® 20. Incubate a minimum of 1 hour at room temperature with gentle shaking.
5. Wash membrane 4 times for 5 minutes in a generous amount of PBS plus 0.1% Tween 20 with gentle shaking.
6. Dilute the IRDye® secondary antibody (1:20,000; consult the IRDye secondary antibody pack insert) in Odyssey Blocking Buffer plus 0.1 - 0.2% Tween 20 (also add 0.01% - 0.02% SDS for PVDF) and incubate, protected from light, 1 hour with gentle shaking.
7. Wash membrane as in step 5; protect from light.
8. Rinse membrane in 1X PBS.
9. Scan on an Odyssey Imaging System.

Two-Color Detection

Note: Two-color detection requires careful selection of primary and secondary antibodies. For details, refer to *Western Blot Analysis and Western Blot and In-Cell Western™ Assay Detection Using IRDye Subclass Specific Antibodies*.

- In step 4, combine two primary antibodies and incubate simultaneously with the membrane.
- In step 6, combine the IRDye 680LT and IRDye 800CW secondary antibodies that recognize the primary antibodies and incubate simultaneously with membrane.

For Best Results

- Let membrane dry after transfer for 1 hour or overnight.
- Clean forceps, incubation trays, and Odyssey scanning surface (if applicable) before use with 100% methanol to remove any residual dye signal from previous use; rinse with a small volume of distilled water followed by isopropanol. Dry with a lint-free wipe.
- Handle membranes by the edges only. Be sure to use clean forceps with smooth edges.
- Avoid pouring antibody or wash solutions directly onto membranes; pour onto the bottom of the dish or down the side.
- Do not include detergents during the blocking step.

©2011 LI-COR, Inc. LI-COR, Odyssey, IRDye and In-Cell Western are trademarks and registered trademarks of LI-COR, Inc. Tween is a registered trademark of ICI Americas. All other trademarks belong to their respective owners.

Technical Support: 800-645-4260, or biohelp@licor.com

LI-COR

Biosciences

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