

Product Number

**827-40000**

Quantity: 25 µL

Storage: -20 °C

June 2013

Updates available at:

<http://biosupport.licor.com>

**Limitation of Liability and Limited Use Label License**

LI-COR IRDye® Infrared Dyes and reagent products are offered for research purposes only and are not intended for human therapeutic or diagnostic use. The purchase of this product conveys to the buyer the non-transferable right to use the amount of product purchased and the components of the product in research conducted by the buyer (whether the buyer is a not-for-profit, academic or for-profit entity). The buyer shall not sell or otherwise transfer this product, its components, or materials made there from to any third party. Buyer shall not use this product or its components for commercial purposes. The term "commercial purposes" shall mean any activity by a party for consideration and may include, but is not limited to, use of the product or its components (i) in manufacturing, (ii) to provide a service, information or data, (iii) for therapeutic, diagnostic or prophylactic purposes, or (iv) for resale, whether or not such product or its components are resold for use in research. The use of this product by the buyer constitutes agreement with the terms of this limited use label license for LI-COR IRDye infrared dyes and reagent products. Inquiries regarding the licensing of one or more IRDye infrared dyes should be submitted by e-mail to [busdev@licor.com](mailto:busdev@licor.com).

LI-COR BIOSCIENCES DOES NOT PROVIDE RESEARCH ADVICE OR DETERMINE OR RECOMMEND ANY POTENTIAL USES FOR IRDYE INFRARED DYES AND REAGENT PRODUCTS. LI-COR BIOSCIENCES MAKES NO WARRANTIES OF ANY KIND, EITHER EXPRESS OR IMPLIED, AS TO ANY MATTER INCLUDING, BUT NOT LIMITED TO, WARRANTY OF FITNESS FOR PURPOSE, OR MERCHANTABILITY OR RESULTS OBTAINED FROM USE OF IRDYE INFRARED DYES. IN NO EVENT SHALL LI-COR BE LIABLE FOR LOST PROFITS, CONSEQUENTIAL, EXEMPLARY, SPECIAL, DIRECT, INCIDENTAL, OR PUNITIVE DAMAGES, OR ATTORNEY FEES, EVEN IF LI-COR HAD BEEN ADVISED OF, KNEW, OR SHOULD HAVE KNOWN, OF THE POSSIBILITIES THEREOF. NO EMPLOYEE, AGENT OR REPRESENTATIVE OF LI-COR HAS THE AUTHORITY TO BIND LI-COR TO ANY ORAL REPRESENTATION OR WARRANTY EXCEPT AS SPECIFICALLY SET FORTH HEREIN.

© 2013 LI-COR, Inc. LI-COR is an ISO 9001 registered company. LI-COR and Odyssey are registered trademarks of LI-COR, Inc., in the United States and other countries. All other trademarks belong to their respective owners. For patent information, visit [www.licor.com/patents](http://www.licor.com/patents)

Doc #988-13756

**LI-COR®**

4647 Superior Street • P.O. Box 4000  
Lincoln, Nebraska 68504 USA  
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)

## Odyssey® Protein Molecular Weight Marker – Sample Pack

Odyssey Protein Molecular Weight Marker is a mixture of ten blue-stained, recombinant proteins (10 - 250 kDa), including three reference bands (25, 50, and 75 kDa). The Marker is designed to provide a ladder of convenient, consistent sizes for SDS-PAGE applications and is recommended for monitoring electrophoretic separation, molecular weight sizing, and monitoring blot transfer efficiency. The Marker also provides fluorescence detection in the 700 channel of all Odyssey Family Imaging Systems.

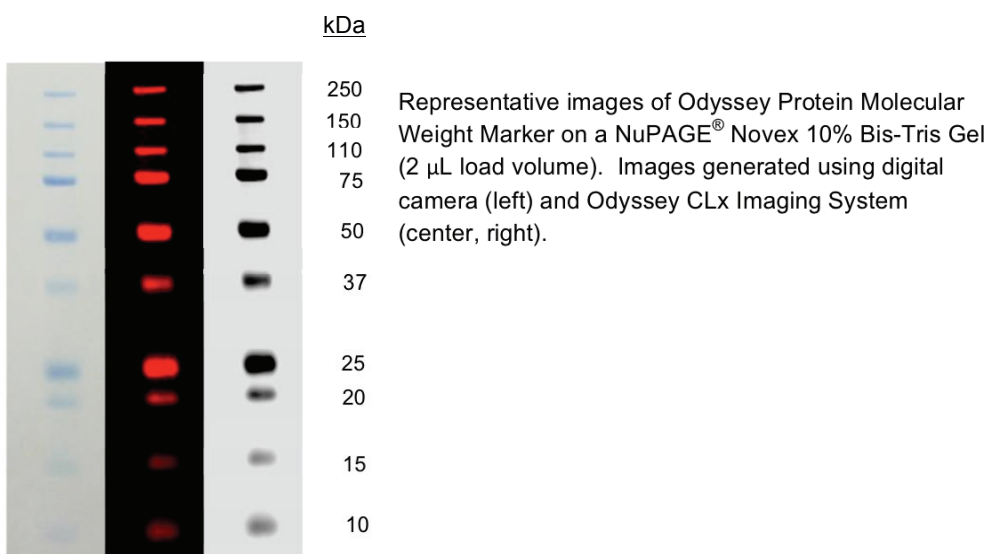
### Specifications

Odyssey Protein Molecular Weight Marker contains ten recombinant protein bands of 10, 15, 20, 25, 37, 50, 75, 100, 150, and 250 kDa. Store at -20 °C upon arrival.

**Volume:** 25 µL

**Loading Buffer:** 30% (w/v) glycerol, 2% SDS, 62.5 mM Tris, pH 6.8, 50 mM DTT, 5 mM EDTA, 0.02% NaN<sub>3</sub>

**Shelf Life:** 1 year at -20 °C



### Instructions

Odyssey Protein Molecular Weight Marker is provided in loading buffer, ready to load, with no dilution required. Allow the tube to reach room temperature and thoroughly mix before use to dissolve any solids that may have precipitated at -20 °C.

Do not heat above room temperature. Recommended loading volumes:

Load Volume	PAGE Application
1 - 3 µL	Mini gels
1 - 3 µL	Blotting, to monitor transfer efficiency (mini gel)
3 - 5 µL	Large gels

### NOTICE TO PURCHASER: LIMITED LICENSE

*This product is 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 re-sell this product or its components.*