

## Components

**928-90000** Chameleon Kit Pre-stained Protein Ladder, 2 x 250 µL

## Specifications

- Channel: 700 and 800 nm
- Size: 2 x 250 µL
- Storage: Store in small aliquots at -15 to -25 °C and always protect from light. Do not subject the ladder to more than 25 cycles of freezing and thawing.
- Shelf Life: 6 months from date of receipt
- Compatible Imagers Include: Odyssey® M Imager, Odyssey DLx Imager, Odyssey CLx Imager, Odyssey XF Imager, Odyssey Fc Imager

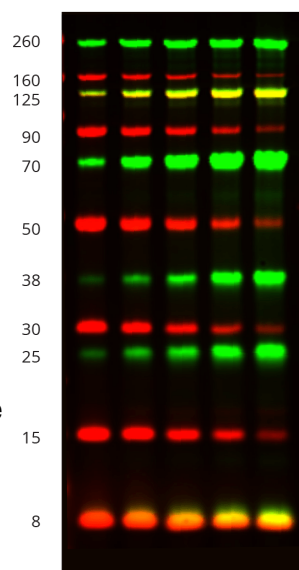
## Product Description

The Chameleon Kit Pre-stained Protein Ladder provides multicolored, pre-stained, and near-infrared fluorescent bands of convenient and consistent protein sizes.

In gels, the ladder can be used to visualize progress of the protein separation during electrophoresis and to estimate the molecular weight of unknown proteins based on their relative mobility. In blotting applications, the ladder can be used to monitor protein transfer and as a reference to estimate the molecular weight of proteins of interest.

The example shows a two-color, near-infrared image of custom mixtures prepared using ratios suggested below. Samples were resolved on a 4 - 12% Bis-Tris Gel and transferred to nitrocellulose via wet tank transfer. The blot was scanned on an Odyssey® CLx Imager. Red and green colors represent the 700 and 800 nm channels, respectively.

The ladder is optimized for use with Bis-Tris and Tris-Glycine systems. To see a complete portfolio of LI-COR molecular weight markers, visit [licor.com/proteinmarkers](http://licor.com/proteinmarkers).



Chameleon Kit  
LI-COR P/N 928-90000

**Signal Optimization for Chameleon Kit:** The kit is designed to independently control the fluorescence signal intensity of the marker in the 700 and 800 nm channels on the Odyssey family of imagers. Customer optimization is required. Test samples composed of mixtures prepared at the ratios listed below as shown.

Channel	Amount to Add (µL)				
700 nm	5	4	3	2	1
800 nm	1	2	3	4	5

**Note:** These ratios are for use with the Chameleon Kit components only. Mixing the Chameleon 700 and Chameleon 800 products (P/Ns 928-70000 and 928-80000) will not result in the same appearance.

## Loading Volumes

Application	Recommended Loading Volume
Protein Gel	10 µL
Western Blot	1 - 5 µL*

\*Ladder volume may require optimization to obtain uniform signal compared to your sample.

## Protocol

Chameleon Pre-stained Protein Ladders arrive ready to use in loading buffer. Bring the product to room temperature while protecting from light. Mix the product thoroughly before use to dissolve any solids that may have precipitated during storage. **Do not heat above room temperature.**

During electrophoresis, the ladder will separate into 11 visible and near-infrared bands in the 8 - 260 kDa range. The ladder can be used for easy visualization or two-color near-infrared detection. For individual molecular weights, please refer to the image above.

**Note:** *The Chameleon Pre-stained Protein Ladders are not standards, and protein band migration may appear different depending on the gel type used.*

## Analysis

### Empiria Studio® Software

Molecular weights for Chameleon Pre-Stained Protein ladders are preset in Empiria Studio Software, allowing you to estimate band size without entering the molecular weights. For more information, see the instructions for the Experiment workflow you are using at [licor.com/WesternWorkflowOverview](http://licor.com/WesternWorkflowOverview).

### Image Studio™ Software

For instructions on entering marker molecular weights and sizing bands in Image Studio Software, visit [licor.com/chameleon-sizing](http://licor.com/chameleon-sizing). Although Image Studio Software is no longer in development, we continue to sell and support it.

### Limitation of Liability and Limited Use Label License

For Research Use Only. Not intended for human therapeutic or diagnostic use.

By using this product, you agree to the Limitation of Liability and the Limited Use Label License available online at [licor.com/packinserts](http://licor.com/packinserts) under the **Pack Inserts** heading (the "License"). The License is incorporated herein by this reference, and the License may be updated from time to time.



© 2022 LI-COR, Inc. LI-COR, Odyssey, Chameleon, Empiria Studio, and Image Studio 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 Biosciences  
4647 Superior Street  
Lincoln, NE 68504  
Phone: +1-402-467-0700  
Toll free: 800-645-4267  
biosales@licor.com  
licor.com/bio

#### Regional Offices

LI-COR Biosciences GmbH  
Siemensstraße 25A  
61352 Bad Homburg  
Germany  
Phone: +49 (0) 6172 17 17 771  
licor.eu@licor.com

LI-COR Biosciences UK Ltd.  
St. John's Innovation Centre  
Cowley Road • Cambridge  
CB4 0WS • United Kingdom  
Phone: +44 (0) 1223 422104  
bio-eu@licor.com