

E-PAGE™ MagicMark™ Unstained Protein Standard

Catalog no. LC5701 **Size** 250 µL **Store at** -30°C to -10°C

Pub. Part no. LC5701.pps **MAN0001522** **Rev. Date** 17 January 2012

Description

The E-PAGE™ MagicMark™ Unstained Protein Standard is suitable for estimating the molecular weight of proteins on E-PAGE™ pre-cast gels after staining or western blotting. The E-PAGE™ MagicMark™ Unstained Protein Standard allows you to directly visualize protein standard bands on a western blot without the need for modifying the proteins or using special detection reagents. Each standard protein contains an IgG binding site.

The important features of the standard are:

- Consists of 5 recombinant protein bands in the range of 20–220 kDa
- Suitable for western blotting and estimating molecular weight
- Designed for E-PAGE™ pre-cast gels
- Visualized with alkaline phosphatase or peroxidase conjugated antibody using chromogenic or chemiluminescent substrates
- Visualized also with SimplyBlue™ SafeStain, silver stain, or fluorescent stains on E-PAGE™ pre-cast gels

Specifications

Contents: 250 µL E-PAGE™ MagicMark™ Protein Standard

Storage Buffer: 125 mM Tris-HCl, pH 6.8; 10 mM DTT; 1% (w/v) SDS;
17.4% (w/v) glycerol; 0.025% Bromophenol Blue

Storage: Store at -30°C to -10°C. To avoid repeated freezing and thawing, aliquot in small volumes and store.

Stability: 6 months at -20°C

For research use only. Not for human or animal therapeutic or diagnostic use.

Directions

The E-PAGE™ MagicMark™ Unstained Protein Standard is supplied in a ready-to-use format. You do not need to heat or reduce the standard.

1. Load the standard into the marker well of an E-PAGE™ 96 gel using the loading volumes listed below to obtain the best results.

| Gel | Electrophoresis | Blotting |
|----------------|-----------------|----------|
| E-PAGE™ 96 gel | 10 µL | 5–10 µL* |

***Note:** We recommend testing different amounts of the standard to determine the optimal amount of standard to use with your western blotting conditions. The amount of standard will depend on the binding affinity of MagicMark™ for your antibody species (next page), sensitivity of your detection system, and exposure time.

2. Perform electrophoresis as described in the E-PAGE™ manual.

Blotting

1. After transferring proteins to a suitable membrane, perform the blocking step, primary antibody incubation step, and (if necessary) secondary antibody incubation step with the blot using your method of choice.
2. Visualize proteins using a colorimetric, chemiluminescent, or fluorescent detection system using the manufacturer's recommendations. After detection, you should observe 5 protein standard bands as shown on page 3.

Staining

Stain the E-PAGE™ gel with SimplyBlue™ SafeStain, Coomassie R-250 stain, SilverXpress® Silver Stain, or fluorescent stain as described in the E-PAGE™ manual.

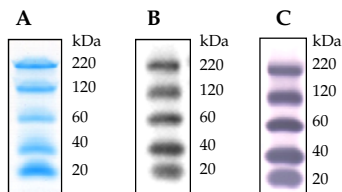
Affinity of E-PAGE™ MagicMark™ Proteins to Antibodies

| Species | Affinity of MagicMark™ |
|--|------------------------|
| Human, Horse, Cow | ++++ |
| Pig, Rabbit | +++ |
| Goat, Sheep, Hamster, Guinea Pig, Rat, Mouse | ++ |
| Chicken | + |

Example

E-PAGE™ MagicMark™ Unstained Protein Standard (10 μ L) was electrophoresed on a 6% E-PAGE™ 96 Gel and stained with SimplyBlue™ SafeStain (A) as described in the E-PAGE™ manual. The standard (5 μ L) was also blotted onto a 0.45- μ m nitrocellulose membrane, and detected with Anti-V5 Antibody (1:5000 dilution) from Life Technologies using the WesternBreeze® Anti-mouse Chemiluminescent Kit (B) or WesternBreeze® Anti-mouse Chromogenic Kit (C).

Note: For troubleshooting, refer to the MagicMark™ Western Protein Standard manual available at www.lifetechnologies.com/manuals.



Product Qualification

The Certificate of Analysis (CofA) provides detailed quality control information for each product. The CofA is available on our website at www.lifetechnologies.com/support, and is searchable by product lot number, which is printed on each box.

Accessory Products

The following products are available separately. For more information, visit www.lifetechnologies.com or contact Technical Support.

| Product | Quantity | Catalog no. |
|-------------------------------------|----------|--------------|
| E-PAGE™ 96 Gels | 8 gels | EP096-06 |
| Mother E-Base™ | 1 unit | EB-M03 |
| Daughter E-Base™ | 1 unit | EB-D03 |
| WesternBreeze® Immunodetection Kits | 1 kit | See web site |
| Nitrocellulose Membrane, 0.45-μm | 20 blots | LC2001 |
| Invitrolon™ PVDF Membranes, 0.45-μm | 20 blots | LC2005 |
| SimplyBlue™ SafeStain | 1 L | LC6060 |
| SilverXpress® Silver Staining Kit | 1 kit | LC6100 |

Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

©2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

For support visit www.lifetechnologies.com/support or email techsupport@lifetech.com

www.lifetechnologies.com

