

E-PAGE™ SeeBlue® Pre-Stained Standard

Cat. no. LC5700 Size 500 µL Store at 2°C to 8°C

Pub. Part no. LC5700.pps MAN0005692 **Rev. Date:** 17 January 2012

Description

The E-PAGE[™] SeeBlue[®] Pre-Stained Standard is specifically designed for use with E-PAGE[™] pre-cast gels. It allows you to visualize protein molecular weight ranges during electrophoresis and evaluate western transfer efficiency.

Important features of the standard are:

 Consists of 5 pre-stained protein bands (3 blue and 2 contrasting colors) ranging from 20–260 kDa

Designed for E-PAGE[™] pre-cast gels

Supplied in a ready-to-use format

Specifications

Contents: 500 µL E-PAGE™ SeeBlue® Pre-Stained Standard

Storage Buffer: Tris-HCl, Formamide, and SDS

Storage: Store at 2°C to 8°C

Stability: 4 months at 4°C

Directions

The E-PAGE[™] SeeBlue[®] Pre-Stained Standard is supplied in a ready-to-use format. **Do not** heat or add reducing agent.

Load the standard into the marker well of an E-PAGE^T 96 gel using the loading volumes listed in the following table to obtain the best results:

Gel	Electrophoresis	Blotting
E-PAGE™ 96 Gel	10 μL	5 μL

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

Example

The apparent molecular weights of protein bands in the E-PAGE™ SeeBlue® Pre-Stained Standard are shown below. Ten microliters of E-PAGE™ SeeBlue® Pre-Stained Standard was separated on a 6% E-PAGE™ 96 gel as described in the E-PAGE™ manual.

k Da	Protein
261— —— 173— ——	Myosin Phosphorylase B
97-	Bovine serum albumin
42-	Alcohol dehydrogenase
21-	Myoglobin

Certificate of Analysis

The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to www.lifetechnologies.com/support and search for the Certificate of Analysis by product lot number, which is printed on the box.

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