

Product Number
926-85000

WesternSure™ ULTRA Chemiluminescent Substrate

Kit Contents:

WesternSure ULTRA
Luminol Enhancer Solution,
50 mL

WesternSure ULTRA Stable
Peroxide Solution, 50 mL

Storage: Room Temperature

February 2013

Updates available at:

<http://biosupport.licor.com>

Limitation of Liability

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. For patent information, visit www.licor.com/patents. LI-COR, Odyssey, WesternSure, and C-DiGit are trademarks or registered trademarks of LI-COR, Inc. in the United States and other countries. All other trademarks belong to their respective owners.

Doc #988-13466



4647 Superior Street • P.O. Box 4000
Lincoln, Nebraska 68504 USA
Biotechnology Support U.S.: 800-645-4260
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

Introduction

WesternSure™ ULTRA Chemiluminescent Substrate is a highly sensitive enhanced substrate for detecting horseradish peroxidase (HRP) on immunoblots and is optimized for use with the LI-COR® C-DiGit™ Blot Scanner and the Odyssey® Fc Imaging System. This substrate's intense signal output enables detection of picogram amounts of antigen. The sensitivity and intensity allow for easy detection of HRP using imaging systems or film.

Protocol

NOTE: Optimize the antigen and antibody concentrations. WesternSure HRP-conjugated secondary antibodies (LI-COR P/N 926-80010 or 926-80011) are optimized for use with this substrate. HRP-conjugated secondary antibodies from other vendors may be suitable; optimization may be required.

1. Dilute primary antibody as recommended by vendor.
2. Dilute WesternSure HRP Secondary Antibody 1:20,000 - 1:100,000. (LI-COR P/N 926-80010 or 926-80011).
3. Prepare Working Solution by mixing equal parts of the Stable Peroxide Solution and the Luminol Enhancer Solution. Use 0.1 mL Working Solution per cm² of membrane. The Working Solution is stable for 8 hours at room temperature.
4. Incubate blot 5 minutes in substrate Working Solution.
5. Drain excess reagent.
NOTE: Light emission is most intense during the first 5 - 30 minutes after substrate incubation. Light emission will continue for several hours, but will decrease with time.
6. For imaging on the C-DiGit Blot Scanner, place the wet blot directly on the scanning surface, protein side facing down. The blot will not dry out during imaging. The blot can be placed in a clear plastic sleeve, if desired. Select 'Standard' or 'High' sensitivity setting, then select 'Start'.
7. For imaging on the Odyssey Fc Imaging System, cover blot with a clear plastic sleeve. Place the protected blot in the tray, protein side facing up. The blot needs to be covered to avoid drying during imaging. Select the acquisition time, then select 'Start'.
8. For film, cover the blot with clear plastic wrap or a clear plastic sleeve. Place the protected blot in a film cassette with the protein side facing up. Place film on top of blot and expose for desired length of time.