

Texas Red®-X Protein Labeling Kit

Catalog no. T10244

Pub. no. MAN0006894

Rev. date 29 August 2012

Detailed protocol is available online at www.lifetechnologies.com/manuals.

WARNING! For every chemical, read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from www.lifetechnologies.com/support.

Chemical	CAS	Warning
DMSO	67-68-5	Components of the product may be absorbed into the body through the skin.

1. Add 1 mL of deionized water to Component B to prepare a 1 M sodium bicarbonate solution.
2. Add 50 μ L of the 1 M sodium bicarbonate solution (from step 1) to 0.5 mL of a 2 mg/mL protein solution.
3. Add 10 μ L DMSO to Texas Red®-X Reactive Dye (Component A) that has equilibrated to room temperature.

For research use only. Not for use in diagnostic procedures.

4. Transfer the protein solution (from step 2) to the vial of reactive dye with DMSO (from step 3). Stir the reaction mixture for 1 hour at room temperature.
5. Assemble the Purification Column.
6. Stir the Purification Resin (Component C) and pipet the resin into the column until it is ~3 cm from the top.
7. Allow excess buffer to drain into the column bed and load the labeled protein (from step 4) onto the column resin.
8. Slowly add 1X elution buffer (10-fold dilution of Component D) and collect the first fluorescent band, which contains the labeled protein.

Approximate fluorescence excitation and emission maxima of Texas Red®-X Reactive Dye: 595/615 nm.

Limited Product Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

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

DISCLAIMER — LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

For support visit www.lifetechnologies.com/support
www.lifetechnologies.com