

Glycoprotein Western Detection Kit Protocol-at-a-Glance

(PT4056-2)

The **Glycoprotein Western Detection Kit** (Cat. No. 635648) is designed for detection of glycoprotein sugar moieties on Western blotting membranes. The kit contains sufficient materials to stain 20 Western blots. The blots may be generated by performing gel electrophoresis and Western blotting to nitrocellulose or PVDF membranes using your standard laboratory methods, according to the instructions provided by the respective manufacturer of your reagents.

A. Kit Components

- Oxidizing Reagent (3 x 85 ml)
- Staining Reagent (3 x 85 ml)
- HRP Positive Control (250 µl; apply 5–10 ul per lane)

B. Additional Reagents Required

- 3% acetic acid

Mix 15 ml of glacial acetic acid with 485 ml of ultrapure water. Store at room temperature.

C. Detection Procedure

Perform this procedure in a fume hood, since the Staining Reagent is toxic.

1. Wash membrane with 20 ml of 3% acetic acid for 5 min on an orbital shaker.
Repeat this step once.
2. Transfer membrane to a container holding 10 ml of Oxidizing Reagent and gently agitate on an orbital shaker for 20 min.
3. Wash membrane with 10 ml of 3% acetic acid for 5 min on an orbital shaker.
Repeat this step twice.
4. Transfer membrane to a container holding 10 ml of Staining Reagent and gently agitate on an orbital shaker for 15 min.
5. Wash membrane extensively with 3% acetic acid and then with ultrapure water.
Glycoproteins will appear as magenta bands.

NOTE: Oxidizing Reagent may be stored at room temperature. If a precipitate is present after warming to room temperature, then spin briefly in a centrifuge and use supernatant for detection.



Clontech

United States/Canada
800.662.2566

Asia Pacific
+1.650.919.7300

Europe
+33.(0)1.3904.6880

Japan
+81.(0)77.543.6116

Clontech Laboratories, Inc.
A Takara Bio Company
1290 Terra Bella Ave.
Mountain View, CA 94043
Technical Support (US)
E-mail: tech@clontech.com
www.clontech.com

Notice to Purchaser

Clontech products are to be used for research purposes only. They may not be used for any other purpose, including, but not limited to, use in drugs, *in vitro* diagnostic purposes, therapeutics, or in humans. Clontech products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without written approval of Clontech Laboratories, Inc.

Clontech, the Clontech logo and all other trademarks are the property of Clontech Laboratories, Inc. unless otherwise noted. Clontech is a Takara Bio Company. ©2008 Clontech Laboratories, Inc.

(PR862560; published 18 June 2008)