

ab308244 – Glycoprotein Staining Kit

Detection of glycoproteins on a PVDF or Nitrocellulose membrane.

For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit: [abcam website](#)

Storage and Stability

Store the kit at 4°C, protected from light. Briefly centrifuge all small vials prior to opening. Read the entire protocol before performing the assay procedure.

Materials Supplied

Item	Quantity	Storage Condition after preparation
Oxidizing Reagent (10X) (Avoid light)	20 ml	4 °C
Staining Reagent A (6X) (Avoid light)	40 ml	-
Staining Reagent B (6X) (Avoid light)	40 ml	-
Glycoprotein Positive Control	1 vial	-20°C
Glycoprotein Negative Control	1 vial	-20°C

Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

- Equipment's to run SDS-PAGE and transfer proteins from the gel to a PVDF or nitrocellulose membrane.
- 3% glacial acetic acid (i.e., 15 ml of glacial acetic acid in 485 ml of water). Store at room temperature (RT).

Reagent Preparation

Oxidizing Reagent Working Solution: Prepare the Oxidizing Reagent Working Solution by mixing one part of Oxidizing Reagent (10X) with nine parts of 3% glacial acetic acid and mix well. Store at 4 °C.

Staining Reagent Working Solution: Prepare fresh Staining Reagent Working Solution by mixing one part of Staining Reagent A (6X), one part of Staining Reagent B (6X) with four parts of water.

Glycoprotein Positive & Negative Controls: Reconstitute each vial in 250 µl of water to make 1 mg/ml solution. Divide into aliquots and store at -20 °C.

Glycoprotein Staining Protocol:

SDS-PAGE:

Run an SDS-PAGE to separate the glycoproteins, following the standard SDS-PAGE procedure.

Gel Transfer:

Transfer the proteins from the gel onto a PVDF or Nitrocellulose membrane using standard gel transfer procedure.

Staining of Glycoproteins on membrane:

1. Wash the membrane with 20 ml of 3% glacial acetic acid for 5 min with gentle shaking. Repeat this step once.
2. Add 10 ml of Oxidizing Reagent Working Solution and gently shake for 20 min.

3. Wash the membrane with 10 ml of 3% glacial acetic acid for 5 min with gentle shaking. Repeat this step twice.
4. Add 12 ml of fresh Staining Reagent Working Solution to the container with the membrane and gently shake for 15 min.
5. Wash the membrane briefly with 10 ml of 3% glacial acetic acid and water.
6. Take the image immediately. Glycoproteins will appear as magenta bands.

Note: The background will become darker over time.

Technical Support

Copyright © 2023 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)