INSTRUCTIONS

Glycoprotein Isolation Kit, WGA



Storage: Upon receipt store kit at 4°C. Kit is shipped at ambient temperature.

Introduction

The Thermo Scientific[™] Glycoprotein Isolation Kit, WGA isolates glycoproteins from complex protein mixtures including serum and tissue and cultured cell lysates using the lectin wheat germ agglutinin (WGA) immobilized on agarose. Lectins are proteins that have a selective affinity for carbohydrate moieties. The WGA lectin preferentially binds N-acetyl glucosamine (GlcNAC) and terminal GlcNAC structures that are commonly present in many serum and membrane glycoproteins. WGA also has affinity for sialic acid.

Protein glycosylation is a common post-translational modification. Asparagine (N-linked) and serine/threonine residues (Olinked) are glycosylated during passage through the endoplasmic reticulum and golgi apparatus in eukaryotic and prokaryotic (i.e., Archaea and Eubacteria) systems. Glycoconjugates are important for immune regulation, inflammation, cell-to-cell adhesion and contact inhibition, cell signaling, protection against proteolytic degradation, and other biological processes.

This kit is easy to use and contains all the necessary components for isolating sialic acid/GlcNAC-containing glycans. A sample containing up to 1.5mg of total protein is first diluted with the Binding/Wash Buffer and applied to the WGA resin bed. Following incubation, the resin is washed and the bound glycoproteins are eluted. Glycoproteins have been successfully isolated from serum as well as HeLa and CHO cell lysates in approximately 50 minutes.

Important Product Information

- If needed, add protease inhibitors to samples; however, avoid cocktails containing EDTA or other metal chelators.
- Samples purified with this kit are compatible with 1D gel electrophoresis and the Thermo ScientificTM Coomassie PlusTM (Bradford) Assay Kit (Product No. 23236). Many other downstream applications require sample processing to remove incompatible substances in the Elution Buffer. To quantify protein using the Thermo Scientific BCA Protein Assay (Product No. 23225), desalt sample using a 5mL Thermo ScientificTM ZebaTM Spin Desalting Column (Product No. 89891). For 2D gel electrophoresis, remove interfering substances by precipitation or dialysis.
- For serum samples, isolation of less-abundant glycoproteins is improved by first removing the albumin and IgG. This can be accomplished with an Albumin/IgG Removal Kit (see Related Products Section).







Procedure Summary Flow Chart



Procedure for Glycoprotein Isolation using WGA

A. Preparation of sample

- 1. Equilibrate the Binding/Wash and Elution Buffers to room temperature.
- 2. Dilute sample containing 1 to 1.5mg of total protein 4:1 with 5X Binding/Wash Buffer stock solution (e.g., mix 400µL sample with 100µL 5X Binding/Wash Buffer). The total volume, including dilution, must not exceed 800µL.

B. Isolation of glycoproteins

- 1. To prepare the 1X Binding/Wash Buffer, dilute 460µL 5X Binding/Wash Buffer with 1840µL of ultrapure water, which is sufficient volume to process one sample.
- 2. Insert a column into a collection tube.
- 3. Gently swirl the bottle of WGA Lectin Resin to obtain a homogeneous suspension. Use a wide-bore or cut pipette tip to transfer 200μL of 50% resin slurry to the column.
- 4. Centrifuge 1 minute at $1000 \times g$ and discard the storage buffer. Reuse the collection tube through Step B12.
- 5. Place column in collection tube and add 200 μ L of 1X Binding/Wash Buffer to the resin. Close the top cap and centrifuge column for 1 minute at 1000 × g and discard rinse. Repeat this step two times.
- 6. Place bottom cap on column and add sample, from Step A1, to the resin. Close the top cap.
- 7. Incubate column for 10 minutes at room temperature with end-over-end mixing using a rotator (e.g., Labquake[™] Shaker by Thermolyne). Alternatively, rock back and forth on a rocking platform.
- 8. Remove top cap and then bottom cap from column. Place column in the collection tube, and replace top.

Note: Remove top cap before bottom cap to prevent sample from leaking from the bottom of the column.

9. Centrifuge column for 1 minute at $1000 \times g$ and discard flow-through.

Note: If desired, save the flow-through for SDS-PAGE or protein assay analysis.

- 10. Reinsert column and add 400 μ L 1X Binding/Wash Buffer to the resin. Cap column and centrifuge for 1 minute at $1000 \times g$ and discard flow-through. Repeat this step.
- 11. Place bottom cap on column and add 400µL of 1X Binding/Wash Buffer to the resin. Cap column and incubate for 5 minutes at room temperature with end-over-end mixing using a rotator.
- 12. Remove top cap and then bottom cap from column. Place column in the collection tube, and replace top cap. Centrifuge column for 1 minute at $1000 \times g$ and discard flow-through.
- 13. Repeat Steps B11-B12.
- 14. Replace bottom cap on column. Add 200µL of Elution Buffer to resin and cap column. Incubate column for 10 minutes at room temperature with end-over-end mixing using a rotator.
- 15. Remove top cap and then bottom cap from column. Place column in a new collection tube. Replace top cap and centrifuge column for 1 minute at $1000 \times g$.
- 16. Carefully set aside the collection tube and remove top cap.
- 17. Repeat Steps B14-B16. Collect eluate in the same collection tube that contains eluate from the first elution. Store eluted glycoproteins on ice for immediate use or freeze for later analysis.



Troubleshooting

Problem	Possible Cause	Solution
Low glycoprotein recovery in elution fraction	Some glycoproteins have high affinity for the immobilized WGA and will not elute with Elution Buffer	Increase incubation with elution buffer to 15 minutes, or boil resin in 200µL of SDS-PAGE sample buffer for 5 minutes and then centrifuge column in a 2mL tube for 1 minute at $1000 \times g$ to collect eluate Note: Boiling the resin results in detachment of some lectin and also may release nonspecifically bound proteins.
Glycoprotein is not binding to the resin	Sample contains metal chelator(s)	Confirm EDTA or other metal chelators are not present in the sample

Related Thermo Scientific Products

89804	Glycoprotein Isolation Kit, ConA
89875	Pierce Albumin/IgG Removal Kit
89161	Top 2 Abundant Protein Depletion Columns (Human Albumin and IgG)
89891	Zeba Spin Desalting Columns
23225	BCA Protein Assay Kit
23236	Coomassie Plus (Bradford) Assay Kit

General References

Cooper, C.A., *et al.* (2001) GlycoSuiteDB: A new curated relational database of glycoprotein glycan structures and their biological sources. *Nucl Acid Res* **20**(1):332-5.

Cummings, R.D. (1997). Affinity chromatography of oligosaccharides and glycopeptides. Affinity Separations: A Practical Approach (Matejschuk, P., Ed.), Oxford Univ. Press, London, pp. 123-39.

Ghosh, D., et al. (2004). Lectin affinity as an approach to the proteomic analysis of membrane glycoproteins. J Proteome Res 3:841-50.

Young, N.M., et al. (2002). Structure of the N-linked glycan present on multiple glycoproteins in the gram-negative bacterium, Campylobacter jejuni. J Biol Chem 277:42530-9.

Products are warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation"). No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

NO OTHER WARRANTIES, EXPRESS OR IMPLIED, ARE GRANTED, INCLUDING WITHOUT LIMITATION, IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR ANY PARTICULAR PURPOSE, OR NON INFRINGEMENT. BUYER'S EXCLUSIVE REMEDY FOR NON-CONFORMING PRODUCTS DURING THE WARRANTY PERIOD IS LIMITED TO REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER'S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR PRODUCTS AS THE RESULT OF (I) ACCIDENT, DISASTER OR EVENT OF FORCE MAJEURE, (II) MISUSE, FAULT OR NEGLIGENCE OF OR BY BUYER, (III) USE OF THE PRODUCTS IN A MANNER FOR WHICH THEY WERE NOT DESIGNED, OR (IV) IMPROPER STORAGE AND HANDLING OF THE PRODUCTS.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

Current product instructions are available at <u>www.thermoscientific.com/pierce</u>. For a faxed copy, call 800-874-3723 or contact your local distributor. © 2013 Thermo Fisher Scientific Inc. All rights reserved. All (other) trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.