



# **Enzymatic Deglycosylation Kit**

## **Instruction Manual**

**Catalog Number**  
**170-6500**

**BIO-RAD**

# Table of Contents

<b>Section 1</b>	<b>Introduction .....</b>	<b>1</b>
<b>Section 2</b>	<b>Kit Components and Specifications ....</b>	<b>2</b>
<b>Section 3</b>	<b>Protocols .....</b>	<b>4</b>
3.1	Sample Preparation .....	4
3.2	Non-Denaturing Protocol.....	5
3.3	Denaturing Protocol.....	5
3.4	Experimental Control.....	8
<b>Section 4</b>	<b>Product Information.....</b>	<b>8</b>
<b>Section 5</b>	<b>Technical Assistance.....</b>	<b>9</b>

# Section 1

## Introduction

The Enzymatic Deglycosylation kit enzymatically cleaves all N and most O-linked oligosaccharides from glycoproteins or glycopeptides. PNGase F removes all Asn-linked oligosaccharides while the combination of NANase II and O-Glycosidase DS removes all Ser/Thr linked Gal( $\beta$ 1,3)GalNAc( $\alpha$ 1) and all sialic acid substituted Gal( $\beta$ 1,3)GalNAc( $\alpha$ 1). Modifications of the Ser/Thr linked Gal( $\beta$ 1,3)GalNAc core other than by neuraminic acid (such as galactose, GlcNAc or fucose substitutions) will inhibit O-Glycosidase cleavage. To insure complete deglycosylation of such modified cores, the use of additional enzymes may be required.

## Section 2

### Kit Components and Specifications

Component	Specificity	Concentration	Volume	Storage
NANase II	Releases $\alpha$ 2-3 and $\alpha$ 2-6 linked N-acetylneuraminic acid from complex oligosaccharides	10 U/ml (in 20 mM Tris-HCl pH 7.5, 25 mM NaCl)	40 $\mu$ l	4 °C
O-Glycosidase DS	Releases unsubstituted Gal( $\beta$ 1,3)GalNAc( $\alpha$ 1) disaccharide attached to serine or threonine.	1 U/ml (in 20 mM Tris-HCl pH 7.5, 25 mM NaCl)	40 $\mu$ l	4 °C
PNGase F	Releases all Asn-linked oligosaccharides from glycoproteins.	2.5 U/ml (in 20 mM Tris-HCl pH 7.5, 50 mM NaCl, 1 mM EDTA)	40 $\mu$ l	4 °C
Bovine Fetuin Control	N/A	500 $\mu$ g	N/A	4 °C
5x Reaction Buffer	N/A	250 mM sodium phosphate, pH 6.0	200 $\mu$ l	RT
pH Adjustment Buffer	N/A	0.5 M sodium phosphate dibasic	200 $\mu$ l	RT
Denaturing Solution	N/A	2% SDS, 1 M $\beta$ -Mercaptoethanol	200 $\mu$ l	RT
NP-40	N/A	>99 %	100 $\mu$ l	RT

RT = Room Temperature

## Section 3 Protocols

Two deglycosylation protocols are provided: one with and one without a denaturing step. The choice of protocols depends on the glycoprotein. Some proteins require denaturation prior to PNGase F digestion. Initially, carry out both protocols with your protein. If denaturation releases more oligosaccharides than without, use the denaturing protocol for all subsequent studies.

### 3.1 Sample Preparation

Isolate glycoprotein according to your usual procedures.

**Solid Sample:** Dissolve up to 100 µg glycoprotein in 12 µl of distilled water. Add 4 µl of 5x Reaction Buffer.

**Liquid Sample:** Add 4 µl of 5x Reaction Buffer to 12 µl of glycoprotein solution (up to 100 µg glycoprotein).

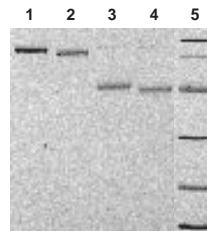
### 3.2 Non-Denaturing Protocol

1. Add 2 µl of NANase II to the reaction vial.
2. Add 2 µl of O-Glycosidase to the reaction vial.
3. Incubate at 37 °C for 1 hour.
4. Add 10 µl of distilled water and 10 µl of pH Adjustment Buffer.
5. Add 2.0 µl of PNGase F to the reaction vial.
6. Incubate at 37 °C for 24 hours.
7. Check the efficiency of the deglycosylation reaction by running samples (before and after deglycosylation) on a SDS-PAGE gel. Stain with Coomassie Blue or Silver Stain to observe the shift in mobility of your sample. Alternatively, use the Immun-Blot® Kit for Glycoprotein Detection to determine the deglycosylation efficiency.

### 3.3 Denaturing Protocol

1. Add 2 µl of NANase II to the reaction vial.
2. Add 2 µl of O-Glycosidase to the reaction vial.
3. Incubate at 37 °C for 1 hour.

4. Add 10  $\mu$ l of deionized water and 10  $\mu$ l of pH Adjustment Buffer.
5. Add 2.5  $\mu$ l of Denaturing Solution and heat for 5 minutes at 100  $^{\circ}$ C.
6. Cool on ice for 5 minutes.
7. Add 2.5  $\mu$ l of NP-40 to the reaction vial and mix.
8. Add 2.0  $\mu$ l of PNGase F to the reaction vial.
9. Incubate at 37  $^{\circ}$ C for 3 hours.
10. Check the efficiency of the deglycosylation reaction by running samples (before and after deglycosylation) on a SDS-PAGE gel. Stain with Coomassie Blue or Silver Stain to observe the shift in mobility of your sample. Alternatively, use the Immun-Blot Kit for Glycoprotein Detection to determine the deglycosylation efficiency.



A. 12 % Ready Gel stained with Coomassie<sup>®</sup> blue. Mobility shift indicates proteins were deglycosylated.



B. A parallel gel blotted onto nitrocellulose. Glycosylation determined with the Immun-Blot Kit for Glycoprotein Detection. Deglycosylated proteins do not react and are not detected.

**Fig. 1. Human transferrin and Ovalbumin were deglycosylated with the Enzymatic Deglycosylation Kit and detected with the Immun-Blot Kit for Glycoprotein Detection.** Lane 1: Human transferrin; Lane 2: Deglycosylated Human transferrin; Lane 3: Ovalbumin; Lane 4: Deglycosylated Ovalbumin; Lane 5: Biotinylated SDS-PAGE Standards, low range.

### 3.4 Experimental Control

Bovine Fetuin is included in the kit for use as a positive control. Use this control to verify that the deglycosylation reaction was carried out correctly. Dilute the Bovine Fetuin Control (500 µg) in 50 µl of distilled water to yield a 10 mg/ml working solution and perform the deglycosylation reaction protocol. Store reconstituted control at -20 °C.

## Section 4 Product Information

<b>Catalog Number</b>	<b>Product Description</b>
170-6500	<b>Enzymatic Deglycosylation Kit</b>
170-6490	<b>Immun-Blot Kit for Glycoprotein Detection</b>
170-6501	<b>N-Linked Oligosaccharide Profiling Kit</b>
170-6510	<b>N-Linked Oligosaccharide Sequencing Kit</b>

## Section 5 Technical Support

If you require additional technical assistance contact your local Bio-Rad representative or in the U.S. dial 1-800-4BIORAD and press 2 for the technical service department.

Coomassie is a trademark of ICI.