



Deglycosylation Enhancement Kit

Instruction Manual

Catalog Number
170-6508

BIO-RAD

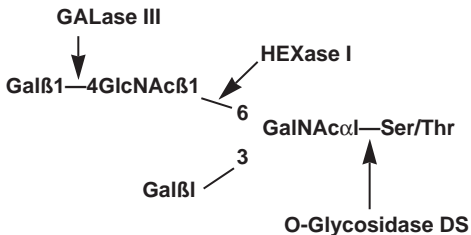
Table of Contents

Section 1	Introduction	1
Section 2	Kit Components and Specifications	3
Section 3	Protocols	4
3.1	Sample Preparation	5
3.2	Non-Denaturing Protocol.....	5
3.3	Denaturing Protocol.....	6
Section 4	Product Information.....	8
Section 5	Technical Assistance.....	10

Section 1

Introduction

The Deglycosylation Enhancement Kit (catalog number 170-6508) enhances deglycosylation by the additional release of β -galactose and β -GlcNAc modified cores. The kit includes HEXase I which releases non-reducing terminal β -linked N-acetyl-glucosamine and GALase III which releases non-reducing terminal (β 1,4) linked galactose. These enzymes together with the enzymes provided in the Enzymatic Deglycosylation Kit (170-6500) will release Gal(β 1,3)GalNAc(α 1) with β -galactose or β -GlcNAc substitutions. For example:



Additional modifications such as α -fucose or α -galactose to these structures require additional enzymes for complete O-linked deglycosylation. To insure complete deglycosylation of such modified cores, the use of additional enzymes may be required.

Note: The Deglycosylation Enhancement Kit protocol differs from the one contained in the Enzymatic Deglycosylation Kit. To use the Deglycosylation Enhancement Kit effectively, follow the deglycosylation protocol contained in this instruction manual.

Section 2 Kit Components and Specifications

	HEXase I	GALase III	2x Reaction Buffer pH 7.5
Specificity	Non-reducing terminal β -linked N-acetylglucosamine	Non-reducing terminal (β 1,4) linked galactose	N/A
Concentration	8 U/ml	1.5 U/ml	100 mM sodium phosphate, pH 7.5
Volume	40 μ l	40 μ l	
Storage	4 °C	4 °C	4 °C

Section 3 Protocols

The HEXase I and GALase III supplied in this kit will degrade N-linked oligosaccharides present in your sample. To minimize N-linked oligosaccharide degradation, release N-linked oligosaccharides with PNGase F prior to O-linked oligosaccharides release.

Two deglycosylation protocols are provided. One protocol includes a denaturing step and the other omits the denaturing step. Some glycoproteins require denaturation prior to PNGase F digestion. Denaturation of these glycoproteins with SDS/ β -mercaptoethanol will increase the efficiency of N-linked oligosaccharide release 10 fold. The choice of protocols depends on the glycoprotein. 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 usual procedures.

Solid Sample: Dissolve up to 100 μ g of glycoprotein in 20 μ l of distilled water. Add 25 μ l of 2x Reaction Buffer, pH 7.5.

Liquid Sample: Dilute 20 μ l of glycoprotein solution (containing up to 100 μ g) with 25 μ l of 2x Reaction Buffer, pH 7.5.

3.2 Non-Denaturing Protocol

1. Aliquot 45 μ l of prepared sample (Section 3.1).
2. Add 2 μ l of PNGase F and incubate for 24 hours at 37 °C.
3. Add 3 volumes of cold ethanol. Chill on ice for 5 minutes.
4. Centrifuge the sample to pellet the glycoprotein and decant the supernatant.
5. Wash the pellet with cold 70% ethanol and decant the supernatant.

6. Dry the pellet in a centrifugal vacuum evaporator.
7. Add 30 μl of distilled water to re-suspend the glycoprotein pellet.
8. Add 10 μl of 5x Reaction Buffer 6.
9. Add 2 μl of each:
 - NANase II
 - GALase III
 - HEXase I
 - O-Glycosidase DS
10. Incubate for 2 hours at 37 °C.
11. Run the treated and untreated glycoprotein in separate lanes in a SDS-PAGE gel. Deglycosylated proteins will exhibit an increase in mobility due to the reduction in molecular weight. Alternatively use the Immun-Blot® Kit for glycoprotein detection (catalog number 170-6490) to determine the efficiency of the deglycosylation reaction.

3.3 Denaturing Protocol

1. Aliquot 45 μl of prepared sample (Section 3.1).

2. Add 2.5 μl of Denaturing Solution to sample. Mix gently and heat at 100 °C for 5 minutes.
3. Chill on ice and add 2.5 μl of NP-40 solution.
4. Add 2 μl of PNGase F and incubate for 3 hours at 37 °C.
5. Add 3 volumes of cold ethanol. Chill on ice for 5 minutes.
6. Centrifuge the sample to pellet the glycoprotein and decant the supernatant.
7. Wash the pellet with cold 70% ethanol and decant supernatant.
8. Dry the pellet in a centrifugal vacuum evaporator..
9. Add 30 μl of distilled water to re-suspend the glycoprotein pellet.
10. Add 10 μl of 5x Reaction Buffer.
11. Add 2 μl of each:
 - NANase II
 - GALase III
 - HEXase I
 - O-Glycosidase DS
12. Incubate for 2 hours at 37 °C.

13. Run the treated and untreated glycoprotein in separate lanes in a SDS-PAGE gel. Deglycosylated proteins will exhibit an increase in mobility due to the reduction in molecular weight. Alternatively use the Immun-Blot[®] kit for glycoprotein detection (catalog number 170-6490) to determine if deglycosylation is complete.

Section 4 Product Information

Catalog Number	Product Description
<i>Carbohydrate Analysis Kits</i>	
170-6490	Immun-Blot Kit for Glycoprotein Detection
170-6500	Enzymatic Deglycosylation Kit
170-6508	Deglycosylation Enhancement Kit
170-6513	GALase III , 1.5 U/ml, 0.04 ml
170-6880	HEXase I , 42 U/ml, 0.04 ml
170-6881	O-Glycosidase DS , 1 U/ml, 0.04 ml
170-6882	NANase II , 5 U/ml, 0.04 ml
170-6883	PNGase F , 2.5 U/ml, 0.04 ml

Catalog Number	Product Description
170-6501	N-Linked Oligosaccharide Profiling Kit
170-6510	N-Linked Oligosaccharide Sequencing Kit
170-6502	N-Linked Oligosaccharide Gel Refill , 6
170-6514	N-Linked Oligosaccharide Gel and Buffer Refill
170-6815	O-Linked Oligosaccharide Profiling Kit
170-6816	O-Linked Oligosaccharide Gel and Buffer Refill Pack
170-6817	O-Linked Oligosaccharide Gel Refill Pack
170-6503	Oligosaccharide Electrophoresis Buffer Refill
170-6811	Monosaccharide Composition Analysis Kit
170-6812	Monosaccharide Gel and Buffer Refill Pack
170-6813	Monosaccharide Gel Refill Pack
170-6814	Monosaccharide Buffer Refill Pack
<i>Carbohydrate Analysis Instruments</i>	
170-6555	Glyco Doc™ Imager , 100/120 V
170-6557	Glyco Doc Imaging System , 100/120 V
170-6559	Glyco Doc Analytical Software

Section 5 Technical Support

If you require additional technical assistance contact your local Bio-Rad representative or in the US dial 1-800 4BIORAD.