

Enzymatic Deglycosylation of a Protein Containing Core 1 O-glycans with O-Glycosidase

Paula Magnelli, Dave Landry and Alicia Bielik

For many cancers, such as colon, ovary, uterus and bladder (mucosas), tumor progression and its poor prognosis strongly correlates with alterations in the patterns of mucin O-glycosylation. For instance, β 3Gn-T6 (the enzyme responsible for Core 3 O-glycan synthesis), is abundant in normal colon tissue while its expression is strongly downregulated in adenocarcinoma (1,2,3). As a result, mucin glycosylation switches from common Core 3 O-glycan structures to short Core 1 structures. These Core 1 structures, T and Tn, are hallmark epitopes of cell malignancy (4).

The study of glycan aberrations in cancer opens new avenues in the development of novel therapies. In this regard, the endo-O-glycosidase is proving useful to characterize and quantitate a variety of commonly found O-glycan forms.

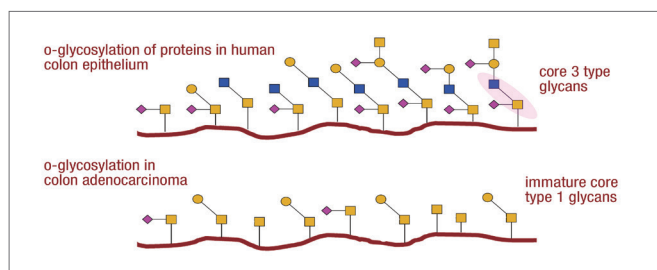
This application note describes the use of a recombinant O-Glycosidase that can cleave core 1 O-linked disaccharides (5) and immature Tn core (GalNAc). Conditions have been determined that allow this O-Glycosidase to be used under the same denaturing conditions used for PNGase F digestion.

Materials

O-Glycosidase & Neuraminidase Bundle
(NEB #E0540)

Fetuin (NEB #P6042)

Core 1 disaccharide (Galb1,3GalNAc;
Accurate Chemical #BCR20/06)



General Protocol

1. Set up the reaction in 200 μ l tubes as follows:

COMPONENTS	DENATURING CONDITIONS	NATIVE CONDITIONS
Fetuin (10 mg/ml) in water	18 μ l	18 μ l
10X Denaturing Buffer	2 μ l	2 μ l
Incubate 95°C for 5 minutes. Cool on ice, spin		
10X NP-40	4 μ l	–
10X GlycoBuffer 2	4 μ l	4 μ l
Milli-Q Water	18 μ l	24 μ l
Neuraminidase	2 μ l	2 μ l
O-Glycosidase	2 μ l	2 μ l

2. Incubate at 37°C for 1 to 4 hours. After incubation, add 1 μ l of 4 M KCl followed by 150 μ l of methanol. Chill overnight at 4°C to precipitate proteins. After the overnight precipitation, spin the sample at 14,000 rpm for 30 minutes, and reserve the supernatant.

- Concentrate supernatant to dryness with a Speed Vac set at medium heat (Savant; equipped with a high vacuum pump and finger trap immersed in a Dewar containing isopropanol and dry ice). Reconstitute with 400 μ l Milli-Q® water.
- De-ionize the sample from step 4 by gently rocking in 200 μ l of prepared mixed bed ion exchange resin AG 501-X8 (Bio-Rad; #142-6424) for 5 minutes. Collect the supernatant with a 1 ml syringe using a 23 gauge needle. **Note:** before use, the resin must be converted to the acetate form by soaking in an equal volume of 1 M acetic acid followed by washing ten times with equal volumes of water.
- Remove the needle and load the entire sample (400 μ l) from Step 5 to an activated Sep-Pak® cartridge (Waters; #WAT051910). Collect the entire flow through (400 μ l). Wash the Sep-Pak 2 times with 400 μ l of Milli-Q water and pool the washes with the flow through. Concentrate to 70 μ l using a Speed Vac. **Note:** before use, the Sep-Paks are activated by washing two times with 400 μ l methanol followed by 4 times with 400 μ l Milli-Q water.

- Detect free sugars by HPAEC-PAD Chromatography using the following conditions:

Column: CarboPac 20 with Amino Guard.

Elution: 40 mM NaOH isocratic for 12 minutes, 150 mM regeneration for 10 minutes, flow rate: 0.5 μ l/min.

Detection: Pulse electrochemical, Au electrode, quadruple potential.

Injection sample: 30 μ l, with or without internal Gal β 1,3GalNAc standard (30 nanograms).

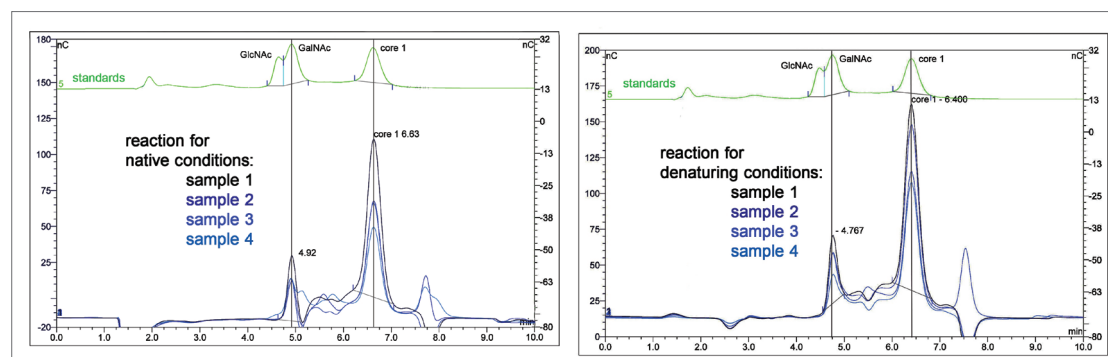
References:

- Iwai, T. et al., (2005) *Proceedings of the National Academy of Sciences U.S.A.* 102, 4572-4577.
- Iwai, T. et al., (2002) *Journal of Biological Chemistry* 277, 12808-12809.
- Capon, C. et al., (2001) *Biochemical Journal* 358, 657-664.
- Springer, G.F. et al., (1997) *Journal of Molecular Medicine* 75, 594-602.
- Koutsouliis, D. et al., (2008) *Glycobiology* 18, 799-805.

Results



Superimposed chromatograms of released sugars with decreasing amounts of enzyme shown in table. The chromatograms represent a 5-fold increase of all materials in Step 2.



Sample #	units	O-GLYCOSIDASE ON NATIVE FETUIN		O-GLYCOSIDASE ON DENATURED FETUIN	
		GalNAc Peak Area nC*min	Core 1 Peak Area nC*min	GalNAc Peak Area nC*min	Core 1 Peak Area nC*min
1	200,000	9.55	36.92	9.84	43.29
2	10,000	6.23	21.85	8.02	39.91
3	50,000	6.74	23.01	8.07	27.12
4	12,500	4.51	18.08	5.09	25.11

One or more of these products are covered by patents, trademarks and/or copyrights owned or controlled by New England Biolabs, Inc. For more information, please email us at gbd@neb.com. The use of these products may require you to obtain additional third party intellectual property rights for certain applications.

Your purchase, acceptance, and/or payment of and for NEB's products is pursuant to NEB's Terms of Sale at www.neb.com/support/terms-of-sale. NEB does not agree to and is not bound by any other terms or conditions, unless those terms and conditions have been expressly agreed to in writing by a duly authorized officer of NEB.

SEP-PAK® is a registered trademark of Waters.

MILLI-Q® is a registered trademark of Millipore Corporation.

© Copyright 2018, New England Biolabs, Inc.; all rights reserved.



www.neb.com

