

# Removal of terminal galactose from a glycoprotein containing tri- and tetra-antennary *N*-linked sugars with $\alpha$ 1-3, 6 Galactosidase

Paula Magnelli, Alicia Bielik and Dave Landry

With advances in transplantation and stem cell research, there has been a renewed interest in the study of glycoforms carrying the Gal $\alpha$ 1-3Gal epitope.

This motif is widely present in non-primate mammalian cells, while absent in Old World monkeys and humans (1). Naturally occurring high levels of anti-Gal antibodies cause xenotransplantations to fail within a few hours (2). This ability to ablate Gal-exposing cells has been exploited to develop safer human tissue grafts (3).

Specific glycosidases are required to characterize these kinds of systems. This application note describes the use of an  $\alpha$ 1-3,6 Galactosidase from *Xanthomonas manihotis* (recombinant expressed in *E. coli*) to remove terminal galactose residues from the tri- and tetra- antennary *N*-glycoprotein Bovine Thyroglobulin (4).

## Materials

$\alpha$ 1-3,6-Galactosidase  
(NEB #P0731)

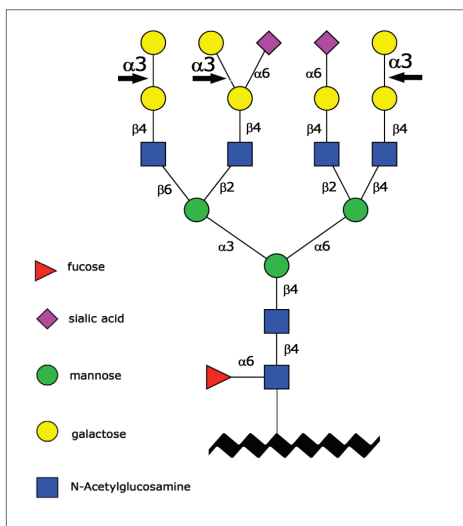
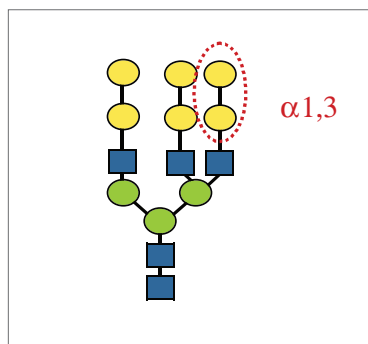
Galactose standard  
(Sigma #G0750)

Bovine Thyroglobulin (Calbiochem; #609310)

10X G6 buffer  
(supplied with enzyme)



**Structure of the Bovine Thyroglobulin tetra-antennary carbohydrate moiety. Arrows denote the  $\alpha$ 1-3,6 Galactosidase cleavage sites.**



## General Protocol

1. Preparation of Glycoprotein substrate: Dialyze 1  $\mu$ l of a 10 mg/ml solution of Bovine Thyroglobulin in water against 100 volumes of G6 buffer, for 4 hours at 4°C. The dialyzed solution can be stored in aliquots of 100  $\mu$ l.

Glycoprotein Substrate 10 mg/ $\mu$ l	85 $\mu$ l
G6 Buffer (10X)	10 $\mu$ l
$\alpha$ 1-3,6 Galactosidase	5 $\mu$ l (20 units)
Total volume	100 $\mu$ l

2. Incubate at 37°C for 4 hours. Add 200  $\mu$ l water followed by 600  $\mu$ l methanol (1)\*. Chill overnight at 4°C to precipitate proteins. After the overnight precipitation, spin the sample at 14 K 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  $\mu$ l Milli-Q™ water.
- De-ionize the sample from step 4 by gently rocking in 200  $\mu$ l of prepared mixed bed ion exchange resin AG 501-X8 for 5 minutes (Bio-Rad; #142-6424). Collect the supernatant with a 1ml 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  $\mu$ l) from Step 5 to an activated Sep-Pak® cartridge (Waters; #WAT051910). Collect the entire flow through (400  $\mu$ l). Wash the Sep-Pak 2 times with 400  $\mu$ l of Milli-Q water and pool the washes with the flow through. Concentrate to 70  $\mu$ l using a Speed Vac. Note: before use, the Sep-Paks are activated by washing two times with 400  $\mu$ l methanol followed by 4 times with 400  $\mu$ l Milli-Q water.
- Detect free galactose by HPAEC-PAD Chromatography using the following conditions:  
Column: CarboPac 20 with Amino Guard. Elution: 20mM NaOH isocratic for 12 minutes, 150 mM regeneration for 10 minutes, flow rate: 0.5  $\mu$ l/min. Detection: Pulse electrochemical, Au electrode, quadruple potential. Injection sample: 30  $\mu$ l, with or without internal Galactose standard (30 nanograms).

#### References

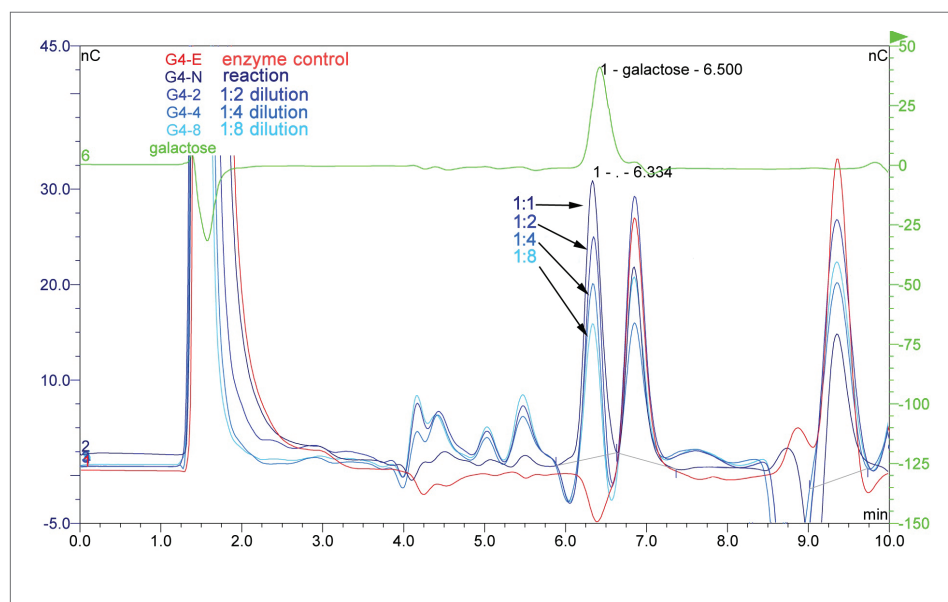
- Koike, C. et al., (2002) *The Journal of Biological Chemistry* 277, 10114–10120.
- Cooper, D.K., et al., (1993) *Lancet* 342, 682–683.
- Hewitt, Z., et al., (2007) *Stem Cells* 25 10–18.
- Wong-Madden, T., et al., (1995) *Glycobiology* 5, 19–28.

## Results:



### FIGURE 1: Superimposed chromatograph of released sugars.

Chromatogram showing galactose peak released by serial decreasing amounts of  $\alpha$ 1-3,6 Galactosidase for the same amount of substrate. The superimposed peaks are designated 1:1 (20 units); 1:2 (10 units), 1:4 (5 units) and 1:8 (0.5 units).



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](mailto: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](http://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.

PRO-Q® is a registered trademark of Invitrogen, Inc.

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



[www.neb.com](http://www.neb.com)

