

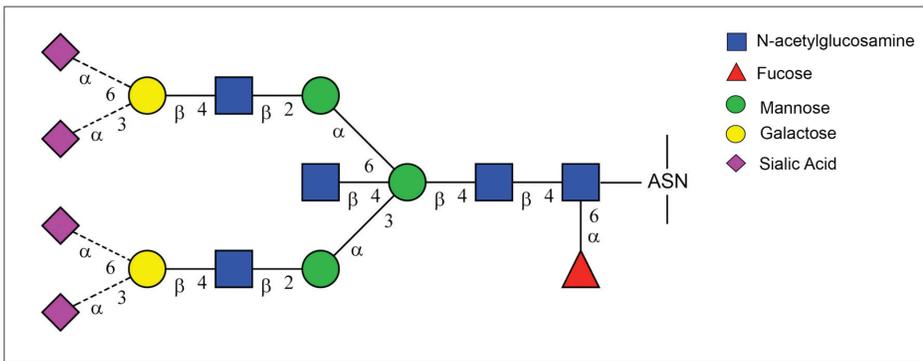
# $\beta$ -N-Acetylglucosaminidase

## Application Note

Paula Magnelli, Alicia Bielik and Dave Landry

### Removal of terminal N-acetylglucosamine from the biantennary N-linked sugars of IgG.

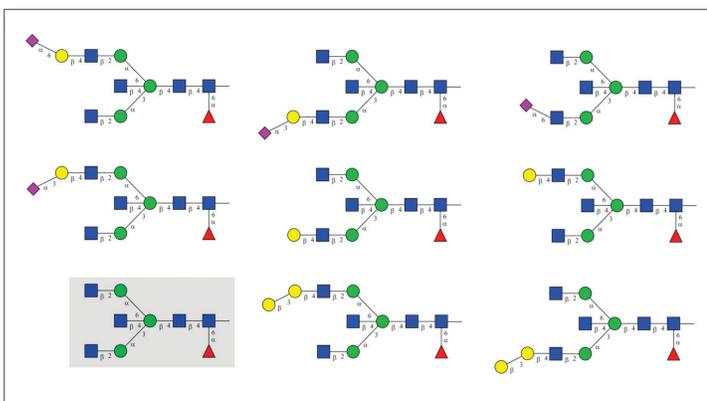
IgG contains one asparagine-linked biantennary carbohydrate in the Fc region of the heavy chain (1). Even though the heterogeneous biantennary structure of different molecules can be characterized by the variability of the presence of terminal galactose, the overall oligosaccharide pattern is constant for healthy individuals (2).



Mature biantennary carbohydrate structure for IgG.

Changes in the extent of galactosylation in the N-glycan of IgG are related to several diseases such as cancer (3), arthritis (4), Castleman's disease (5), periodontitis (6), and coeliac disease (7). Other emerging developments are in the use of recombinantly produced glycoproteins for therapeutic applications. Functional consequences can result from improper glycosylation. In the production of monoclonal antibodies it has been shown that varying dissolved oxygen concentrations in murine hybridoma cell lines during growth can influence the extent of galactosylation (8).

Specific glycosidases can be used to characterize the diversity of the carbohydrate structures in IgG. This application note describes the use of a  $\beta$ -N-Acetylglucosaminidase (recombinant expressed in *E. coli*) to remove terminal N-acetylglucosamine from the biantennary N-glycoprotein IgG. The amount of N-acetylglucosamine that is released is a direct result of the extent of galactosylation.



Nine of the detected thirty immature biantennary N-glycan structures have cleavable terminal N-acetylglucosamines. The only bisecting terminal N-acetylglucosamine cleaved by the N-Acetylglucosaminidase is the N-glycan framed in the highlighted box.

(see other side)

- DNA CLONING
- DNA AMPLIFICATION & PCR
- EPIGENETICS
- RNA ANALYSIS
- LIBRARY PREP FOR NEXT GEN SEQUENCING
- PROTEIN EXPRESSION & ANALYSIS**
- CELLULAR ANALYSIS

### Materials

- $\beta$ -N-Acetylglucosaminidase (NEB #P0732)
- N-acetylglucosamine standard (Sigma #A8625)
- IgG from human serum (Sigma; #4506)
- 10X G1 buffer (supplied with enzyme)

## General Protocol

- Prepare 10 mg/ml solution of IgG in water (store in aliquots of 100  $\mu$ l @ -20°C).
- Set up the exo-glycosidase digestion reaction:
 

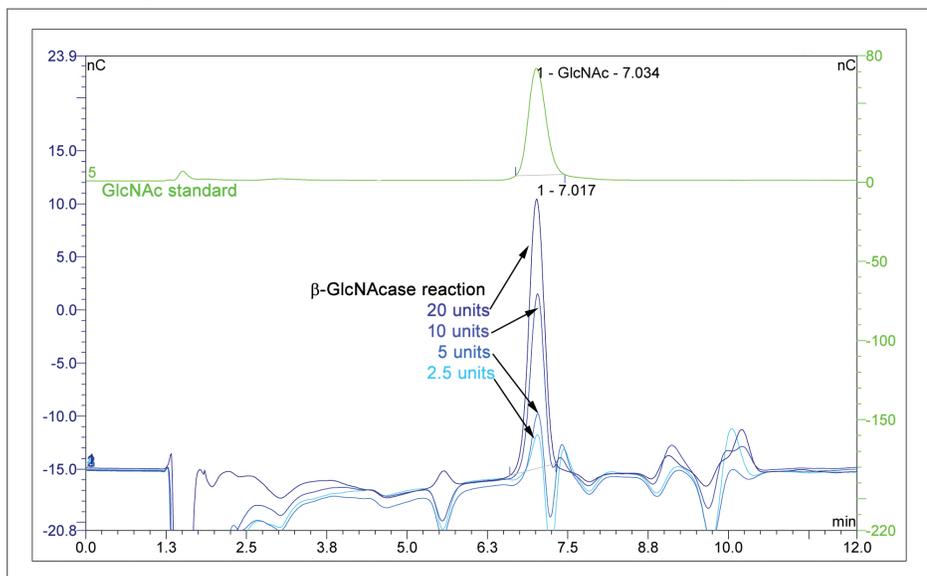
Glycoprotein Substrate 10 mg/ $\mu$ l	85 $\mu$ l
G1 buffer (10x)	10 $\mu$ l
$\beta$ -N-Acetylglucosaminidase	<u>5 <math>\mu</math>l (20 units)</u>
Total volume	100 $\mu$ l
- Incubate at 37°C for 4 hours. Add 200  $\mu$ l water followed by 600  $\mu$ l methanol. Chill overnight at 4°C to precipitate proteins. After the overnight precipitation, spin the sample at 14K 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 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 were activated by washing two times with 400  $\mu$ l methanol followed by 4 times with 400  $\mu$ l Milli-Q water.
- Detect free N-acetylglucosamine (GlcNAc) by HPAEC-PAD Chromatography using the following conditions:
 

*Column:* CarboPac 20 with Amino Guard.

*Elution:* 20mM NaOH isocratic for 12 minutes, 150mM 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 GlcNAc standard (30 nanograms).



Superimposed chromatograms of released N-acetylglucosamine for serial decreasing amounts of  $\beta$ -N-Acetylglucosaminidase.

## References

- Kornfeld, R. et al., (1971) *Journal of Biological Chemistry* 246, 3259-3268.
- Parekh, R. B. et al., (1988) *Lancet* 1, 966-969.
- Fleming, S. C. et al., (1998) *Journal of Clinical Pathology* 51, 825-830.
- Watson, M. et al., (2001) *Arthritis and Rheumatism* 42, 1682-1690.
- Kanayam, Y. et al., (1991) *Clinica Chimica Acta* 197, 00221-00228.
- Gordana, S. et al., (2006) *Journal of Periodontology* 77, 1887-1893.
- Cremata, J. A. et al., (2003) *Clinical and Experimental Immunology* 133, 422-429.
- Kunkel, J.P. et al., (1998) *Journal of Biotechnology* 62, 55-71.

One or more of these products are covered by one or more patents, trademarks and/or copyrights owned or controlled by New England Biolabs, Inc. For more information, please contact NEB's Global Business Development team at [gbd@neb.com](mailto:gbd@neb.com).

Your purchase, acceptance, and/or payment of and for NEB's products is pursuant to NEB's Terms of Sale at <https://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. MILLI-Q® is a registered trademark of Millipore Corporation. SEP-PAK® is a registered trademark of Waters.



v1.0

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



New England Biolabs, Inc., 240 County Road, Ipswich, MA 01938-2723 Telephone: (978) 927-5054  
Toll Free: (USA Orders) 1-800-632-5227 (USA Tech) 1-800-632-7799 Fax: (978) 921-1350 e-mail: [info@neb.com](mailto:info@neb.com)