

Glycan Analysis of Murine IgG by Enzymatic Digestion with Remove-iT™ Endo S and PNGase F, Followed by Mass Spectrometric Analysis

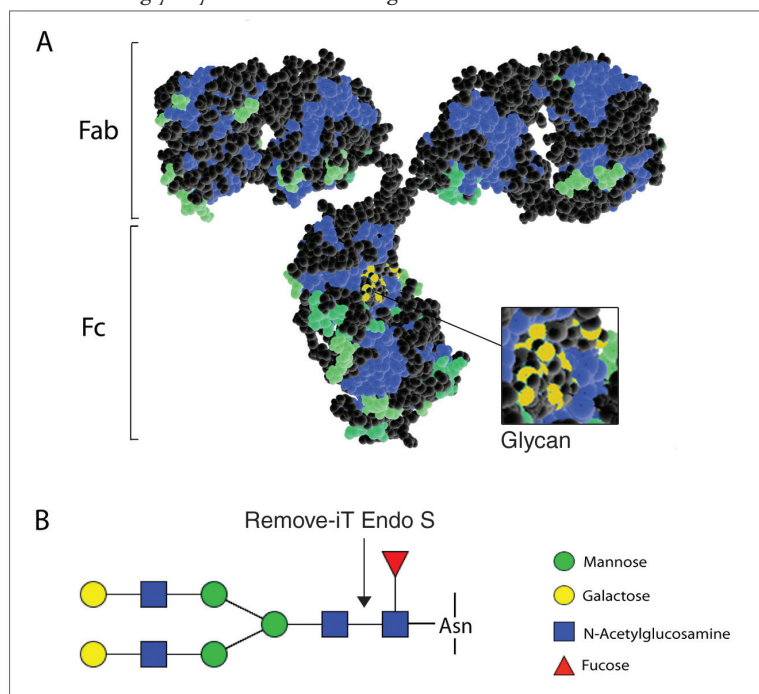
Immunoglobulin Gs (IgGs) are antibody molecules that are composed of four peptide chains — two heavy chains and two light chains. There are four IgG subclasses (IgG1, 2a, 2b, and 3) in mice. The heavy chains are known to be glycosylated, while the light chains are not. The *N*-glycan moiety attached to the asparagine 297 residue in the Fc domain of the antibody is critical for the structure and biological activity of the molecule (1). A growing number of monoclonal IgG antibodies are currently being developed and used as therapeutic agents and it is apparent that there are many variables in cell culture systems that can greatly influence the heterogeneity of the glycans on IgGs. Therefore, it has become increasingly important to monitor the glycosylation profiles of these molecules in the production process.

Endoglycosidase S (Endo S) isolated from *Streptococcus pyogenes* has been shown to specifically and completely cleave the biantennary complex *N*-glycan at asparagine 297 of IgG under native conditions (2). This enzyme removes the *N*-glycan moiety after the first *N*-acetylglucosamine (GlcNAc) residue on the chitobiose core, leaving only a GlcNAc with or without a core fucose residue on the protein. In contrast, PNGase F cleaves between the innermost GlcNAc and asparagine residue of high mannose, hybrid and complex oligosaccharides from *N*-linked glycoproteins.

Remove-iT Endo S is cloned and expressed in *E. coli* as a fusion to the chitin binding domain (CBD). The specificity of Remove-iT Endo S is identical to Endo S with the added benefit of the CBD tag to remove the enzyme following a deglycosylation reaction if needed.

This application compares the enzymatic removal of glycans on murine IgG using Remove-iT Endo S and PNGase F Glycerol free under native conditions. Remove-iT Endo S is a more robust enzyme for this purpose, completely removing the sugar residues from monoclonal mouse IgG. Conversely, the PNGase F digest does not result in a complete digestion under native conditions.

Structure and glycosylation of a murine IgG



(A) Structural model of murine IgG. In the IgG, beta-sheets are colored blue, loops are colored black and the helices are colored green. The brackets indicate the antigen-binding Fab portion and the Fc effector portion of IgG. The inset highlights the two conserved glycans (yellow) attached to Asn-297 of the heavy chains. The model was generated using JMOl 12.2.23 from a model deposited in the Protein Data Bank by L. Harris (University of California, Riverside).
(B) Schematic representation of the fully substituted IgG heavy-chain glycan and the location of the Remove-iT Endo S cleavage.

DNA CLONING

DNA AMPLIFICATION & PCR

EPIGENETICS

RNA ANALYSIS

SAMPLE PREP FOR NEXT GEN SEQUENCING

PROTEIN EXPRESSION & ANALYSIS

CELLULAR ANALYSIS

Materials

- Remove-iT Endo S (NEB #P0741)
- PNGase F (Glycerol Free) (NEB #P0705)
- Anti-MBP Monoclonal Antibody (Murine IgG2a) (NEB #E8032)
- G7 Reaction Buffer (10X, supplied with PNGase F (Glycerol free))
- G6 Reaction Buffer (10X, supplied with Remove-iT Endo S)
- 3K Millipore Amicon Ultra Filter Unit (cat. #UFC500324)
- Dilution Buffer: 20mM Tris HCl pH 7.5, 50 mM NaCl, 1mM EDTA
- Protein Ladder (NEB #P7703)
- Agilent 6210 TOF MS with both 1200 Series Capillary and Nano pumps with a ChipCube. Custom PLRP-S Chip (75 μm x 150 mm with 40 nl trap) or equivalent TOF or Q-TOF and nanoLC system.

General Protocols

Deglycosylation of IgG using Remove-iT Endo S occurs optimally under native conditions.

Deglycosylation using PNGase F occurs optimally under denaturing conditions (using SDS, DTT and heat); however, this is not optimal for downstream mass spectrometry analysis and thus the reaction needs to be performed under native conditions without the addition of detergents.

Removal of Glycerol from Murine IgG Substrate

1. Dilute 50 μ l of 1 mg/ml Anti-MBP Monoclonal Antibody (Murine IgG2a, NEB #E8032S) with 450 μ l of the Dilution Buffer: 20mM Tris HCl pH 7.5, 50 mM NaCl, 1mM EDTA.
2. Apply to a 0.5 ml 3K Millipore Amicon Ultra Filter Unit (cat. # UFC500324) and spin in a microcentrifuge for 30 minutes at 12,000 rpm.
3. Discard flow-through and add an additional 450 μ l of Dilution Buffer to the sample. Spin in a microcentrifuge for 30 minutes at 12,000 rpm.
4. Place the Amicon filter device upside-down in a clean microcentrifuge tube and spin for 2 minutes at 1,000 rpm to transfer glycerol-free murine IgG to the tube.

Deglycosylation of IgG with Remove-iT Endo S

Keep all enzyme solutions on ice.

1. Add 10 μ l prepared glycerol free Anti-MBP Monoclonal Antibody (Murine IgG2a, NEB #E8032) at 1 mg/ml (10 μ g total) to a 200 μ l tube.
2. Add 5 μ l of 10X G6 Reaction Buffer. (500 mM NaAcetate pH 5.5, 50 mM CaCl₂)
3. Add 34 μ l of water.
4. Add 1 μ l (200 units) of Remove-iT Endo S.
5. Mix with pipette and incubate at 37°C for 1 hour.
6. Reserve 10 μ l for SDS-PAGE gel analysis, if desired.
7. Analyze remainder by nanoLC-TOF MS.

Deglycosylation of IgG with PNGase F (Glycerol Free)

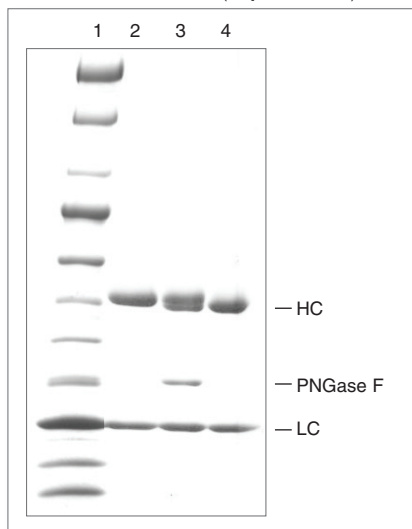
1. Add 10 μ l prepared glycerol free Anti-MBP Monoclonal Antibody (Murine IgG2a, NEB #E8032) at 1 mg/ml (10 μ g total) to a 200 μ l tube.
2. Add 5 μ l of 10X G7 Reaction Buffer. (500 mM Sodium Phosphate pH 7.5)
3. Add 34 μ l of water.
4. Add 1 μ l (500 units) of PNGase F (Glycerol Free).
5. Mix with pipette and incubate at 37°C for 1 hour.
6. Reserve 10 μ l for SDS-PAGE gel analysis if desired
7. Analyze remainder by nanoLC-TOF MS.

Results:

N-glycan moiety removal under native conditions

Remove-iT Endo S has a high specificity for removing the *N*-glycan moiety of IgG under native conditions. Digestion of IgG with PNGase F (Glycerol Free) under native conditions (Lane 3) yields a partial deglycosylation as seen by the presence of a doublet band of the heavy chain (HC). Remove-iT Endo S (Lane 4) yields a complete deglycosylation of IgG under native conditions as shown by a complete shift of the band.

SDS-PAGE gel shift analysis of the deglycosylation of IgG under native conditions using Remove-iT Endo S and PNGase F (Glycerol Free)

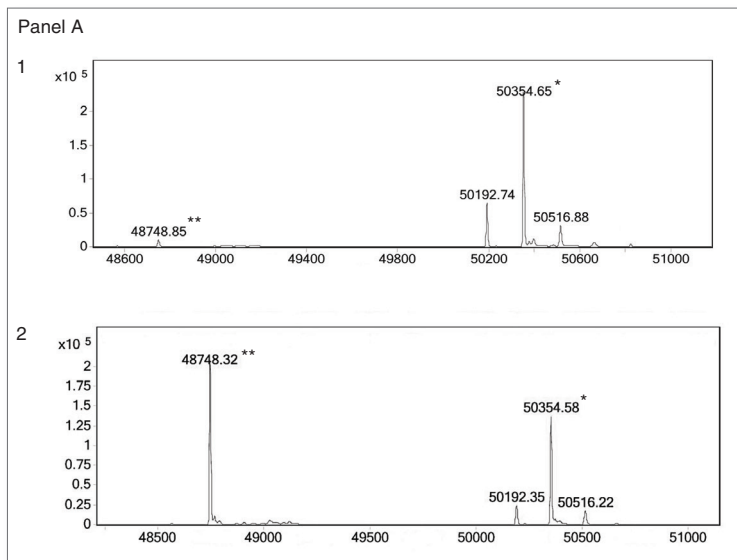


Deglycosylation was assessed by SDS-PAGE: following IgG deglycosylation for 1 hour at 37°C with 500 units of PNGase F (Glycerol Free), in lane 3; or 200 units of Remove-iT Endo S, in lane 4. Samples were standardized against an IgG control, with no deglycosylation enzymes, in lane 2. For reference, Protein Ladder (NEB #P7703) was run in lane 1.

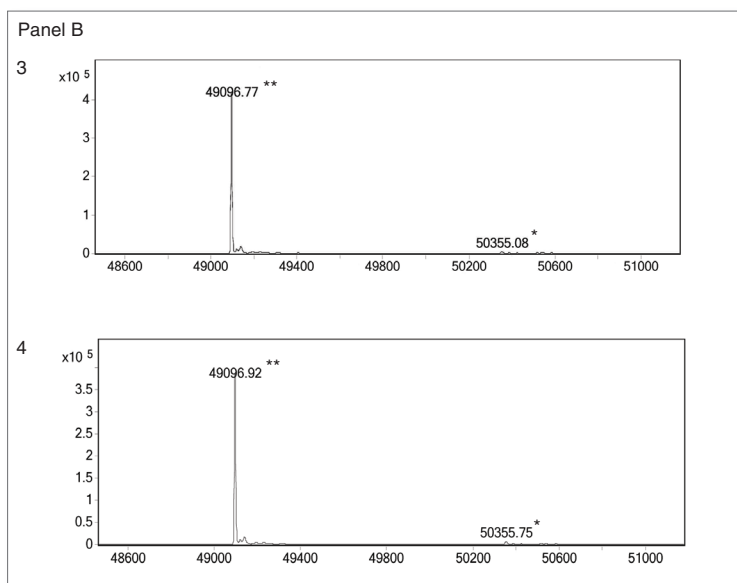
Mass spectrometric analysis

Deglycosylated Murine IgG samples were denatured using 10 mM DTT for 30 minutes at room temperature, and diluted to 0.1% formic acid with water. Samples were analyzed by reverse phase liquid chromatography (LC) and electrospray ionization time-of-flight mass spectrometry (ESI-TOF MS). A custom reverse-phase chip, containing an integrated trapping column (40 nl capacity), separation column and nano-ESI emitter (75 μ m x 150 mm both packed with PLRP-S, 5 μ m particles, 1000 Å pore size) was used for the separation of proteins (3). The chip trap column was loaded at 2 μ l/min and the separation column developed at a flow rate of 500 nl/min using an Agilent 1200 series nano LC connected directly to an Agilent 6210 series ESI-TOF MS. The column was equilibrated with 0.1% formic acid in water containing 5% acetonitrile. 1 μ l of deglycosylated murine IgG sample was injected onto the column and was developed after two minutes with a fifteen-minute linear gradient from 5% to 95% acetonitrile and then held at 95% acetonitrile for five minutes. Protein was found to elute at approximately 10 minutes after injection. The mass spectra were acquired from 150 to 3200 m/z, one cycle/sec and 10,000 transients per scan using an ionization energy of 1800 V, fragmentor of 215 V and drying gas of 325° C at 4.0 l/min. The acquired spectra were extracted and the protein spectra were deconvoluted.

ESI-TOF MS analysis of monoclonal mouse IgG digested with PNGase F (Glycerol Free) (Panel A) or Remove-iT Endo S (Panel B) under native conditions.



ESI-TOF MS : PNGase F digested monoclonal mouse IgG incubated 1 hour at 37 °C (figure 1) or 24 hrs at 37°C (figure 2). * indicates IgG with intact glycan MW 50,354, ** indicates deglycosylated heavy chain of IgG MW 48,748. PNGase F does not completely remove glycans from the IgG substrate under native conditions after a 24 hr digest.



ESI-TOF MS : Remove-iT Endo S digested monoclonal mouse IgG incubated 1 hour (figure 3) or 24 hrs (figure 4). * indicates IgG with intact glycan MW 50,355, ** indicates deglycosylated heavy chain of IgG MW 49,096. Remove-iT Endo S completely removes glycans from IgG after 1 hour at 37°C. Differences in MW of deglycosylated IgG between PNGase F and Remove-iT Endo S are the result of the enzyme specificity. Remove-iT Endo S leaves a GlcNAc residue on the glycoprotein, a feature which can be used to determine occupancy on a given glycoprotein.

Conclusion

Comparison of the deglycosylation activities of PNGase F and Remove-iT Endo S demonstrates that Remove-iT Endo S completely removes glycans from IgG after a 1 hour digestion at 37°C, whereas PNGase F is unable to completely remove glycans from IgG, under native conditions, after a 24 hour digestion at 37°C. Remove-iT Endo S is a superior choice for glycobiology applications that demand rapid and reliable deglycosylation of IgG under native conditions.

References:

1. Arnold, J.N. et al. (2007) *Annual Rev. Immunol.* 25, 21-50.
2. Colin, M. and A. Olsén (2001) *EMBO Journal* 20, 3046-3056.
3. Vollmer, M and van de Goor (2009) *Methods in Molecular Biology* 544, 3-15.