

**NOTICE:** ProZyme was purchased by Agilent in July 2018. Documents for products and product lots manufactured before August 2019 will contain references to ProZyme. For more information about these products and support, go to: [www.agilent.com/en/contact-us](http://www.agilent.com/en/contact-us).



**PRO**zyme

## TECHNICAL NOTE

### Gly-Q

### Gly-X

GlykoPrep

### Glyko Enzymes

Glyko Standards

### InstantPC

InstantAB

### InstantQ

2-AB

APTS Express

PhycoLink

PhycoPro

RPE & APC Conjugates

Streptavidins

#### Keywords

Biotherapeutic

Exoglycosidase

N-Glycans

InstantPC

InstantQ

Capillary Electrophoresis

HILIC

# An Intact Protein Approach to Exoglycosidase Sequencing Followed by Released N-Glycan Analysis from Biotherapeutic Glycoproteins

Andres Guerrero, Anna Fong, John Yan, Tom Rice, Justin Hyché, Aled Jones  
ProZyme, Hayward, CA

## SUMMARY

- Analysis of N-glycans released from biotherapeutics frequently relies on derivatization with a fluorescent label.
- Exoglycosidases cleave specific terminal monosaccharides allowing for sequential digestion thus aiding in the identification of complex N-glycans.
- Glycoprotein samples can be first treated in parallel by a panel of specific exoglycosidases followed by N-glycan sample preparation using Gly-X with InstantPC or InstantQ.

## INTRODUCTION

The structure of N-linked glycans can play a critical role in the pharmacology of therapeutic proteins, potentially affecting immunogenicity, pharmacokinetics and pharmacodynamics, making the characterization of N-glycans an essential part of the biotherapeutic development process (1). A common analytical approach for the analysis of N-glycans includes the enzymatic release of the glycans with PNGase F, derivatization with a fluorophore (InstantPC, 2-AB, InstantQ) and analysis by liquid

chromatography (LC) with fluorescence detection (FLD) with potential hyphenation to mass spectrometry (MS), or capillary electrophoresis (CE) with LED-induced fluorescence (LEDIF). In addition, the use of exoglycosidase enzymes (Figure 1) offers a specific set of tools for the sequential digestion and release of common monosaccharides (sialic acid, galactose and N-acetylglucosamine) from the non-reducing terminus of the N-glycan, aiding in structural characterization of biotherapeutic proteins containing these important post-translational modifications.

Here we present a workflow for the sequential digestion of terminal monosaccharides from N-glycans found on biotherapeutic proteins. In this unique approach, exoglycosidases are used on the intact glycoprotein prior to PNGase F digestion and labeling of the released N-glycans. This approach allows for a simple workflow that can be applied for different labeling chemistries while providing flexibility in fluorescent dye selection.

## METHODS

### Materials

MabThera Lot # H0102803, Enbrel Lot # R180844

Sialidase A [GK80040]

$\beta$ (1-3,4)-Galactosidase [GKX-5013]

$\beta$ (1-2,3,4,6) N-Acetylhexosaminidase, Sequencing-grade [GKX-5023]

### N-Glycan sample preparation kits

Gly-X N-Glycan Rapid Release and Labeling with InstantPC Kit (96-ct) [GX96-IPC]

Gly-X N-Glycan Rapid Release and Labeling with InstantQ Kit (96-ct) [GX96-IQ]

### Exoglycosidase digestion and N-glycan sample preparation conditions

Glycoprotein samples were treated with exoglycosidase enzymes under the following conditions prior to release and labeling of N-glycans (Figure 2).

- 1. Denaturation:** 20  $\mu$ L of 2 mg/mL (40  $\mu$ g) glycoprotein stock along with 2  $\mu$ L Gly-X denaturant was incubated at 90°C for 3 minutes.
- 2. Exoglycosidase digestion:** After denaturation, separate glycoprotein samples were incubated overnight at 42°C with the following sequences of exoglycosidase enzymes (2  $\mu$ L of each enzyme) after the addition of 6  $\mu$ L of 0.5 mM Ammonium Acetate pH 5.5.

- a. Sialidase A [GK80040]
- b. Sialidase A [GK80040] +  $\beta$ -Galactosidase [GKX-5013]
- c. Sialidase A [GK80040] +  $\beta$ -Galactosidase [GKX-5013] +  $\beta$ -N-Acetylhexosaminidase [GKX-5023]

- 3. N-Glycan sample preparation:** Following exoglycosidase digestion N-glycans were prepared following standard Gly-X (with InstantPC and InstantQ) protocols starting at the 5 minute deglycosylation step.

### N-Glycan separation and analysis

#### HILIC separation conditions

60-minute UHPLC high resolution method

Column: 2.1 x 150 mm, 1.7  $\mu$ m

Flow rate: 0.4 mL/minute

Gradient: 25–38% 50 mM ammonium formate pH 4.4, between 2.5–50 minutes

Column temperature: 60°C

#### Gly-Q CE separation

InstantQ labeled N-glycans were analyzed with Gly-Q, a capillary electrophoresis-based 2-minute separation along with LED induced fluorescence detection (LEDIF)

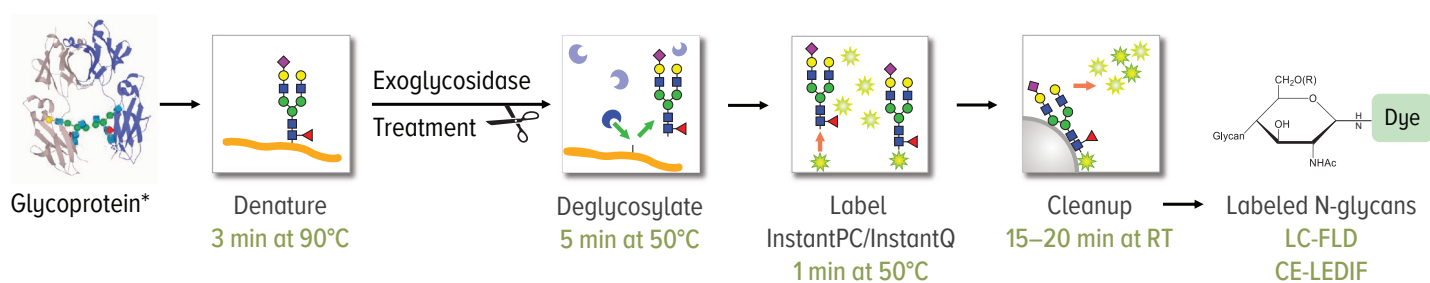
Instrument method: N-glycans ProZyme

Processing methods: Rituxan\_Method/Enbrel\_Method

Exoglycosidase [Product Code]	Short Name	Example Glycan	Specificity
Sialidase A (Arthrobacter) [GK80040]	ABS		$\alpha$ (2–3, 6, 8, 9) N-Acetylneuraminic acid linkages
$\beta$ -Galactosidase (Bovine testis) [GKX-5013]	BTG		$\beta$ (1–3, 4) Galactose linkages
$\beta$ -N-Acetylhexosaminidase (Jack bean) [GKX-5023]	JBH		B (1–2, 3, 4, 6) GlcNAc and GalNAc linkages

◆ N-Acetylneuraminic acid (Neu5Ac)   ● Galactose   ■ N-Acetylglucosamine (GlcNAc)   ● Mannose   ▼ Fucose

**Figure 1:** Exoglycosidase enzymes along with example glycans and substrate specificity.

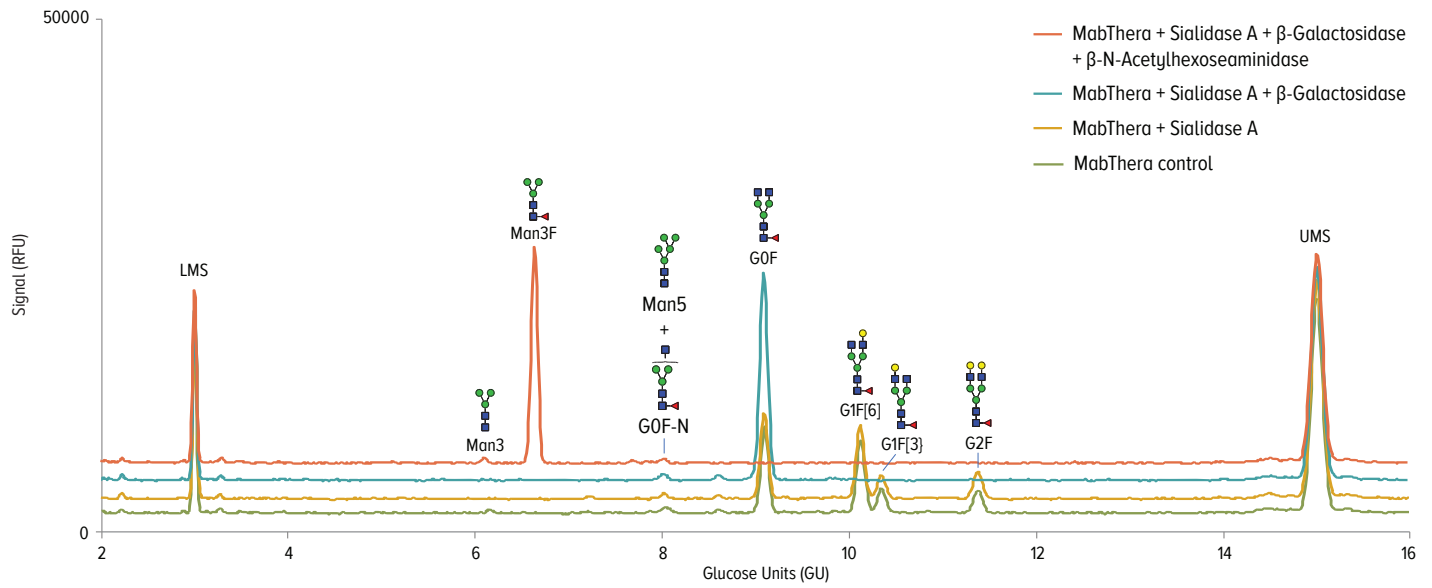


**Figure 2:** General workflow for exoglycosidase treatment, PNGase F digestion, InstantDye labeling and cleanup prior to LC-FLD and CE-LEDIF analysis.

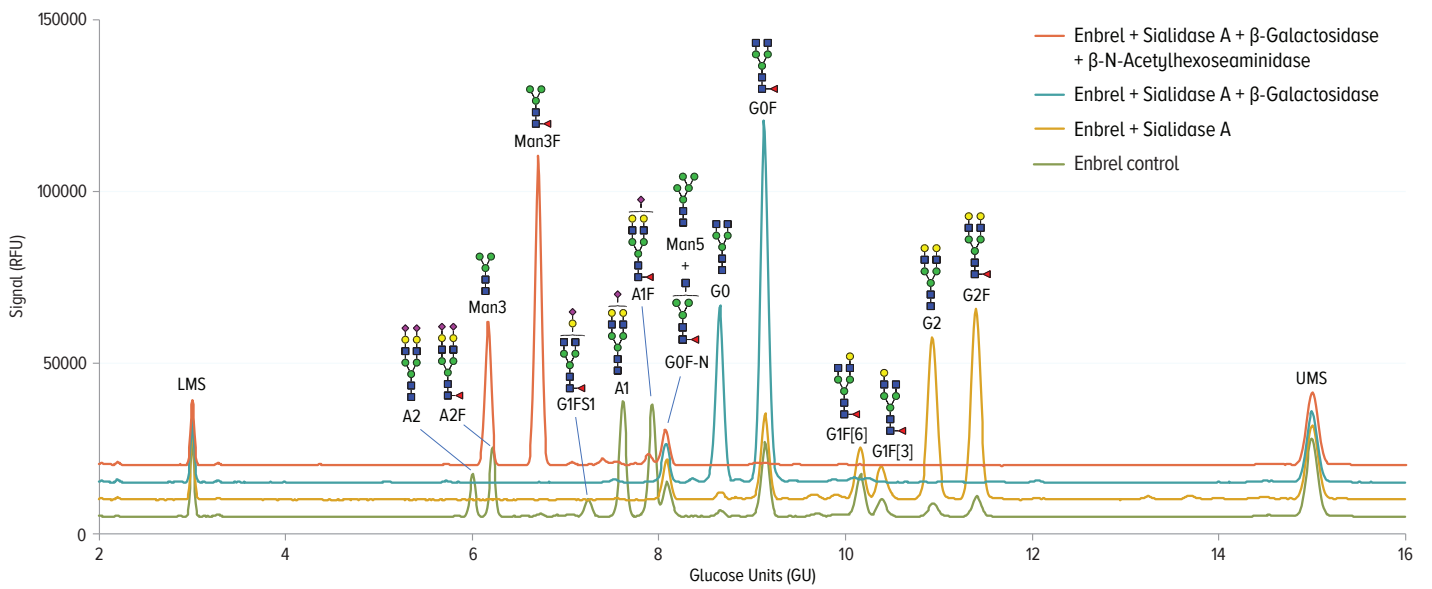
\* Glycoprotein structure adopted from PDB entry DOI: 10.2210/pdb5VGP/pdb  
 PDB structure modified and image generated with Lite Mol: <https://webchemdev.ncbr.muni.cz/LiteMol/>

## RESULTS

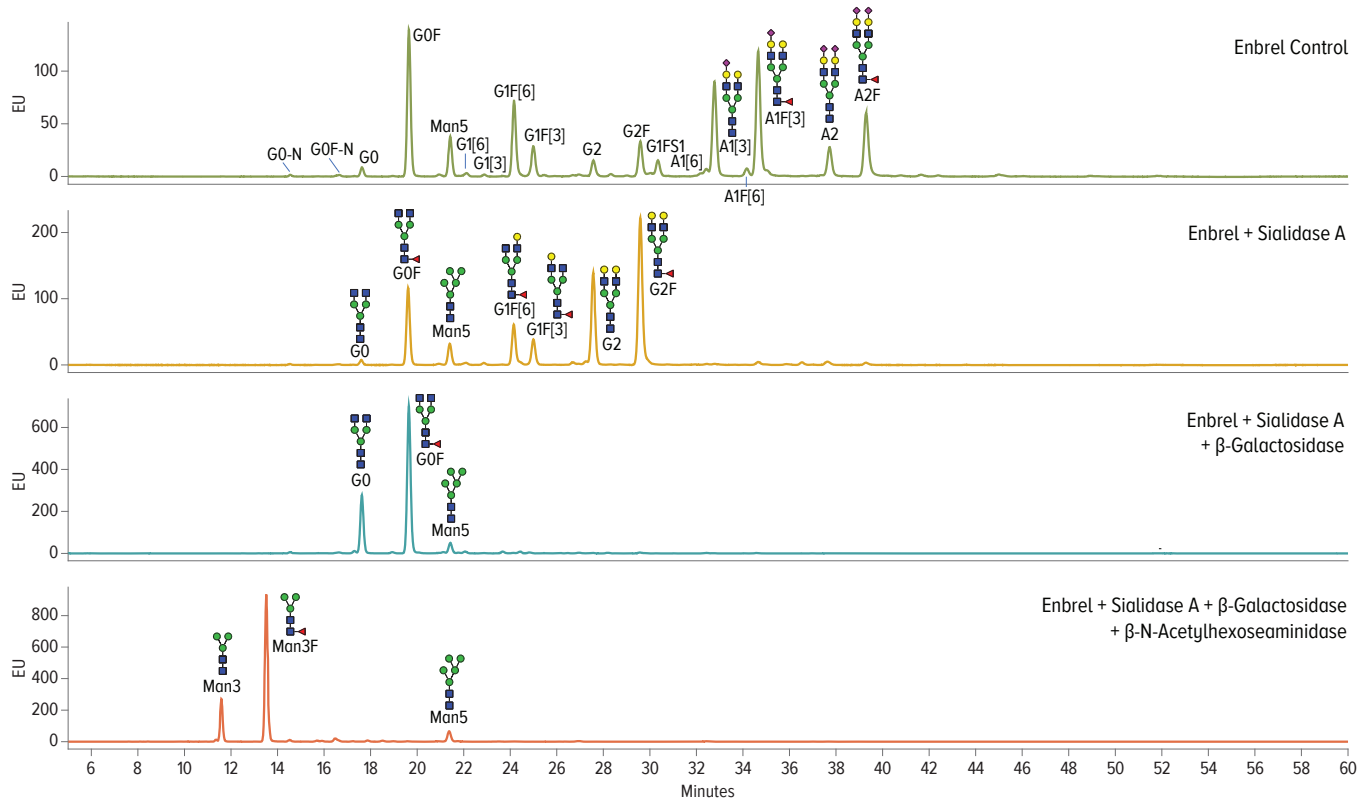
- In parallel digestion of glycoprotein samples with Sialidase A, Sialidase A +  $\beta$ -Galactosidase and Sialidase A +  $\beta$ -Galactosidase +  $\beta$ -N-Acetylhexosaminidase results in the sequential release of terminal sialic acid followed by galactose and N-acetylglucosamine residues.
- Exoglycosidase digestions of MabThera and Enbrel were prepared as described above and N-glycans released and labeled with InstantQ and InstantPC. InstantQ labeled samples from MabThera were analyzed by Gly-Q (Figure 3) and samples from Enbrel were analyzed by both InstantQ and InstantPC (Figures 4 and 5).
- Removal of terminal sialic acid residues results in the shift of sialylated glycans A1, A1F, A2, A2F and G1FS1 to asialylated species G1F, G2 and G2F.
- After the removal of terminal sialic acid, removal of the subsequent terminal galactose residues with  $\beta$ -Galactosidase further shifts the glycan profile towards G0 and G0F.
- Finally, the treatment of glycoprotein samples with all three exoglycosidase enzymes results in the trimming of N-glycans down to the core structures Man3 and Man3F.
- InstantQ labeling and analysis on Gly-Q results in comigrations of Man5 and G0F-N as seen in MabThera (Figure 3). Digestion of MabThera N-glycans with the combination of exoglycosidase enzymes removes the N-Acetylglucosamine from G0F-N resulting in its shift to Man3F.



**Figure 3:** Overlaid Gly-Q electropherograms for sequential digestions of MabThera LMS - lower migration standard (DP3), UMS - upper migration standard (DP15).



**Figure 4:** Overlaid Gly-Q electropherograms for in parallel digestions of Enbrel.



**Figure 5:** Stacked UHPLC chromatograms for Enbrel N-glycans. A) Total N-glycan profile, B) treatment with Sialidase A, C) Sialidase A + β-Galactosidase, D) Sialidase A + β-Galactosidase + β-N-Acetylhexosaminidase.

## CONCLUSIONS

1. Exoglycosidases may be used on intact and denatured glycoprotein samples prior to the release and labeling of N-glycans.
2. The use of exoglycosidases on intact glycoprotein samples allows for a standard digestion protocol prior to release and labeling allowing for flexibility in fluorescent dye selection.
3. Exoglycosidase digestion can aid in resolving peaks which would otherwise coelute or comigrate.

## References

1. Antibody glycosylation and its impact on the pharmacokinetics and pharmacodynamics of monoclonal antibodies and Fc-fusion proteins. Liu L. J. Pharm. Sci. 2015;104(6):1866–1884





**Advancing Glycosciences**

---

Hayward, California **Toll Free** +1 (800) 457-9444 **Phone** +1 (510) 638-6900 **Fax** +1 (510) 638-6919 **Website** [www.prozyme.com](http://www.prozyme.com)  
**Technical** [info@prozyme.com](mailto:info@prozyme.com) **Orders** [orders@prozyme.com](mailto:orders@prozyme.com) **Virtual Patent Marking** [www.prozyme.com/patents](http://www.prozyme.com/patents)

*ProZyme, GlykoPrep, N-Glycanase, InstantQ and Gly-X are trademarks of ProZyme, Inc. in the United States and other countries. All other trademarks are the property of their respective owners.*