

Poster Reprint

**ASMS 2019 ThP-697**

# Streamlined Workflows for N-Glycan Analysis of Biotherapeutics Using InstantPC and 2-AB with LC-FLD-MS

John Yan, Andres Guerrero, Ace G. Galermo,  
Ted Haxo, Sergey Vlasenko, Justin Hyche,  
Tom Rice, Aled Jones

ProZyme, A part of Agilent, Hayward CA

## Introduction

The characterization of N-glycans is an essential part of the biotherapeutic development process. The structure of N-linked glycans can influence the function of glycosylated biotherapeutics<sup>1</sup>, frequently making glycosylation a critical quality attribute (CQA). N-Glycan analysis often involves the labeling of released glycans with a tag to allow for detection by fluorescence (FLD) and mass spectrometry (MS). Many of the frequently used fluorescent tags such as 2-AB<sup>2</sup> are limited with regard to MS sensitivity compared with recently introduced dyes such as InstantPC, and pre-existing N-glycan sample preparation workflows can be time-consuming.<sup>3</sup>

Here we present streamlined workflows for preparation of InstantPC and 2-AB labeled N-glycans coupled with analysis using Agilent LC/FLD/MS instrumentation.

## Experimental

### Sample Preparation

Gly-X with InstantPC and 2-AB workflows (Figure 1) were used to prepare labeled N-glycans from monoclonal antibody rituximab (Rituxan, lot # M190170) and Fc fusion protein etanercept (Enbrel, lot # 1092537), 40 µg protein per preparation.

InstantPC labeled samples were prepared using a developmental protocol on an Agilent Bravo liquid handling system with an adapted Gly-X in-solution deglycosylation protocol followed by instant glycosylamine labeling of released N-glycans (Figure 2), followed by vacuum-driven cleanup of free dye using HILIC SPE.

2-AB labeled samples were prepared per standard Gly-X 2-AB Express manual method with reductive amination chemistry. Released N-glycans were desolvated by vacuum filtration on a solid-state matrix followed by an on-matrix 2-AB labeling step, eliminating the need for glycan drying prior to the labeling step, thereby reducing total sample preparation time. Four replicates of each sample were analyzed with fluorescence/MS detection and relative percent peak areas calculated.

### N-Glycan Analysis

InstantPC and 2-AB labeled N-glycans were separated by hydrophilic interaction liquid chromatography (HILIC) using an Agilent AdvanceBio Glycan Mapping column along with an Agilent 1290 Infinity II UHPLC system with in-line fluorescence detection (Table 1) coupled to an Agilent 6545 LC/Q-TOF mass spectrometer (Table 2).

## Experimental

All HILIC separations were conducted under the conditions described in Table 1. A fixed flow splitter was utilized post-FLD, diverting approximately 50% of the flow to waste and 50% to the MS. MassHunter BioConfirm software was used for data processing, with a personal compound database (PCD).

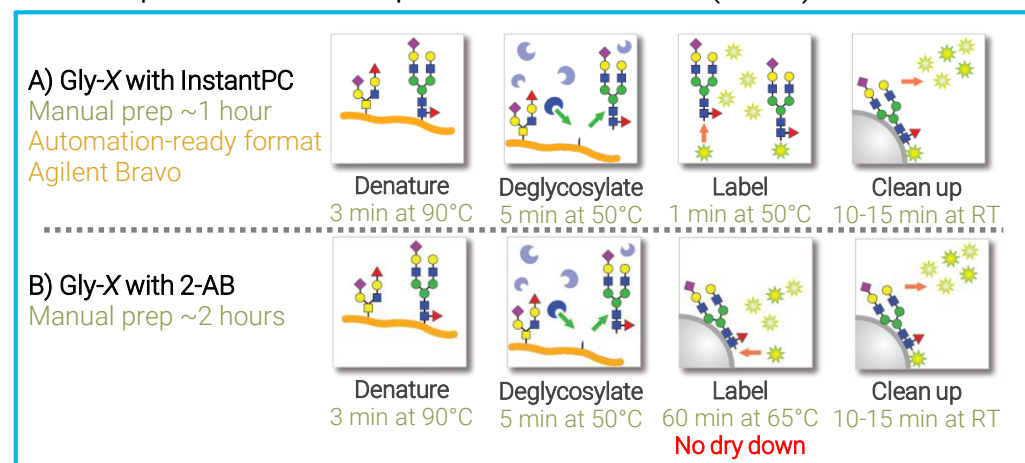


Figure 1. Gly-X N-glycan sample prep. A) InstantPC workflow with in solution deglycosylation and labeling followed by on-matrix cleanup; B) 2-AB workflow with deglycosylation in solution, followed by on-matrix labeling and cleanup.

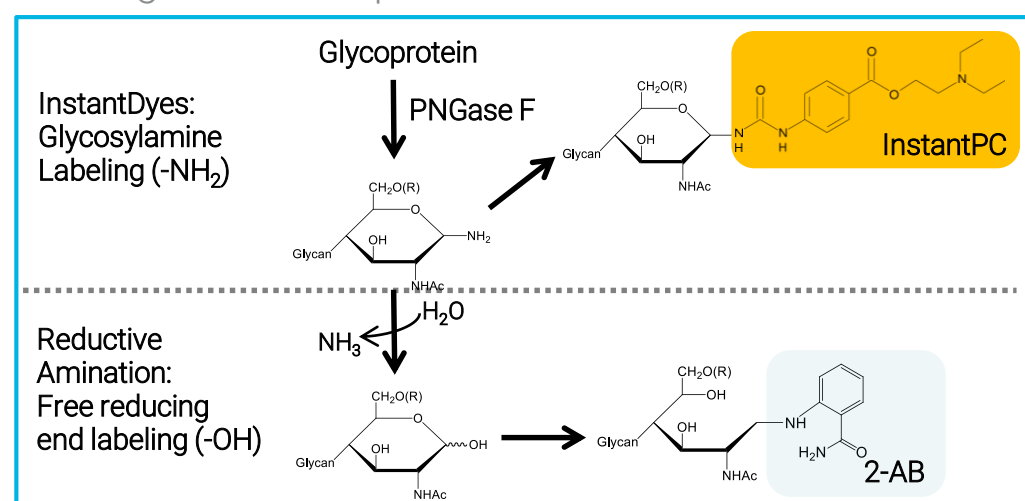


Figure 2. Comparison of InstantPC glycosylamine labeling and traditional reductive amination with 2-AB.

Table 1. UHPLC HILIC/FLD conditions (Agilent 1290).

Parameter	Value																																																																														
Column	Agilent AdvanceBio Glycan Mapping, 2.1 x 150 mm, 1.8 μm (p/n 859700-913)																																																																														
Column Temp	40 °C																																																																														
Mobile Phase	A) 50 mM ammonium formate, pH 4.5 B) Acetonitrile																																																																														
Gradient Program	<table><tr><th colspan="3">InstantPC</th><th colspan="3">2-AB</th></tr><tr><th>Time (min)</th><th>%B</th><th>Flow rate (mL/min)</th><th>Time (min)</th><th>%B</th><th>Flow rate (mL/min)</th></tr><tr><td>0</td><td>80</td><td>0.5</td><td>0</td><td>82</td><td>0.4</td></tr><tr><td>2</td><td>75</td><td>0.5</td><td>2</td><td>82</td><td>0.4</td></tr><tr><td>48</td><td>62</td><td>0.5</td><td>2.5</td><td>77</td><td>0.4</td></tr><tr><td>49</td><td>40</td><td>0.5</td><td>48</td><td>62</td><td>0.4</td></tr><tr><td>51.5</td><td>80</td><td>0.5</td><td>49</td><td>40</td><td>0.4</td></tr><tr><td>52</td><td>80</td><td>0.5</td><td>51.5</td><td>40</td><td>0.4</td></tr><tr><td>60</td><td>80</td><td>0.5</td><td>52</td><td>82</td><td>0.4</td></tr><tr><td></td><td></td><td></td><td>54</td><td>82</td><td>0.4</td></tr><tr><td></td><td></td><td></td><td>58</td><td>82</td><td>0.6</td></tr><tr><td></td><td></td><td></td><td>58.5</td><td>82</td><td>0.6</td></tr><tr><td></td><td></td><td></td><td>60</td><td>82</td><td>0.4</td></tr></table>	InstantPC			2-AB			Time (min)	%B	Flow rate (mL/min)	Time (min)	%B	Flow rate (mL/min)	0	80	0.5	0	82	0.4	2	75	0.5	2	82	0.4	48	62	0.5	2.5	77	0.4	49	40	0.5	48	62	0.4	51.5	80	0.5	49	40	0.4	52	80	0.5	51.5	40	0.4	60	80	0.5	52	82	0.4				54	82	0.4				58	82	0.6				58.5	82	0.6				60	82	0.4
InstantPC			2-AB																																																																												
Time (min)	%B	Flow rate (mL/min)	Time (min)	%B	Flow rate (mL/min)																																																																										
0	80	0.5	0	82	0.4																																																																										
2	75	0.5	2	82	0.4																																																																										
48	62	0.5	2.5	77	0.4																																																																										
49	40	0.5	48	62	0.4																																																																										
51.5	80	0.5	49	40	0.4																																																																										
52	80	0.5	51.5	40	0.4																																																																										
60	80	0.5	52	82	0.4																																																																										
			54	82	0.4																																																																										
			58	82	0.6																																																																										
			58.5	82	0.6																																																																										
			60	82	0.4																																																																										
Injection Volume	1 μL (equivalent to glycans from 0.4 μg protein)																																																																														
Detection	Agilent 1260 Infinity II FLD InstantPC: λ <sub>Ex</sub> 285 nm, λ <sub>Em</sub> 345 nm 2-AB: λ <sub>Ex</sub> 260 nm, λ <sub>Em</sub> 430 nm																																																																														

Table 2. 6545XT Q-TOF

6545XT Q-TOF	
Source	Dual AJS ESI
Gas Temperature	150 °C
Drying Gas Flow	9 L/min
Nebulizer	35 psi
Sheath Gas Temperature	300 °C
Sheath Gas Flow	10 L/min
Vcap	3000 V
Nozzle Voltage	500 V
Fragmentor	120 V
Skimmer	65 V
Mass Range	m/z 600-3000
Scan Rate	1 spectra/sec
Acquisition Mode	High resolution (4 GHz)

HILIC Separation of InstantPC and 2-AB N-Glycans

HILIC separation with fluorescence detection of InstantPC and 2-AB labeled N-glycans from Rituxan and Enbrel results in well resolved peaks for major glycan species from Rituxan and Enbrel (Figures 3 & 4). The HILIC retention time of 2-AB N-glycans is shorter than InstantPC N-glycans, although elution order of N-glycan species is comparable. Critical pairs such as G0F/Man5 and G1F[6]/[3], which are often monitored during the development process of biotherapeutics are well separated with both InstantPC and 2-AB labels, lending to confident determination of relative percentage composition determination. An added benefit of InstantPC is the separation of isoforms G2S1[6]/[3] and G2FS1[6] from Enbrel (Figure 3) compared to 2-AB (Figure 4) using above described chromatography conditions. Analysis with fluorescence detection of InstantPC and 2-AB labeled N-glycans from biotherapeutics Rituxan and Enbrel results in comparable relative percent areas for major glycoforms G0F, G1F[6]/[3], G2F, G2S2 and G2FS2.

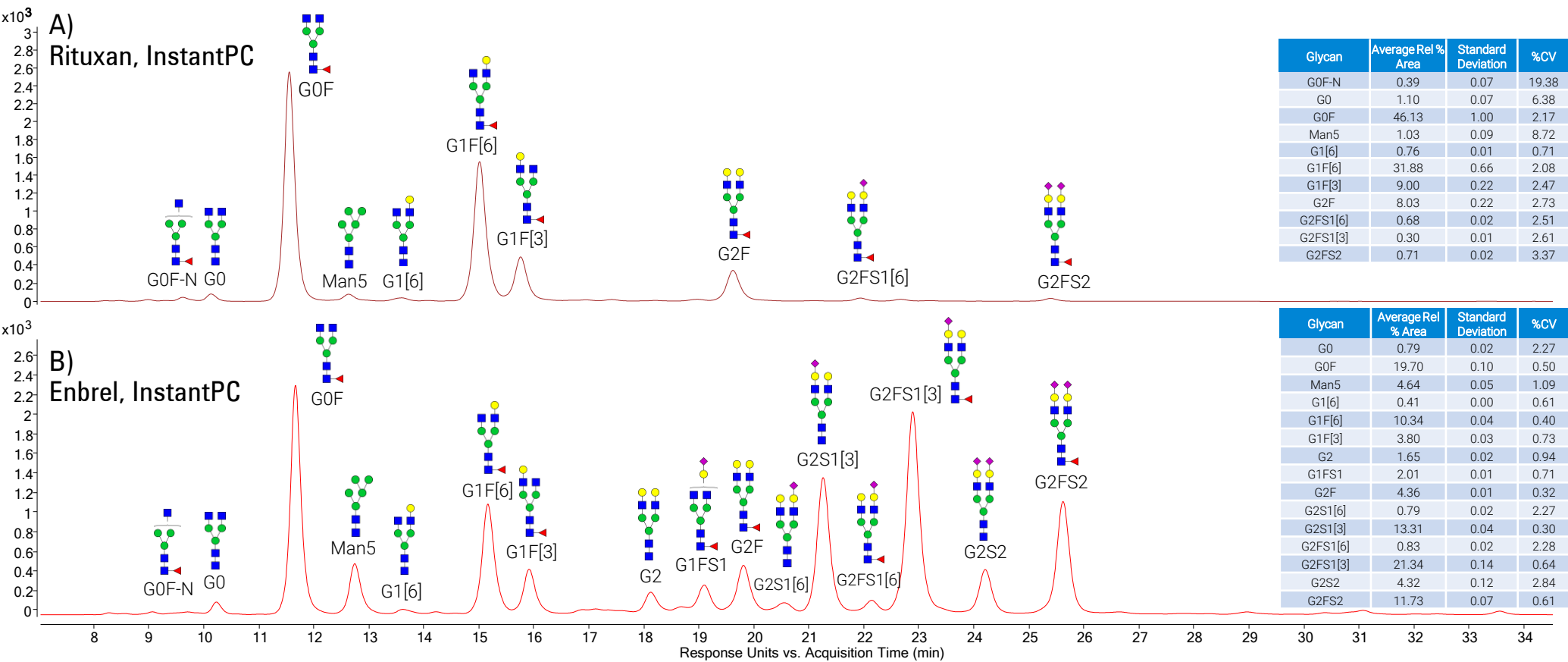


Figure 3. HILIC-UHPLC fluorescence profile of A) Rituxan and B) Enbrel N-glycans labeled with InstantPC. N-Glycan relative percent areas are shown in the inset tables, n = 4.

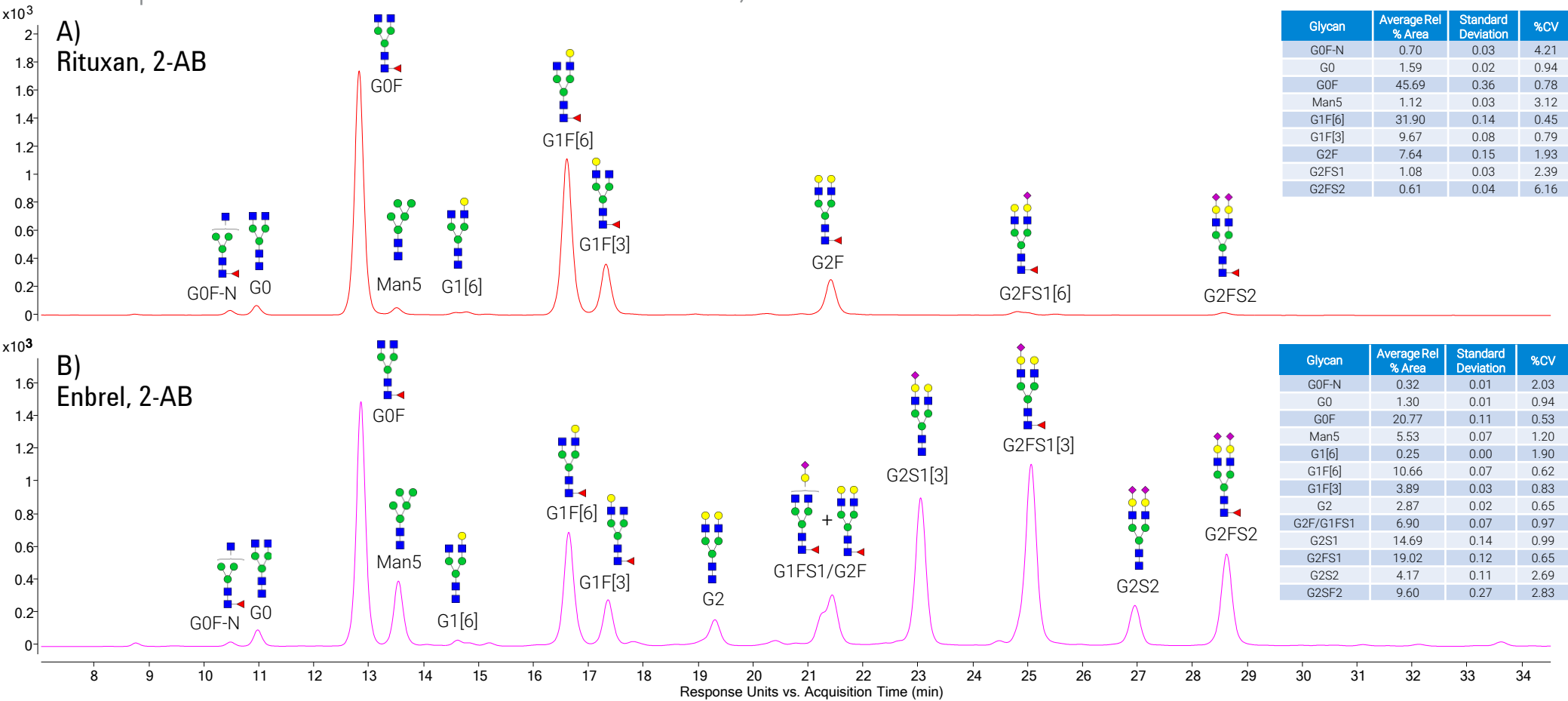


Figure 4. HILIC-UHPLC fluorescence profile of A) Rituxan and B) Enbrel N-glycans labeled with 2-AB. N-Glycan relative percent areas are shown in the inset tables, n = 4.

## FLD and MS detection of InstantPC and 2-AB N-Glycans

InstantPC displays higher fluorescence and MS signal compared to 2-AB (Figure 5). Individual spectra for InstantPC and 2-AB labeled Man5 also show higher MS signal of InstantPC (Figure 6).

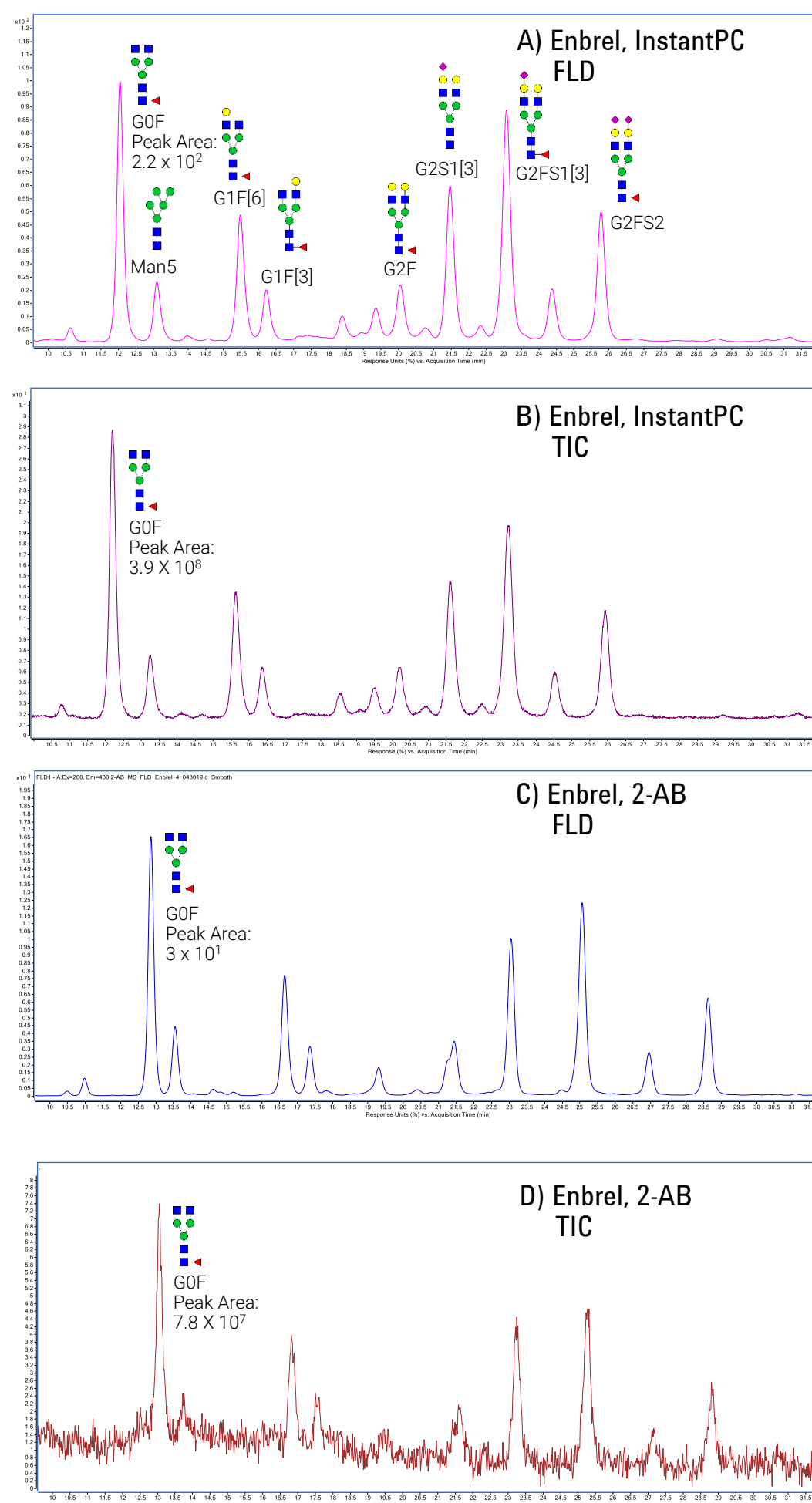


Figure 5. FLD and MS of InstantPC and 2-AB labeled N-glycans from Enbrel. A) InstantPC FLD; B) InstantPC TIC (total ion chromatogram); C) 2-AB FLD; D) 2-AB TIC.

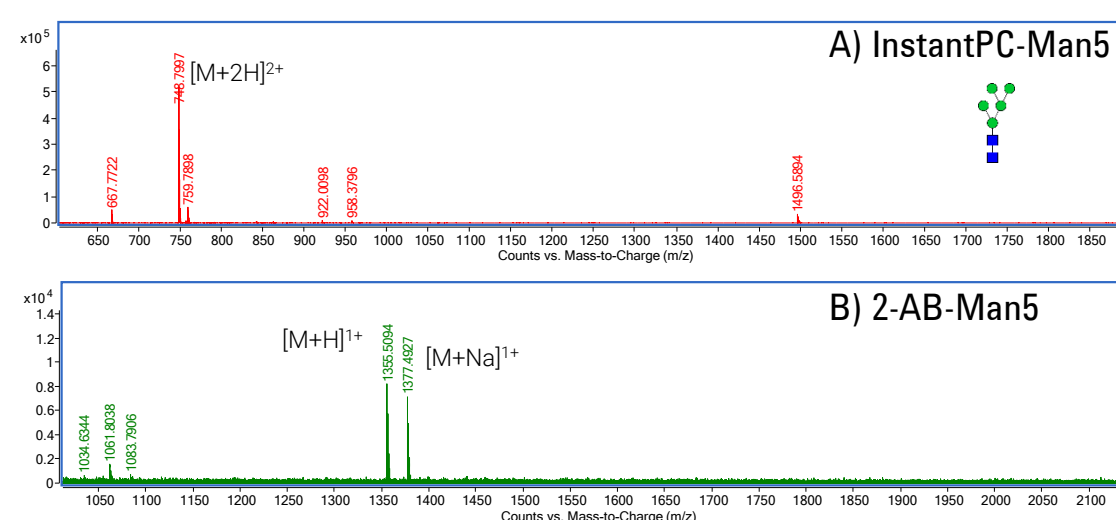


Figure 6. Mass spectrum comparison of Man5 from Enbrel, labeled with A) InstantPC and B) 2-AB.

## Conclusions

- Gly-X sample prep enables 5 minute release of N-linked glycans suitable for labeling by both glycosylamine labeling with InstantPC and reductive amination chemistry with 2-AB.
- On-matrix 2-AB labeling eliminates dry down step.
- Glycan species were profiled by relative fluorescence peak area % and structurally assigned using high resolution tandem mass spectrometry. Reproducibility between sample preparation replicates is high.
- Compared with 2-AB, InstantPC glycans display higher FLD signal and greater MS response in positive mode, allowing for confident detection of low abundance glycan species.

## References

- Liu, L. Antibody Glycosylation and its Impact on the Pharmacokinetics and Pharmacodynamics of Monoclonal Antibodies and Fc-Fusion Proteins. *J. Pharm. Sci.* **2015**, *104*(6), 1866–1884.
- Anumula, K. R. Advances in fluorescence derivatization methods for high-performance liquid chromatographic analysis of glycoprotein carbohydrates. *Anal. Biochem.* **2006**, *350*, 1-23.
- Kimzey, M.; et al. Development of a 5-Minute Deglycosylation Method for High Throughput N-Glycan Analysis by Mass Spectrometry. *ProZyme Technical Note*, Bulletin 4001, Rev E.

For Research Use Only. Not for use in diagnostic procedures.