# UHPLC Analysis of 2-Aminobenzamide-Labeled Glycans with the Vanquish Flex System

Alexandra Manka and Mauro De Pra Thermo Fisher Scientific, Germering, Germany

#### **Key Words**

Biosimilars, Biotherapeutics, Bovine Fetuin, Fluorescence Detector, Glycan Analysis, IgG

#### Goal

To demonstrate successful separation and detection of 2-aminobenzamide (2-AB)-labeled complex glycan libraries on a Thermo Scientific<sup>™</sup> Vanquish<sup>™</sup> Flex UHPLC system equipped with a Vanquish Fluorescence Detector F. Retention time precision for these glycan libraries as well as for a 2-AB-labeled single glycan standard are reported. Area precision and linearity are evaluated for three labeled glycan standards. Further, recommendations for the selection of the right detector flow cell are given.

#### Introduction

Glycans are involved in many biological and physiological processes, such as cell and pathogen recognition, gene expression, or cellular immunity, communication, growth, and development. In biological systems, glycans can be found in a "free state" or in conjugated forms like glycoproteins, glycolipids, or proteoglycans. N-linked glycans are considered an important species in therapeutic protein drug development because there is strong evidence that bioactivity and efficiency of proteins depend on glycosylation. Examples for glycosylated protein-based drugs are recombinant proteins and monoclonal antibodies. The function of these proteins often depends on the structure and type of glycans attached to them.<sup>1</sup> Further, glycans can serve as biomarkers for severe diseases such as cancer.<sup>2,3,4</sup> Thus, glycan analysis is of tremendous interest, especially in the biopharmaceutical industry.

In protein glycosylation studies, typically the glycans are analyzed qualitatively and quantitatively after enzymatical or chemical release from the protein. This approach often leads to quite complex and heterogeneous mixtures of oligosaccharides that require high-efficiency separation techniques prior to characterization, usually done by mass spectrometry (MS). Characterization cannot be accomplished by ultraviolet (UV) detection because glycans lack chromophores; this requires a modification to the detection strategy.



A common approach is to label the glycans with fluorescent tags, separate them with high performance liquid chromatography (HPLC), and detect them with fluorescence detectors (FLD). Even though the tagging procedure and sample cleanup are time consuming, no further sample treatment prior to MS-analysis is needed. Consequently, an HPLC system with a fluorescence detector can be coupled directly to an MS. In addition, fluorescently labeled glycans generally provide better and more MS/MS fragmentation peaks.<sup>5</sup> Therefore, HPLC with fluorescence detection is a commonly used approach for glycan analysis, especially when sensitivity is critical and analysis by MS is also required.



This technical note shows the separation of two 2-AB-labeled glycan libraries and linearity data for three different 2-AB-labeled glycan standards on a Thermo Scientific Vanquish Flex system equipped with a Vanquish Fluorescence Detector F. The system, including two different flow cells, is fully biocompatible. Measurements were conducted with two different columns, the Thermo Scientific<sup>™</sup> GlycanPac<sup>™</sup> AXR-1 column and the Thermo Scientific<sup>™</sup> Accucore<sup>™</sup> 150 Amide HILIC column (hydrophilic interaction chromatography).

### **Experimental**

## Instrumentation

Vanquish Flex UHPLC system (Figure 1), equipped with:

- System Base (P/N VF-S01-A)
- Quarternary Pump F (P/N VF-P20-A)
- Split Sampler FT with 25 µL sample loop (P/N VF-A10-A)
- Column Compartment H (P/N VH-C10-A) with active preheater (P/N 6732.0110)
- Fluorescence Detector F (P/N VH-D51-A) equipped with either 2 μL flow cell (P/N 6079.4330) or 8 μL flow cell (P/N 6079.4230)

#### **Reagents and Materials**

- Deionized (DI) water 18 MΩ×cm resistivity or higher
- Acetonitrile, Optima<sup>™</sup> LC/MS grade (Fisher Scientific P/N 10001334)
- Ammonium formate, Optima LC/MS grade (Fisher Scientific P/N 11377490)
- Formic acid, Optima LC/MS grade (Fisher Scientific P/N 10596814)

#### Standards

- 2-AB-labeled N-glycan library from bovine fetuin (Prozyme P/N GKSB-002)
- 2-AB-labeled N-glycan Immunoglobulin G (IgG) library (Ludger P/N CAB-IGG-01)
- 2-AB labeled NA2 glycan (Ludger P/N CAB-NA2-01)
- 2-AB labeled A1 glycan (Ludger P/N CAB-A1-01)
- 2-AB labeled A2 glycan (Ludger P/N CAB-A2-01)

#### List of Abbreviations

2-AB-NA2:	2-AB-labeled asialo, galactosylated biantennary glycan
2-AB-A1:	2-AB-labeled mono-sialylated, galactosylated biantennary glycan
2-AB-A2:	2-AB-labeled di-sialylated, galactosylated biantennary glycan
ACN:	Acetonitrile
FLD:	Fluorescence detector
PMT:	Photomultiplier tube
RSD:	Relative standard deviation

#### Preparation of Ammonium Formate Buffers

Dissolve 6.30 g (100 mM) or 3.15 g (50 mM) of ammonium formate in 1 L of DI water, respectively. Adjust pH to 4.4 by carefully adding formic acid.

#### Settings for the Vanquish Fluorescence Detector F

The Vanquish Fluorescence Detector F was used with two different excitation/emission wavelength pairs. Sensitivity was set to 8 unless stated otherwise. The lamp mode was set to HighPower during data acquisition and to LongLife during re-equilibration.

The emission signal stray light can be suppressed by setting the appropriate cut-off filter-wheel. Several settings are available for the emission filter in the Vanquish Fluorescence Detector F: Auto, 280, 370, 435, and 530 nm. With the Auto setting for the filter wheel, the most suitable cut-off filter is selected automatically, which was 370 nm for the experiments described here.



Figure 1. Vanquish Flex UHPLC System with Fluorescence detector.

Chromatographi	c Conditions I			Chrom
Column:	GlycanPac AXR-1, 2.1 × 250 mm, 1.9 µm (P/N 088135)			Column
Mixer volume:	350 μL + 50 μl			Mixer v
Mobile phase A:	Acetonitrile			Mobile
Mobile phase B:	DI water			Mobile
Mobile phase C:	Ammonium formate, 100 mM, pH 4.4			Gradien
Gradient:	Time (min)	%В	%C	_
	-5.0	93	7	_
	0	93	7	_
	60	31	50	_
	60.5	0	100	
	65.5	0	100	_
	75	93	7	Flow ra
	100	93	7	Temper
Flow rate:	0.40 mL/min			
Temperature:	Column compartment:30 °CDetector flow cell:35 °C		Detection parame	
Detection	Excitation wave	length:	250 nm	
parameters:	Emission wavelength:		430 nm	—
	Lamp mode:		HighPower	—
	Sensitivity:		7	—
	Data collection rate:		10 Hz	Sample
	Response time: 0.2 s			
Sample:	1 nmol/mL 2-AB-labeled N-glycan library from bovine fetuin Sample was dissolved in DI water			_

hromatographic Conditions II			
olumn:	Accucore 150-Amide-HILIC, 2.1 × 150 mm, 2.6 μm (P/N 16726-152130)		
lixer volume:	350 μL + 50 μL		
obile phase A:	Acetonitrile		
obile phase B:	Ammonium formate, 50 mM, pH 4.4		
radient:	Time (min)	%В	
	0	20	
	30	42	
	30.5	50	
	32	50	
	32.5	20	
	60	20	
ow rate:	0.40 mL/min		
emperature:	Column compartment: Detector flow cell:	50 °C 45 °C	
etection	Excitation wavelength:	320 nm	
arameters:	Emission wavelength:	420 nm	
	Lamp mode:	HighPower	
	Sensitivity:	8	
	Data collection rate:	5 Hz	
	Response time:	1 s	
amples:	1 nmol/mL 2-AB labeled <i>N</i> -glycan lgG library 1 nmol/mL 2-AB labeled NA2 glycan 1 nmol/mL 2-AB labeled A1 glycan 1 nmol/mL 2-AB labeled A2 glycan Samples were dissolved with appropriate amounts of 82/18 acetonitrile/ammonium formate 50 mM pH 4.2.		

## **Data Processing**

Thermo Scientific<sup>™</sup> Dionex<sup>™</sup> Chromeleon<sup>™</sup> Chromatography Data System software, version 7.2 SR 3 **Results and Discussion** 

## Separation of 2-AB-Labeled Glycan Library Samples

The *N*-linked glycans from bovine fetuin are heterogeneous in terms of charge, as their sialic acid content can vary between 0 and 5. Here, we show that these rather diverse glycans samples can be successfully separated and detected with the Vanquish Flex system equipped with a Vanquish Fluorescence Detector F. Over 60 peaks were detected in the sample of 2-AB-labeled *N*-fetuin glycans. All peaks eluted within 50 min in a ternary gradient separation (Figure 2).



Figure 2. Separation of 50 pmol bovine fetuin 2-AB-N-glycans measured with the Vanquish FLD F using a 2  $\mu$ L flow cell. Chromatographic conditions correspond to those described in the table "Chromatographic Conditions I".

The IgG sample is less heterogeneous, consisting primarily of neutral and mono-sialylated glycans. The IgG 2-AB-*N*-glycans could be separated on an Accucore 150-Amide-HILIC column, separating the sample by size, delivering over 20 peaks within 30 minutes (Figure 3).



Figure 3. Separation of 5 pmol IgG 2-AB-*N*-glycans measured with the Vanquish FLD F using an 8 µL flow cell. Chromatographic conditions correspond to those described in the table "Chromatographic Conditions II".

#### **Retention Time Precision**

Retention time precision (RSD) is essential in labeled glycan analysis for peak identification. This applies particularly to cases where the relative retention times are expressed based on the retention time of glucose oligomers, with the purpose of structure assignment based on a database.<sup>6</sup> A method with repeatable retention times allows the injection of the unknown sample and the glucose ladder standard in separated runs, and subsequent peak identification by comparison of the chromatograms. This avoids the practice of spiking the sample with the glucose standard, hence eliminating the risk of co-elution between glucose oligomers and analyte peaks.

Retention time precision was measured for repeated injections of the bovine fetuin 2-AB-*N*-glycans and the IgG 2-AB-*N*-glycans, as well as for a single standard of 2-AB-NA2 glycan



For the bovine fetuin 2-AB-*N*-glycan separation with a ternary gradient shown in Figure 2, 10 arbitrarily selected peaks were evaluated. The retention time precision for these peaks is shown in Figure 4. The RSD for this separation is generally  $\leq 0.1\%$ , which is excellent for a ternary gradient separation.

For the IgG 2-AB-*N*-glycan separation shown in Figure 3, 10 arbitrarily selected peaks were evaluated. The retention time precision for these peaks is shown in Figure 5. For this separation, the retention time RSD is even better and < 0.04% for all peaks.

For all peaks shown in this technical note, the RSD was far below the value 0.3% that is commonly reported.<sup>7</sup> In addition, retention time precision was evaluated for six consecutive injections of the single standard 2-AB-NA2; an RSD of 0.008% was measured in this case. Retention time precision data can be found in Table 1.



Figure 4. Retention time precision for 10 arbitrarily selected peaks from the bovine fetuin 2-AB-*N*-glycan separation shown in Figure 2.

Figure 5. Retention time precision for 10 arbitrarily selected peaks from the IgG 2-AB-*N*-glycan separation shown in Figure 3.

Table 1. Retention time precision (n=6). Fetuin glycans from Figure 2, RSD% represents average of 10 peaks; lgG glycans from Figure 3, RSD% represents average of 10 peaks; 2-AB-NA2 standard, measured with "Chromatographic Conditions II".

	Fetuin 2-AB- <i>N</i> -glycans	lgG 2-AB- <i>N</i> -glycans	2-AB-NA2
Retention time precision (RSD%)	0.079	0.023	0.008

## **Area Precision and Linearity**

Area precision was evaluated for five consecutive injections of 1 pmol 2-AB-NA2, and the measured RSD was 0.92%. Linearity was evaluated for three different glycan standards (2-AB-NA2, 2-AB-A1, and 2-AB-A2) and two different flow cells (2  $\mu$ L and 8  $\mu$ L), respectively. Linearity was excellent in all cases and interpolation resulted in r<sup>2</sup> well above 0.99 (Table 2).

Table 2. Linearity obtained for 3 different glycan standards with different sample amounts ( $n \ge 5$ ) in two different flow cells, respectively. Same chromatographic conditions as in Figure 3.

Sample	Range (pmol)	r² in 2 µL flow cell	r² in 8 µL flow cell
2-AB-NA2	0.05 - 5.00	0.9998	0.9999
2-AB-A1	0.50 – 10.00	0.9976	0.9976
2-AB-A2	0.25 - 5.00	0.9996	0.9976



## Conclusions

The Vanquish Flex system equipped with the Vanquish Fluorescence Detector F is a powerful tool for the analysis of 2-AB-labeled glycans. The retention time precision of the system is excellent. Hence, this system can be effectively used for separation of labeled glycans and peak recognition by customized or public databases.

Linearity of the method was excellent for detection of samples in the range spanning from low femtomole to low picomole, independent of the flow cell. Most likely the linearity range stretches to much higher concentrations, however this aspect is outside the purposes of this study and was not investigated.

The Vanquish Fluorescence Detector F can be equipped with two different flow cells, an 8  $\mu$ L standard flow cell and a 2  $\mu$ L micro flow cell for minimized extra-column band dispersion, both biocompatible. Figure 6 shows an overlay of measurements obtained by use of these two different flow cells.



Figure 6. Overlay of three peaks from Figure 3, measured using an 8  $\mu$ L flow cell (black), with peaks detected with an identical experiment but a 2  $\mu$ L flow cell (blue).

Although the size of the flow cell is not critical for this application, with the 8  $\mu$ L flow cell peaks were higher and broader, compared to those measured with the 2  $\mu$ L flow cell. The increase in peak height leads to lower limits of detection while the increased peak width impairs resolution. Consequently, the 8  $\mu$ L flow cell should be the preferred choice for methods with very low amounts of sample and very well resolved peaks (as in Figure 3), whereas the 2  $\mu$ L flow cell should be chosen when higher resolving power is required (as in Figure 2).<sup>8</sup>

## References

- Bertozzi, C. R.; Freeze, H. H.; Varki; A.; Esko, J. D. Glycans in Biotechnology and the Pharmaceutical Industry, Essentials of Glycobiology, Second Edition; Cold Spring Harbor Laboratory Press: New York, 2009; Chapter 51.
- Kobata, A.; Amano, J. Altered glycosylation of proteins produced by malignant cells, and application for the diagnosis and immunotherapy of tumours. *Immunol. Cell. Biol.*, 2005, *83*, 429-439.
- 3. Dube, D.H.; Bertozzi, C.R., Glycans in cancer and inflammation--potential for therapeutics and diagnostics. *Nat Rev Drug Discovery* **2005**, *4*, 477-488.
- Arnold J. N.; Saldova R.; Galligan M.C.; Murphy T. B.; Mimura-Kimura Y.; Telford J. E.; Godwin A.K.; Rudd P.M., Novel Glycan Biomarkers for the Detection of Lung Cancer. *J. Proteom. Res.* 2011, *10*(4), 1755-1764.
- 5. Thermo Scientific PN 70513: Novel Glycan Column Technology for the LC-MS Analysis of Labeled and Native N Glycans Released from Proteins and Antibodies, Thermo Fisher Scientific. <u>http://www.thermoscientific.com/content/dam/tfs/ATG/</u> <u>CMD/CMD%20Documents/posters/PN-70513-Novel-</u> <u>Glycan-Column-Technology-PN70513-EN.pdf</u>
- Thermo Scientific TN 147: UHPLC Analysis of 2-ABlabeled Dextran Ladder and Assignment of Glucose Units to Unknown Glycans, Thermo Fisher Scientific. http://www.thermoscientific.com/content/dam/tfs/ ATG/CMD/CMD%20Documents/TN-147-Analysis-2-AB-labeled-Dextran-Ladder-Glucose-Units-TN70926-E.pdf
- Domann P. J.; Pardos-Pardos A. C.; Ludger Fernandes D.; Spencer D. I. R.; Radcliffe C. M.; Royle L. Dwek R. A.; Rudd P. M., Separation-based glycoprofiling approaches using fluorescent labels. *Practical Proteomics*, 2007, 1/2007, 70-76.
- Dionex Technical Note 92: Combining Fluorescence Detection with UHPLC: An Overview of the Technical Requirements; Dionex Corp.

#### www.thermofisher.com/chromatography

©2016 Thermo Fisher Scientific Inc. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

Africa +43 1 333 50 34 0 Australia +61 3 9757 4300 Australia +43 810 282 206 Belgium +32 53 73 42 41 Brazil +55 11 3731 5140 Canada +1 800 530 8447 China 800 810 5118 (tree call domestic) 400 650 5118 TN71673-EN 09/16S

Japan +81 6 6885 1213 Korea +82 2 3420 8600 Latin America +1 561 688 8700 Middle East +43 1 333 50 34 0 Netherlands +31 76 579 55 55 New Zealand +64 980 6700 Norway +46 8 556 468 00

Russia/CIS +43 1 333 50 34 0 Singapore +65 6289 1190 Sweden +46 8 556 468 00 Switzerland +41 61 716 77 00 Taiwan +886 2 8751 6655 UK/Ireland +44 1442 233555 USA +1 800 532 4752

