

Towards a More Complete Glycome: Advances in Ion Chromatography-Mass Spectrometry (IC-MS) for Improved Separation and Analysis of Carbohydrates

Neil G. Rumachik,* Tian Tian,* Yuanxue Hou, Charanjit Saini, Zoltan Szabo, Yan Liu, Chris Pohl, Thermo Fisher Scientific, 1228 Titan Way, Sunnyvale, CA, 94086; *contributed equally to this work

ABSTRACT

Purpose: Demonstrate the coupling of high performance anion exchange (HPAE) chromatography with Orbitrap mass spectrometry to enable more complete glycome characterization

Methods: Developed novel Thermo Scientific™ Dionex™ CarboPac™ PA300 column to use with a Thermo Scientific™ Dionex™ ICS-6000 HPIC™ coupled to a Thermo Scientific™ Q Exactive™ Hybrid Quadrupole Orbitrap™ mass spectrometer for structural characterization of neutral and charged glycans

Results: Novel anion exchange column technology improves resolution of neutral glycans while maintaining capabilities to separate charged species. Improved platform performance is demonstrated via analysis of complex oligosaccharides from biologics and food samples

INTRODUCTION

Glycosylation is one of the most prevalent and diverse types of post-translational modification (PTM), and it is known to be important for myriad cellular and developmental processes, including cell adhesion, molecular trafficking and clearance, receptor activation, signal transduction, and endocytosis. This PTM is often very challenging to study, as glycosylation is non-template driven, heterogeneous, and substoichiometric. O-linked glycan sample preparation is particularly challenging, as there is no known enzyme for the universal release of O-glycans from biomolecules. Additionally, glycans are often complex both in structure and composition, often creating high-density, heterogenous clusters of glycans on regions of biomolecules like mucins. Therefore, high resolution chromatographic separations become important for subsequent mass spectrometric characterization.

To date, few tools are available for the analysis of the glycome without derivatization, a process which is known to introduce issues such as differential loss of sialic acid and incomplete labeling. We have previously reported the use of ion chromatography-mass spectrometry (IC-MS) to analyze native sialylated and sulfated glycans [1]. Here, we expand upon our previous reports, introducing the Dionex CarboPac PA300 column, a new introduction into the Dionex CarboPac suite of columns that uses novel column technology to enable the simultaneous separation of neutral and charged glycans without the need of derivatization. Using the column in an IC-MS workflow takes advantage of the improved chromatographic separation as well as the high resolution and mass accuracy of the Orbitrap mass analyzer. The MS data and tandem MS/MS spectra provide diagnostic fragment ions, allowing for the reliable structural annotation of heterogenous glycans. The unique HPAE-PAD/MS workflow presented here provides complementary often orthogonal information to that of other standards for O-glycan analysis.

MATERIALS AND METHODS

Sample Preparation

For glycan release, samples were incubated in 50 mM NaOH containing 1 M NaBH₄ at 50 °C for 16 h. After incubation, samples were kept on ice and neutralized by careful addition of formic acid. The neutralized samples were evaporated to dryness in a speedvac. The samples were then washed three times with 0.5 mL methanol. Afterwards, samples were dissolved in 0.5 mL of 0.1 % TFA solution (in water). Free and released glycan samples were passed through porous graphitic carbon (Thermo Scientific™ HyperSep™ filter plates, 40 µL). After washing the resin twice with 0.5 mL of 0.1 % TFA solution, oligosaccharides were eluted in 400 µL of 40 % acetonitrile containing 0.1 % TFA and evaporated to dryness.

High performance anion exchange chromatography (HPAEC)

A Dionex ICS-6000 HPIC dual channel IC system coupled with a Thermo Scientific™ Dionex™ AS-AP Autosampler was used for glycan separations. The system was equipped with a Thermo Scientific™ Dionex™ BorateTrap Inline Trap Column between the gradient mixer and injection valve. A Dionex CarboPac PA300 analytical column (2 x 250 mm) and guard column (2 x 50 mm) were used at 0.25 mL/min flow rate for separations. The column effluent was sent either to the PAD or MS for detection of glycans. In the system configuration with MS, a Thermo Scientific™ Dionex™ ERD 500 electrolytically regenerated desalter (anion, 2 mm) was placed prior to MS ionization.

Mass spectrometry

All MS experiments were performed on a Q Exactive Hybrid Quadrupole Orbitrap mass spectrometer in negative mode ESI with Labile Compound User License. The mass spectrometer was operated with full mass scan: m/z 400-2,000 at a resolution of 60,000 (FWHM at m/z 200). The AGC target was set to 3×10^6 with maximum IT: 120 ms. Data dependent MS² were acquired using the Top 10 method. MS² spectra were interpreted using the UniCarb-DB database.

RESULTS

Figure 1 shows the improved separation of neutral glycan species with the new Dionex CarboPac PA300 column. While maintaining the ability separate charged species towards the latter half of the run, the new column technology affords improved separation of early eluting glycans. This introduces the ability to tailor HPAE-PAD/MS workflows based on sample composition, expanding the toolbox of technologies to study and characterize the glycome.

Figure 1. Chromatograms of released O-glycans from a porcine gastric mucin type III sample for the Dionex CarboPac PA200 (bottom trace) and Dionex CarboPac PA300 (top trace). The window in which neutral glycans elute from the columns is shown with dashed lines.

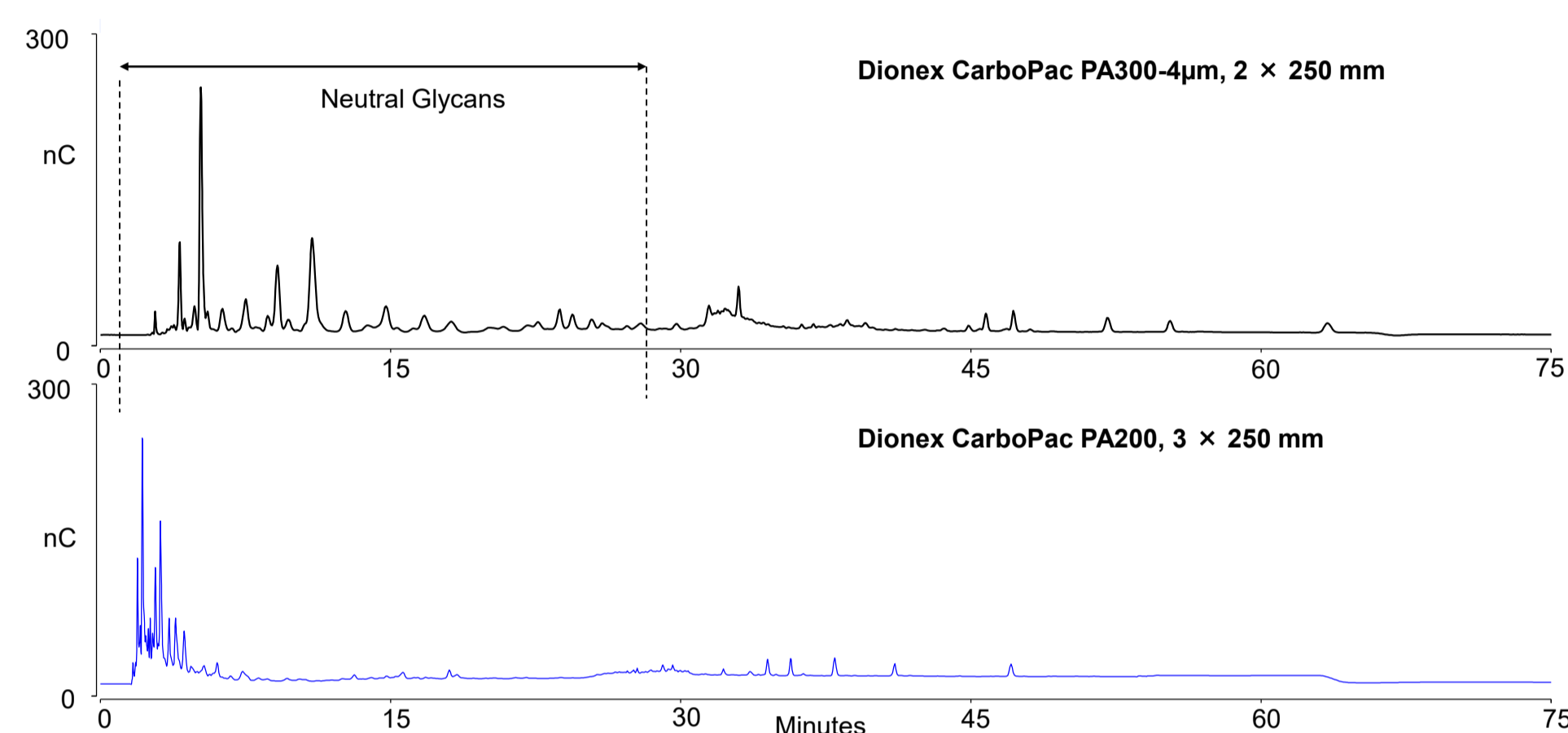


Figure 2. Base peak chromatogram of porcine gastric mucin type III O-glycans injected into a Q Exactive HF-X Hybrid Quadrupole-Orbitrap Mass Spectrometer.

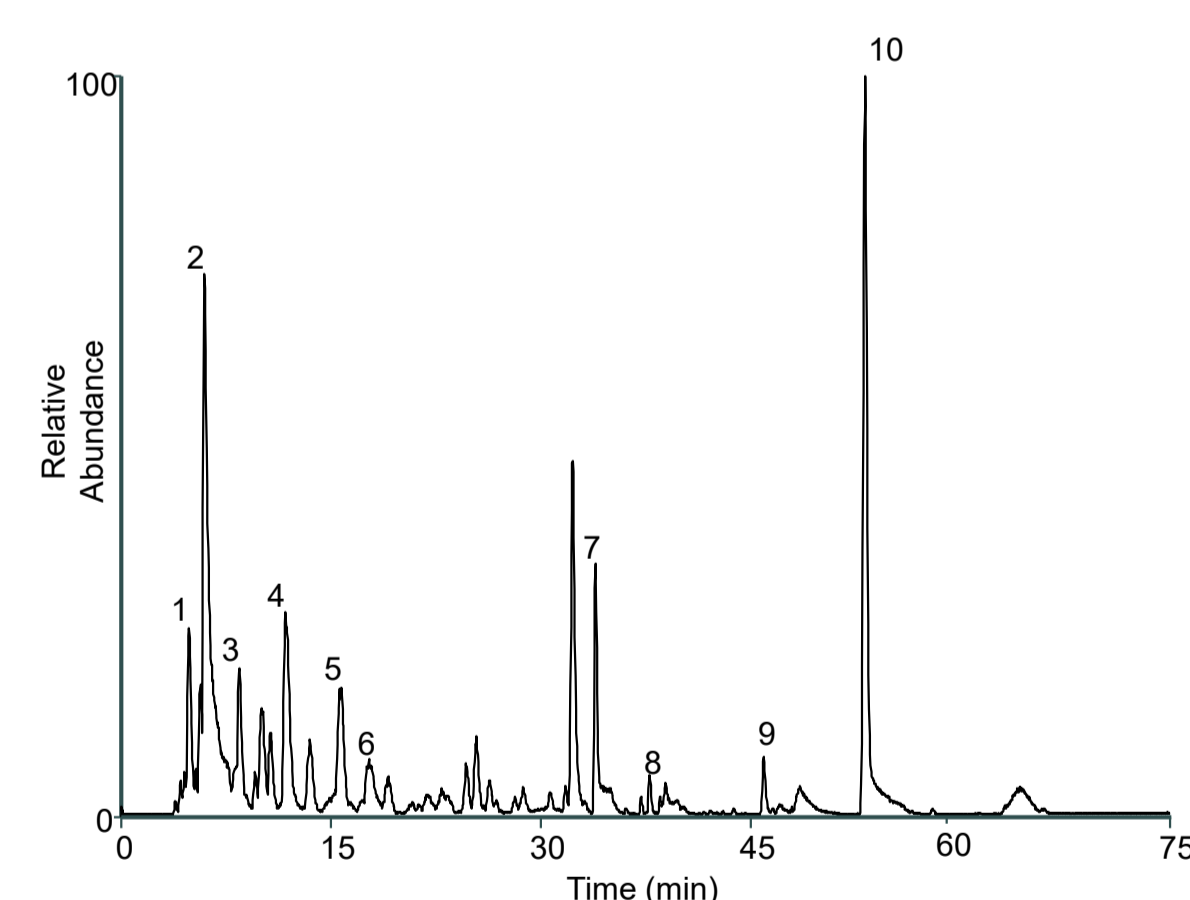
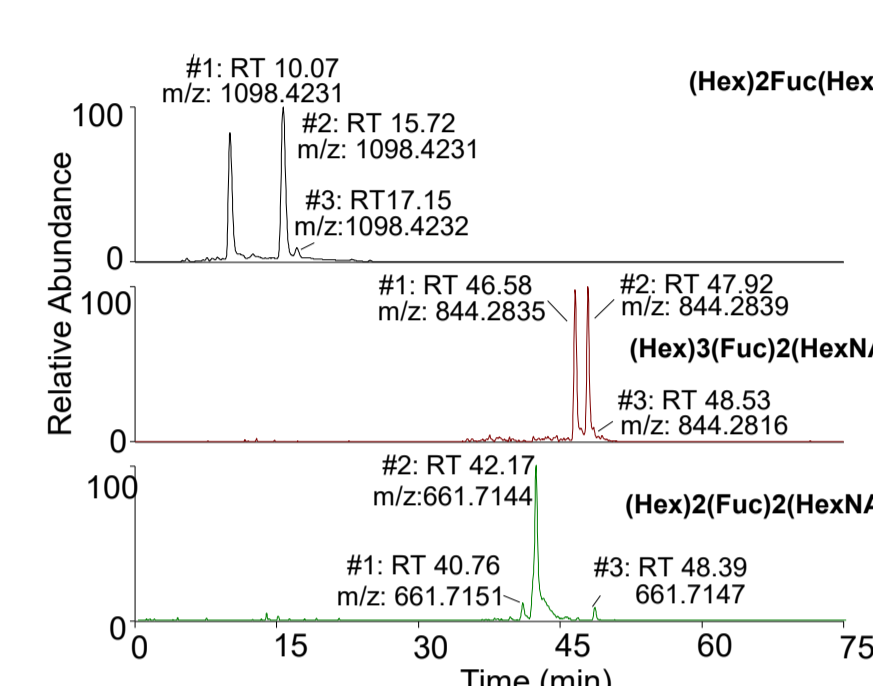


Table 1. Glycan identities of the labeled peaks in Figure 3.

Peak	RT (min)	Observed m/z	Theoretical m/z	Mass accuracy (ppm)	Composition	Ion
1	4.85	733.2896	733.2884	1.6465	HexFuc(HexNAc)2	[M-H] ⁻
2	6.08	530.2080	530.2090	1.8860	HexFucHexNAc	[M-H] ⁻
3	8.46	733.2892	733.2884	1.0910	HexFuc(HexNAc)2	[M-H] ⁻
4	11.97	1041.4006	1041.3992	1.4403	(Hex)2(Fuc)2(HexNAc)2	[M-H] ⁻
5	15.58	1098.4233	1098.4206	2.4581	(Hex)2Fuc(HexNAc)3	[M-H] ⁻
6	17.66	1203.4506	1203.4520	1.1633	(Hex)3(Fuc)2(HexNAc)2	[M-H] ⁻
7	34.01	821.3063	821.3045	2.1916	NeuAcHexFucHexNAc	[M-H] ⁻
8	38.19	852.2811	852.2848	4.3413	N-glycan, hybrid, sulfated	[M-2H] ²⁻
9	45.92	1121.3582	1121.3560	1.9619	(Hex)2(Fuc)2(HexNAc)2-S	[M-H] ⁻
10	53.19	813.2474	813.2452	2.7052	HexFuc(HexNAc)2-S	[M-H] ⁻

Hex: Hexose; Fuc: Fucose; HexNAc: N-Acetylhexosamine; NeuAc: N-Acetylneuraminic acid; S: Sulfate

Figure 3. The Dionex CarboPac PA300 column provides isomer resolution of both neutral and sulfated O-glycans from a porcine gastric mucin type III sample.



The excellent resolution achieved by the Dionex CarboPac PA300 column is especially important for compositionally identical but chemically distinct isomeric species. Coupled with the high mass accuracy of the Orbitrap mass analyzer, separation by HPAE chromatography affords the identification of otherwise coeluting isomers, as in the three examples shown in Figure 3 for (Hex)₂Fuc(HexNAc)₃, (Hex)₃(Fuc)₂(HexNAc)₄-S, and (Hex)₂(Fuc)₂(HexNAc)₃-S.

CONCLUSIONS

- High performance anion exchange (HPAE) offers highly efficient chromatographic O-glycan separations
- HPAE shows outstanding separation of sulfated O-Glycan alditols and their isomers
- The coupling of HPAE to an Orbitrap mass spectrometer enables the collection of highly informative MS² spectra, allowing reliable structural annotation
- The Dionex CarboPac suite of columns affords workflow solutions to tailor analysis based on specific sample composition
- IC-MS provides access to the acidic subset of the glycome and offers complementary data to RPLC-MS-based approaches

REFERENCES

- Anal. Bioanal. Chem. 2017;409(12):3089-3101.
- Mol. Cell Proteomics. 2017;16(5):743-758.

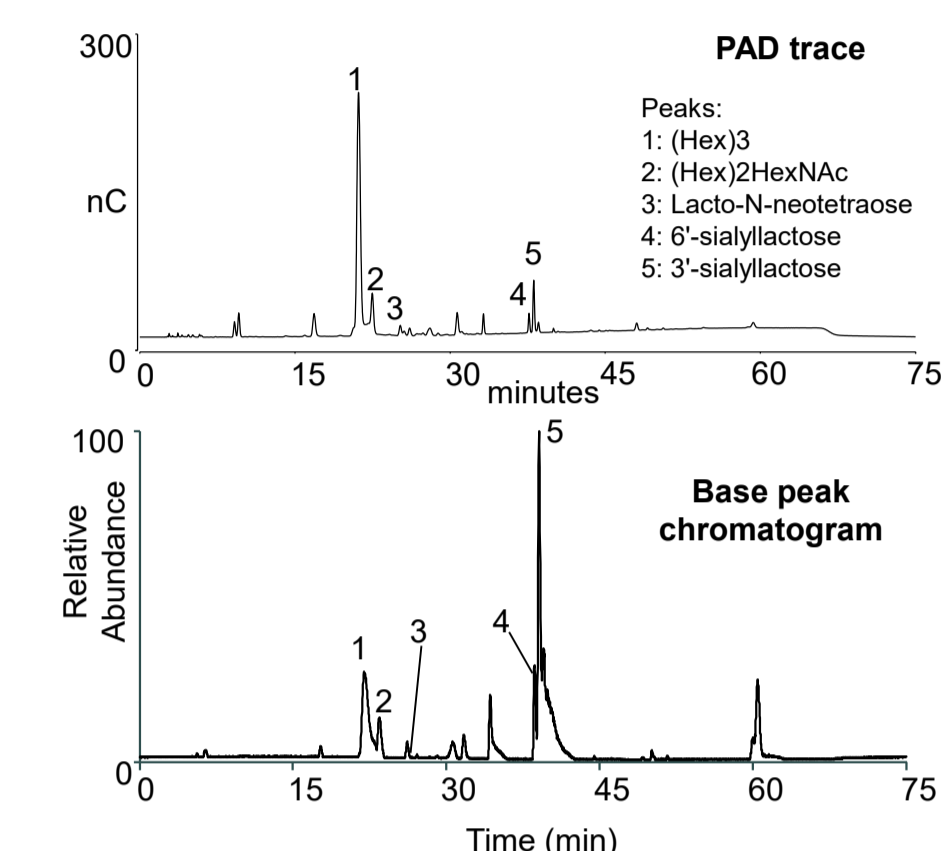
ACKNOWLEDGEMENTS

We would like to thank Jeff Rohrer, Sachin P. Patil, and Pranathi R. Perati at Thermo Fisher Scientific for helpful discussions and critical review.

TRADEMARKS/LICENSING

© 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.

Figure 4. Free oligosaccharides from commercial lactose-free milk with the Dionex CarboPac PA300 column.



Structural annotation of oligosaccharides in foods has become increasingly important, as these glycans are often used as prebiotics to promote healthy gut microbiomes. Figure 4 shows the ability of CarboPac™ PA300 column to separate neutral, non-sialylated glycans from charged, sialylated structures. A selection of peaks are numbered, and the corresponding peak identity is shown in the legend of the upper panel.