

Glycopeptide Analysis Using Electron Transfer Dissociation and Porous Graphite Chromatography

Rosa Viner, Terry Zhang, Vlad Zabrouskov

Thermo Fisher Scientific, San Jose, CA

Abstract

Purpose: N-glycopeptide analysis by nano-LC/MS² using Electron Transfer Dissociation (ETD).

Methods: Glycoprotein digests were analyzed by nano-LC ESI using different types of HPLC columns and a Thermo Scientific LTQ XLTM mass spectrometer equipped with an ETD source option.

Results: A graphic carbon column (HypercarbTM, Thermo Scientific) demonstrated excellent capabilities for glycopeptide analysis especially in the case of short hydrophilic peptides containing bi- or tri-antennary glycan chains without any enrichment. Formation of metal adducts on the Hypercarb column promotes higher charged species and, as a result, promotes ETD fragmentation of glycopeptides. The combination of porous graphite chromatography and ETD-MS/MS is demonstrated to be a useful and flexible tool for studying glycosylation and identifying PTM sites.

Introduction

Glycosylation plays a key role in controlling numerous biological processes. It is one of the most diverse and complex post-translational modifications (PTMs) found on proteins and its characterization remains a great analytical challenge. LC/MS/MS is the most powerful and versatile technique for glycopeptide structure elucidation. However, commonly used collisional-induced dissociation (CID) has limitations for determining the modification site due to the labile nature of glycan modifications. As a new dissociation technique, Electron Transfer Dissociation (ETD) preserves labile PTMs and provides a new and powerful tool that makes the identification of modification sites possible. Since glycosylated proteins and their resulting peptides are most often highly heterogeneous, high quality liquid chromatography is critical for glycopeptide analysis.

In this study, two reasonably well-characterized glycoproteins, bovine α -acid glycoprotein and human α -acid glycoprotein were analyzed using nano-LC MS/MS with ETD. Liquid chromatography separation conditions were systematically optimized for glycopeptide analysis with mass spectrometry using different stationary phase columns. Two reversed phase columns, C18 and C18, and a graphic carbon column were evaluated.

Materials & Methods

Sample:

Reduced and alkylated enzymatic digests of bovine and human α -acid glycoproteins. Glycoproteins were purchased from Sigma.

LC/MS:

HPLC System: SurveyorTM MS pump with a flow splitter
Column: AgilentTM ZORBAX[®] 300SB C8 column (75 μ m x 5 cm)
Microvial: C18 column (150 μ m x 150 cm)
Thermo Scientific Hypercarb column (75 μ m x 5 cm)
Mobile Phase: A: water, 0.1% formic acid; B: Acetonitrile, 0.1% formic acid
Gradient: 5-50% B in 30 minutes
Mass Spectrometer: Thermo Scientific LTQ XL linear ion trap mass spectrometer equipped with ETD and nano-ESI source

Spray Voltage: 2 kV
Capillary Temp: 160 °C
Capillary Voltage: 25 V
Tube Length: 125 V
MS1 Target: 1e4
Mass range: 50-2000 m/z or 100-4000 m/z
Ion Reagent: Fluorobenzene
Ion Reagent Isolation: On
Ion Target: 2e5
Max Ion Injection Time: 50 ms
ETD Reaction Time: 100 ms
Mass Spectrometer: Thermo Scientific LTQ Orbitrap XLTM
Mass range: 600-2000 m/z, resolution 100000
Full MS Scan: 5e6
Survey Scan: Source CID at 65 V for m/z 200

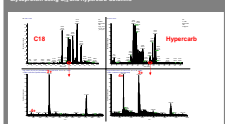
Data Processing:

Data were processed using BioWorksTM 3.3.1 with SEQUEST[®]. The Xtract program (Thermo Scientific) was used for deconvolution of multiply charged precursors. The GlycoMod tool from the Swiss-Prot website was used to assign possible oligosaccharide structures and compositions.

Results

Figure 1 shows the results of the LC/MS analysis of bovine α -acid glycoprotein on C₁₈ and Hypercarb columns. The top panels are the base peak chromatograms and the bottom panels are the bi-antennary glycopeptide 91-99 MS/MS profiles. One pmol of protein digest was injected into the C₁₈ column versus 500 fmol on the Hypercarb column resulting in a C₁₈ base peak intensity three times higher than the one from the Hypercarb column. However, as shown in the bottom panel of Figure 1, chromatography on the Hypercarb column promotes additional higher charge state precursor ions than C₁₈ which is instrumental in obtaining high quality ETD MS/MS. Overall intensity of glycopeptide precursor ions with the Hypercarb column is the same as, or higher than with the C₁₈ column.

FIGURE 1. Comparison of Base Peak Chromatograms of Bovine α -acid Glycoprotein using C₁₈ and Hypercarb Columns.



In addition, as shown in Figure 2, chromatography on the Hypercarb column promotes metal adduct formation. Figure 2 shows a high resolution deconvoluted spectrum of 91-99 bovine glycopeptides acquired on an LTQ Orbitrap XLTM. At least one metal adduct was observed for each glycopeptide. As demonstrated in Figure 1, formation of metal adducts is likely responsible for producing high charge state ions: the precursor of the 4+ charge state (567.3) was the dominant peak for glycopeptide 5 while the 3+ charge state (1134.6) was the dominant peak for glycopeptide 2. These results can be explained by partial neutralization of the negative charge of the sialic acid by metal cations⁽⁶⁾ and because the metal stabilizes the labile glycans and promotes formation of a higher charge state precursor⁽⁶⁾.

FIGURE 2. Deconvoluted Full MS Spectrum of Bovine α -acid Bi-antennary Glycopeptide *CVY*CSFQK.

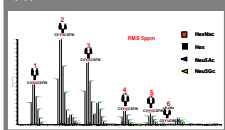


FIGURE 3. ETD spectrum of CVY*CSFQK (charge 4+, potassium adduct ion, m/z 663).

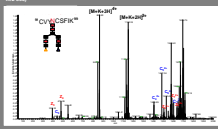
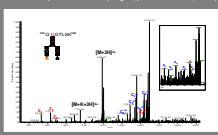


FIGURE 4. ETD spectrum of Q*GTSLK (charge 4+, potassium adduct ion, m/z 752).

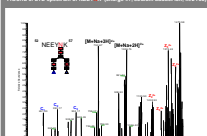


Figures 3 and 4 show ETD spectra of a potassium adduct of bovine α -acid glycoprotein peptides 91-99 (4+ charge) and peptides 103-109 (4+ charge) are detected as 4 on Figure 2. Glycosylation sites and peptide sequences were identified based on an almost complete series of c/z ions. No significant loss of carbohydrate was detected and, as expected, all observed glycan-containing fragments retained the metal ion. The overall difference in the performance of reversed phase chromatography and porous graphite carbon columns can be explained by higher trapping efficiency of the graphite column for hydrophilic peptides. Bovine α -acid glycoprotein contains five N-glycosylation sites with complex-type glycan structures⁽¹⁾.

TABLE 1. Bovine α -acid glycopeptides detected by nano-LC/MS/MS

Peptide/Type of LC column	Graphite	C ₁₈	C ₁
103*GVGSLK ⁽¹⁾	+	+	+
103*GVGSLK ⁽²⁾	+	+	+
103*GVGSLK ⁽³⁾	+	+	+
103*GVGSLK ⁽⁴⁾	+	+	+
103*GVGSLK ⁽⁵⁾	+	+	+
103*GVGSLK ⁽⁶⁾	+	+	+
103*GVGSLK ⁽⁷⁾	+	+	+
103*GVGSLK ⁽⁸⁾	+	+	+
103*GVGSLK ⁽⁹⁾	+	+	+
103*GVGSLK ⁽¹⁰⁾	+	+	+
103*GVGSLK ⁽¹¹⁾	+	+	+
103*GVGSLK ⁽¹²⁾	+	+	+
103*GVGSLK ⁽¹³⁾	+	+	+
103*GVGSLK ⁽¹⁴⁾	+	+	+
103*GVGSLK ⁽¹⁵⁾	+	+	+
103*GVGSLK ⁽¹⁶⁾	+	+	+
103*GVGSLK ⁽¹⁷⁾	+	+	+
103*GVGSLK ⁽¹⁸⁾	+	+	+
103*GVGSLK ⁽¹⁹⁾	+	+	+
103*GVGSLK ⁽²⁰⁾	+	+	+
103*GVGSLK ⁽²¹⁾	+	+	+
103*GVGSLK ⁽²²⁾	+	+	+
103*GVGSLK ⁽²³⁾	+	+	+
103*GVGSLK ⁽²⁴⁾	+	+	+
103*GVGSLK ⁽²⁵⁾	+	+	+
103*GVGSLK ⁽²⁶⁾	+	+	+
103*GVGSLK ⁽²⁷⁾	+	+	+
103*GVGSLK ⁽²⁸⁾	+	+	+
103*GVGSLK ⁽²⁹⁾	+	+	+
103*GVGSLK ⁽³⁰⁾	+	+	+
103*GVGSLK ⁽³¹⁾	+	+	+
103*GVGSLK ⁽³²⁾	+	+	+
103*GVGSLK ⁽³³⁾	+	+	+
103*GVGSLK ⁽³⁴⁾	+	+	+
103*GVGSLK ⁽³⁵⁾	+	+	+
103*GVGSLK ⁽³⁶⁾	+	+	+
103*GVGSLK ⁽³⁷⁾	+	+	+
103*GVGSLK ⁽³⁸⁾	+	+	+
103*GVGSLK ⁽³⁹⁾	+	+	+
103*GVGSLK ⁽⁴⁰⁾	+	+	+
103*GVGSLK ⁽⁴¹⁾	+	+	+
103*GVGSLK ⁽⁴²⁾	+	+	+
103*GVGSLK ⁽⁴³⁾	+	+	+
103*GVGSLK ⁽⁴⁴⁾	+	+	+
103*GVGSLK ⁽⁴⁵⁾	+	+	+
103*GVGSLK ⁽⁴⁶⁾	+	+	+
103*GVGSLK ⁽⁴⁷⁾	+	+	+
103*GVGSLK ⁽⁴⁸⁾	+	+	+
103*GVGSLK ⁽⁴⁹⁾	+	+	+
103*GVGSLK ⁽⁵⁰⁾	+	+	+
103*GVGSLK ⁽⁵¹⁾	+	+	+
103*GVGSLK ⁽⁵²⁾	+	+	+
103*GVGSLK ⁽⁵³⁾	+	+	+
103*GVGSLK ⁽⁵⁴⁾	+	+	+
103*GVGSLK ⁽⁵⁵⁾	+	+	+
103*GVGSLK ⁽⁵⁶⁾	+	+	+
103*GVGSLK ⁽⁵⁷⁾	+	+	+
103*GVGSLK ⁽⁵⁸⁾	+	+	+
103*GVGSLK ⁽⁵⁹⁾	+	+	+
103*GVGSLK ⁽⁶⁰⁾	+	+	+
103*GVGSLK ⁽⁶¹⁾	+	+	+
103*GVGSLK ⁽⁶²⁾	+	+	+
103*GVGSLK ⁽⁶³⁾	+	+	+
103*GVGSLK ⁽⁶⁴⁾	+	+	+
103*GVGSLK ⁽⁶⁵⁾	+	+	+
103*GVGSLK ⁽⁶⁶⁾	+	+	+
103*GVGSLK ⁽⁶⁷⁾	+	+	+
103*GVGSLK ⁽⁶⁸⁾	+	+	+
103*GVGSLK ⁽⁶⁹⁾	+	+	+
103*GVGSLK ⁽⁷⁰⁾	+	+	+
103*GVGSLK ⁽⁷¹⁾	+	+	+
103*GVGSLK ⁽⁷²⁾	+	+	+
103*GVGSLK ⁽⁷³⁾	+	+	+
103*GVGSLK ⁽⁷⁴⁾	+	+	+
103*GVGSLK ⁽⁷⁵⁾	+	+	+
103*GVGSLK ⁽⁷⁶⁾	+	+	+
103*GVGSLK ⁽⁷⁷⁾	+	+	+
103*GVGSLK ⁽⁷⁸⁾	+	+	+
103*GVGSLK ⁽⁷⁹⁾	+	+	+
103*GVGSLK ⁽⁸⁰⁾	+	+	+
103*GVGSLK ⁽⁸¹⁾	+	+	+
103*GVGSLK ⁽⁸²⁾	+	+	+
103*GVGSLK ⁽⁸³⁾	+	+	+
103*GVGSLK ⁽⁸⁴⁾	+	+	+
103*GVGSLK ⁽⁸⁵⁾	+	+	+
103*GVGSLK ⁽⁸⁶⁾	+	+	+
103*GVGSLK ⁽⁸⁷⁾	+	+	+
103*GVGSLK ⁽⁸⁸⁾	+	+	+
103*GVGSLK ⁽⁸⁹⁾	+	+	+
103*GVGSLK ⁽⁹⁰⁾	+	+	+
103*GVGSLK ⁽⁹¹⁾	+	+	+
103*GVGSLK ⁽⁹²⁾	+	+	+
103*GVGSLK ⁽⁹³⁾	+	+	+
103*GVGSLK ⁽⁹⁴⁾	+	+	+
103*GVGSLK ⁽⁹⁵⁾	+	+	+
103*GVGSLK ⁽⁹⁶⁾	+	+	+
103*GVGSLK ⁽⁹⁷⁾	+	+	+
103*GVGSLK ⁽⁹⁸⁾	+	+	+
103*GVGSLK ⁽⁹⁹⁾	+	+	+
103*GVGSLK ⁽¹⁰⁰⁾	+	+	+

FIGURE 5. ETD spectrum of NEEY*AK (charge 4+, sodium adduct ion, m/z 755).



Four out of five glycopeptides were detected and identified using a Hypercarb column compared to two on C₁₈ and three on C₁ columns (Table 1) without any enrichment. Only the largest and most hydrophilic peptide was not detected.

Similar results were obtained for human α -acid glycoprotein which also contains five N-glycosylation sites⁽¹⁾. All its glycopeptides can be detected after C₁₈ or C₁ chromatography but mostly as charge 2+ or 3+ species and as a result, MS² ETD spectra did not generate enough information for unambiguous identification. On the other hand, multiple alkali metal on acid glycopeptides were observed after separation on the Hypercarb column at the higher charge states. Figure 5 shows an example of an ETD spectrum of a 4+ charge bi-antennary peptide 52-57 (sodium adduct), which was successfully identified by BioWorks 3.3.1. This peptide contains two asparagine residues but the site of glycosylation was still easily identified as Asn 56.

Conclusions

- A graphic carbon column (Hypercarb, Thermo Scientific) demonstrated excellent capabilities for glycopeptide analysis especially for short hydrophilic peptides containing bi- or tri-antennary glycan chains without any enrichment.
- Formation of metal adducts on Hypercarb columns promote higher charged species and as a result, promotes ETD fragmentation of glycopeptides.
- Optimization is required for successful glycoprotein nano-LC-ETD MS analysis.
- As a gentle fragmentation technique, ETD preserves labile glycans, facilitating the identification of both the peptide of interest and its site of modification.

References

1. S. I. Smolke, V. C. Chen, O. Krohlin, and H. Pernas, *Anal. Chem.*, **2006**, *78*, 6556-63.
2. S. M. Peterson and J. J. Mulholland, *J. Am. Soc. Mass Spectrom.*, **2006**, *17*(2), 168-79.
3. K. A. Neman, R. Amunugama, and S. A. Mulcahy, *J. Phys. Chem. A*, **2005**, *109*(16), 3638-16.
4. K. F. Medvedev, S. Guan, D. A. Malby, and A. L. Burlingame, *J. Am. Soc. Mass Spectrom.*, **2007**, *18*(6), 1617-24.
5. M. J. Treuhart, C. E. Costello, and H. B. Nalati, *Biochem. J.*, **1992**, *283*, 105-112.

Agilent and ZORBAX are registered trademarks of Agilent Technologies, Inc. SEQUEST is a registered trademark of the University of Washington. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

Thermo
SCIENTIFIC