POSTER NOTE

Complete Characterization of a Cysteinelinked Antibody-Drug Conjugate Performed on a Hybrid Quadrupole-Orbitrap Mass Spectrometer with High Mass Range

Aaron O. Bailey¹, Eugen Damoc², Stephane Houel¹, Kai Scheffler³, Jonathan L. Josephs¹; Thermo Fisher Scientific, ¹San Jose, CA, USA; ²Bremen, Germany; ³Dreieich, Germany

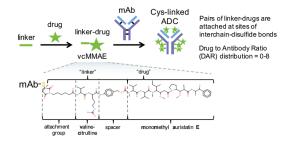
ABSTRACT

We have modified the instrument control software of a benchtop quadrupole-Orbitrap mass spectrometer to add native MS capability. In this study we demonstrate complete characterization of Brentuximab vedotin, a cysteine-linked ADC, which requires native MS conditions for intact analysis. We demonstrate preservation of non-covalent bonding of antibody subunits during electrospray ionization. HMR mode can be turned off for peptide mapping. We use trypsin peptide mapping approach with HCD fragmentation to achieve 99% coverage of the *Brentuximab vedotin* sequence using a single LC-MS analysis of a 90 min reverse phase gradient. Finally, we demonstrate that signature ions specific for HCD fragmentation of Brentuximab vedotin can be utilized to increase MS/MS assignment confidence

INTRODUCTION

The complexity of modern therapeutic proteins, such as antibody-drug conjugates (ADCs), present a great analytical challenge which requires high resolution chromatography combined with high resolution mass spectrometry. Complementary MS approaches such as peptide mapping and intact mass analysis are needed for complete characterization of therapeutic proteins. Cysteine-linked ADCs present a unique challenge for characterization as proper intact analysis requires native MS conditions to preserve structurally-critical non-covalent binding between antibody chains. We have modified commercially-available Thermo Scientific[™]Q Exactive[™] Plus and Q Exactive[™] HF Orbitrap[™] mass spectrometers to perform native LC-MS experiments. In the present study, we demonstrate this capability with intact analysis of *Brentuximab vedotin*, a cysteine-linked ADC (Figure 1). Additionally, we have performed denaturing LC-MS and peptide mapping on these same instruments to generate complementary datasets for complete characterization.

Figure 1. Schematic for Constructing Cysteine-Linked ADC Brentuximab vedotin is a cysteine-linked ADC which is constructed by modifying an antibody with vcMMAE, a preformed linker-drug comprised of a valine-cirtuline-based linker and a monomethyl auristatin E toxic drug. Saturaturated (8 drugs) cys-linked ADCs are held intact with only non-covalent binding.



MATERIALS AND METHODS

Brentuximab vedotin was prepared for peptide mapping (reduction, alkylation, and trypsin digestion) or intact analysis (no treatment). For denaturing LC-MS intact analysis 1 µg of protein samples were separated using a 10 min gradient of 10-90% ACN in H₂O and 0.1% formic acid (Thermo ScientifeTM MAbPacTM column RP; flow rate 250 µL/min). For native LC-MS intact analysis 10 µg of sample was desalted online using size exclusion chromatography (WatersTM BEH SEC 4.6x150mm; 50 mM NH₄OAc isocratic elution, flow rate 300 µL/min) and directly presented to the mass spectrometer via electrospray ionization. Peptide mapping was performed using 2.5 µg of sample separated using a 90 min gradient of 2-90% ACN in H₂O and 0.1% formic acid (Acclaim RSLC 120 C18: flow rate 250 uL/min). Commercially-available Orbitrap mass spectrometers (Q Exactive HF and Q E20 C19, how rate 250 µL7min), commercially-available Orbitral mass spectrometers (0L Exactive PL) with PL and 0L Exactive PL) were modified to include High Mass Range (HMR) mode to allow improved high mass transmission and scanning up to m/2 8000. Native intact and denaturing MS data were acquired in HMR mode at setting of R=15k or 17.5k and deconvolved using the ReSpectTM algorithm and Silding Window integration in Thermo ScientificTM BioPharma FinderTM 1.0 SP1 software. Deconvolution species were identified automatically using the publicly available FASTA sequence for Brentwirmab vodotin, a mass tolerance of 50 ppm, and a statie modification of GinPyrvo-Cill for the heavy chain. Peptide mapping data were acquired by data dependent selection with R=60k or 70k for Full MS and R=15k or 17.5k for MS/MS. Peptide mapping data were searched prior the Mach Apping reaching the prior Biopers Endocomposition and a static termore the mach apping data were searched prior the Mach Apping the prior the prior the prior to the prior to fee for point and the prior to the prior to the prior to the prior to the perior to the prior to the perior to the termoson of fee point and the perior to the termoson of the prior to the perior to the perior to the termoson of fee point and the perior to the termoson of the termoson of the period to the termoson of the perior to the termoson of the termoson of the termoson of the perior to the termoson of the t using the MassAnalyzer algorithm in BioPharma Finder software with a tolerance of 5 ppm

Figure 2. LC-MS Instrumentation for Complete ADC Characterization All experiments were performed using a Thermo Scientific™ Vanquish™ UHPLC connected to a Exactive HF or Q Exactive Plus instrument with High Mass Range (HVR) mode.



RESULTS

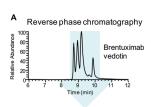
DENATURING LC-MS, CYSTEINE-LINKED ADC

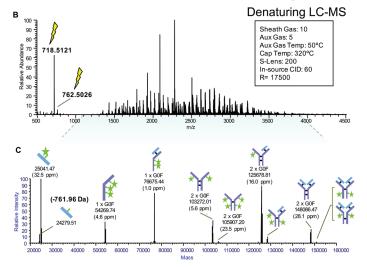
Intact protein LC/MS analysis conventionally involves using mobile phases which are comprised of organic and acidic/basic pH, often suited specifically for reverse phase chromatography. This strategy can be useful for achieving high resolution protein separations. Conditions such as these, however, are not compatible for performing intact analysis on certain classes of compounds which require preservation of non-covalent bonds to maintain structural integrity, such as cysteine-linked ADCs. We demonstrate this phenomenon using the cysteine-linked ADC Brentuximab vendotin. Denaturing (reverse phase) LC/MS analysis of *Brentuximab vedotin* results in detection of roughly six unraveled forms (Figure 3A-C). We observed a previously-reported¹ collisionally-induced m/z 718 fragment of the fragile vcMMAE linker-drug (Figure 3B). Upon deconvolution we also observed a mass corresponding to light chain with addition of one vcMMAE and a loss of approximately 762 Da.



Figure 3. Denaturing LC-MS analysis

(A) Unmodified sample (Jug) was analyzed by reverse phase chromatography coupled to a Q Exactive Plus Orbitrap MS operating in HMR mode and produced several peaks. (B) The resulting averaged MS spectrum is a complex mixture of charge state envelopes as well as a previously described volMAE specific reporter fragment ion at miz 7 18. (C) Data analysis with ReSpect deconvolution and Sliding Window integration show roughly six covalently-structured forms of unraveled cysteine-linked ADC. We detect a protein species which corresponds to a light chain with addition of one linker drug and a loss of 762 Da, which is also present in the raw spectrum





NATIVE INTACT LC-MS, CYSTEINE-LINKED ADC

Native MS intact protein analysis allows direct observation of molecules which rely on noncovalent interactions to preserve critical structural features, such as maintaining interchain associations which hold together cysteine-linked ADCs. The use of 100% aqueous physiological pH buffers in native MS analysis produces decreased charge states (increased m/z) and improves mass separation of heterogeneous mixtures. We performed native size exclusion LC-MS and observed 5 distinct species corresponding to intact Brentuximab vedotin with 0, 2, 4, 6, or 8 vcMMAE linker-drugs (Figure 4). We measured an average drug-to-antibody ratio of 4.07, which is consistent with a previously published studies reporting 3.9-4.2 drugs per antibody²

RESULTS

DENATURING LC-MS, CYSTEINE-LINKED ADC

Intact protein LC/MS analysis conventionally involves using mobile phases which are comprised of organic and acidic/basic pH, often suited specifically for reverse phase chromatography. This strategy can be useful for achieving high resolution protein separations. Conditions such as these, however, are not compatible for performing intact analysis on certain classes of compounds which require preservation of non-covalent bonds to maintain structural integrity, such as cysteine-linked ADCs. We demonstrate this phenomenon using the cysteine-linked ADC Brentuximab vendotin. Denaturing (reverse phase) LC/MS analysis of Brentuximab vedotin results in detection of roughly six unraveled forms (Figure 3A-C). We observed a previously-reported¹ collisionally-induced m/z 718 fragment of the fragile vcMMAE linker-drug (Figure 3B). Upon deconvolution we also observed a mass corresponding to light chain with addition of one vcMMAE and a loss of approximately 762 Da.

Figure 4. Native intact LC-MS analysis Α Size exclusion chromatography (A) Unmodified sample (10 ug) was analyzed using size exclusion chromatography coupled to a Q Exactive Plus Orbitrap MS operating in HMR mode. Buffer exchange occurs online as ADC forms elute as a single peak, Brentuximab 50 mM vedotin NH₄OAc followed by a second peak corresponding separated buffer salts. (B) Averaging 2 min chromatographic time produces a native intact MS spectrum which includes all DAR forms (DAR 0-8). (C) ReSpect deconvolution and 60isocratic elution 40buffe salts 20 Silding Window integration can accommodate peak tailing to report quantitatively accurate abundances for the mixture of DAR forms which have diverse elution profiles. 07 4 Time (mi A pattern of lower abundance species were detected corresponding to a low abundance loss of 762 Da from each glycoform at each DAR value (green arrows). **(D)** Native LC-MS Based on the individual deconvolved abundances of the Sheath Gas: 40 Aux Gas: 5 Aux Gas Temp: 125°C Cap Temp: 275°C G0F/G0F glycoform, we calculated an average DAR value of 4.07, which is consistent with previous reports². в 100 S-Lens: 200 90 In-source CID: 100 R= 17500 80 Abundance 00 20 Relative / 20 10 5400 7400 DAR4 С 脊靴 DAR6 DAR2 2 x G0F 153355.40 100 2 x G0F DAR8 DAR0 80 衸 Ŷ 60 (-762.25 Da 40 2 x G0F 158632.20 2 x G0 20 íц (1 0 150000 152000 151000 154000 155000 157000 158000 Mass D GOF/GOF Mass Accu (ppm) Relative Abundance Average DARC 11.7 6.77 Drug-to-Antibody DAR2 23.1 69.23 Ratio (DAR) DAR4 22.4 100.00 4.07 DAR6 40.5 69.75 DAR8 17.6 10.61

PEPTIDE MAPPING, CYSTEINE-LINKED ADC

A fundamental component of biotherapeutic protein characterization is peptide mapping. Whereas intact mass analysis aims to detect the abundances and distributions of mass deviation combinations, peptide mapping allows highly sensitive analysis of site-specific sequence features. The vcMMAE linker-drug on Brentuximab vedotin poses particular challenges when attempting to identify drug conjugation sites. We prepared a sample for peptide mapping using reduction and alkylation to block non-drug-conjugated cysteines, followed by trypsin digestion. In one 90 min LC-MS gradient we were able to achieve 99% sequence coverage for both light and heavy chains and detect peptides spanning all four drug conjugation sites. HCD fragmentation allowed detection of a peptide in the hinge region of the heavy chain that is differentially modified with 0-2 vcMMAE drugs. As a result efficient elution requires sustained delivery of high organic mobile phase.

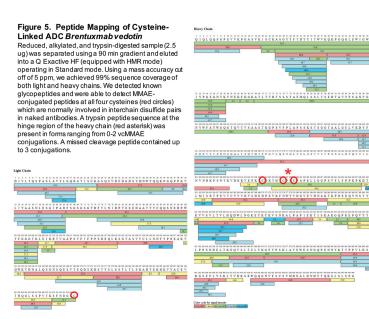


Figure 6. Hinge Region Peptide of Cysteine-linked ADC is Site of Multiple Conjugations (A) Our data analysis in BioPhama Finder software resulted in detection of peptides which covered the hinge region of the heavy chain (red asterisk). A faithfully-trypsin-cleaved THT-KPK peptide was detected with 0-2 vcMMAE conjugations at cysteines (red circles) normally involved in interchain disulfide pairs. (B) Addition of vcMMAE to peptides dramatically increases hyphobicity which results in poor elution and increased retention time. MS/MS analysis of the (C) 1 linker-drug (both positional isomers) and (D) two linker-drug forms in BioPharma Finder al clear sequencing of y-ions in the hinge peptides, and thus facilitated automatic detection. na Finder allowed

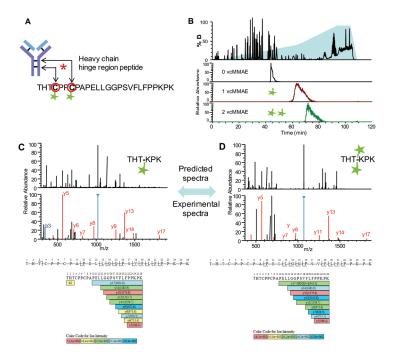
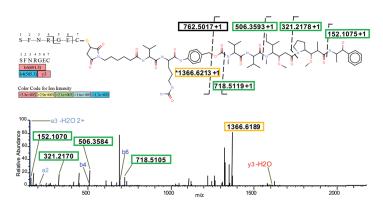


Figure 7. HCD Signature Fragment lons for vcMMAE Linker-Drug The light chain C-terminal peptide SFN-GEC is a conjugation site for vcMMAE. This modified peptide was automatically identified by BioPhama Finder (left side top panel). Further manual inspection produced additional fragment assignments for vcMMAE signature ions (right side top parel). Theoretical masses (top panel) were calculated manually and matched to experimental masses (bottom panel) within 5ppm (green boxes). A cleavage site for the loss of 762 Da is shown (black box; theoretical monoisotopic mass = 762.5017). We observed a high abundance ion at m/z 1366.6189 (orange box, asterisk) which corresponds to the peptide-retaining fragment pair of a 762 Da loss with an additional loss of 2 protons, presumably due to formation of a seven-membered aromatic ring



CONCLUSIONS

•We have modified the control software in Q Exactive Plus and Q Exactive HF mass spectrometers to add native MS capability.

•Native LC/MS intact analysis of Brentuximab vedotin resulted in detection of intact ADC forms DAR0-8. ReSpect deconvolution and Sliding Window integration showed an average DAR of 4.07, consistent with previous studies.

•Acquisition of MS/MS spectra with HCD fragmentation on Q Exactive Plus and Q Exactive HF Orbitrap mass spectrometers followed by data analysis with BioPharma Finder resulted in 99% sequence coverage from a single 90 min gracient using 5 ppm mass tolerance.

•Addition of vcMMAE linker drug dramatically increases peptide hydrophobicity and retention time.

•Signature HCD fragment ions of linker-drug may allow additional means for identifying drugconjugated peptides.

REFERENCES

1. Janin-Bussat MC, Dillenbourg M, Corvaia N, Beck A, Klinguer-Hamour C. Characterization of antibody drug conjugate positional isomers at cysteine residues by peptide mapping LC-MS analysis. J Chromatogr B Analyt Technol Biomed Life Sci. 2015 Feb 15;981-982:9-13.

2. Debaene F, Boeuf A, Wagner-Rousset E, Colas O, Ayoub D, Corvaïa N, Van Dorsselaer A, Beck A, Cianférani S. Innovative native MS methodologies for antibody drug conjugate characterization: High resolution native MS and IM-MS for average DAR and DAR distribution assessment. Anal Chem. 2014 Nov 4;86(21):10674-83.

For Research Use Only. Not to be used in diagnostic procedures.

Find out more at thermofisher.com

© 2016 Thermo Fisher Scientific Inc. All rights reserved. All rights reserved. Waters is a trademark of Waters Corporation. ReSpect™ is atrademark of Positive Probability Ltd. All otherAll trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. PN64858-EN 08/16S

