Poster ID: P-II-0565

Cross-linking Mass Spectrometry

Thermo Fisher SCIENTIFIC

Optimized XL-MS workflows for heterobifunctional crosslinkers SDA and DizSEC

Yi He¹, Erum Raja², Leigh Foster², Anthony Ciancone³, Anneliese Faustino⁴, Piyoosh Sharma⁴, Stephen D Fried^{4,5}, Francis O'Reilly³, Ryan Bomgarden², Rosa Viner¹

¹Thermo Fisher Scientific, San Jose, CA, USA; ²Thermo Fisher Scientific, Rockford, Illinois, USA; ³Center for Structural Biology, National Cancer Institute, NIH, Frederick, MD, USA; ⁴Department of Chemistry, Johns Hopkins University, Baltimore, MD, USA; ⁵ Department of Biophysics, Johns Hopkins University, Baltimore, MD, USA

Abstract

Purpose: To optimize XL-MS workflows for heterobifunctional photoactivatable -crosslinkers SDA and DizSEC.

Methods: Heterobifunctional photoactivatable -crosslinkers SDA (succinimidyl 4,4'-azipentanoate), sulfo-SDA (sulfosuccinimidyl 4,4'-azipentanoate) and DizSEC (2,5-Dioxopyrrolidin-1-yl (2-(3methyl-3H-diazirin-3-yl)ethyl)carbamate) were used to crosslink standard proteins. Separation was achieved using Thermo Scientific[™] Vanquish[™] Neo LC system with a 60 min gradient using Thermo Scientific[™] EASY-Spray[™] PepMap[™] Neo column. Following separation, the peptides were detected on Thermo Scientific[™] Orbitrap Ascend[™] or Exploris 480 [™] mass spectrometers, and data were analyzed using XlinkX node in Thermo Scientific[™] Proteome Discoverer[™] v3.2 software and xiSEARCH/xiFDR. Identified XLs were visualized using the XMAS plug-in for ChimeraX.

Results

 Table 1. Orbitrap Exploris and Ascend MS acquisition parameter
settings for SDA and DizSEC MS2 methods.

Settings	Exploris	Ascend
MS1 scan	Orbitrap	Orbitrap
Resolution	120K	240K
Normalized AGC Target (%)	300%	150%

Results

Methods for DizSEC crosslinked samples

DisSEC is a novel MS-cleavable heterobifunctional crosslinker that connects Lys with any amino acid in proteins and forms Ureas. Ureas are doubly fissile under HCD/CID (across each N-carbonyl bond) and generate two distinct sets of fragment ions (as shown in Figure 5). Collisional energy is crucial for cleavable crosslinkers. To improve accuracy in identification, we developed a MS2-MS3 method for DizSEC crosslinked peptides to compared with the MS2 method. Optimized parameters are listed below.

Results

Figure 7. (A) HSA DizSEC crosslinks and looplinks identified using XlinkX MS2 search in different MS methods. (B) Venn diagram showing the overlap of identified HSA crosslinks on OT Ascend.



Results: With optimized MS2 and MS2-MS3 methods, SDA and DizSEC crosslinked peptides can be identified with high confidence.

Introduction

Cross-linking mass spectrometry (XL-MS) has grown dramatically as a key workflow for elucidating protein higher-order structure and mapping protein-protein or protein-nucleic acid interaction networks on a proteome-wide scale. Photoactivatable, heterobifunctional crosslinkers like SDA can link lysine residues with any residue, providing detailed information for protein structures, but also pose challenges due to lack of MS-cleavability and increased search space during data analysis. Recently, a new cleavable crosslinker, DizSEC, was developed to overcome this issue. DizSEC links the same residues as SDA, but DizSEC crosslinked peptides can generate unique peptide pairs in MS/MS to facilitate data analysis.

In this work, we optimized XL-MS workflows for SDA and DizSEC crosslinked samples and developed a new MS2-MS3 method to improve DizSEC XL identifications.

Materials and methods

Sample preparation

Heterobifunctional photoactivatable -crosslinkers SDA (succinimidyl 4,4'-azipentanoate), sulfo-SDA (sulfosuccinimidyl 4,4'-azipentanoate) and DizSEC (2,5-Dioxopyrrolidin-1-yl (2-(3methyl-3H-diazirin-3-yl)ethyl)carbamate) were used to crosslink standard proteins.

Max. injection time (ms)	20	100	
Charge state	3-8	3-8	
MS mass range, m/z	375-1600	380-1400	
RF lens	50	60	
MS2 scan	Orbitrap	Orbitrap	
Resolution	45k	60K	
Normalized AGC Target (%)	200%	500%	
Max. injection time (ms)	105	123	
Charge states	3-8	3-6 (prioritizing +4 to +6)	
First mass (m/z)	150	150	
Intensity threshold	1e⁴	2.5e⁴	
Isolation width	1.6	1.4	
NCE (%)	SCE 20,25, 30	SCE 20, 25, 30	
Top speed		5s	

Figure 2. Orbitrap Ascend MS acquisition methods for SDA and DizSEC crosslinked peptides.



Figure 5. Structure and MS fragmentation of DizSEC crosslinked peptides.



Table 3. Orbitrap Ascend MS acquisition parameter settings for **DizSEC MS2-MS3 method.**

Settings	MS2	MS3	
Detector	Orbitrap	lon trap	
Resolution/Scan Rate	30K	Rapid	
Normalized AGC Target (%)	100%	200%	
Max. injection time (ms)	59	80	
Charge states	3-8	2-6	
Intensity threshold	1e⁴	NA	
Isolation width	1.6	2.5	
NCE (%)	CID 25	HCD 30	
Top speed/scans	5s	4 scans	
Delta Mass (Da)	25.9942		

Conclusions

- Three different SDA crosslinkers provided a similar number of crosslinked peptides for monomeric proteins.
- SDA and sulfo-SDA outperformed DizSEC for multimeric proteins probably due to the length of the linker.

Liquid chromatography and mass spectrometry

Samples were separated by reverse phase-HPLC on a Vanquish Neo system using Thermo Scientific[™] EASY-Spray[™] PepMap[™] RSLC C18 column. Peptides were eluted over a 60 min 6-50% gradient (A: water, 0.1% formic acid; B: 80% acetonitrile, 0.1% formic acid) at 300 nL/min flow rate. The crosslinked samples were analyzed on Thermo Scientific[™] Orbitrap[™] Ascend[™] or Thermo Scientific[™] Exploris[™] 480 mass spectrometers in DDA modes. Specific LC and MS settings are shown in Tables 1, Figure 2, and Table 3.

Data analysis

The acquired spectra were analyzed using Proteome Discoverer 3.2 software: the XlinkX node with the noncleavable, MS2 or MS2-MS3 search option for crosslinked peptides and SEQUEST™ HT search engine for unmodified, looplinks and monolinks peptides. Detailed settings for DizSEC is shown in Figure 5. Data were also analyzed using xiSEARCH/xiFDR. Carbamidomethylation (+57.021 Da) was used as a static modification for cysteine. Different crosslinked mass modifications for lysine were used as variable modifications for lysine or Nterminus in addition to methionine oxidation (+15.996 Da). Data were searched against a protein database containing 10 proteins including human serum albumin (P02768) or yeast enolase (P00924). The false discovery rate (FDR) was set to 1% at CSM and cross-link levels.

Results

Comparison of three crosslinkers

We compared the performance of three heterobifunctional photoactivatable -crosslinkers SDA, sulfo-SDA and DizSEC with standard proteins. The structures of the crosslinkers are shown in Figure 1.

Table 2. Summary of XlinkX non-cleavable search results from three crosslinkers on OT Exploris 480.

		Looplinks	CSMs	Crosslinks
HSA	SDA	148	200	113
	Sulfo-SDA	128	171	99
	DizSEC	117	164	108
Enolase Dimer	SDA	75	107	53
	Sulfo-SDA	75	93	60
	DizSEC	51	52	17

Figure 3. Venn diagram showing the overlap of identified enolase crosslinks on OT Exploris 480.



Figure 4. Intra (yellow) and inter (red) crosslinks identified in the enolase dimer structure with three crosslinkers and their distance distributions.

Figure 6. Examples of HSA DizSEC crosslinkers identified by (A) MS2 method and (B) MS2- MS3 method.

2024021_HSA_DizSEC_SCE202530_nofaims.raw #18853_RT: 45.5608 min FTMS: 764.1809@hcd25.00, z=+5, Mono m/z=763.58026 Da, MH+=3813.87220 Da, Match Tol.=0.5 Da Α α: RHPYFYAPELLFFAKR 763.97943 YKAAFTECCQAADK 763.77991 5 300 764.18030 $\alpha b_6^{2+}, \alpha b_6^{2+}, \alpha b_6^{2+}, \alpha b_6^{2+}$ 64.37860 432.71124 ab7+,ab7+,ab7+,ab7+ 1217.70386 935.45245 718.3889 100 -539.04340 1662.72583 1076.57861 1688.70288 2000 В a: SLHTLFGDKLCTVATLR LVNEVTEFAK Crosslink Peptides 20240430_HSA_DizSEC_MS2MS3_26Da_4scans.raw #30739 RT: 51.8337 min FTMS, 795.4291@cid25.00, z=+4, Mono m/z=795.18054 Da, MH+=3177.70034 Da, Match Tol.=0.5 Da



- Increased numbers of SDA and DizSEC crosslinked peptides were identified using the instrument method prioritizing high charge states 4-6.
- An optimized MS2-MS3 method for DizSEC crosslinker was developed to reduce false positives in XL identifications.

References

- 1. Tanaka Y, Kohler JJ. Photoactivatable crosslinking sugars for capturing glycoprotein interactions. J Am Chem Soc. 2008 Mar 19;130(11):3278-9.
- 2. Mendes ML, Fischer L, Chen ZA, Barbon M, O'Reilly FJ, Giese SH, Bohlke-Schneider M, Belsom A, Dau T, Combe CW, Graham M, Eisele MR, Baumeister W, Speck C, Rappsilber J. An integrated workflow for crosslinking mass spectrometry. Mol Syst Biol. 2019 Sep;15(9):e8994.
- 3. Fischer L, Rappsilber J. Quirks of Error Estimation in Cross-Linking/Mass Spectrometry. Anal Chem. 2017 Apr 4;89(7):3829-3833.
- 4. Faustino AM, Sharma P, Manriquez-Sandoval E, Yadav D, Fried SD. Progress toward Proteome-Wide Photo-Cross-Linking to Enable Residue-Level Visualization of Protein Structures and Networks In Vivo. Anal Chem. 2023 Jul 18;95(28):10670-10685.
- 5. Lagerwaard IM, Albanese P, Jankevics A, Scheltema RA. Xlink Mapping and AnalySis (XMAS)—Smooth Integrative Modeling in ChimeraX. bioRxiv: 2022.04.21.489026.

Figure 1. Structure and reactions of three heterobifunctional photo-cross-linkers SDA, LCSDA and DizSEC.



Conflict of interest

YH, ER, LF, RB and RV are employees of Thermo Fisher Scientific.

Trademarks/licensing

© 2024 Thermo Fisher Scientific Inc. All rights reserved. SEQUEST is a trademark of the University of Washington. Aurora Ultimate is a trademark of IonOpticks Pty Ltd. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.

Science at a scan Scan the QR code on the right with your mobile device to download this and many more scientific posters.



Learn more at thermofisher.com/us/en/home/life-science/structural-biology/integrative-structural-biology.html