Thermo Fisher S C I E N T I F I C

Optimized XL-MS workflows for protein-protein and proteinnucleic acid interactions

Yi He¹, Erum Raja², Leigh Foster², Olexandr Dybkov³, Sergey Moshkovskiy³, Henning Urlaub^{3,4}, Ryan Bomgarden², Rosa Viner¹ ¹Thermo Fisher Scientific, San Jose, CA, USA; ²Thermo Fisher Scientific, Rockford, Illinois, USA; ³Bioanalytical Mass Spectrometry Group, Max Planck Institute for Multidisciplinary Sciences, 37077 Göttingen, Germany; ⁴Bioanalytics Group, Institute of Clinical Chemistry, University Medical Center Göttingen, 37075 Göttingen, Germany

Abstract

Purpose: To optimize XL-MS workflows for protein-protein and protein-nucleic acid interactions.

Methods: For protein-protein interaction, heterobifunctional photoactivatable crosslinker DizSEC (2,5-Dioxopyrrolidin-1-yl (2-(3-methyl-3H-diazirin-3-yl)ethyl)carbamate) were used to crosslink standard proteins. For protein-nucleic acid interaction, *E. coli* ribosomes were exposed to UV or nitrogen mustard and enriched using TiO₂. Separation was achieved using Thermo Scientific[™] Vanquish[™] Neo LC system with a 60 min gradient using EASY-Spray[™] PepMap[™] Neo column. Following separation, the peptides were detected on Thermo Scientific[™] Orbitrap[™] Ascend[™] or Astral[™] mass spectrometers, and data were analyzed using the XlinkX node or the NuXL node in Thermo Scientific[™] Proteome Discoverer[™].

Results

Methods for DizSEC crosslinked samples

Collisional energy is crucial for cleavable crosslinkers. We compared three MS methods: MS2 SCE, MS2 SCE prioritizing charge states +4 to +6, and a new MS2-MS3 method for DizSEC. Optimized parameters are listed below.

Figure 2. Orbitrap Ascend MS acquisition methods for DizSEC crosslinked peptides prioritizing charge states +4 to +6.



Figure 4. (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.**



Figure 7. Example of Annotated MS2 spectrum of a UV induced RNA-peptide crosslink from OT-Astral. The MS2 spectrum view in Proteome Discoverer of the

peptide ATLGEVGNAEHmLR crosslinked with U



Results: With optimized MS methods, DizSEC crosslinked peptides can be identified with high confidence. Moreover, we identified two to three times the number of unique RNA crosslinks on Orbitrap Astral MS.

Introduction

Crosslinking 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. UV light is one of the common ways to create crosslinks to investigate both types of interactions. However, challenges remain in sample preparation, instrument methods and data analysis. In this work, we optimized XL-MS workflows for photoactivatable crosslinkers SDA and DizSEC (cleavable) crosslinked samples and protein-RNA crosslinks. A new MS2-MS3 method was developed to improve DizSEC XL identifications for protein-protein interaction. Importantly, we further optimized protein-nucleic acid workflows on the Thermo Fisher Scientific[™] Orbitrap[™] Astral[™] mass spectrometer and achieved two-to-three times identifications for XL-MS analysis.

Materials and methods

Sample preparation

Heterobifunctional photoactivatable DizSEC (2,5-Dioxopyrrolidin-1-yl (2-(3-methyl-3H-diazirin-3-yl)ethyl)carbamate) was used to crosslink standard proteins. E. coli ribosomes were exposed to UV or chemical crosslinkers. Crosslinked samples were digested, purified using TiO_2 .

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 4-50% 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 Figure 2, Tables 1 and 2.

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	

Protein-nucleic acid interactions

We developed DDA methods for NuXL identification and quantitation on OT Astral. E. coli ribosomes were crosslinked with UV light (215 nm) or nitrogen mustard .Optimized parameters are listed below.

Table 2. Orbitrap Astral MS crosslinker specific LC-MS acquisition parameter settings.

Settings	OT-OT	OT-Astral	
LC gradient	4-50% B in 60 min	4-50% B in 60 min	



Conclusions

- Increased numbers of 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.
- NuXL workflow was benchmarked on the Orbitrap Astral[™] mass spectrometer with UV- or chemically induced *E. coli* ribosome crosslinked samples.
- Using the Astral analyzer's high sensitivity and fast scan rate (up to 200 Hz), we identified two to three times the number of RNA crosslinks and spectrum matches for protein-nucleic acid interactions.

Availability

The NuXL node is released for Proteome Discoverer 3.0 and 3.1. The binary installer, documentation, and example data are available at https://openms.de/applications/nuxl/.

References

- 1. 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.
- 2. Röst HL, Sachsenberg T, Aiche S, Bielow C, Weisser H, ... & Kohlbacher O. OpenMS: a flexible open-source software platform for mass spectrometry data analysis. Nat Methods.

Data analysis

Protein-protein crosslinking data 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. Data were also analyzed using xiSEARCH/xiFDR. Carbamidomethylation (+57.021 Da) was used as a static modification for cysteine and methionine oxidation (+15.996 Da) as variable modification. 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.

Protein-nucleic acid crosslinking data was analyzed using the NuXL node in PD 3.1. We calculate match-odds and subscores and employ semi-supervised score calibration using Percolator. Moreover, entrapment experiments have been employed using manually curated data ensuring proper false discovery rate (FDR) control.

Results

Protein-protein interactions

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 Ncarbonyl bond) and generate two distinct sets of fragment ions . The structure of the crosslinker is shown in Figure 1.

Figure 3. Examples of HSA DizSEC crosslinkers identified by (A) MS2 method and (B) MS2- MS3 method.





MS1 scan	Orbitrap	Orbitrap
Resolution	120K	180K
Normalized AGC Target (%)	300%	500%
Max. injection time (ms)	20	5
Charge state	2-6	2-6
MS mass range, m/z	400-1400	400-1400
Intensity threshold	1e⁴	5e ³
RF lens	50	50
MS2 scan	Orbitrap	Astral
Resolution	45K	Standard
Normalized AGC Target (%)	200%	100%
Max. injection time (ms)	105	20
Charge states	2-6	2-6
First mass (m/z)	150	150
Intensity threshold	1e⁴	5e ³
Isolation width	1.6	1.6
NCE (%)	30	30
Top speed	2s	1s

Figure 5. Comparison of unique identified RNA crosslinks and spectrum matches from NM duplicate samples in OTOT and OT Astral DDA runs.



2016 Aug 30;13(9):741-8.

- 3. Stützer, A., Welp, L. M., Raabe, M., Sachsenberg, T., Kappert, C., Wulf, A., ... & Urlaub, H. (2020). Analysis of protein-DNA interactions in chromatin by UV induced cross-linking and mass spectrometry. Nature Communications, 11(1), 1-12
- 4.Welp,L.M., Sachsenberg,T., Wulf, A., Chernev, A., Horokhovskyi, Y., Neumann, P., Pašen, M., Siraj, A., Raabe, M., Johannsson, S., et al. (2024). Chemical crosslinking extends and complements UV crosslinking in analysis of RNA/DNA nucleic acid-protein interaction sites by mass spectrometry. bioRxiv 2024.08.29.610268

Acknowledgements

This project has received funding from the European Union's Horizon 2020 research and innovation program under the Marie Skłodowska-Curie grant agreement Nº 956148.

Trademarks/licensing

© 2025 Thermo Fisher Scientific Inc. All rights reserved. SEQUEST is a trademark of the University of Washington. All other 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. **PO003796 0325**







8

Figure 1. Structure and reactions of MS-cleavable heterobifunctional photo-cross-linkers DizSEC.



Peptide A (CO) m/z = 979.5129	Peptide A (Pep1) m/z = 966.5229 (Not Used for Verification)
#30746 RT: 51.8440 min ITMS, 795.4291@cid25.00 979.5129@hcd30.00 $\sum_{100}^{5} 60 - \frac{1}{40} \frac{\alpha b_4^+}{439.23} \frac{\alpha y_{15}^{2+}}{879.76} \frac{\alpha y_{11}}{\alpha y_9} \frac{\alpha y_{12}}{\alpha y_{9}} \frac{\alpha y_{14}}{\alpha y_{9}} \frac{\alpha y_{14}}{\alpha y_{14}} \frac{\alpha y_{14}}{\alpha y_{14}} \frac{\alpha y_{14}}{\alpha y_{14}} \frac{\alpha y_{14}}{\alpha y_{14}}$	#30747 RT: 51.8453 min ITMS, 795.4291@cid25.00 966.5229@hcd30.00 $\sum_{150}^{150} \frac{1}{4} \alpha y_{15}^{2+} \alpha y_{12}^{+} \alpha y_{12}^{+}$ 338.20 866.62 αy_8 1380.80 αy_{14} 0 1000 1500 m/z
Peptide B (Amine) m/z = 610.8481	Peptide B (Isocyanate) m/z = 623.8391 (Not Used for Verifica
#30744 RT: 51.8400 min ITMS, 795.4291@cid25.00 610.8481@hcd30.00 200 $\frac{1}{\beta a_2^+}$ $\frac{\beta y_8^{2+}}{504.97}$ $\frac{\beta y_9^{2+}}{554.52}$ $\frac{\beta y_8^+}{1008.60}$ 100 $\frac{1}{200}$ $\frac{1}{400}$ $\frac{1}{600}$ 800 1000 1200 m/z	#30745 RT: 51.8428 min ITMS, 795.4291@cid25.00 623.8391@hcd30.00 $\sum_{strong}^{10} \frac{\beta b_2^+}{213.22} \beta y_5 \beta a_5 \frac{\beta y_7^+}{920.55} 986.81$ $\frac{\beta y_3}{200} \frac{\beta y_3}{400} \frac{\beta y_5}{600} \frac{\beta a_5}{1000} \frac{920.55}{1200} \frac{986.81}{1200}$ m/z

■ OTOT ■ OTAS

Figure 6. Venn diagram showing the overlap of unique identified RNA crosslinks in duplicates from (A) UV- or (B) NM samples in OT Astral DDA runs







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/xlms