

Iman Mohtashemi, Jie Qian, and David M. Horn  
Thermo Fisher Scientific, San Jose, CA, USA

Overview

**Purpose:** Disulfide mapping with minimal sample preparation and data interpretation

**Methods:** Following limited proteolysis, beta lactoglobulin peptides were analyzed using electron transfer dissociation (ETD)-triggered MS<sup>3</sup> experimental paradigm and disulfide bonds were identified by multiple software platforms

**Results:** We have demonstrated a workflow for discovery of disulfide bonds using ETD triggered MS<sup>n</sup> techniques.

Introduction

The biological function of proteins is a result of a number of factors including conformation, alternative splicing and a plethora of known and unknown modifications. Among the important contributors to tertiary and quaternary structure is disulfide linkages. Disulfide linkages influence not only the structural integrity of proteins but also their biological functions. Characterization of these linkages is important in many biological assay protocols. Protein structural characterization continues to present major challenges for the biological mass spectrometry community. Traditional collision induced dissociation (CID) is inefficient at cleaving disulfide bonds and produces difficult-to-interpret heterogeneous MS/MS ion populations. We present a complimentary approach leveraging multiple software platforms along with a novel ETD triggered MS<sup>3</sup> experimental paradigm.

Methods

Sample Preparation

Beta lactoglobulin was used as a model protein to study disulfide linkages in these experiments. The native protein was partially digested using trypsin for limited proteolysis. This method obviates the need for reduction/alkylation and extensive sample manipulation. The UniProt entry for beta lactoglobulin is shown below. The disulfide bonds from the UniProt entry were taken from two literature references.<sup>1, 2</sup>

UniProt ID P02754:

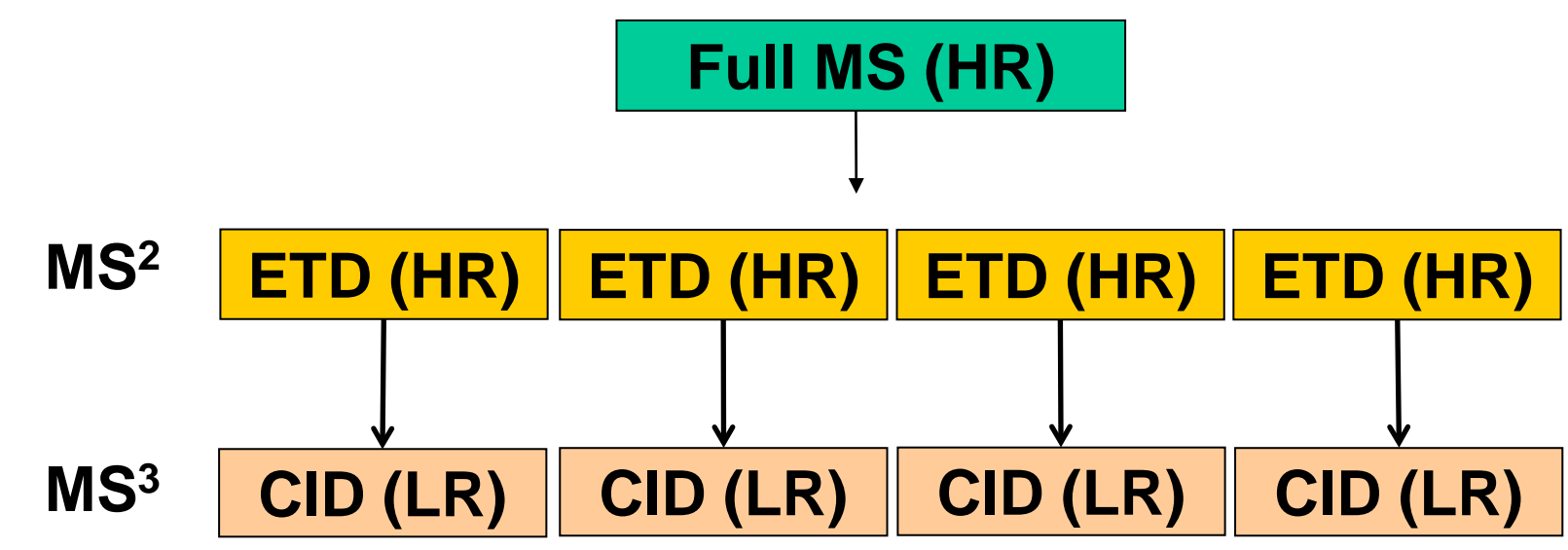
10	20	30	40	50	60
MKCLLLAL	TCGAQALIVT	QTHKGLDIQK	VAGTWYSLAM	AASDISLDA	QSAPLRVYVE
70	80	90	100	110	120
ELKFTPEGDL	EILQKWENG	ECAQKIIAE	KTQIPAVFKI	DALNENKVLV	LDTVYKKYLL
130	140	150	160	170	
FCMENSAPFE	QSLACQLVLR	TPEVDDEALE	KFDKALKALP	MHIRLSFNPT	QLEEQCHI

Disulfide bond	82 ↔ 176	(Ref.13) (Ref.19)
Disulfide bond	122 ↔ 137	Alternate (Ref.13)
Disulfide bond	122 ↔ 135	(Ref.13) (Ref.19)

Data Acquisition

An aliquot of the digest was loaded onto a microcapillary analytical column. Peptides were gradient-eluted and analyzed using a hybrid ion trap- Orbitrap™ mass spectrometer for ETD triggered CID MS<sup>3</sup> experiment. In this method, the mass spectrometer was utilized to acquire high-resolution precursor ion spectra with top 4 high-resolution tandem mass spectra (MS/MS). Also, each MS/MS spectrum was followed by a CID MS<sup>3</sup> spectrum of the most intense ion from the corresponding MS/MS scan.

**FIGURE 1. Schematic view of an ETD triggered MS<sup>3</sup> experimental paradigm. HR represents high resolution data detected by the Orbitrap mass spectrometer while LR represents low resolution data acquired in the ion trap.**



Data Analysis

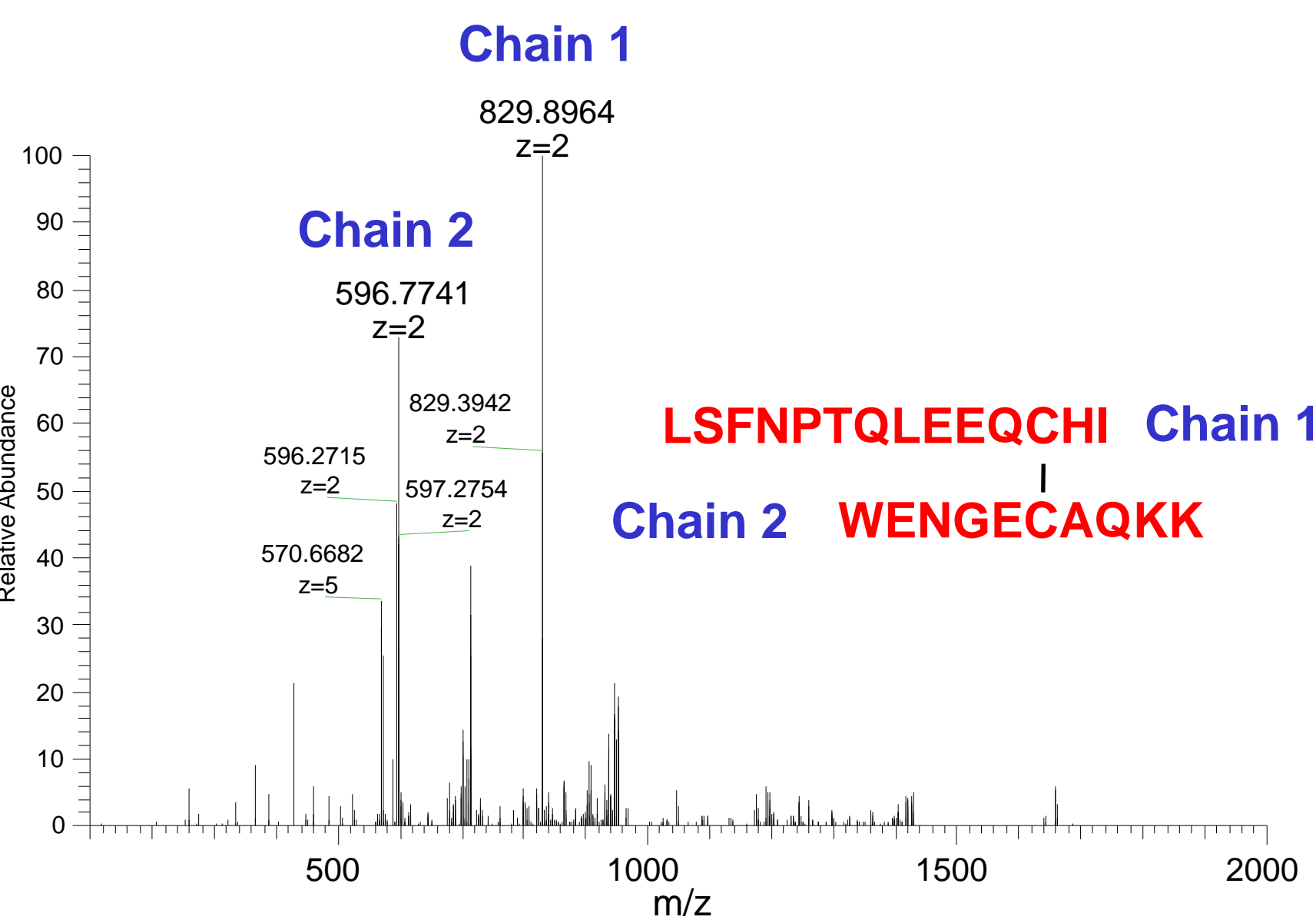
For disulfide mapping, the data were first analyzed using Thermo Scientific Proteome Discoverer software version 1.3. A workflow was created to search the CID MS<sup>3</sup> spectra using SEQUEST, with the precursor mass for the spectrum chosen to be the MS2 fragment that triggered the MS3 event. Thermo Scientific ProSightPC software version 2.0 was subsequently used to confirm the peptides identified by the Proteome Discoverer™ software by searching the high resolution ETD MS/MS spectra using the “Delta M” error tolerant search mode with a precursor mass tolerance of 3000 Da and a fragment tolerance of 10 ppm.

Results

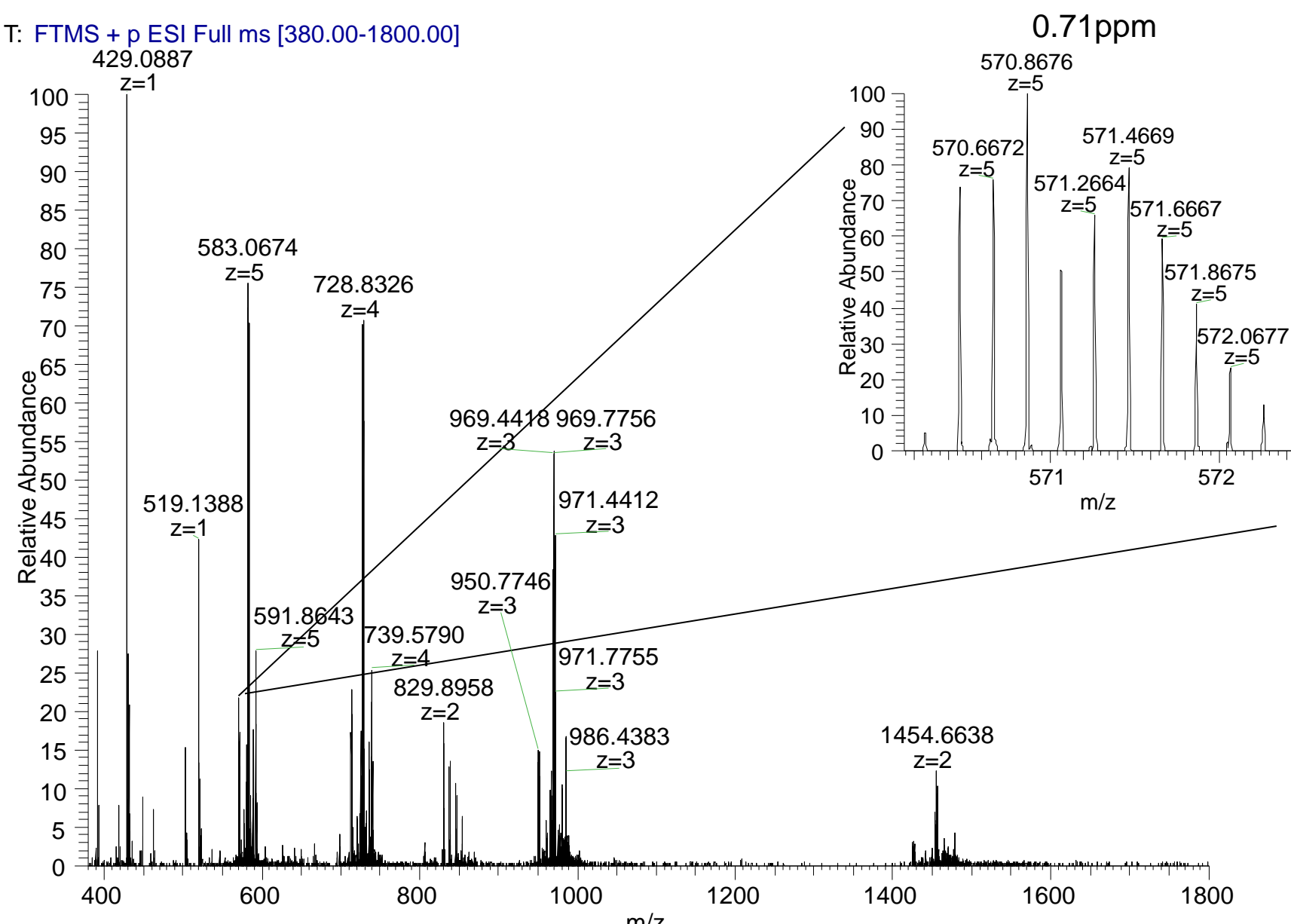
Unlike collision-induced dissociation (CID), ETD favors fragmentation of disulfide linkages.<sup>3</sup> For disulfide bound peptide, we expected that ETD fragmentation will favor dissociation of the two peptides at the disulfide bond over backbone cleavage, producing abundant fragments corresponding to the unlinked peptides. Figure 2 shows one such case, where the two major fragments in the MS/MS spectrum correspond to two tryptic fragments from beta-lactoglobulin that are known to be disulfide bound.

These peptides could then be isolated and fragmented using CID in the ion trap for subsequent identification by database search. The high mass measurement accuracy of the precursors (Figure 3) and the high sequence coverage of the CID MS<sup>3</sup> data lead to very confidently identified disulfide bound peptides.

**FIGURE 2. ETD MS<sup>2</sup> spectrum of disulfide linked peptides. ETD dissociated the two peptides linked by a disulfide bond, producing intense fragments corresponding to the unlinked peptides. Masses corresponding to the dissociated chains were among the most abundant species since ETD favors fragmenting disulfide bonds.**



**FIGURE 3. Detection of disulfide linked peptides LSFNPTQLEEQCHI and WENGECQKK in full MS scan.**



Disulfide Mapping Using Proteome Discoverer Software

When searching only the ETD MS/MS data by SEQUEST, no cysteine-containing peptides were identified as shown in the sequence map of beta-lactoglobulin shown in Figure 4.

**FIGURE 4. Protein coverage by ETD MS<sup>2</sup> spectra. Green highlights represent the portions of the sequence where peptides were identified with high confidence.**

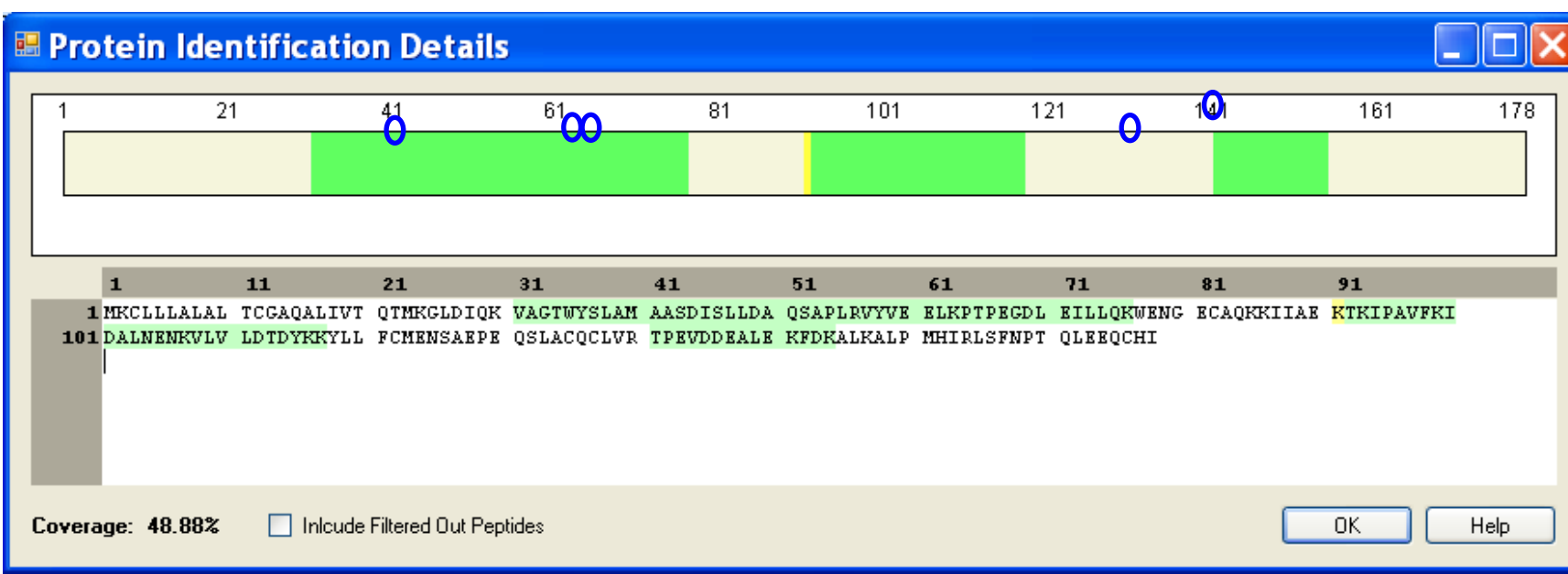
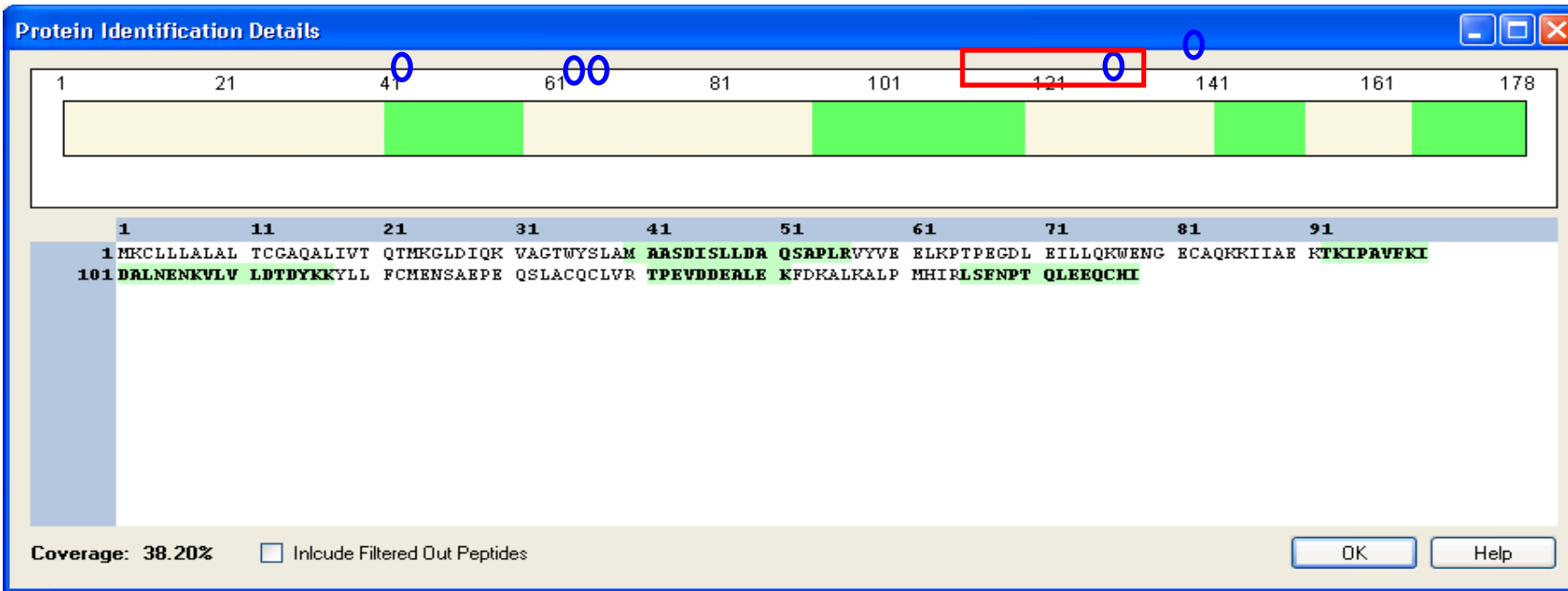
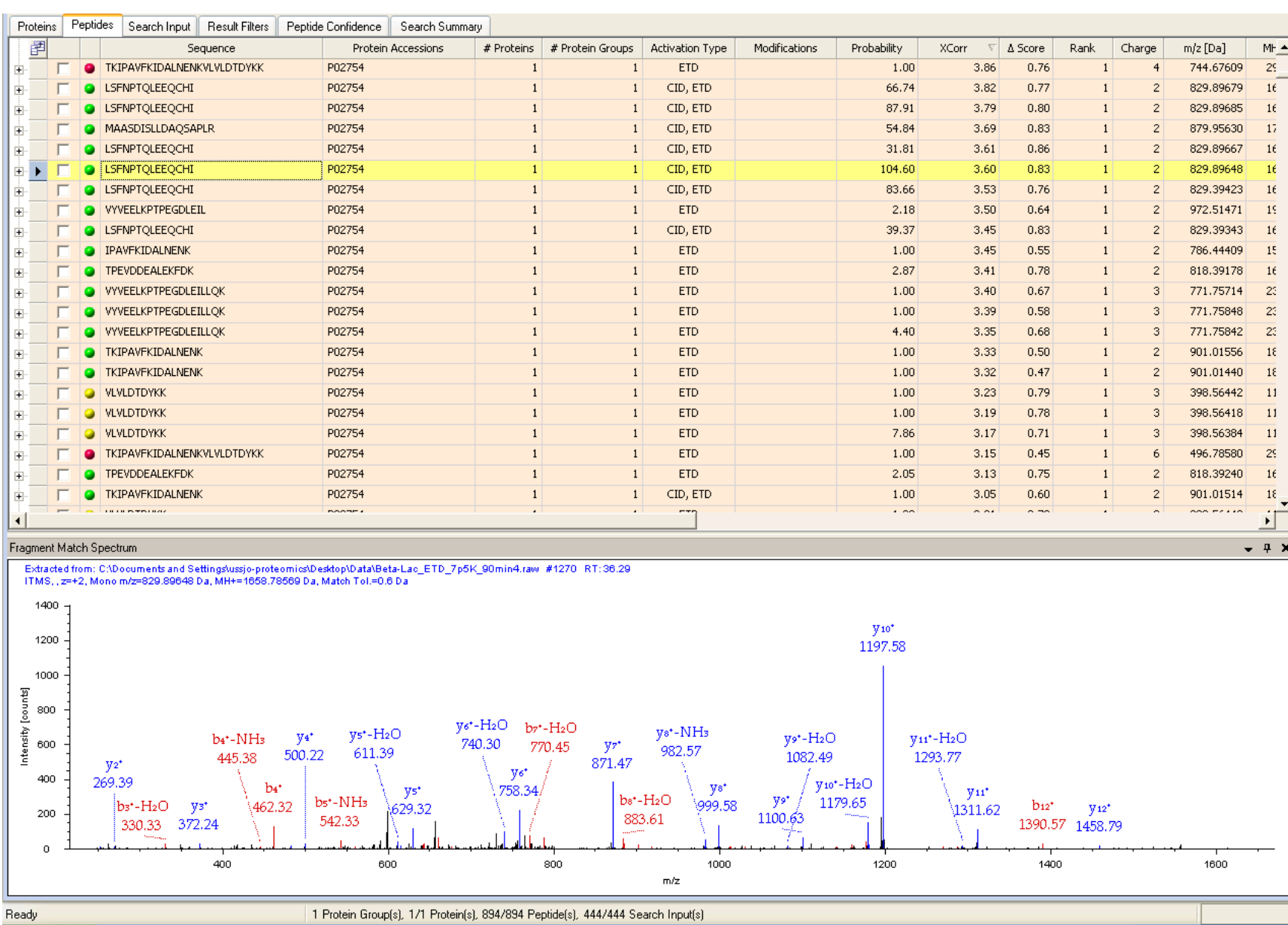


Figure 5 shows the sequence coverage when the CID MS<sup>3</sup> are included, where an additional cysteine-containing peptide at the C-terminus was identified. Figure 6 shows the Proteome Discoverer software results highlighting the C-terminal peptide detected by MS<sup>3</sup>.

**FIGURE 5. Protein coverage by combined ETD MS/MS and CID MS<sup>3</sup> spectra.**



**FIGURE 6. From CID MS<sup>3</sup>, identification of the cysteine-containing peptide LSFNPTQLEEQCHI using SEQUEST in Proteome Discoverer software. The upper panel shows a list of peptides identified by CID MS<sup>3</sup> spectra including LSFNPTQLEEQCHI; bottom panel demonstrates a CID MS<sup>3</sup> spectrum for LSFNPTQLEEQCHI with fragment ions annotated.**



A dynamic chemical modification for the mass of the peptide LSFNPTQLEEQCHI that included loss of two hydrogen [M-2H] was created in Proteome Discoverer software and the ETD MS/MS data was searched again using this modification with SEQUEST. As seen in Figure 7, the complimentary chain for the disulfide linked peptides was confidently identified as WENGECQKK.

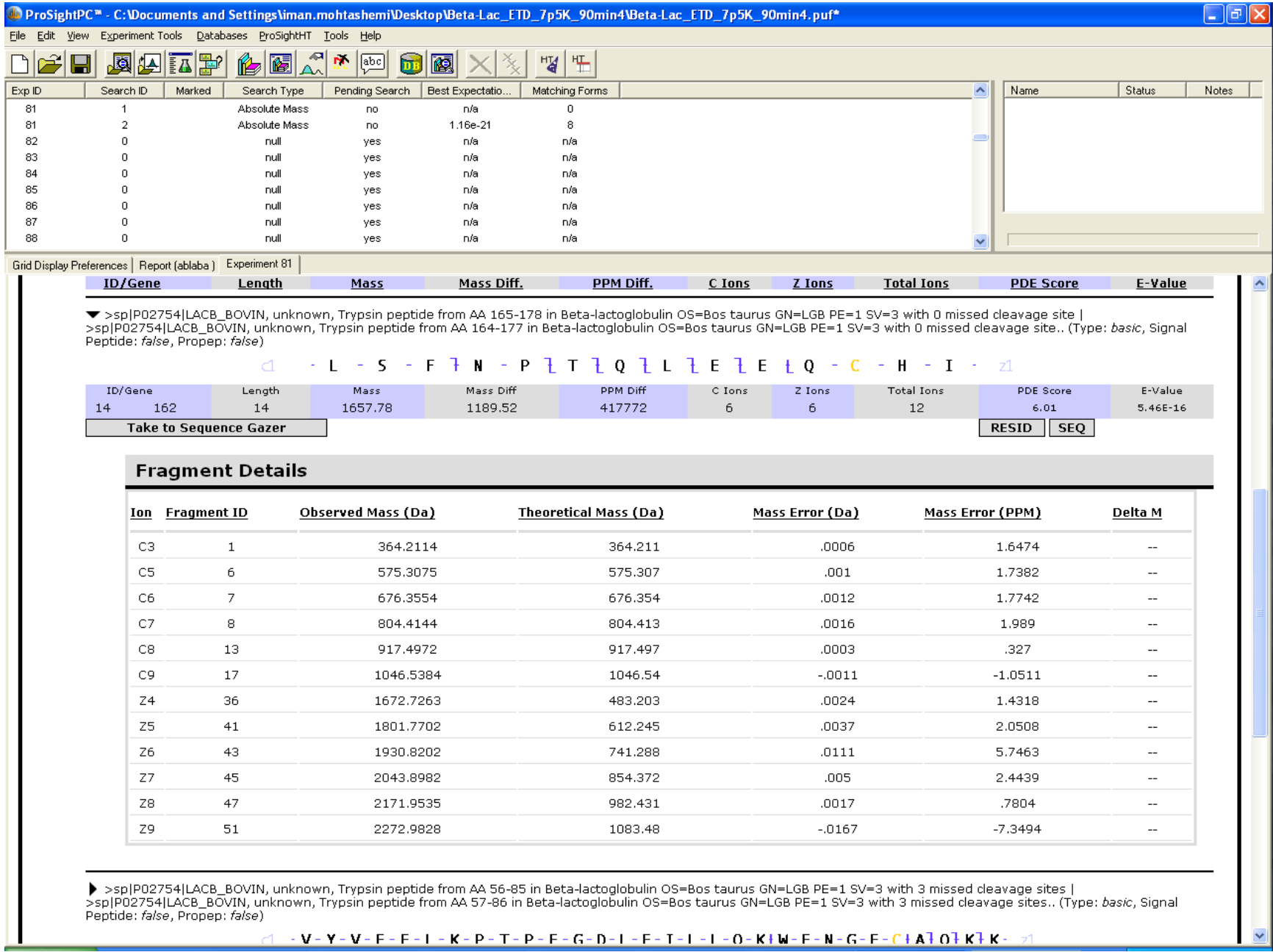
**FIGURE 7. From ETD MS<sup>2</sup>, WENGECQKK was identified as the other chain of the disulfide linked peptides using LSFNPTQLEEQCHI with loss of two hydrogen [M-2H] as a dynamic chemical modification.**

Protein	Peptides	Search Input	Filters	Peptide Confidence	Search Summary
WENGECQKK	WENGECQKK	ETD	CALSFNPTQLEEQCHI	5	2846.30488
WENGECQKK	WENGECQKK	ETD	CALSFNPTQLEEQCHI	4	2846.30556
WENGECQKK	WENGECQKK	ETD	CALSFNPTQLEEQCHI	4	2846.30556
WENGECQKK	WENGECQKK	ETD	CALSFNPTQLEEQCHI	3	2846.30556
WENGECQKK	WENGECQKK	ETD	CALSFNPTQLEEQCHI	3	2846.30556

Disulfide Mapping Using ProSightPC Software

ProSightPC software does not have a “variable” modification search mode for unknown post translational modification identification. Instead, it can searches for any single modification of any mass using a “delta-M” mode. In delta-M mode, another set of fragments are matched to the sequence: b + “delta-M” and y + “delta-M”. The delta-M is calculated as the mass difference between the measured precursor and the matched sequence. In this case, to each chain of the disulfide linked peptides, mass of the other chain with loss of two hydrogen would be the calculated delta mass. In the example shown in Figure 8, the peptide LSFNPTQLEEQCHI was very confidently identified by ProSightPC software with an E-value of 5.46e-16 from the ETD MS/MS data with a delta mass calculated to be 1189.52, which is 2 Da smaller than the theoretical mass of WENGECQKK. This further confirms the disulfide bond detected by the CID MS<sup>3</sup> method and Proteome Discoverer software.

**FIGURE 8. Identification of peptide LSFNPTQLEEQCHI by ProSightPC software using the delta-M mode. E-value was 5.46e-16 and the delta mass was 1189.52 Da. This confirmed the disulfide bond detected by CID MS<sup>3</sup> and Proteome Discoverer software.**



Conclusion

Here we present a comprehensive approach leveraging multiple software platforms along with a novel ETD triggered MS<sup>3</sup> experimental paradigm.

- ETD fragmentation dissociated the two peptides linked by a disulfide bond, producing intense fragments corresponding to the unlinked peptides which could then be isolated and fragmented by CID for identification.
- Without knowledge a priori of the disulfide linkages, we could detect inter-peptide disulfide linkages using this method. In this case, disulfide linkage between Cys82 and Cys176 was identified unambiguously.
- Analysis of the data using Proteome Discoverer software allowed the unlinked peptide in the CID MS3 spectra to be identified and then to be used as a modification to identify the linked peptide in the ETD MS<sup>2</sup> spectrum.
- Analysis of the data using delta-M mode in ProSightPC software also allowed identification of the interpeptide disulfide linkage.

References

- McKenzie HA, Ralston GB, Shaw DC. Location of sulphydryl and disulfide groups in bovine - lactoglobulins and effects of urea. *Biochemistry*. 1972 Nov 21;11(24):4539-47.
- Kuwata K, Hoshino M, Forge V, Era S, Batt CA, Goto Y. Solution structure and dynamics of bovine beta-lactoglobulin A.. *Protein Sci*. 1999 Nov;8(11):2541-5.
- Zubarev RA, Kruger, NA, Fridriksson EK, Lewis MA, Horn DM, Carpenter BK, McLafferty FW. Electron capture dissociation of gaseous multiply-charged proteins is favored at disulfide bonds and other sites of high hydrogen atom affinity. *J Am. Chem. Soc*. 1999; 121 (12):

SEQUEST is a registered 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 manners that might infringe the intellectual property rights of others.