

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 MS³ 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 MSⁿ 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 MS³ 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.^{1, 2}

UniProt ID P02754:

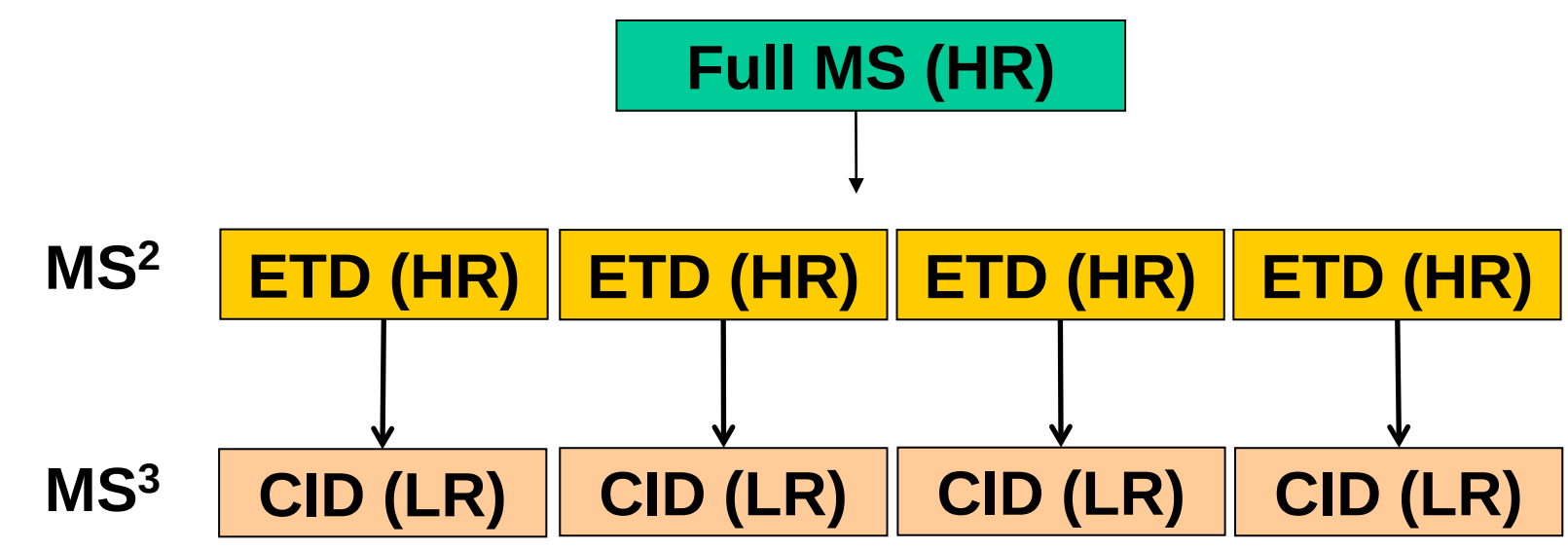
10	20	30	40	50	60
MKCLLLAL	TCGAQALIVT	QTHKGLDIQK	VAGTWYSLAM	AASDISLDA	QSAPLRVYVE
70	80	90	100	110	120
ELKFTPEGL	EILQKWMNG	ECAQKIIAE	KTQIPAVFKI	DALMENKVLV	LDTVYKKYLL
130	140	150	160	170	
FCMENSAEPE	QSLACQLVLR	TPEVDDEALE	KFDKALKALP	MHIRLSFNPT	QLEEQCHI

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 MS³ 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 MS³ spectrum of the most intense ion from the corresponding MS/MS scan.

FIGURE 1. Schematic view of an ETD triggered MS³ 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 MS³ 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.³ 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 MS³ data lead to very confidently identified disulfide bound peptides.

FIGURE 2. ETD MS² 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.

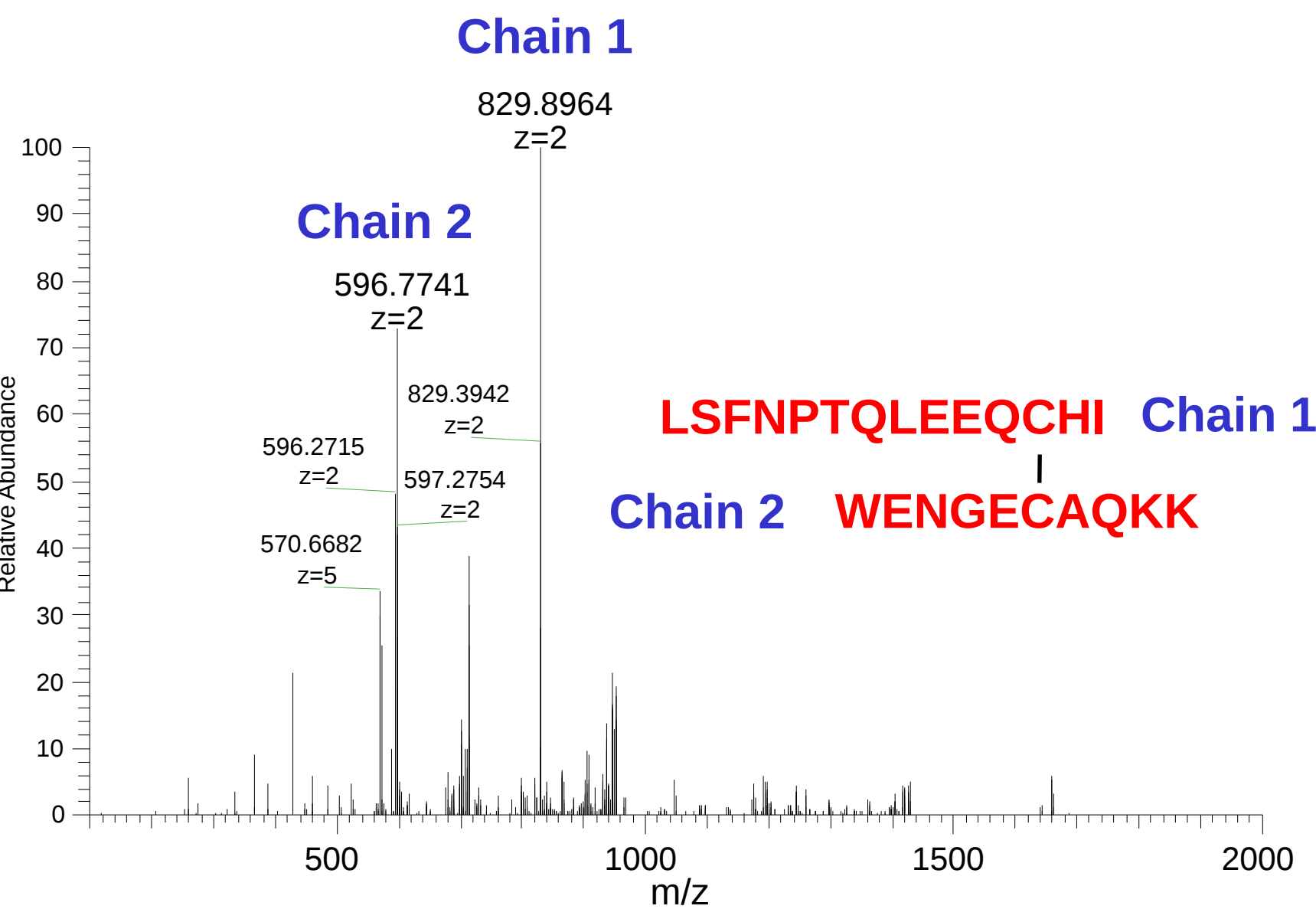
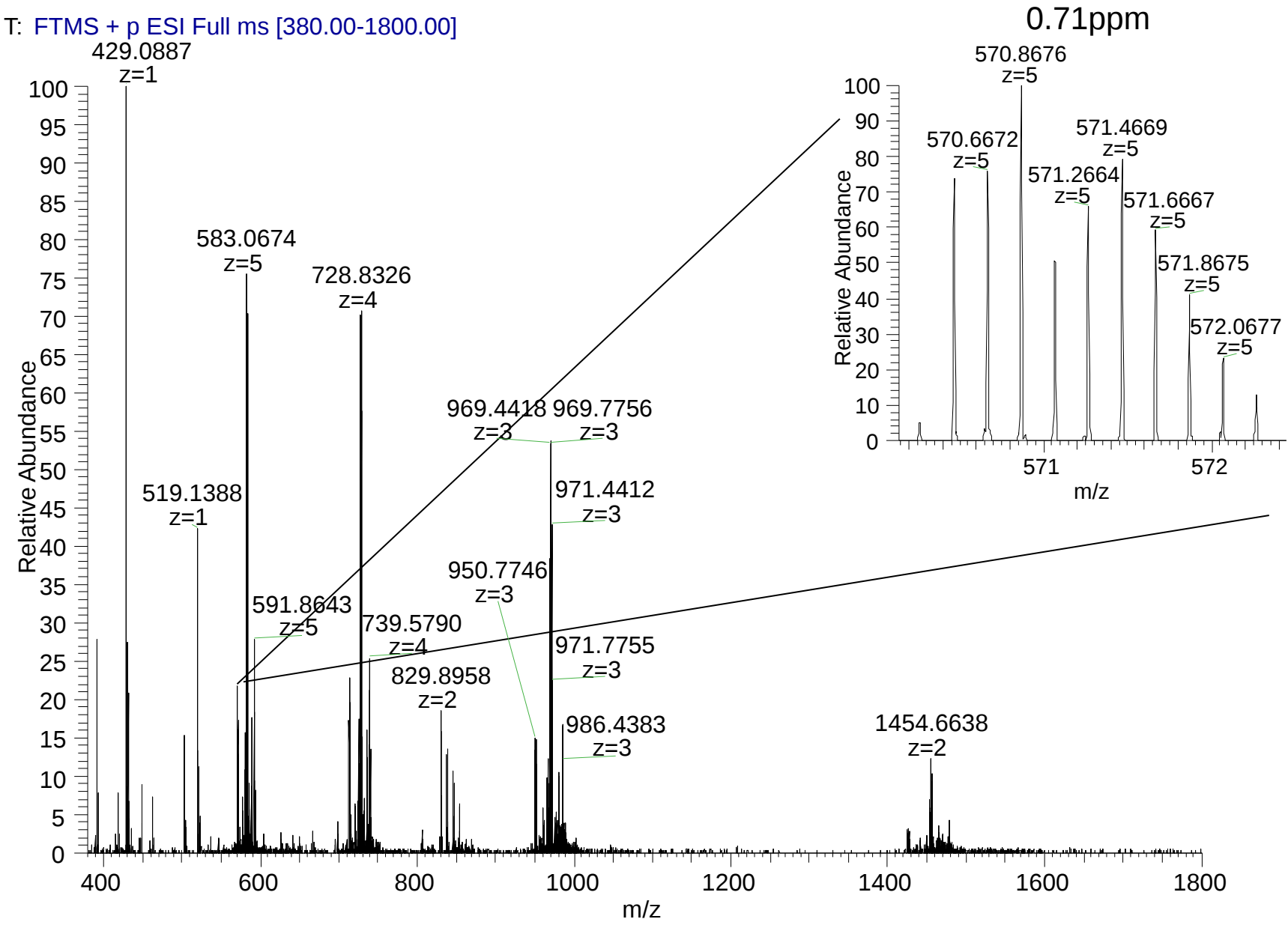


FIGURE 3. Detection of disulfide linked peptides LSFNPTQLEEQCHI and WENGCAQKK in full MS scan.



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 MS² spectra. Green highlights represent the portions of the sequence where peptides were identified with high confidence.

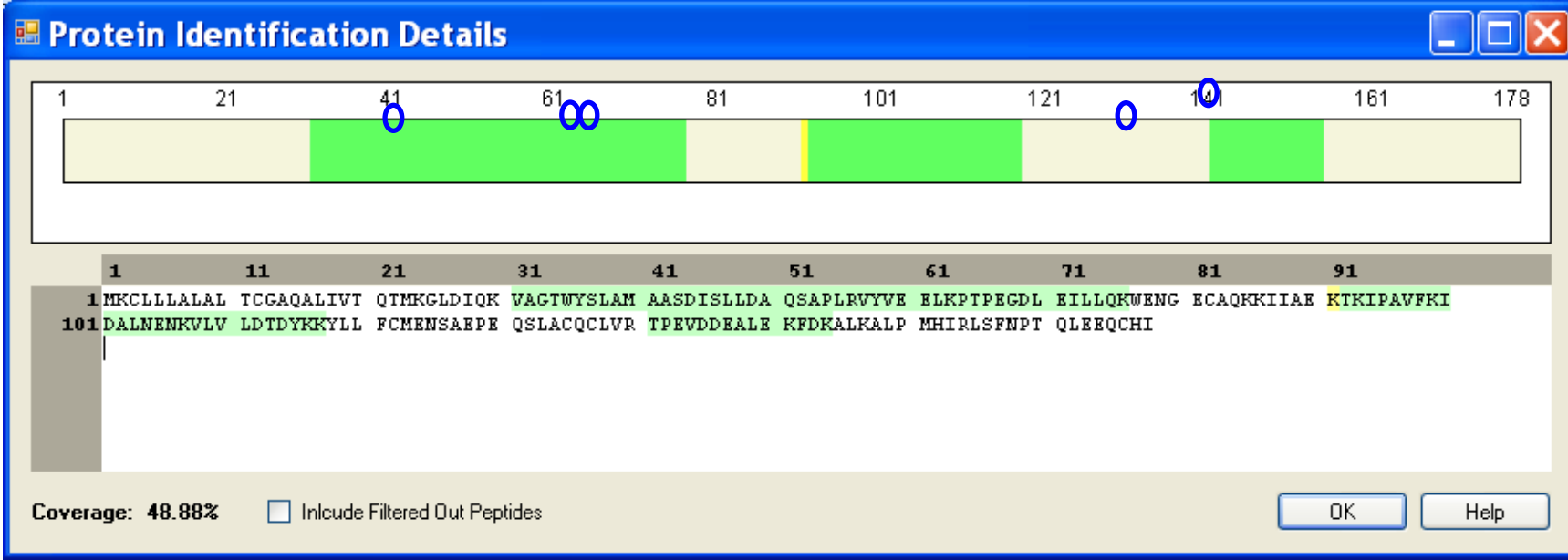


Figure 5 shows the sequence coverage when the CID MS³ are included, where an additional cysteine-containing peptide at the C-terminus was identified. Figure 6 shows the Proteome Discoverer software results highlighting the C-terminal peptide detected by MS³.

FIGURE 5. Protein coverage by combined ETD MS/MS and CID MS³ spectra.

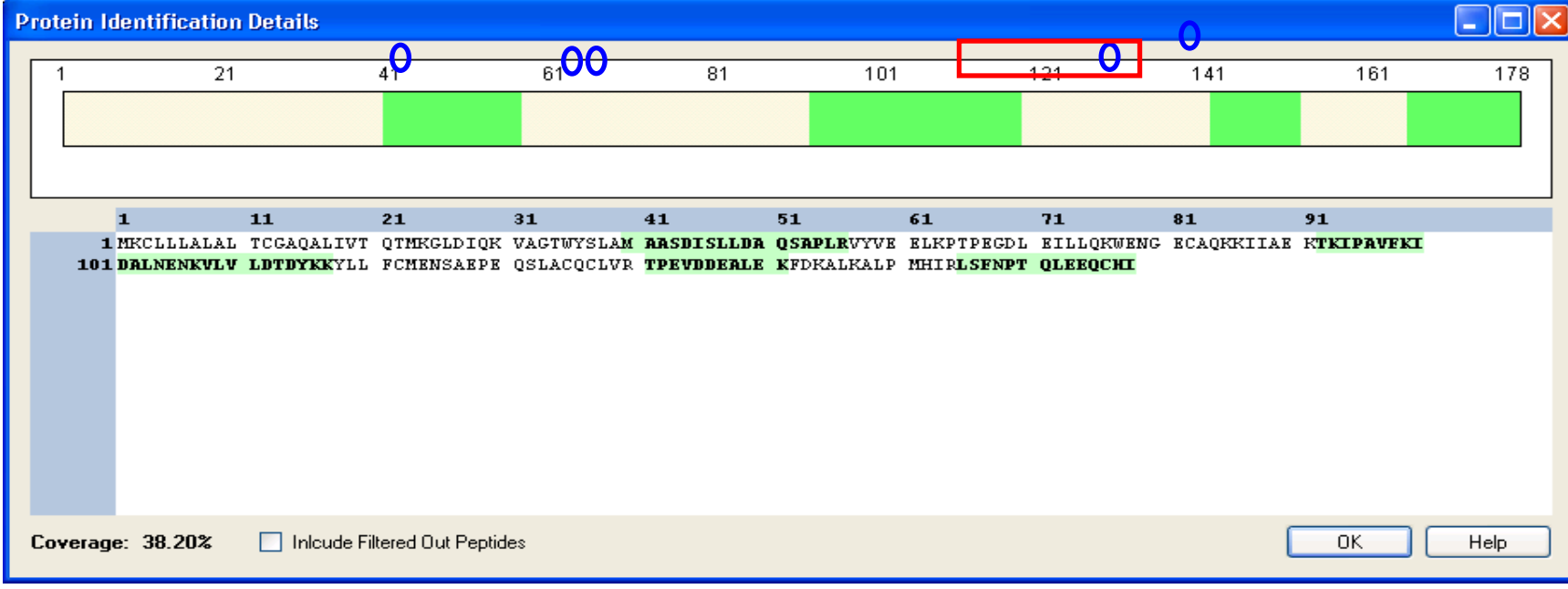
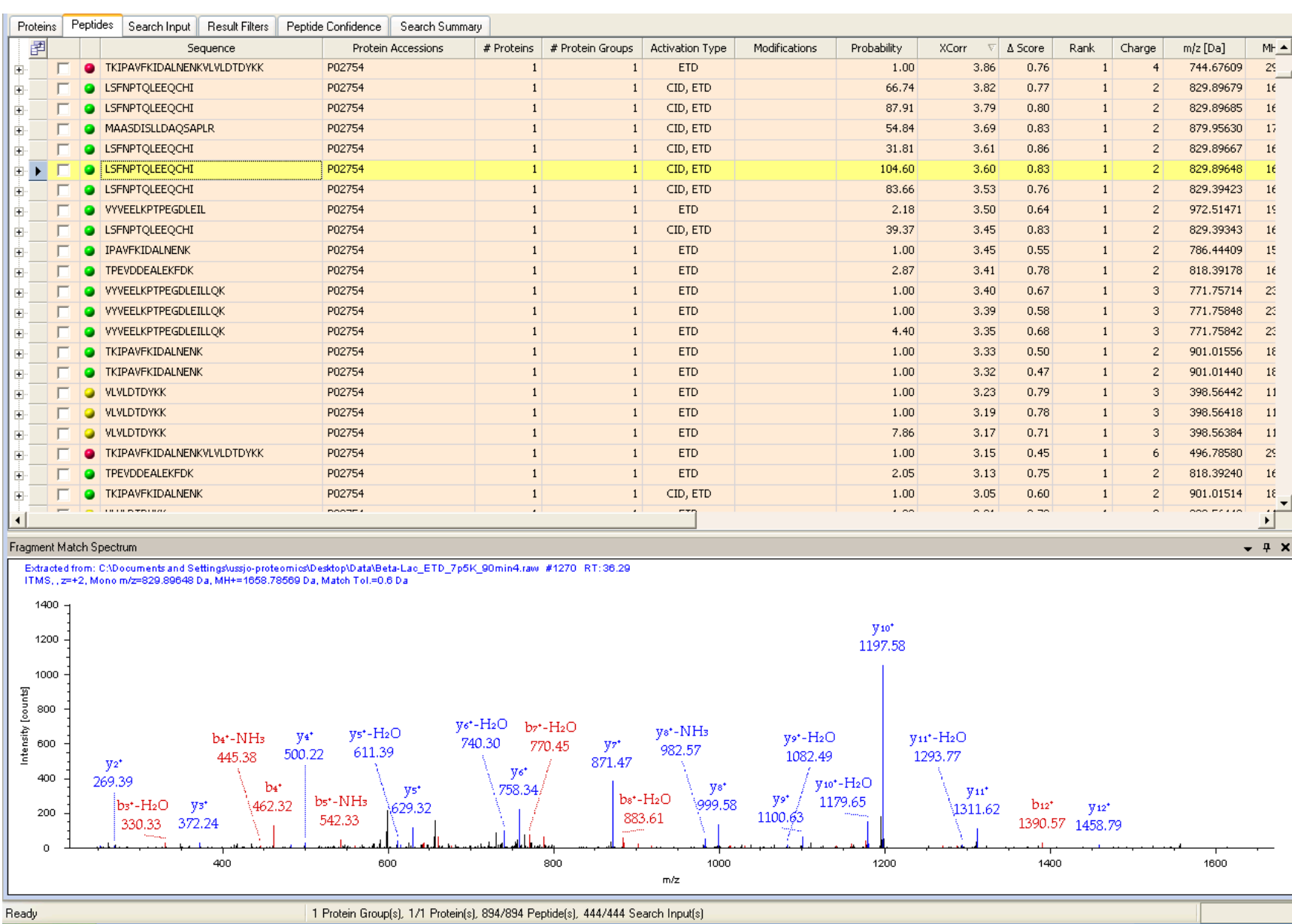


FIGURE 6. From CID MS³, identification of the cysteine-containing peptide LSFNPTQLEEQCHI using SEQUEST in Proteome Discoverer software. The upper panel shows a list of peptides identified by CID MS³ spectra including LSFNPTQLEEQCHI; bottom panel demonstrates a CID MS³ spectrum for LSFNPTQLEEQCHI with 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 WENGCAQKK.

FIGURE 7. From ETD MS², WENGCAQKK 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	Peptide	Search Input	Filters	Peptide Confidence	Search Summary					
	Sequence	Activation Type			Modifications	Y	Charge	MMH [Da]	ΔM [ppm]	First Scan
	☒ WENGCAQKK	ETD			CGSLNPTQLEEQCHI	5	2846.30488	1.15	871	
	☒ WENGCAQKK	ETD			CGSLNPTQLEEQCHI	4	2846.30556	1.39	875	
	☐ IDALNENKVLDTDWKK	ETD			4	2091.1540	1.00	887		
	☐ IDALNENKVLDTDWKK	ETD			3	2091.13474	0.69	889		
	☐ TRPEVDALFKFK	ETD			3	1635.77561	0.43	841		