

# Detection and Identification of New Protein Features using Multi-Attribute Monitoring (MAM) as Part of Mass Spectrometry-based Quality Control (QC)

Michael Blank, Stephane Houel, Kevin McCowen and Jonathan Josephs  
Thermo Fisher Scientific, 355 River Oaks Pkwy, San Jose, California 95134



CONTACT INFORMATION: Kevin McCowen  
Kevin.mccowen@thermofisher.com

## PURPOSE

The GMP compliant detection and quantitation of low level proteins in a drug substance modeling host cell proteins.

## OBJECTIVE

In addition to the expression of the desired drug molecule, transgenic host cells such as the widely employed Chinese Hamster Ovary cell (CHO) manufacture a number of native proteins associated with cell life cycle and maintenance. While these proteins are often excluded from the final drug product through a series of purification and chromatographic polishing steps, some are retained and present immunogenic effects that can adversely affect the safety of the therapeutic drug. Consequently the detection and quantification of these proteins are of great interest, but are challenging as many of these host cell proteins (HCPs) are at extremely low stoichiometry (on the order of parts per million) compared to the drug molecule. This work incorporates targeted ultra low level protein quantitation and non-targeted detection using a GMP-compliant multi attribute (MAM) quality control method using mass spectrometry.

## METHODS

### Sample Preparation

The NIST mAb standard commercially available at 10 mg/mL was adjusted to 1 mg/mL using 90 µL of 7M guanidine HCl with TRIS buffer. To this solution, 500 mM DTT was added as a reducing agent and reacted for 30 minutes at RT. Iodoacetic acid at 500 mM was added to alkylate the sample for 20 minutes. The sample was thoroughly buffer exchanged to remove the guanidine, following digestion with Pierce™ LCMs Grade trypsin in a 1:10 ratio for 30 minutes at 37 C. Digestion was quenched using 10% formic acid. The Pierce™ 6 Protein digest was spiked in at a calculated 1% molar ratio to the NIST mAb. A serial dilution was performed to reduce the concentration of the 6 proteins to 0.1% and ultimately to 0.001% (10 ppm concentration).

### Test Method(s)

Proteins and peptides present in the product and spiked samples were detected using data dependent acquisition (DDA). On the Thermo Scientific™ Q Exactive™ HF mass spectrometer under control of Thermo Scientific™ Chromleon™ CDS software, a 120,000 FWHM @ 200 m/z MS scan was followed by the HCD fragmentation of the top 5 most abundant precursors and detection of fragment ions at 30,000 resolution. For the MAM method, the data were acquired in triplicate using MS1 only from 300 to 1800 m/z at 120,000 resolution at an AGC target of 36k.

Chromatographic separation and delivery to the mass spectrometer was performed with a Thermo Scientific™ Dionex™ Vanquish™ Horizon uHPLC system coupled to a Thermo Scientific™ Accucore™ C18 uHPLC, 1.5 µm, 2.1 x 150 mm column maintained at 60 C. The binary solvent system consisted of 0.1% formic acid in water (A) and 0.1% formic acid in acetonitrile (B) flowed at a fixed rate of 0.20 mL/min. A multistage gradient was employed beginning with an equilibration phase for 5 minutes at 1% B before increasing abruptly to 10% at 6 minutes and eventually attaining 35% after 70 minutes. A column regeneration cycle at 90% B followed from 72 to 77 minutes before returning to 1% B at 79 minutes. A second "blanking gradient" see below brought the total run time to 115 min.

3 µL containing approximately 2.5 µg (15 pmol) of NISTmAb and as little as 0.15 fmol (10 ppm) of Pierce 6 Protein digest were loaded onto column for each injection which was performed in triplicate for each concentration down to 10 ppm. The eluent from the LC was diverted to waste for 2 minutes prior to being switched in line with the mass spectrometer. It was again diverted to waste at 70 minutes concurrent with the beginning of the column regeneration step.

### Data Analysis

The MS/MS data of the neat injection of the 6 protein digest were searched against a combined FASTA file of their respective sequences in Thermo Scientific™ BioPharma Finder™ software with carboxymethylation (+58.005 Da) as a fixed modification and without dynamic modifications, as these were expected to be of extremely low abundance and unsuitable for protein quantitation. The resultant list of peptides with a confidence of greater than 0.90 were exported as a target list for monitoring in a soon to be released version of Chromleon 7 software based on 7.2 SRS.

The MS spectral raw data acquired by Chromleon software were extracted using 5 ppm extraction tolerance for the targeted analysis with confirmation of peptides based upon their isotopic correlation being greater than 0.90 and their mass accuracy all confirming isotopes being less than 5 ppm. A weak moving average smoothing was applied.

Figure 1: Multi Attribute Method (MAM) workflow summary. A peptide mapping experiment followed by analysis in BioPharma Finder comprises the discovery phase to identify salient peptides. A targeted peptide list is then handed off to Chromleon to routinely and GMP compliantly monitor these targets, detect new features, and report out.



## RESULTS

### Peptide Mapping

A search of the peptide mapping data for the neat injection of the Pierce 6 Protein Digest revealed a total 384 components identified from their MS/MS spectra. Identification took into account not only mass accuracies of both the precursor and the fragment ions but also the correlation between the theoretically predicted fragmentation spectra and the observed spectra. For many of the detected peptides, multiple charge states were observed. This list was filtered only for those of the very highest confidence and then exported into Chromleon software for method optimization.

The imported list was substantially larger than any that would be typically used for a MAM method but was able to be processed handily by this version of Chromleon software without issue. This enabled an opportunity to execute further method development that sought to eliminate peptides with poor signal to noise, interference from the NIST mAb, low ionization efficiency, or unfavorable chromatographic behavior leaving a refined optimized method. Additionally, in the cases of multiple charge states usually at least one would be eliminated. While the inclusion of multiple charges is highly beneficial for drug product attribute monitoring as they can serve as additional confirmatory evidence and be used as part of the scoring algorithm, for the detection and quantitation of low level foreign proteins, some charge states were simply unsuitable as they did not have sufficient signal to noise at lower concentrations.

Figure 2: BioPharma Finder results showing the identification of a peptide based on the correlation between the observed MS/MS spectra and the theoretically predicted fragmentation pattern. High correlation and accuracy less than 3 ppm lead to a confident assignment. This peptide and many others can then be directly exported to Chromleon for monitoring as part of MAM.



### Targeted Quantification

As part of a system suitability evaluation, several critical criteria were evaluated prior to commencing primary data analysis, including retention time, peak area reproducibility, and mass accuracy. Several representative peptides from the NIST mAb sequence were used to evaluate this over three replicate runs. Average retention time variation was 0.1%, average peak area variation was 5%, and the median mass accuracy was 1 ppm, which was deemed acceptable for analysis to continue.

With the optimized list, it was possible to detect at least two peptides from each of the six proteins spiked into the NIST mAb digest to quantitatively model the behavior of HCPs (Table 1). In many cases all four confirming isotopes could clearly be observed at 5 ppm mass accuracy (Figures 3, 4, and 5) and could be automatically detected and quantified, especially when matched against the imported retention time values, lending a high degree of confidence to the assignment of these peptides and their resulting integration. This is coupled together with the fact that these peptides have already been identified previously using MS/MS.

Not unexpectedly, the two smallest proteins (Bovine Cytocrome C and Chicken Lysozyme) were the most difficult to detect as the pool of tryptic peptides to monitor was much smaller than the remaining four, not to mention that as these were by far the lowest in concentration (around 1 ppm). Overall, each of the peptides in Table 1 could be detected in each of the replicates, even at 10 ppm concentration, but not every targeted peptide was able to be effectively quantified according to ICH guidelines.

Table 1. Proteins and their peptides detected and quantified at 10 ppm concentration relative to the NIST mAb drug substance. At least two peptides for each protein could be detected and many quantified based on their linearity and coefficients of variation between technical replicates

Protein/Peptide	Peptide Name	Observed	Theoretical	Peak Area (mV*min)	Retention Time (min)	Mass Accuracy (ppm)	CV (%)
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.52	2.52	10.0	2.52	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.53	2.53	10.0	2.53	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.54	2.54	10.0	2.54	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.55	2.55	10.0	2.55	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.56	2.56	10.0	2.56	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.57	2.57	10.0	2.57	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.58	2.58	10.0	2.58	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.59	2.59	10.0	2.59	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.60	2.60	10.0	2.60	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.61	2.61	10.0	2.61	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.62	2.62	10.0	2.62	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.63	2.63	10.0	2.63	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.64	2.64	10.0	2.64	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.65	2.65	10.0	2.65	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.66	2.66	10.0	2.66	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.67	2.67	10.0	2.67	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.68	2.68	10.0	2.68	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.69	2.69	10.0	2.69	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.70	2.70	10.0	2.70	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.71	2.71	10.0	2.71	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.72	2.72	10.0	2.72	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.73	2.73	10.0	2.73	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.74	2.74	10.0	2.74	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.75	2.75	10.0	2.75	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.76	2.76	10.0	2.76	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.77	2.77	10.0	2.77	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.78	2.78	10.0	2.78	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.79	2.79	10.0	2.79	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.80	2.80	10.0	2.80	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.81	2.81	10.0	2.81	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.82	2.82	10.0	2.82	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.83	2.83	10.0	2.83	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.84	2.84	10.0	2.84	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.85	2.85	10.0	2.85	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.86	2.86	10.0	2.86	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.87	2.87	10.0	2.87	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.88	2.88	10.0	2.88	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.89	2.89	10.0	2.89	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.90	2.90	10.0	2.90	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.91	2.91	10.0	2.91	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.92	2.92	10.0	2.92	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.93	2.93	10.0	2.93	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.94	2.94	10.0	2.94	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.95	2.95	10.0	2.95	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.96	2.96	10.0	2.96	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.97	2.97	10.0	2.97	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.98	2.98	10.0	2.98	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	2.99	2.99	10.0	2.99	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.00	3.00	10.0	3.00	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.01	3.01	10.0	3.01	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.02	3.02	10.0	3.02	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.03	3.03	10.0	3.03	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.04	3.04	10.0	3.04	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.05	3.05	10.0	3.05	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.06	3.06	10.0	3.06	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.07	3.07	10.0	3.07	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.08	3.08	10.0	3.08	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.09	3.09	10.0	3.09	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.10	3.10	10.0	3.10	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.11	3.11	10.0	3.11	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.12	3.12	10.0	3.12	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.13	3.13	10.0	3.13	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.14	3.14	10.0	3.14	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.15	3.15	10.0	3.15	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.16	3.16	10.0	3.16	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.17	3.17	10.0	3.17	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.18	3.18	10.0	3.18	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.19	3.19	10.0	3.19	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.20	3.20	10.0	3.20	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.21	3.21	10.0	3.21	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.22	3.22	10.0	3.22	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.23	3.23	10.0	3.23	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.24	3.24	10.0	3.24	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.25	3.25	10.0	3.25	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.26	3.26	10.0	3.26	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.27	3.27	10.0	3.27	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.28	3.28	10.0	3.28	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.29	3.29	10.0	3.29	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.30	3.30	10.0	3.30	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.31	3.31	10.0	3.31	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.32	3.32	10.0	3.32	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.33	3.33	10.0	3.33	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.34	3.34	10.0	3.34	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.35	3.35	10.0	3.35	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.36	3.36	10.0	3.36	1.0	1.0
Bovine Cytocrome C	YVYVYVYVYVYVYVYVYVYV	3.37	3.37	10.0	3.37	1.0	