

# High Throughput Quantitation of 46 Histone PTMs through Unscheduled SRM-based Method Development on a Nano-HPLC Triple Quadrupole Platform

Jenny Chen,<sup>1</sup> Tommy Cheung,<sup>2</sup> David Arnott,<sup>2</sup> Yan Chen,<sup>1</sup> Keith Waddell,<sup>1</sup> Cindy Lai<sup>1</sup>

<sup>1</sup>Thermo Fisher Scientific, San Jose, CA; <sup>2</sup>Genentech, South San Francisco, CA

## Overview

Histone posttranslational modifications (PTM) was analyzed on a high throughput, unscheduled quantification platform using nano HPLC coupled with a Thermo Scientific™ TSQ™ Quantiva™ mass spectrometer. Forty-six modified histone peptides were selected for quantitation using 186 transitions in 35 and 60 minute program. The sensitivity and reproducibility of these research methods are evaluated. Results here establish a baseline to guide MS-based quantification method development for research.

## Introduction

Histone posttranslational modifications (PTM) are important aspects of epigenetic gene regulation, and are linked to many diseases. The core histones (H2A, H2B, H3, and H4) are densely populated with numerous PTMs. Most histone PTM analysis has involved antibody-based techniques, but mass spectrometry (MS) has emerged as an attractive alternative, overcoming limitations of antibodies such as poor specificity and epitope occlusion by nearby PTMs. The most common MS strategy is to analyze enzymatically-digested histones, with chemical modification of lysine residues for better detection. The complexity of histone peptide digests, with many closely related peptides and large dynamic ranges make fast and accurate quantitation challenging. We therefore developed a workflow for high throughput quantitation of histone peptides using nano-HPLC triple quadrupole instrumentation.

## Methods

### Sample Preparation

Core histones were acid-extracted from cultured HEK293T cells, column-purified by ion exchange, and desalted by perchloric acid precipitation. The purified histones were prepared for MS using a hybrid chemical derivatization strategy, whereby an initial conversion of free lysines to their propionylated forms under mild aqueous conditions was followed by trypsin digestion and labeling of new peptide N-termini with phenyl isocyanate. Labeled histone samples were desalted using C18-stage-tips. Approximately 100 ng of peptides were loaded onto a Thermo Scientific™ PepMap™ C18, 75µm x15cm column and separated by a linear gradient of 35 or 60 minutes. MS data were acquired in selected reaction monitoring (SRM) mode using "Scheduled" and "Unscheduled" SRM lists on TSQ Quantiva triple quadrupole MS.

### Mass Spectrometry

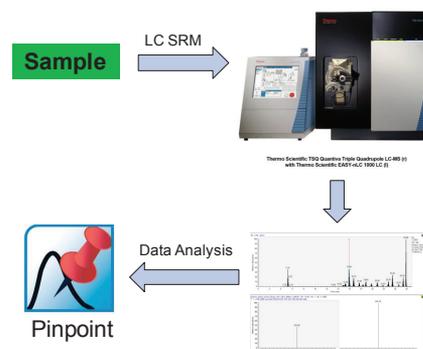
A mixture of peptides from human histones H3 and H4 was quantified on a Thermo Scientific™ Easy1000™ nano™ HPLC coupled with TSQ Quantiva triple quadrupole MS. These samples had been previously analyzed by data-dependent tandem mass spectrometry on a Thermo Scientific™ Orbitrap Elite™ mass spectrometer where a 90-minute HPLC program was required to resolve closely related and isobaric peptides. As a more rapid and targeted alternative for quantitation, 186 SRM transitions for "light" and "heavy" isotopically labeled forms of 46 peptides were established, first in "Unscheduled", and then "Scheduled" modes. In "Unscheduled" mode, all SRM transitions were performed in each duty cycle throughout the experiment, using 0.7 unit mass resolution in both Q1 and Q3, and chrome filter at 7 sec. Cycle times and collision gas pressures in the 60 minute program were investigated. All 186 SRMs were quantified by using targeted SRM methods at both 35 and 60 minute gradients.

A "Scheduled" mode was then implemented to improve sensitivity. Eight time segments in a 60-minute program were created. An even higher throughput was investigated using a 35min program in "Scheduled" mode.

### Data Analysis

Data was acquired using Thermo Scientific™ Xcalibur™ software version 2.2. Thermo Scientific™ Pinpoint™ software version 1.4 was used for simultaneous quantitative data processing.

FIGURE 1. Peptide Quantitation Workflow using nano-HPLC coupled with Quantiva MS



## Histone Peptide Information

A total of 46 histone peptides from 5 backbone peptides (the list below) were quantified with their isotopic <sup>13</sup>C-labeled heavy pairs using 186 SRMs. The peptide backbones are:

TKQTAR  
KSTGGKAPR  
KQLATKAAR  
KSAPATGGVKKPHR  
GKGGKGLGKGGAKR

## Results

### Quantitation Optimization using "Unscheduled" Mode

The quantitation of histone peptides are optimized using the heavy histone peptides in "unscheduled" mode. All the heavy peptide transitions were subjected to detection through the 60min LC run. Optimization of the cycle time and collision gas pressure found best signal for most peptides at a cycle time of 800ms with 1mTorr of collision gas. For each peptide, its collision energy was optimized based on its predicted collision energy by Pinpoint. In addition, with 500ms cycle time and 1ms interscan delay, all 186 SRMs were detected.

Table 1. Target List for "Unscheduled" SRM at 60min Run Time

SRM Table							
Compound	Start Time (min)	End Time (min)	Polarity	Precursor (m/z)	Product (m/z)	Collision Energy (V)	
1	K(Me2)SAPATGGV(KMe2)K(un)P(HR) light	5	55	Positive	416.996	434.24	24
2	K(Me2)SAPATGGV(KMe2)K(un)P(HR) heavy	5	55	Positive	416.996	532.322	24
3	K(Me2)SAPATGGV(KMe2)K(un)P(HR) heavy	5	55	Positive	418.493	440.231	24
4	K(Me2)SAPATGGV(KMe2)K(un)P(HR) heavy	5	55	Positive	418.493	532.322	24
5	K(Me2)SAPATGGV(KMe2)K(un)P(HR) light	5	55	Positive	420.499	434.24	22
6	K(Me2)SAPATGGV(KMe2)K(un)P(HR) light	5	55	Positive	420.499	539.33	22
7	K(Me2)SAPATGGV(KMe2)K(un)P(HR) light	5	55	Positive	420.501	448.256	22
8	K(Me2)SAPATGGV(KMe2)K(un)P(HR) light	5	55	Positive	420.501	532.322	22
9	K(Me2)SAPATGGV(KMe2)K(un)P(HR) heavy	5	55	Positive	421.997	440.231	22
10	K(Me2)SAPATGGV(KMe2)K(un)P(HR) heavy	5	55	Positive	421.997	539.33	22
11	K(Me2)SAPATGGV(KMe2)K(un)P(HR) heavy	5	55	Positive	421.999	454.247	22
12	K(Me2)SAPATGGV(KMe2)K(un)P(HR) heavy	5	55	Positive	421.999	532.322	22
13	TK(Me2)Q(TAR) light	5	55	Positive	426.241	475.262	28
14	TK(Me2)Q(TAR) light	5	55	Positive	426.241	631.388	28
15	TK(Me2)Q(TAR) heavy	5	55	Positive	429.236	475.262	28
16	TK(Me2)Q(TAR) heavy	5	55	Positive	429.236	631.388	28
17	TK(Me2)Q(TAR) light	5	55	Positive	433.23	475.262	23
18	TK(Me2)Q(TAR) light	5	55	Positive	433.23	645.367	23
19	TK(Me2)Q(TAR) light	5	55	Positive	433.249	373.729	28
20	TK(Me2)Q(TAR) light	5	55	Positive	433.249	620.341	28
21	TK(Me2)Q(TAR) heavy	5	55	Positive	436.226	475.262	23
22	TK(Me2)Q(TAR) heavy	5	55	Positive	436.226	645.367	23
23	TK(Me2)Q(TAR) heavy	5	55	Positive	436.244	373.73	28
24	TK(Me2)Q(TAR) heavy	5	55	Positive	436.244	626.36	28

Figure 2. Base Peak Chromatogram of 60 min "Unscheduled" Analysis of 186 SRMs

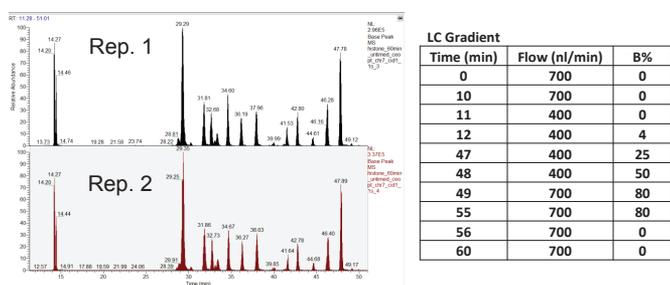


Figure 3. Optimization of Collision Energy (CE) for Heavy Peptide K(Me3)SAPATGGV(KMe2)K(un)P(HR)

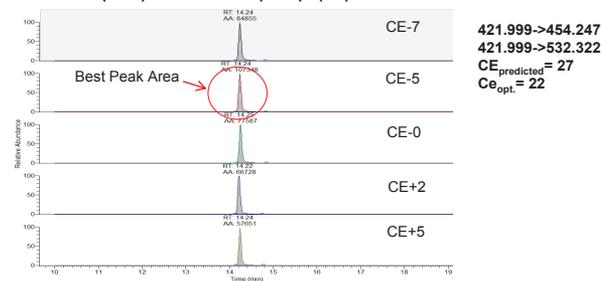


Figure 4. Cycle Time Optimization of Low Abundant Peptide [N-light]K[tri-Methyl]SAPATGGV(KMe2)K(un)P(HR) using Best Peak Area

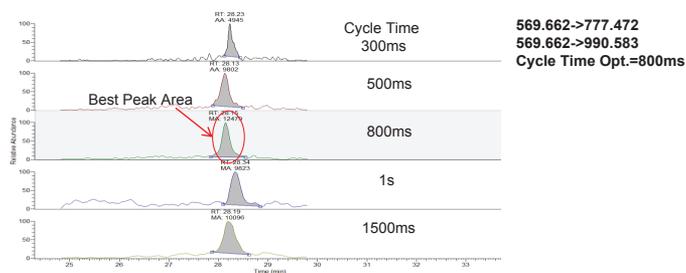
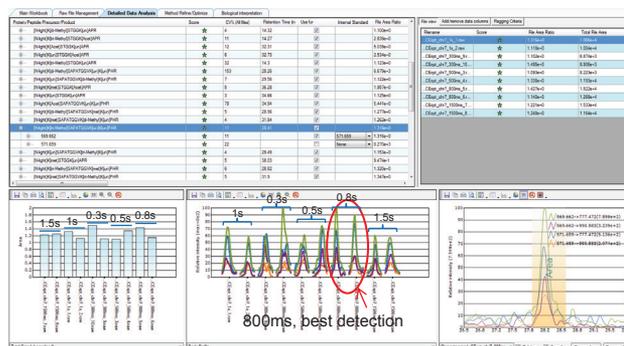


Figure 5. Pinpoint Comparison of Histone Peptide Detection at Different Cycle Time



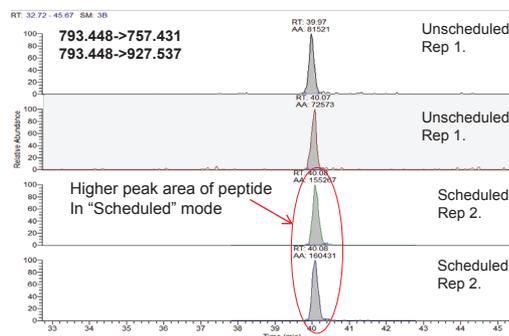
### Quantitation using "Scheduled" Mode

A "Scheduled" mode was then developed to improve sensitivity. Eight time segments in a 60-minute program were created, with less than 30 SRMs in each time segment. Quantitation of the 46 histone peptides via the Pinpoint program found high reproducibility, with CV < 15% (n=4). To achieve higher throughput, a 35min program in "Scheduled" mode was developed.

Table 3. Example of SRM List in "Scheduled" Mode

Compound	Start Time (min)	End Time (min)	Polarity	Precursor (m/z)	Product (m/z)	Collision Energy (V)	
1	TK(Me2)Q(TAR) light	11.7	16.7	Positive	433.249	373.729	28
2	TK(Me2)Q(TAR) light	11.7	16.7	Positive	433.249	620.341	28
3	TK(Me2)Q(TAR) heavy	11.7	16.7	Positive	436.244	373.73	28
4	TK(Me2)Q(TAR) heavy	11.7	16.7	Positive	436.244	626.36	28
5	K(Me2)SAPATGGV(KMe2)K(un)P(HR) light	11.8	16.8	Positive	416.996	434.24	24
6	K(Me2)SAPATGGV(KMe2)K(un)P(HR) heavy	11.8	16.8	Positive	416.996	532.322	24
7	K(Me2)SAPATGGV(KMe2)K(un)P(HR) heavy	11.8	16.8	Positive	418.493	440.231	24
8	K(Me2)SAPATGGV(KMe2)K(un)P(HR) heavy	11.8	16.8	Positive	418.493	532.322	24
9	K(Me2)SAPATGGV(KMe2)K(un)P(HR) light	11.8	16.8	Positive	420.499	434.24	22
10	K(Me2)SAPATGGV(KMe2)K(un)P(HR) light	11.8	16.8	Positive	420.499	539.33	22
11	K(Me2)SAPATGGV(KMe2)K(un)P(HR) light	11.8	16.8	Positive	420.501	448.256	22
12	K(Me2)SAPATGGV(KMe2)K(un)P(HR) light	11.8	16.8	Positive	420.501	532.322	22
13	K(Me2)SAPATGGV(KMe2)K(un)P(HR) heavy	11.8	16.8	Positive	421.997	440.231	22
14	K(Me2)SAPATGGV(KMe2)K(un)P(HR) heavy	11.8	16.8	Positive	421.997	539.33	22
15	K(Me2)SAPATGGV(KMe2)K(un)P(HR) heavy	11.8	16.8	Positive	421.999	454.247	22
16	K(Me2)SAPATGGV(KMe2)K(un)P(HR) heavy	11.8	16.8	Positive	421.999	532.322	22

Figure 6. Detection Comparison of Peptide "GK[un]GGK[un]GLGK[Acet]GGAK[Acet]R Using "Scheduled" vs "Unscheduled" Mode



The cycle time for "unscheduled" mode was 800ms for all 186 transitions. The cycle time for "Scheduled" mode was also 800ms for less than 30 transitions in each of the eight time segments. The detection sensitivity of "scheduled" mode was higher, because MS now spent longer time on each transition. In addition, quantitation of the 46 histone peptides by "scheduled" mode found higher reproducibility with tighter CV (CV<15%, n=4, Figure 7.). However, "Unscheduled" mode still achieved reasonable sensitivity performance for all transitions and served as a quick and straightforward starting point for method optimization.

### Histone Peptide Quantitation using 35 min HPLC run

A even higher throughput was achieved using a 35min program in "Scheduled" mode, maintaining sensitivity and quantitative performance. Figure 7. shows the base peak chromatogram of a 35min run for "Scheduled" quantitation of 46 histone peptides using 186 SRMs using also eight time segments. With shorter gradient, all the 186 transitions were still detected with good sensitivity and reproducibility (CV<15%, n=3). For example, Pinpoint analysis found that in 35min runs (Figure 9.) the histone peptide "[N-light]K[di-methyl]STGGK[un]APR was well detected compared to 60min runs (Figure 8.).

Figure 7. Base Peak Chromatogram of 35 min "Scheduled" Analysis of 186 SRMs

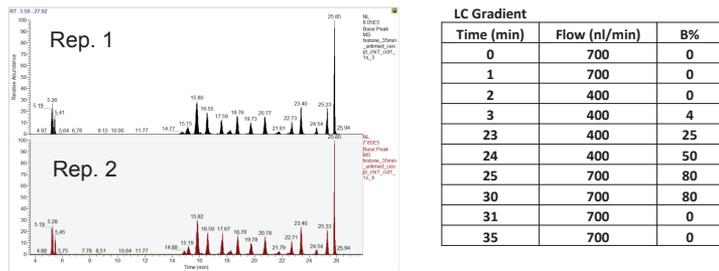


Figure 8. Pinpoint Analysis of Histone Peptide Detection at "Scheduled" Mode at 60min Run Found Higher Reproducibility with Tighter CV

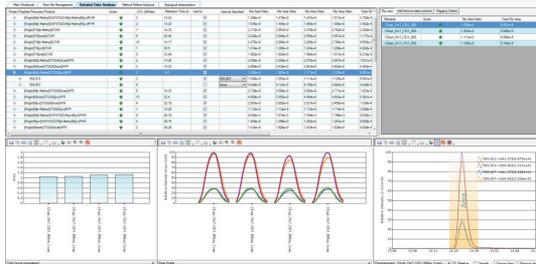
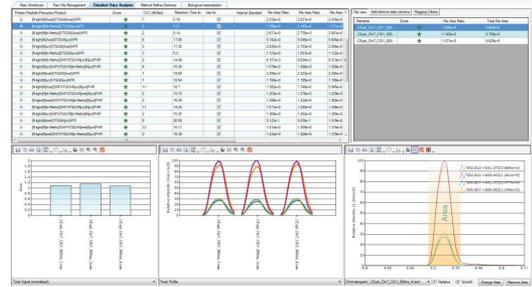


Figure 9. Pinpoint Analysis of Histone Peptide Detection at "Scheduled" Mode at 35min Run



## Conclusion

We developed a workflow for sensitive and high throughput quantitation of 46 modified histone peptides through 186 transitions in 60min runs using nano-HPLC couple with triple quadrupole MS. The high scanning speed of TSQ Quantiva triple quadrupole MS allows MS detection sensitivity optimization through "Unscheduled" mode. Further improvement of detection sensitivity was achieved using "Scheduled" time segments. In addition, we demonstrated quantitation of the 46 histone peptides using 35min runs with better throughput.

The nano-HPLC and TSQ Quantiva MS platform allowed histone peptides with abundances spanning three orders of magnitude to be quantified in a single experiment, separating closely related peptides with identical precursor masses. The simplicity and convenience of this workflow on the TSQ Quantiva MS allows triple quadrupole mass spectrometry to contribute powerfully to epigenetic research programs.

**For research use only. Not for use in diagnostic procedures.**

[www.thermoscientific.com](http://www.thermoscientific.com)

©2015 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

**Africa** +43 1 333 50 34 0  
**Australia** +61 3 9757 4300  
**Austria** +43 810 282 206  
**Belgium** +32 53 73 42 41  
**Canada** +1 800 530 8447  
**China** 800 810 5118 (free call domestic)  
 400 650 5118  
 PNE64449-EN 0615S

**Denmark** +45 70 23 62 60  
**Europe-Other** +43 1 333 50 34 0  
**Finland** +358 10 3292 200  
**France** +33 1 60 92 48 00  
**Germany** +49 6103 408 1014  
**India** +91 22 6742 9494  
**Italy** +39 02 950 591

**Japan** +81 45 453 9100  
**Korea** +82 2 3420 8600  
**Latin America** +1 561 688 8700  
**Middle East** +43 1 333 50 34 0  
**Netherlands** +31 76 579 55 55  
**New Zealand** +64 9 980 6700  
**Norway** +46 8 556 468 00

**Russia/CIS** +43 1 333 50 34 0  
**Singapore** +65 6289 1190  
**Spain** +34 914 845 965  
**Sweden** +46 8 556 468 00  
**Switzerland** +41 61 716 77 00  
**UK** +44 1442 233555  
**USA** +1 800 532 4752

**Thermo**  
 SCIENTIFIC  
 A Thermo Fisher Scientific Brand