Demonstration of SRM Method Visualization and Automated On-the-fly Retention Time Updating for Targeted Peptide Quantitation

Vane D. Sanghvi¹, Susan E. Abbatiello², Mary Blackburn¹, Qingyu Song¹, Jacek W. Sikora³, Yupeng Zheng³, Paul M. Thomas³, Neil Kelleher³ Thermo Fisher Scientific, ¹San Jose, CA; ²Cambridge, MA; ³Kelleher Research Group, Northwestern University, Evanston IL

ABSTRACT

It is a challenge to target hundreds of peptides in a single LC run by SRM-MS. Here we look at two software features that assist in better method definition and improved data acquisition:

- SRM visualization to observe concurrent transitions and approximate dwell times in a given method
- Real-Time RT updating during method acquisition to minimize chances of missing a schedule SRM transition due to chromatographic shifts

Results indicate successful "on-the-fly" updating of scheduled retention time windows as the LC gradient causes peptides to shift, allowing for improved data quality, easier method development, and fewer missed data.

INTRODUCTION

Peptide quantitation using LC-SRM-MS methodologies has progressed over the years to help answer increasingly difficult questions in biology. Targeting hundreds of peptides in a single LC run has resulted in technical challenges that require manual data evaluation to ensure all peptides eluted within their RT windows. Both sample composition and sample load can cause changes in peptide retention times from sample to sample, causing significant RT shifts. Increasing RT windows to ensure peptide detection comes at the cost of either lower dwell times or longer cycle times, either of which can have deleterious effects on LC-SRM-MS precision and sensitivity. These challenges have identified the need for more sophisticated acquisition software. Here we present the application of two software features that assist in visualizing the SRM method design as well as ensuring successful detection of peptide targets in a timed LC-SRM-MS method in the event of chromatographic shifts. The algorithms were evaluated on a set of PRTC peptides in HeLa lysate with several different reversed-phase gradients intended to cause chromatographic shifts. Use of the new features than design and resulted in a higher success rate of peptide detection.

MATERIALS AND METHODS

A sample set consisting of 15 13C/15-labeled peptides (PRTC peptides from Pierce) in HeLa lysate (Pierce) were evaluated on a Thermo Scientific™ TSQ Quantiva™ Triple Quadrupole MS. The LC-SRM-MS method targeted 467 transitions (60 precursors) using timed SRM with variable retention time windows. Out of 467 transitions, 15 were identified as "Retention Time Standards" with Reference Intensity Threshold set as 5.0e4 for 5 precursor *m*/z (493.768, 444.547, 695.832, 558.326, 567.973).

- A Thermo ScientificTM EASY-nLCTM 1000 system was used with a trapping column configuration for sample introduction and gradient delivery.
- Precolumn equilibration was set to 10 uL at 2.5 uL/min or a maximum pressure of 500 bar.
- Analytical column equilibration was set to 10 uL at 0.5 uL/min or a maximum pressure of 500 bar.
- Sample loading was performed at 2.5 uL/min for a total of 10 uL (from a 1 uL sample injection volume).
- Mobile phase A: 2% acetonitrile/0.1% formic acid (v/v); Mobile phase B: 90% acetonitrile/0.1% formic acid (v/v)
- Columns
 - Trap column: ProntoSil C18AQ (Bischoff Chromatography), 3um beads, 150 um x 3 cm
 - Analytical column: Picochip ProntoSil C18AQ (Bischoff Chromatography), 3um beads, 75 um x 10 cm
- Gradients: Seven different gradients were used to artificially shift peptide RTs, and are shown in Figure 3. Flow rate during gradient delivery was 300 nL/min.

To evaluate performance of the instrument control software, LC-SRM-MS data were acquired using three different approaches with each of the 7 chromatographic gradients:

- Unscheduled (all 467 transitions acquired in a 1 second cycle time)
- Scheduled (3 minute RT windows for all transitions with a 1 sec cycle time)
 dRT (1.5 minute RT windows for all transitions except those identified as "RT
- GRT (1.5 minute RT windows for all transitions except those identified as "R Standards", which would have longer RT windows)

Data were analyzed in Thermo Scientific[™] Xcalibur[™] software and Skyline (University of Washington).

SRM Visualization

The use of timed SRM acquisition is often a requirement for the detection and quantitation of a large number of peptides in complex sample matrices to increase the dwell time of transitions and improve analyte detection. Determination of the retention time windows used for each analyte is a balancing game to ensure adequate dwell times per transition without inflating the cycle time, causing reduced sampling across the chromatographic peak. SRM visualization in the method editor allows the user to see the dwell time and retention time windows of all the precursor masses by plotting it in a graph.

Use of SRM visualization allowed for improved detection in the low abundance peptides by pinpointing congested regions of the chromatogram and indicating the average dwell time per transition in those regions. Retention time windows in these regions were easily reduced in duration so that the average dwell time increased.



Figure 1. SRM Visualization Plots Visible in the Method Editor. Two new plots are displayed when "SRM Visualization" is enabled: "Dwell Time Per Transition" (upper right). And "Number of Transitions" (lower right). These plots help illustrate the approximate dwell times per precursor *m*/z and show where in the chromatographic gradient the most concurrent transitions elute, based on the scheduled RT method.





Graphical Representation

The dwell time graph is plotted for each scan event (precursor/product ion) showing the start and end time. The dwell time for each scan event at a given time and precursor is shown in the tooltip as the mouse hovers over each plot (tooltip shown in the lower right corner of each plot). The x axis shows the time scaled at 0.001 minutes and sorted based on the start time, while the y axis is sorted by the precursor mass.

The transitions graph plots the number of transitions at any given time between the experiment start and end time. The x axis shows the time scaled at the cycle time and y axis is the number of transitions.

RESULTS

Gradients for dynamic RT testing 100 90 80 + grad01 70 grad02 60 grad03 %B 50 grad04 40 grad05 30 20 ----- grad06 10 grad07 0 10 15 20 25 30 35 Time (min)

Figure 3. Seven reversed-phase gradients were used to test the ability of the dRT software feature to update scheduled RT windows "on-the-fly" during acquisition.

Samples were acquired on each of the 7 gradients in unscheduled methods to empirically determine the retention time of each of the 60 precursors (data not shown), then each gradient was re-run with a scheduled method. The chromatographic shifts from the different gradients in scheduled SRM mode are shown in Figure 4.



Figure 4. Scheduled SRM-MS chromatograms for each of the 7 gradients used in the study. Peaks are labeled with RT and m/z of the base peak. RT shifts from the "default" gradient (Grad01) were as much as 8 minutes. Four of the 5 peptides used as RT Standards are indicated with colored asterisks (*).

To test the ability of the "Dynamic Retention Time" software feature to adjust RT windows on-the-fly during a scheduled SRM method, the scheduled transition list used for Gradient 1 (GradO1 in Figures 3 and 4) was used for the other six LC gradients. RT windows were reduced to 1 minute except for transitions used as RT Standards, which were set to 5 minutes (see Figure 5).

	COI Global Paramet	scin Para	neters	Summar	· · · · · ·							
Hothed Timeles												
Pathod Duration		5.9		1.1				87	31.5	N	1	Les .
34								_			- 9+	Delete i Clear
AT	Openheed 1											
Name	NOT Falls System 1									5834 Properties		
	Compound	Actentian Time (min)	AT viedos (red	Nets	Present (mid	Product (mild	Collour Drimps (10)	Aehreist Treak *		-	Orongiography Paul 1989	dia.
	E BOATTITT BEAUXIE	10.0	8	Nettra	105.768	176.000	85.9	561			640	n.
Pel ScettO1	2 35A099999 (eevy)(+2	15.0	5	Postilie	41.32	\$72.338	18.9	3.5e4			Use Cycle Time	2
COLUMN TWO IS NOT	3 35449999 (max/y)(+2	2.1	3	Possie	45.58	675.391	2.5	3.044			Cude Tree (sec)	1.
	< 04E14GE000004(+3)	38.27		Pages	+04.547	344,254	24.6	3.564			Use Californial IP Lens	
Product Los Sean	< (ME)/200300(2)	10.27		Parties.	444 242	200.214	14.5	Lint			E CLEMENT FIRM	0.2
-	T THERE AND A AND A	73.47		Party of	000.000	100.00	24.5	1 feet				
Contraction of the local division of the loc	a realization hereit	20.00		inana.	105.812	100.00	25.8	Lies			E GIVENPEN Land	93
Restrict Cost Scat	1 GLUGDER Descrift	27.00	1	Nation	258.326	620.303	26.0	3.0+1			CD-Set (r/for)	1.5
	10 GL1601078 (www.)(+1	27.05	8	Feature	198.526	718.371	21.1	3.6e4			Source Programation-50	e .
LING:	11 012/00/078 (heavy)(+1	27.65	5	Postilier	158.326	832.495	21.4	5.0e4			in the second second	
10102		36.44	4	Castilia.	663.379	141.752	18.3	Lie1			Con consequences	-
	13 AMOLETIDEA(+2)	38.66	5	Pastine	967.975	431.341	18.3	5.044			Die Rotantium filme Rationance	× .
(40)	14 TROPOSAGAQOS (heavy)	22.41	5	Positive	695.832	340,412	25.0	1.044			Digilay Raterban Time	R
	15 APPOLIPTICICA(+2)	38.46	5	Postive	967.973	\$20.51	18.3	1.844			Use Quan Jan	122
	+ 16 STAAMERT Denvil(+2	15.8	1	Position	45.50	175.232	18.9	Natural			Stars Vauderbar	с.
	17 SEAMANNA Deevy (1-2	15.0	8	Pailing	+15.768	741.438	18.9	Netwool			CopyTes	minery free
	13 SEAMININ Deeve)(+2	12.0	1	Posting	413.38	812.405	44.9	Netwed				
	10 SEAMANNA (BENE)(12	18.8		Noise	*15.768	888.497	46.9	Text used				
	20 SSAAV99994(+2)	15.8	1	P00316	486.754	269.224	18.9	Natured				
	21 SGAAPPPWR(+2)	15.0	1	Postive	486.754	466.237	18.9	Netwood				
	22 \$54,0999998(+2)	25.8	4	Positive	401.554	\$65.33	18.9	Netwood				
	23 EEAAMONYM(+2)	21.8	4	Posters	488.754	445.393	18.5	Network				

Figure 5. View of Method Editor including the option for using Retention Time References. Colored asterisks indicate peptide sequence and transitions noted in Figure 4.

Dynamic Retention Time (RT) Adjustment

The software is designed to automatically update the scheduled SRM retention time window during the acquisition, based on the specified reference threshold. The primary goal of this feature is to accommodate the shifts in the retention time due to chromatographic changes that would result in missing the analyte of interest. The reference peak detection thresholds are specified in the acquisition methods (Figure 5). The identification of these reference peaks and corresponding adjustment of retention time windows happens automatically during acquisition. If the reference peaks are not detected, no adjustments are made.

When the RT windows in certain regions are reduced in duration to increase the average dwell time, there is a drawback of ending up with a narrow RT windows, which could result in missing the analyte if the RT shifts outside of the specified window. The use of dynamic RT adjustment allows for increased success of peptide detection, even when small (51 min) RT windows are employed.



Figure 6. Extracted Ion Chromatogram for peptide ISGLIYEETR (+2) to demonstrate the ability of Dynamic RT Adjustment to move the 1 min wide RT windows due to chromatographic shift. Conventional RT scheduling is shown on the left, with 3 min RT windows and manually updated relemtion time data to allow for successful delection of the peptide over different gradients. On the right, Dynamic Retention Time Adjustment was used so that the RT windows could be shortened to 1 min for most transitions, resulting in higher dwell times overall (See Figure 7). Gradients 04 and 05 had RT shifts too large for the RT Standard peptides to be detected with 5 min RT windows, so detection defaulted to the RT defined from Gradient 01.



Figure 7. SRM Visualization comparing conventionally scheduled SRM method (left panes) to Dynamic RT Adjustment scheduling (right panes). Not only does the average dwell time increase when using Dynamic RT Adjustment, but the number of concurrent transitions decreases, due to being able to use smaller RT mindows for scheduling.

Considerations for Dynamic Retention Time (RT) Adjustment

These experiments were conducted as a proof-of-concept to evaluate the Dynamic RT Adjustment feature. The RT windows for the RT Standards were set to 5 minutes, therefore any chromatographic changes causing a RT shift of greater than +/- 2.5 minutes for the RT Standards would result in no peak detected

Additional options for improving flexibility of the method would be to increase the RT windows of the RT Standard peptides, even monitoring for these peptides during the entire chromatographic run

In these experiments, RTs were empirically determined by running an unscheduled method to observe where each peptide eluted. Not only is this time consuming, but lower abundant peptides may not be reliably detected at dwell times near 1 msec. If the same RT Standard peptides are used in all samples, real-time RT updating of scheduled SRMs will ensure a higher success rate of analyte detection with smaller RT windows. The benefit of this approach is increased dwell times (see Figure 7).

CONCLUSIONS

Here we introduced two software features that provide clarity and flexibility to targeted peptide quantitation method design and experiments

SRM Visualization

- · Plots dwell time per transition and concurrent transitions as a function of retention time
- Provides interactivity, providing links between the SRM table and visualization plots
- Improves method design to minimize chances of too many overlapping transitions and/or dwell times that are too low

Dynamic Retention Time Adjustment

- Updates scheduled RT windows during a method if significant chromatographic shifts are observed
- Allows for much narrower RT scheduling windows to be used, which increases dwell times for most transitions
- Minimizes the need for manual data evaluation and re-scheduling of SRM methods
- Minimizes the incidence of "missing data" due to peaks shifting outside of their RT windows

REFERENCES

Escher, C., Reiter, L., MacLean., Ossola, R., Herzog, F., Chilton, J., MacCoss, M.J., Rinner, O. "Using IRT, a normalized retention time for more targeted measurement of peptides". *Proteomics* 2012, (12(8):1111-1121. doi: 10.1002/pmic.201100463.

www.thermofisher.com

©2016 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 Belaium +32 53 73 42 41 Brazil +55 11 2730 3006 Canada +1 800 530 8447 China 800 810 5118 (free call domestic) Italy +39 02 950 591 400 650 5118

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

Japan +81 6 6885 1213 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 Sweden +46 8 556 468 00 Switzerland +41 61 716 77 00 Taiwan +886 2 8751 6655 UK/Ireland +44 1442 233555 USA +1 800 532 4752 PN64750-EN 0616S



A Thermo Fisher Scientific Brand