

Ultra-robust micro-flow LC-MS/MS for targeted high-throughput peptide quantification using the Vanquish Neo UHPLC system

Authors

Stephan Meding, Alexander Boychenko, Runsheng Zheng, Christopher Pynn, Alec Valenta, Martin Samonig
Thermo Fisher Scientific, Germering, Germany

Keywords

micro-flow, cycle time, proteomics, screening, peptide quantification, direct injection, LC-MS, UHPLC, high-throughput

Goal

Demonstrate the robust performance of the Thermo Scientific™ Vanquish™ Neo UHPLC system for micro-flow LC-MS/MS based peptide quantification

Introduction

Peptide analysis with nano-flow gradient separations on long (≥ 25 cm) analytical columns coupled with HRAM (high-resolution accurate-mass) mass spectrometry is the standard for discovery-based “bottom-up” proteomics research. While delivering unsurpassed sensitivity and depth of analysis, nano-flow setups have yet to achieve the throughput and robustness required for large sample cohort analysis. Therefore, nano-flow LC-MS has been of limited use for translational proteomics, in particular for validation studies.

Increasing flow rates in combination with shorter gradients and shorter nano-flow columns (75 μm ID \times 15 cm length) can increase sample throughput.¹ However, higher flow rates result

in reduced electrospray ionization efficiency,² which are difficult to compensate for by larger injection amounts due to the limited sample loading capacity of nano-flow columns. Hence, the uptake of such methods for validation studies is limited.

Micro-flow LC can now overcome this limitation. Its inception dates back to the mid-1970s, but its potential is only now being fully realized. Modern mass spectrometers deliver high detection sensitivity at micro-flow flow rates; improved quality of UHPLC separation columns and packing materials results in better run-to-run and column-to-column reproducibility; and the latest generation of low-flow UHPLC systems, such as the Vanquish Neo UHPLC system, deliver outstanding injection and retention time reproducibility at micro-flow rates and system pressures above 1000 bar. The combination of these technological advances in mass spectrometry and liquid chromatography enables high-throughput proteomic screening of large sample cohorts.

In this Technical Note, a previously published high-throughput micro-flow method^{3,4} has been adapted to the new capabilities of the Vanquish Neo UHPLC system and rigorously tested for reproducibility and robustness. Fast sample loading and accelerated column washing and equilibration result in shorter cycle times without reduced performance or method robustness. With a flow rate of 50 $\mu\text{L}/\text{min}$ during the gradient phase, the method achieves sufficient detection sensitivity and sample throughput while the cycle time is <15 minutes.

Materials and methods

LC-MS grade solvents were used (Table 1). HeLa protein digest spiked with PRTC (Table 1) was diluted to 50 ng/μL and 25 fmol/μL with 200 μL water, 0.1% formic acid.

Table 1. Solvents and additives

Reagent	Grade	Supplier	Part number
HeLa Digest spiked with PRTC, 10 μg and 5 pmol	N/A	Thermo Scientific™ Pierce™	A47996
Acetonitrile with 0.1% formic acid	Optima™ LC-MS	Fisher Chemical	LS120-212
Isopropanol	Optima™ LC-MS	Fisher Chemical	A461-212
Formic acid	Optima™ LC-MS	Fisher Chemical	A117-50
Water with 0.1% formic acid	Optima™ LC-MS	Fisher Chemical	LS118-212
Water	Ultra-Pure, 18.2 MΩ at 25 °C	Thermo Scientific™ Barnstead™ GenPure™ xCAD Plus Ultrapure Water Purification System	

LC-MS analyses were performed on a Vanquish Neo UHPLC system (Table 2) connected to a Thermo Scientific™ TSQ Altis™ triple quadrupole mass spectrometer. For electrospray ionization a Thermo Scientific™ OptaMax™ NG source in HESI configuration was fitted with a 50 μm ID low-flow HESI needle (PN OPTON-30139). For reduced post-column volume the PEEK capillary connecting the grounding union to the HESI needle was replaced by a 50 μm × 150 mm Thermo Scientific™ nanoViper™ capillary (PN 6041.5124).

Table 2. Vanquish Neo UHPLC system with thermostatted column compartment

Module	Part number
Vanquish Neo UHPLC system (comprising Binary Pump N, Split Sampler NT, Solvent Rack, System base with drawer, Ship kit)	VN-S10-A-01
Vanquish Display (required)	6036.1180
Column Compartment N	VN-C10-A-01

The Vanquish Neo UHPLC system was set up in micro-flow direct injection mode with 50 μm ID nanoViper capillaries using the solvents listed in Table 3. A Thermo Scientific™ Acclaim™ PepMap™ 1.0 mm ID × 15 cm column (PN 164711) was placed in the column compartment. LC method details are described in Table 4 and Table 5. MS method details and SRM transitions are described in Table 6 and Table 7 and can be downloaded from Thermo Scientific™ [AppsLab Library of Analytical Applications](#).

The LC-MS system was operated with Thermo Scientific™ Chromeleon™ 7.2.10 MUd. Alternatively, Thermo Scientific™ Xcalibur™ with Thermo Scientific™ SII for Xcalibur™ 1.5.1 software can be used.

Table 3. Solvents of LC-MS method

Solvent	Composition
Eluent A	100% water, 0.1% formic acid
Eluent B	80% acetonitrile, 20% water (v/v), 0.1% formic acid
Weak Wash Liquid of Metering Device	100% water, 0.1% formic acid
Strong Wash Liquid of Metering Device	80% acetonitrile, 20% water (v/v), 0.1% formic acid
Weak (Outer) Needle Wash Liquid	100% water, 0.1% formic acid
Strong (Outer) Needle Wash Liquid	80% acetonitrile, 20% water (v/v), 0.1% formic acid

Table 4. Gradient of LC-MS method

Time (min)	Duration (min)	Flow rate (μL/min)	%B
Gradient separation phase			
0.0	0	100	1
0.2	0.2	100	8
2.2	2.0	50	13
10.7	8.5	50	35
Column wash phase			
10.9	0.2	50	99
11.2	0.3	50	99
11.4	0.2	100	99
12.4	1.0	100	99

Table 5. LC method parameters of LC-MS method

Parameter	Value
Pick-up and loading settings	
Keep Loop Inline	Disabled
Fast Loading	Disabled
Loading Volume	Automatic
Outer Needle Wash Mode	After Draw
Outer Needle Wash Time (Strong)	3.0 s
Outer Needle Wash Speed (Strong)	80.0 μ L/s
Outer Needle Wash Time (Weak)	5.0 s
Outer Needle Wash Speed (Weak)	80.0 μ L/s
Draw Speed	0.2 μ L/s
Draw Delay	2.0 s
Dispense Speed	5.0 μ L/s
Vial Bottom Detection	Enabled
Column equilibration settings	
Fast Equilibration	Disabled
Equilibration Factor	1.0
Temperature settings	
Column Compartment Temperature	50 °C
Autosampler Temperature	4 °C

Table 6. MS method parameters of LC-MS method

Parameter	Value
Source parameters	
Spray Voltage (static, positive)	3500 V
Sheath Gas	25 psig
Auxiliary Gas	5 psig
Sweep Gas	0 psig
Ion Transfer Tube Temperature	325 °C
Vaporizer Temperature	75 °C
Scan parameters	
Acquisition Mode	SRM (selected reaction monitoring)
Polarity	Positive
Cycle Time	0.5 s
Use Calibrated RF Lens	Enabled
Q1 Resolution (FWHM)	0.7 amu
Q3 Resolution (FWHM)	1.2 amu
CID Gas	1.5 mTorr
Source Fragmentation	0 V
Chromatographic Peak Width	6 s
Use Chrome Filter	Enabled

Table 7. Selected reaction monitoring (SRM) transitions of LC-MS method. Transition used for quantitation marked in bold. (Note: PRTC peptides are isotopically labeled. PRTC peptides marked with * were not used for method evaluation in this work due to their low signal response).

PRTC peptide	Start time (min)	End time (min)	Precursor (m/z)	Product (m/z)	Collision energy (V)
SSAAPPPPPR	0	12.4	493.768	476.286	18.9
				573.338	18.9
				670.391	18.9
				741.428	18.9
				812.465	18.9
GISNEGQNASIK	0	12.4	613.316	540.323	23
				725.403	23
				854.446	23
				968.489	23
				1055.521	23
*HVLTSIGEK	0	12.4	496.286	454.275	19
				541.307	19
				642.355	19
				755.439	19
				854.507	19
*DIPVPPPK	0	12.4	451.283	380.275	17.5
				477.327	17.5
				576.396	17.5
				673.449	17.5
				786.533	17.5
IGDYAGIK	0	12.4	422.736	366.194	16.6
				396.270	16.6
				559.333	16.6
				674.360	16.6
				731.381	16.6
*TASEFDSAIAQDK	0	12.4	695.832	582.334	25.8
				740.403	25.8
				855.430	25.8
				1002.498	25.8
				1218.573	25.8
SAAGAFGPPELSR	0	12.4	586.800	611.339	22.1
				668.360	22.1
				815.429	22.1
				886.466	22.1
				943.487	22.1
ELGQSGVDTYLQTK	0	12.4	773.895	660.381	28.5
				761.428	28.5
				876.455	28.5
				1032.545	28.5
				1119.577	28.5

Table 7. (continued)

PRTC peptide	Start time (min)	End time (min)	Precursor (m/z)	Product (m/z)	Collision energy (V)
GLILVGGYGTR	0	12.4	558.325	473.273	21.1
				506.260	21.1
				620.303	21.1
				719.371	21.1
				832.455	21.1
SFANQPLEVVYSK	0	12.4	745.392	603.359	27.5
				942.539	27.5
				1070.597	27.5
				1184.640	27.5
				1255.677	27.5
GILFVSGVSGGEEGAR	0	12.4	801.411	685.314	29.4
				772.346	29.4
				928.436	29.4
				1072.489	29.4
				1171.558	29.4
LTILEELR	0	12.4	498.801	427.254	19.1
				556.296	19.1
				669.381	19.1
				782.465	19.1
				883.512	19.1
NGFILDGFPR	0	12.4	573.302	486.270	21.6
				601.297	21.6
				714.381	21.6
				827.465	21.6
				974.533	21.6
ELASGLSFPVGFK	0	12.4	680.373	555.338	25.3
				702.406	25.3
				789.439	25.3
				959.544	25.3
				1046.576	25.3
LSSEAPALFQFDLK	0	12.4	787.421	658.365	28.9
				805.433	28.9
				918.517	28.9
				1086.607	28.9
				1157.644	28.9

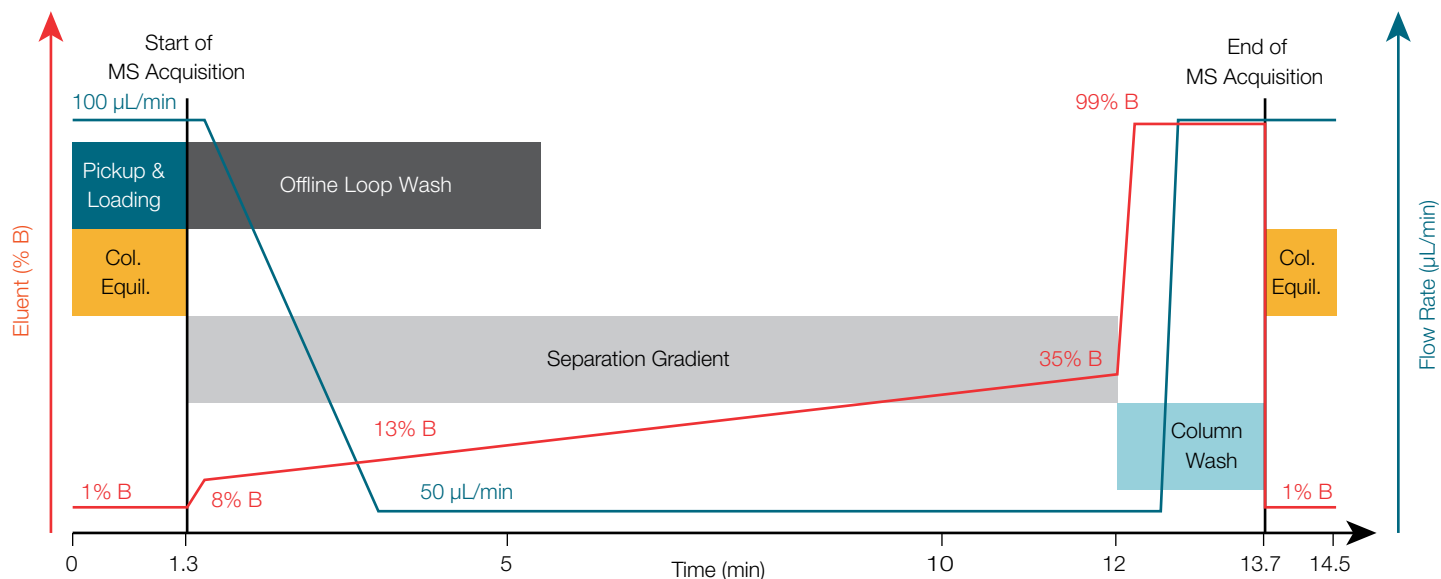


Figure 1. Scheme of the LC method. Both the gradient phase and additional method overhead including sample pickup, loading and column equilibration are depicted. Colored squares indicate the different phases of the run: the injection phase (includes sample aspiration and its loading onto the column), the sample loop wash phase, the column equilibration phase, the column wash phase, and the gradient separation phase. The red trace shows the percentage of eluent B being delivered by the analytical pump. Lowest value is 1% B, highest value is 99% B (Table 4). The blue trace shows the flow rate delivered by the analytical pump. Lowest value is 50 $\mu\text{L}/\text{min}$, highest value is 100 $\mu\text{L}/\text{min}$ (Table 4).

Results and discussion

A high-throughput micro-flow LC-MS method was developed based on previously published work.^{3,4} The goal was to create a highly robust, short method (ca. 14.4 min cycle time enabling 100 samples per day) which delivers good chromatographic separation and sufficient detection sensitivity for proteomic analysis of large sample cohorts. The gradient flow rate of 50 $\mu\text{L}/\text{min}$ was determined as the optimum to deliver both high throughput and high detection sensitivity. For a 1 mm ID column, larger injection amounts are sufficient to compensate for loss in ionization efficiency at micro-flow rate (50 $\mu\text{L}/\text{min}$). Sample loading onto the column, column washing, and column equilibration were performed at 100 $\mu\text{L}/\text{min}$ (Table 4 and 5, Figure 1). After sample loading the sample loop was switched offline to minimize the gradient delay volume. In addition, the flow rate at the start of the gradient was 100 $\mu\text{L}/\text{min}$ and then ramped down to 50 $\mu\text{L}/\text{min}$

within the first 2.2 minutes. This reduces the time until the first peptides are eluted and maximizes the elution window. The total cycle time is ca. 14.4 minutes. The duration of separation gradient phase is 10.7 minutes, or 74% of the cycle time.

To ensure minimal carry-over the sample loop was washed offline using the metering device with both high and low organic containing wash liquids.

The sensitivity of the method was assessed with 1 μL injections corresponding to 50 ng HeLa digest and 25 fmol of each PRTC peptide on column. LC-MS analysis with a TSQ Altis applying default source settings (Table 6) and generic SRM transitions for the PRTC peptides (Table 7) resulted in 12 quantifiable PRTC peptides (Figure 2). The other three PRTC peptides showed low signal intensity and were, therefore, omitted from further analyses.

FWHM \pm SD, sec

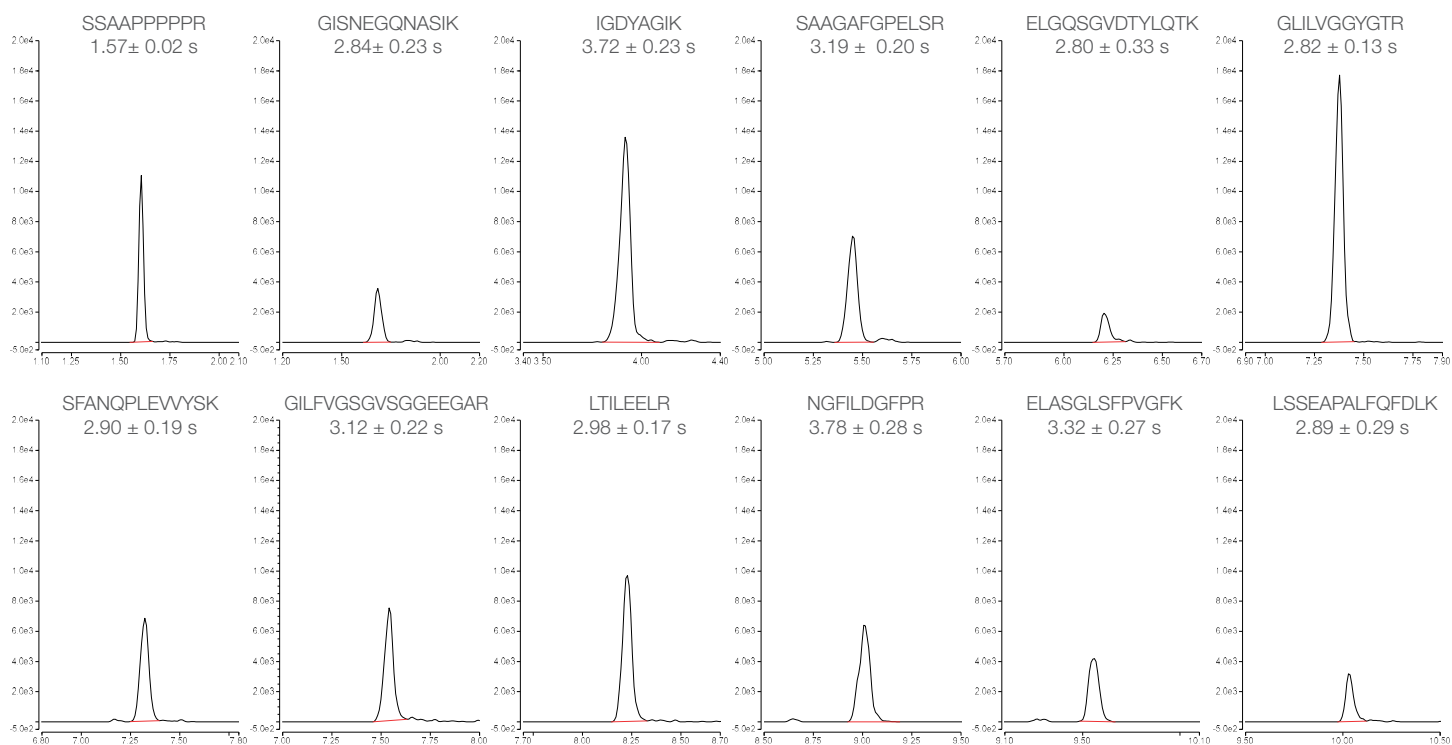


Figure 2: Representative chromatograms of 12 PRTC peptides used for method assessment. The peptide sequence is stated above each chromatogram. The average full width at half maximum (FWHM) and standard deviation over the test sequence is given below.

Since these 12 peptides were detected well above the limit of quantitation (signal-to-noise ratio = 10) they were used for method robustness evaluation. A sequence of 760 injections was run over a period of 7.5 days. The sequence was a repetition of the following injections: 3 sample injections, 3 blank injections, 3 matrix injections resulting in a total of 254 HeLa digest/PRTC injections. After the first 452 injections, sample vials containing fresh matrix and sample were exchanged because the 200 μ L (10 μ g) of matrix and sample were nearly depleted.

Method robustness was evaluated according to the consistency of the pressure traces as well as the retention time stability and peak area reproducibility of the 12 PRTC peptides.

The pressure traces were highly reproducible (Figure 3) with a run-to-run pressure variation of less than 3 bar observed throughout the sequence. It is important to note that the pressure consistency was not only present during the separation gradient phase. The pressure traces were also fully reproducible during the initial flow rate ramp down (0.2 to 2.2 min) and during the column wash phase where the flow rate is ramped up (Figure 3).

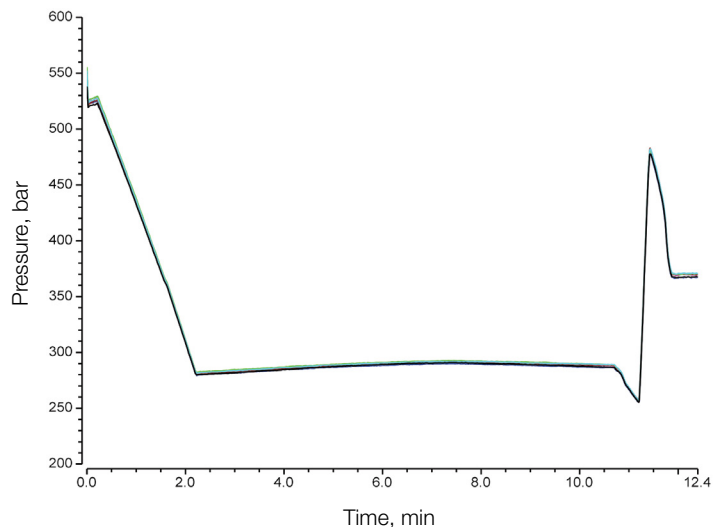


Figure 3. Pressure traces of early (black and purple), intermediate (blue and dark red) and late (green and light blue) runs of the sequence. The pressure traces are highly consistent with a minimal increase of less than 3 bar observed over the entire sequence. Data acquisition starts when the sample loop is switched offline after the sample is loaded onto the column. Therefore, the time scale between Figure 1 where the complete method is outlined and this figure where only the data acquisition during Separation Gradient and Column Wash phases is shown is shifted by 1.3 minutes.

The high congruency of pressure traces was also reflected in the observed retention time stability (Figure 4 and Table 8). The retention times were stable for any of the 12 peptides. Relative standard deviation was well below 0.5% for all peptides during seven days of system operation. Ten peptides had an RSD smaller than 0.1%, one had 0.14% RSD, and one had 0.31% RSD.

There was a minimal retention time reduction (≤ 3 s or $\leq 1\%$) for the first 20 injections (Figure 3). The initial retention time shift is most likely due to stationary phase conditioning. Afterwards, no further retention time changes were observed.

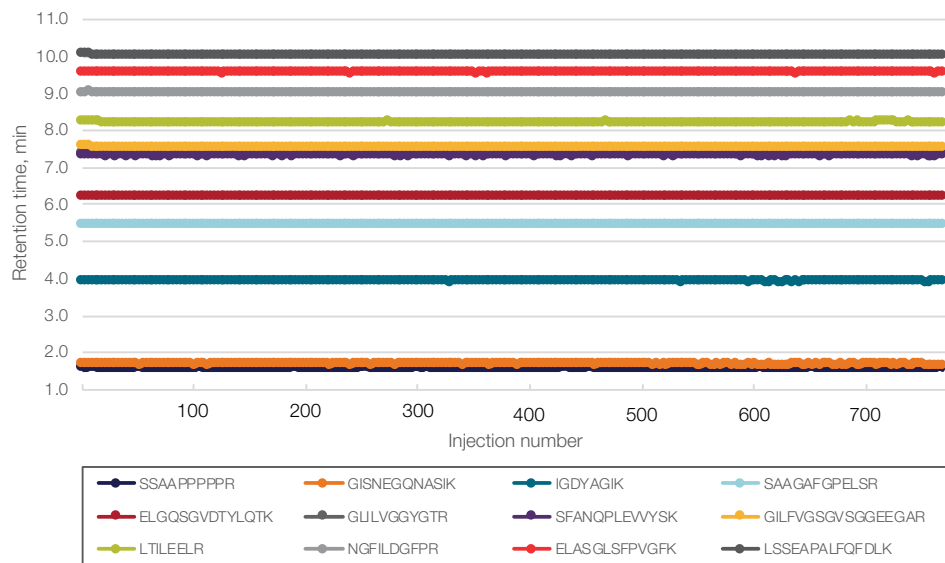


Figure 4. Retention time stability of 12 PRTC peptides for all sample injections of the sequence.

Table 8. Average retention time (RT), standard deviation (SD) and relative standard deviation (RSD, %) for 12 PRTC peptides

Peptide	SSAAPPPIPR	GISNEGQNASIK	IGDYAGIK	SAAGAFGPELSR	ELGQSGVDTYLQTK	GLILVGGYGTR	SFANQPLEVWYSK	GILFVGSGVSGGEEGAR	LTILEELR	NGFILDGFPR	ELASGLSFPVGFK	LSSEAPALFQFDLK
RT (min)	1.604	1.683	3.920	5.449	6.213	7.376	7.322	7.540	8.228	9.010	9.558	10.035
SD (min)	0.000	0.005	0.006	0.004	0.005	0.005	0.006	0.006	0.005	0.007	0.006	0.006
RSD (%)	0.01	0.31	0.14	0.08	0.08	0.07	0.08	0.08	0.06	0.08	0.06	0.06

Peak area reproducibility over the injection sequence was also assessed. As previously stated, a fresh vial of HeLa digest/PRTC was used after the first 452 injections due to sample depletion. Since slight differences in signal intensities between the two sample vials were observed, the peak area stability was considered separately for the first 452 injections and the subsequent 308 injections. A peak area reproducibility of <15% RSD is considered good for quantitative MS analyses.⁵ The observed peak area reproducibility was below 10% for 10 of the peptides. The two peptides with lowest peak areas showed the peak area variabilities of 11.1 and >15%.

In conclusion, a micro-flow LC-MS method on the Vanquish Neo UHPLC system with a cycle time of ca. 14.4 minutes was developed and evaluated. The method delivered highly robust and reproducible results for a sequence of 760 injections over a period of more than one week and did not reveal any LC, MS or column performance declines. Therefore, this method would be suitable for high-throughput proteomic analyses of large sample cohorts needed for validation studies.

Table 9. Average peak area, standard deviation (SD) and relative standard deviation (RSD) for 12 PRTC peptides for injections 1–452 and 453–760

	Peptide	SSAAPPPIPPR	GISNEGQNASIK	IGDYAGIK	SAAGAFPELSR	ELGQSGVDTYLQTK	GLILVGGYTR	SFANQPLEWYSK	GILFVGSVSGGEGAR	LTILEELR	NGFILDGFPR	ELASGLSFPVGFK	LSSEAPALFQFDLK
Injections 1–452	Peak Area (counts * min)	346.8	174.2	925.6	459.5	101.2	881.5	340.8	381.0	478.8	393.3	245.2	136.5
	SD (counts * min)	27.9	15.6	65.6	31.0	11.2	53.9	19.8	27.2	36.9	28.4	22.7	21.2
	RSD (%)	8.0	8.9	7.1	6.7	11.1	6.1	5.8	7.1	7.7	7.2	9.2	15.5
Injections 453–760	Peak Area (counts * min)	412.3	204.8	1,155.4	562.3	119.2	1,148.2	405.0	462.6	605.5	506.5	304.9	168.3
	SD (counts * min)	27.9	15.6	65.6	31.0	11.2	53.9	19.8	27.2	36.9	28.4	22.7	21.2
	RSD (%)	6.8	7.6	5.7	5.5	9.4	4.7	4.9	5.9	6.1	5.6	7.4	12.6

References

1. Boychenko, A., Pynn, C., et al. Technical Note 73208 - Tailored high-throughput low-flow LC-MS methods for large sample cohort analysis. *Thermo Scientific* 2019.
2. Meding, S. and Boychenko, A. Capillary Flow LC-MS Unites Sensitivity and Throughput. *Chromatography Today*, 2016, 3, 43-45.
3. Bian, Y., Zheng, R., et al. Robust, reproducible and quantitative analysis of thousands of proteomes by micro-flow LC-MS/MS. *Nat. Comm.* 2020, 11, 157.
4. Bian, Y., Bayer, F. P. et al. Robust Microflow LC-MS/MS for Proteome Analysis: 38000 Runs and Counting. *Anal. Chem.* 2021, 93, 8, 3686-3690.
5. European Medicines Agency. Guideline on bioanalytical method validation. *EMA/CHMP/EWP/192217/2009 Rev. 1 Corr. 2***, published 21 July 2011.

Learn more at thermofisher.com/vanquishneo