

SOLA μ for reduced sample volume analysis

Jon Bardsley, Ken Meadows, Thermo Fisher Scientific, Runcorn, UK

Key Words

Micro elution, reproducibility, matrix effects, SPE, no dry down, niflumic acid, reduced sample volume

Abstract

This application note demonstrates how the Thermo Scientific™ SOLA μ ™ SPE product can be used to facilitate the scale down of an extraction method for use when sample volume is limited. The use of a Thermo Scientific™ Accucore™ HPLC column provided fast and efficient separation without the need for an ultra high pressure system. MS/MS detection was performed on a Thermo Scientific™ TSQ Vantage™ mass spectrometer.

Introduction

Ethical, analytical and sample availability considerations are a challenge faced by many bioanalytical laboratories and have resulted in a drive to limit sample volume.

In order to achieve the required detection limits many analytical methods utilize dry down and reconstitution steps to remove the dilution effects required by traditional scale SPE when operating with very low sample volumes.

In addition many analytes, such as small volatile molecules or larger biomolecules, suffer from loss of recovery attributed to the evaporation and reconstitution step.

Thermo Scientific™ SOLA μ ™ allows users to directly scale down the volumes used in their analytical methods, allowing for a reduction in sample usage and eliminating issues caused by evaporation without compromising the sensitivity of their assay.

SOLA μ provides reproducibility, robustness and ease of use at low elution volumes by utilizing the revolutionary Thermo Scientific SOLA μ , Solid Phase Extraction (SPE) technology. This removes the need for frits delivering a robust, reproducible format which ensures highly consistent results at low elution volumes.



SOLA μ delivers:

- Lower sample failures due to high reproducibility at low elution volumes
- Increased sensitivity due to lower elution volumes
- The ability to process samples which are limited in volume
- Improved stability of bio-molecules by reduction of adsorption and solvation issues

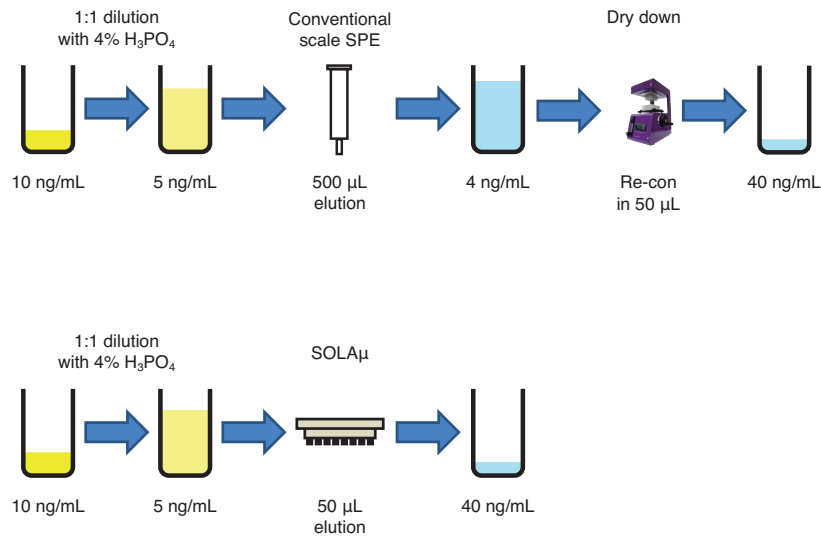


Figure 1: Summary of workflow required to achieve a ten-fold decrease in sample volume without altering the work-flow or compromising the final extracted analyte concentration

Experimental Details

Consumables		Part Number
Fisher Scientific™ LCMS grade water		10777404
Fisher Scientific™ LCMS grade methanol		10653963
Fisher Scientific™ Analytical grade formic acid		10559570
Sample Handling Equipment		Part Number
Liquid handling hardware:		
SPE hardware:	Thermo Scientific™ HyperSep™ 96 vacuum manifold	60103-351
	Vacuum pump, european version	60104-241
Sample handling:	Thermo Scientific™ 96 well square well microplate	60180-P212
	Thermo Scientific™ WebSeal™ mat	60180-M122
Sample Pretreatment		
25 µL of human plasma diluted 1:1 with 4% phosphoric acid		
Sample Preparation		Part Number
Compound(s):	Niflumic acid, niflumic acid d5 (IS)	
Matrix:	Human plasma	
	SOLAµ WAX 2 mg/1 mL 96 well plate	60209-005
Condition:	200 µL methanol	
Equilibrate:	200 µL water	
Load:	Apply sample at 0.5 mL/min	
Wash:	200 µL 25 mM ammonium acetate (pH4)	
	200 µL methanol	
Elute:	2 × 12.5 µL methanol with 2% ammonia	
Direct injection of eluent		

Separation Conditions		Part Number
Instrumentation:	Thermo Scientific™ Dionex™ UltiMate™ 3000 RSLC system	
Column:	Thermo Scientific™ Accucore™ RP-MS HPLC column 50 mm × 2.1 mm 2.6 μm	17626-052130
Guard column:	Thermo Scientific™ Accucore™ RP-MS Defender™ guard cartridge	17626-012105
	Thermo Scientific™ Uniguard™ drop-in guard holder	852-00
Flow rate:	750 μL/min	
Run time:	4 min	
Column temperature:	30 °C	
Injection details:	2 μL full loop injection	
Injection wash solvent 1:	Water	
Injection wash solvent 2:	45:45:10 (v/v/v) propan-2-ol / acetonitrile / acetone (with 5% Ammonia)	
Mobile phase A:	Water with 0.1% formic acid	
Mobile phase B:	Acetonitrile with 0.1% formic acid	

Gradient Conditions

Time (min)	%A	%B
0.0	70	30
2.0	10	90
2.01	70	30
3	70	30

MS Conditions

Instrumentation:	Thermo Scientific™ TSQ Vantage™ triple stage quadruple mass spec
Ionization conditions:	HESI
Polarity:	+ive
Spray voltage (V):	3000
Vaporiser temperature (°C):	475
Sheath gas pressure (Arb):	50
Aux gas pressure (Arb):	60
Capillary temp (°C)	300
Collision pressure (mTorr):	1.5
Scan time (s):	0.02
Q1 (FWHM):	0.7
Q3 (FWHM):	0.7

Compound	Parent (m/z)	S-Lens (V)	Product (m/z)	Collision Energy (V)
Niflumic Acid	283.0	115	265.0	22
Niflumic Acid d5 (IS)	288.8	115	271.1	22

Data processing

Software:	Thermo Scientific™ LC QUAN™ version 2.6 quantitative software
-----------	---

Results

By loading 25 μL of sample onto the SOLA μ plate and eluting in a total of 25 μL a ten-fold decrease in sample volume was achieved when compared to a traditional scale higher bed weight product. The results demonstrate that even with this low elution volume equivalent method performance was achieved.

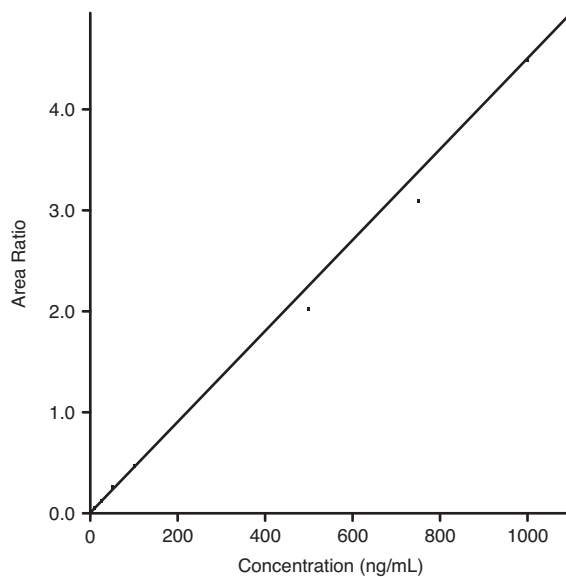


Figure 2: Niflumic acid linearity over the dynamic range 1-1000 ng/mL

Standard	Specified Concentration (ng/mL)	Calculated Concentration (ng/mL)	Accuracy (% difference)	Precision (%RSD n=18)
S1	1.00	0.995	-0.450	-
S2	10.0	10.0	-0.0200	-
S3	25.0	26.4	5.69	-
S4	50.0	55.9	11.9	-
S5	100	102	2.22	-
S6	500	448	-10.4	-
S7	750	685	-8.63	-
S8	1000	997	-0.330	-
QC L	10.0	9.73	2.66	0.355
QC M	500	440	11.9	0.142
QC H	750	671	10.5	0.195

Table 1: Niflumic acid accuracy data for the calibration range 1 to 1000 ng/mL

The assay gave a linear dynamic range from 1 to 1000 ng/mL with an r^2 coefficient of 0.995 (Figure 2, Table 1). The chromatography for the limit of quantitation sample at 1 ng/mL is significantly above the acceptable signal to noise limit (Figure 3).

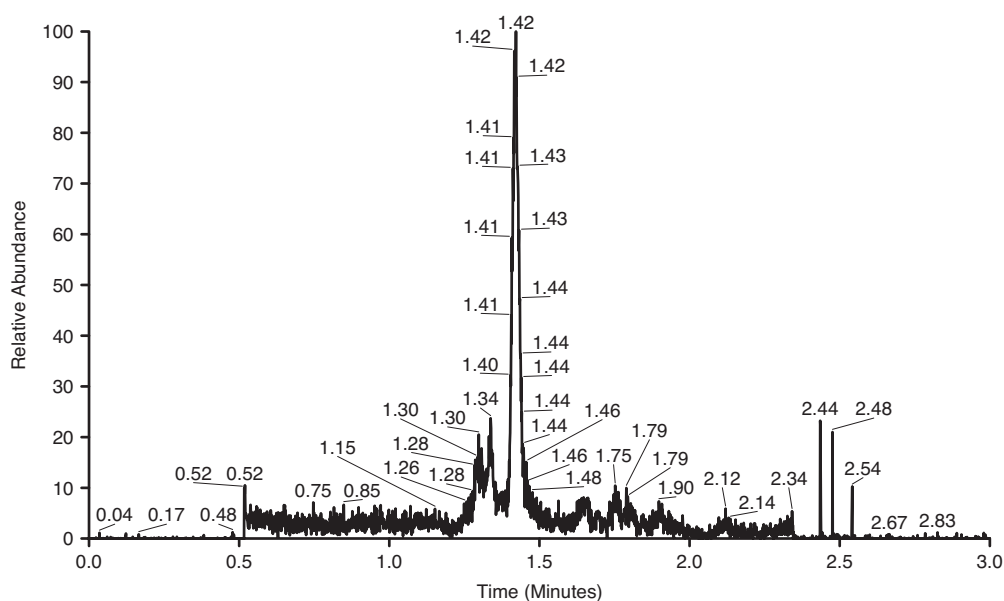


Figure 3: Example chromatogram 1 ng/mL niflumic acid

Low, mid and high QC samples were prepared at concentrations of 10, 500 and 750 ng/mL respectively. Table 1 shows good level of accuracy at all QC levels. Table 2 shows reproducibility data for replicate extractions (n=18) at both high and low QC levels. Relative standard deviation for niflumic acid was less than 8%.

	Precision Data for Niflumic Acid	
	Analyte Peak Area (%RSD)	Peak Area Ratio (%RSD)
Low QC	7.32	0.356
High QC	5.33	0.195

Table 2: Precision data niflumic acid at low QC 10 ng/mL and high QC 750 ng/mL (n=18)

	Recovery of Niflumic Acid (%)
Low QC	101
High QC	98.5

Table 3: Percentage recovery for niflumic acid at low QC 10 ng/mL and high QC 750 ng/mL

	Matrix Effects (%)
Low QC	1.70
High QC	4.78

Table 4: Percentage matrix effects for niflumic acid at low QC 10 ng/mL and high QC 750 ng/mL

Analyte recovery was shown to be greater than 98% by comparison to post extraction fortified blank samples (refer to Table 3). Post extraction fortified blank samples were also compared against pure reference standards to demonstrate matrix effects which were calculated at less than 5% at both high and Low QC levels (refer to Table 4).

A mid QC sample (500 ng/mL) was extracted using SOLA WAX and SOLA_μ WAX. Sample volume was reduced ten fold from 250 μ L to 25 μ L respectively as outlined in Figure 1. A high level of comparison is shown in Figure 4 between the extracts, demonstrating the ability to directly scale volumes with no adverse effect on assay performance.

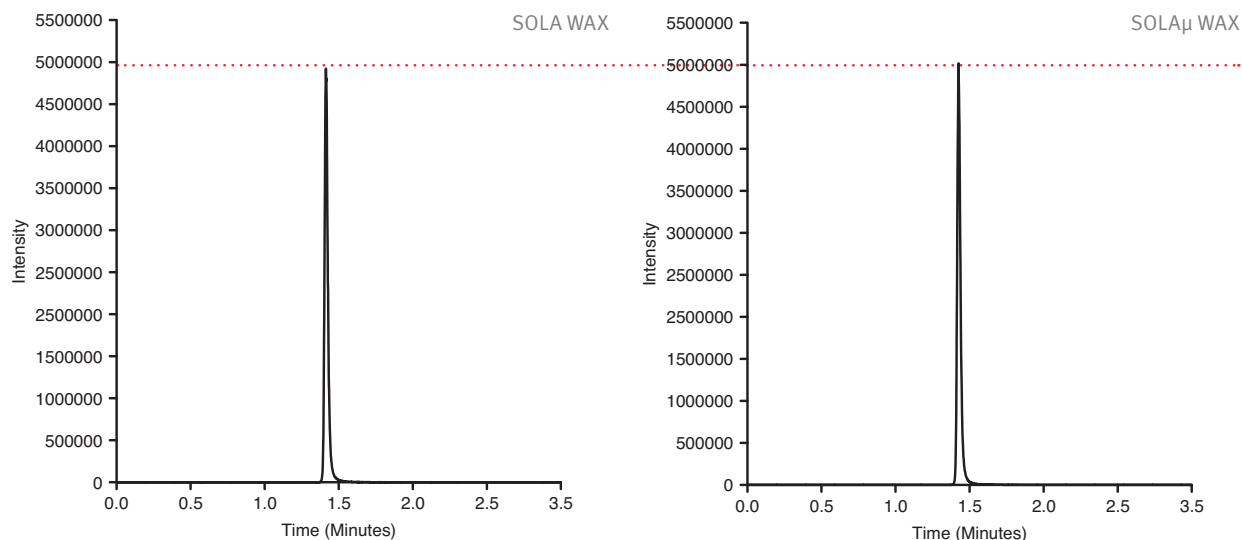


Figure 4: Comparison of niflumic acid extracted with 10 mg SOLA WAX using 250 μ L of sample and from SOLA μ WAX using 25 μ L of sample. Both methods show equivalent assay performance

Conclusion

This application note demonstrates the advantages of SOLA μ for the reduction of sample volume compared to traditional scale SPE while maintaining high levels of precision, accuracy, recovery and sample cleanliness.

By eluting in less solvent it is possible to achieve equivalent limits of detection requiring only a fraction of the sample volume and without relying on lengthy evaporation procedures that may compromise sample integrity. This is particularly important for sample limited applications.

thermoscientific.com/sola-spe

© 2014 Thermo Fisher Scientific Inc. All rights reserved. All 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.

USA and Canada +1 800 332 3331
Australia 1300 735 292 (free call domestic)
China 800 810 5118 (free call domestic)
 400 650 5118
France +33 (0)1 60 92 48 34
Germany +49 (0) 2423 9431 20 or 21

India +91 22 6742 9494
 +91 27 1766 2352
Japan 0120 753 670 (free call domestic)
 0120 753 671 fax
United Kingdom +44 (0) 1928 534 110
New Zealand 0800 933 966 (free call domestic)
Singapore +65 6289 1190
All Other Enquiries +44 (0) 1928 534 050

Technical Support
 For advice and support,
 please visit our website:
www.thermoscientific.com/chromexpert

Thermo
 SCIENTIFIC

A Thermo Fisher Scientific Brand