

LC-MS/MS Quantitative Analysis of the Vitamin K's and their Metabolites in Serum for Research Use

Authors

Rory M Doyle and
Alexander Cherkassky
Thermo Fisher Scientific,
Somerset, New Jersey, USA

ABSTRACT

Purpose: An LC-MS/MS analytical method was developed and verified for the quantitation of the Vitamin K's in serum for research use. Simple sample preparation techniques including protein crash and liquid-liquid extraction were evaluated. A Thermo Scientific™ TSQ Endura triple quadrupole mass spectrometer in positive Electrospray mode with a Thermo Scientific™ Dionex™ Vanquish™ Horizon HPLC system was used.

Methods: 200 μ L (PPX) and 500 μ L (LLE) of serum were used for the analysis of the Vitamin K's and their metabolites. Various columns were evaluated and a Thermo Scientific™ Accucore™ PFP, 100 x 2.1 mm, 2.6 μ m with 0.1% Formic Acid and 5mM Ammonium Formate in water and 0.1% Formic Acid in methanol mobile phases achieved baseline chromatographic separation in approximately 4 minutes run time. Quantitative analysis was performed using scheduled reactive monitoring (SRM) transition pairs for each Vitamin K and internal standard in positive mode and accuracy of the analytical method was verified using pooled reference samples.

Results: Good linearity and reproducibility were obtained across the dynamic range of the Vitamin K's with a coefficient of determination $R^2 > 0.95$ or better for all compounds in the various matrices. The limits of detection and quantitation were determined to range initially from 50 pg/ml to 1 ng/ml with very good reproducibility observed for all compounds.

INTRODUCTION

The Vitamin K's are a group of structurally similar fat soluble compounds that the body requires in order to maintain good physiological function. There are two natural forms - K1 and K2 of which K2 consists of a number of sub types that differ in the length of their carbon side chains and each has a significant influence on the body.

In this case, we evaluated various columns and solvent combinations as well as simple and easy sample preparation techniques in order to develop an LC-MS/MS analytical method that can demonstrate the chromatographic separation, detection and quantification of the Vitamin K's and their metabolites in serum. The Vitamin K's analyzed include Vitamin K1 (Phylloquinone), Vitamin K1-2,3-Epoxyde, Vitamin K2-MK4 (Menaquinone-4), Vitamin K2-MK7 (Menaquinone-7), Vitamin K2-2,3-Epoxyde and Vitamin K3 (Menadiolone). The sample preparation choices were kept simple and included protein crash and a one step liquid-liquid extraction. The methodologies were developed on a TSQ Endura triple quadrupole mass spectrometer in positive Electrospray ionization modes with a Vanquish Horizon HPLC system with a 4 minute analytical gradient.

MATERIALS AND METHODS

Standards

The following analytical reference standards and Internal standards were obtained from -

Isosciences, Inc., King of Prussia, PA

Vitamin K1 (Phylloquinone):	1 mg	Vitamin K1-D7:	1 mg
Vitamin K1-2,3-Epoxyde:	1 mg		
Vitamin K3 (Menadione):	1 mg	Vitamin K3-D8:	1 mg

Cerilliant Corp., Round Rock, TX-

Vitamin K2-MK-4 (Menaquinone-4):	1mg
Vitamin K2-MK-7 (Menaquinone-7):	1 mg

Sigma-Aldrich Corp. St Louis, MO-

Vitamin K2-2,3-Epoxyde:	10 mg
-------------------------	-------

Reagents

The following Fisher Scientific™ acids, reagents and solvents were used

HPLC grade Water	Ethanol
Methanol	Hexane
Ammonium Formate	Formic Acid

Sample Preparation- Protein Crash

- 200 µL of Serum/BSA mixture calibrators, controls and serum sample were added to 1.5 ml eppendorf tubes and 20 µL of Vitamin K ISTD at 1000 ng/mL were added to each tube and vortexed briefly
- 400 µL of Acetonitrile was added to each tube and vortexed for 1 min prior to centrifugation for 10 minutes at 13000 rpm
- The supernatant was transferred to an MS vial and capped.
- All in-house calibrators were prepared in drug-free serum and bovine serum albumin mixture (Golden West Biological, Inc, Temecula, CA)

Sample Preparation- Liquid-Liquid Extraction

- 500 µL of Serum/BSA mixture calibrators, controls and serum samples were added to a test tube and 50 µL of Vitamin K ISTD at 1000 ng/mL were added to each and vortexed briefly
- 1.5 mL of Ethanol was added to each tube and vortexed for 1 min followed by 4 ml of Hexane and again vortexed for 1 min.
- The tubes were centrifugation for 10 minutes at 13000 rpm
- The upper organic layer was transferred to a new test tube and dried down under nitrogen at room temperature
- The extract was reconstituted in 200 µL of 1:3 water and methanol
- The supernatant was transferred to an MS vial and capped.

The calibration curves ranged from 0.1 ng/mL to 1000 ng/mL and various pooled samples were used as control material.

Data Analysis

The software used included for this method included the Thermo Scientific™ Xcalibur™ 3.1 SW, Thermo Scientific™ TSQ Endura Tune™ 2.1 SW, and Thermo Scientific™ Tracefinder™ 4.1 SW

Method

HPLC Conditions-

Vanquish Horizon HPLC binary pump, well plate, thermostatted column compartment

Column:	Accucore PFP, 100 x 2.1 mm, 2.6 µm
Column Temperature:	50 °C
Injection Volume:	20 µL
Sampler Temperature:	4 °C
Needle Wash:	Flush port (50%Methanol:50%Water) 10 seconds
Mobile Phase A:	0.1% Formic Acid + 5mM Ammonium Formate in Water
Mobile Phase B:	0.1% Formic Acid in Methanol
Flow Rate:	0.5 ml/min
Gradient:	0 min- 30%A:70%B 1.0 min- 2%A:98%B 3.0 min- 2%A:98%B 3.1 min- 30%A:70%B
Run time:	40 mins

Method

HPLC Conditions-

Vanquish Horizon HPLC binary pump, well plate, thermostatted column compartment

Column: Accucore PFP, 100 x 2.1 mm, 2.6 μ m
 Column Temperature: 50 °C
 Injection Volume: 20 μ L
 Sampler Temperature: 4 °C
 Needle Wash: Flush port (50%Methanol:50%Water) 10 seconds
 Mobile Phase A: 0.1% Formic Acid + 5mM Ammonium Formate in Water
 Mobile Phase B: 0.1% Formic Acid in Methanol
 Flow Rate: 0.5 ml/min
 Gradient:
 0 min- 30%A:70%B
 1.0 min- 2%A:98%B
 3.0 min- 2%A:98%B
 3.1 min- 30%A:70%B
 Run time: 40 mins

MS and Ion Source Conditions-

TSQ Endura triple quadrupole mass spectrometer

Ion mode: Positive Electrospray (H-ESI) Mode
 Vaporizer Temperature: 325 °C
 Ion Transfer Tube Temperature: 325 °C
 Sheath Gas: 45
 Aux Gas: 25
 Sweep Gas: 15
 Spray Voltage: Positive Ion (V):3500 V
 Q1/Q2 Resolution: 1.2/0.7 (FWHM)
 Cycle time (sec): 0.5
 CID Gas (mTorr): 2
 Chromatographic Peak Width: 6 secs

Table 1. Scan Parameters- SRM Table

Compound	RT (Min)	Polarity	Precursor (m/z)	Product (m/z)	Collision Energies (V)	RF Lens (V)
Vitamin K1	2.59	Positive	451.1	187.1/185.1	23/21	147
Vitamin K1-D7	2.59	Positive	458.3	193.8	23	147
Vitamin K1-2,3-Epoxyde	2.4	Positive	467.2	449.6/161.1	10/17	121
Vitamin K2-MK-4	2.31	Positive	445.1	186.9/227.1	20/15	122
Vitamin K2-MK-7	2.95	Positive	650.3	186.9/94.9	26/31	184
Vitamin K2-MK-4-2,3-Epoxyde	2.19	Positive	461.2	443.1/310.9	10/10	112
Vitamin K3	0.93	Positive	172.9	104.9/77.1	19/33	89
Vitamin K3-D8	0.92	Positive	181.3	108.9/81.3	20/24	99

RESULTS

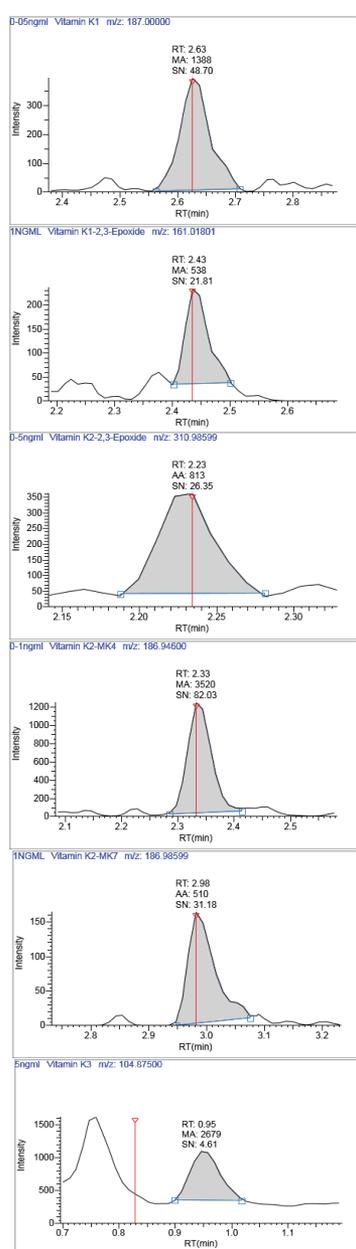
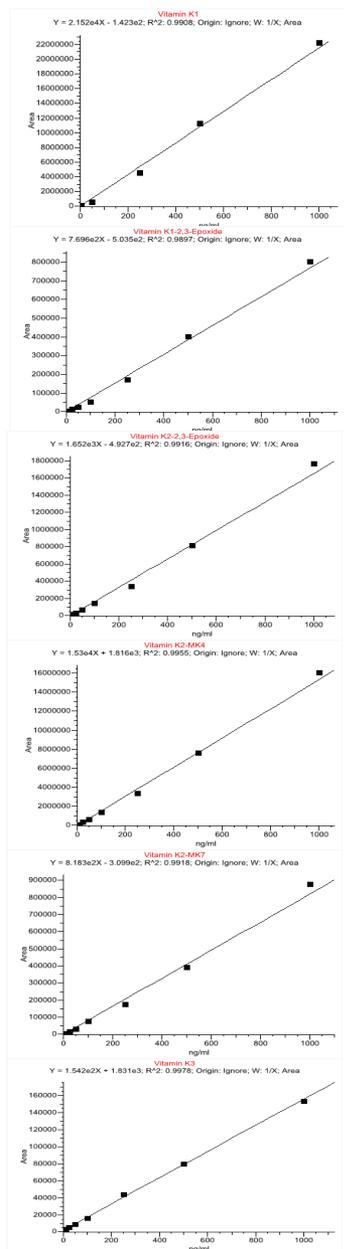
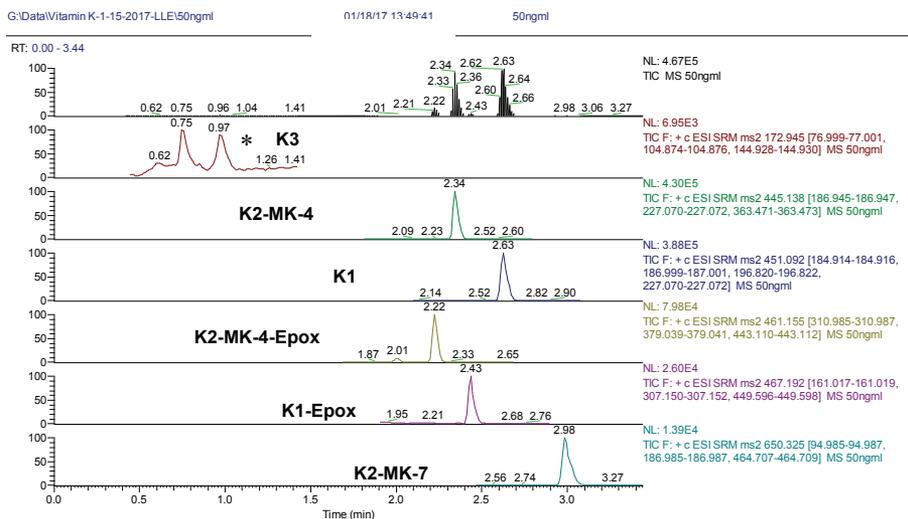
Linearity/Sensitivity

The linear range of the Vitamin K's in serum/BSA matrix was from 100 pg/ml to 1000 ng/ml. The linearity of each extraction was determined in triplicate over 3 days and the results are shown with LOD and LOQ being determined as 3:1 and 10:1 of signal to noise respectively where possible and the mean coefficient of determination (R^2) > 0.98 for each matrix and the %CV for each calibration point were all <10%.

Table 2. Sensitivity.

Compound	Protein Crash LOD/LOQ (pg/ml)	LLE LOD/LOQ (ng/ml)
Vitamin K1	0.25/0.5	0.05/0.1
Vitamin K1-2,3-Epoxyde	2.5/5	1/ 2.5
Vitamin K2-MK-4	0.1/0.25	0.1/0.25
Vitamin K2-MK-7	10/25-Inter	1/ 2.5
Vitamin K2-MK4-2,3-Epoxyde	1/ 2.5	0.5/1
Vitamin K3	100-Inter	5/10

Figure 1: LLE Chromatograms, Curves and LOD



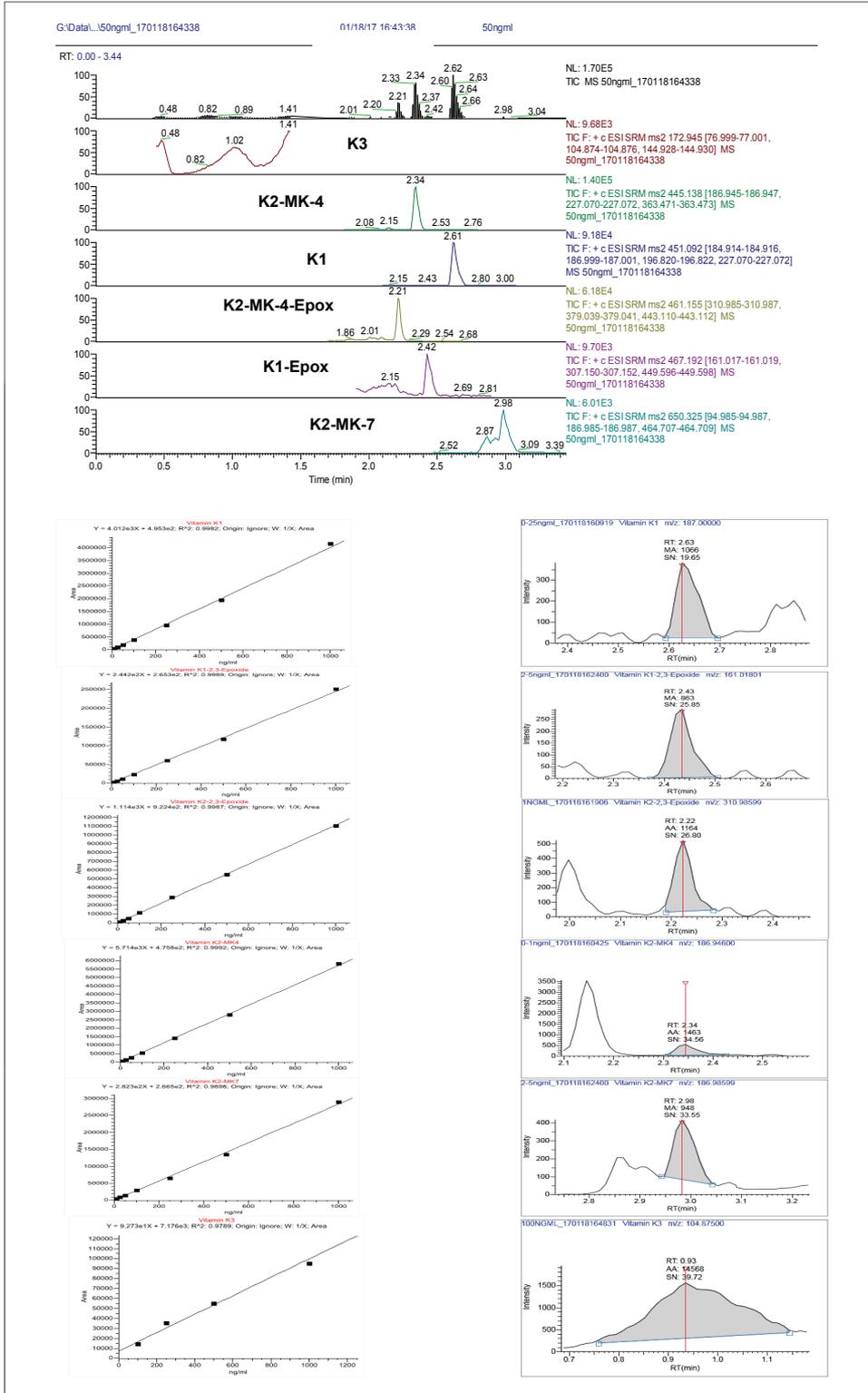
Accuracy

The accuracy was determined by the analysis of pooled sample control material as the percentage deviation from the targeted mean and the results were <10% for all levels. The serum pooled control material concentrations were 25 ng/ml and 250 ng/ml. Therefore, the analytical method can achieve the research laboratory required accuracy for the analysis of the Vitamin K's in serum.

Precision/Specificity

The intra-assay precision (%CV) of the Vitamin K's in serum were determined by extracting and quantifying three replicates of the pooled sample control material. The inter-assay precision was determined over 3 consecutive days and was found to have a %CV <10% for each Vitamin K within their respective linear range for the pooled serum sample control material respectively. Therefore, the analytical method can achieve the required precision for the analysis of the Vitamin K's in serum. Due to the similarity between the Vitamin K's and that it is present in all biological matrices, there were interferences and ion suppression from endogenous lipids which shown in the protein crash chromatogram. Vitamin K's are very sensitive to light and the nitrogen must be passed through steel tubing to avoid loss of compounds.

Figure 2: PPX Chromatograms, Curves and LOD



CONCLUSIONS

- Baseline separation of the Vitamin K's with good LOD/LOQ was achieved in serum particularly by LLE
- Simple sample preparation achieved desirable LOD/LOQ to the relevant levels with further work to be carried out to fine tune these techniques to obtain more sensitive results and to remove the interferences and achieve better sensitivity while maintaining ease of use and low cost
- Good linearity of calibration curves with acceptable accuracy, precision and reproducibility in positive and negative mode was achieved <10% for %CV for each Vitamin K particularly by LLE within their linear range and the sample preparation techniques and analytical methodologies will be further verified and optimized

REFERENCES

1. A Validated HPLC Method for the Determination of Vitamin K in Human Serum – First Application in a Pharmacological Study
The Open Clinical Chemistry Journal, 2011, 4, 17-27
Milka Maranova et al
2. Method for the Determination of Vitamin K Homologues in Human Plasma Using High-Performance Liquid Chromatography-Tandem Mass Spectrometry
Anal. Chem., 2005, 77 (3), pp 757–763
Yoshitomo Suhara, et al

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

Find out more at thermofisher.com

ThermoFisher
SCIENTIFIC