



Fully Automated Sample Preparation for the Determination of Vitamin D on Biomek i7 Hybrid Workstation using Advanced Sample Cleanup with Phree Phospholipid Removal Cartridges

Kerstin Thurow¹, Heidi Fleischer¹, Anna Bach¹, Bhagya Wijayawardena², Miranda Kheradmand²
¹Center for Life Science Automation, University Rostock, Rostock, Germany
²Beckman Coulter Life Sciences, Indianapolis, Indiana, United States

Abstract

Vitamin D is a crucial component of metabolism. Vitamin D deficiency leads to several disease conditions such as rickets and osteomalacia. Therefore, researchers are interested in identification of the role of vitamin D in treating and prevention of metabolic abnormalities. To do so, they often quantify the level of vitamin D in serum using LC-MS based methods. The sample preparation for these analytical methods are labor-heavy and prone to human error. In this application note we evaluate the feasibility of automating the sample prep for vitamin D determination on the Biomek i7 hybrid workstation. We used Phree phospholipid removal cartridges for the protein precipitation prior the sample cleanup by solid phase extraction.

Introduction

The quantitative determination of 25(OH)D₂ and 25(OH)D₃ in blood plasma and blood serum is becoming increasingly important to understand the effects of overdosing and underdosing of vitamins. Due to the complicated matrix, the sample preparation for vitamin D analysis is extensive, including protein precipitation and extraction steps. The protein precipitation and the separation of the metabolites from the vitamin D-binding proteins is usually carried out by adding acetonitrile, methanol, mixtures of acetonitrile and methanol or 2-propanol followed by a centrifugation step [1]. In addition to the proteins, there are other substances in biological matrices, in particular phospholipids (components of the cell membrane), whose highly ionic nature interferes with the ionization in the mass spectrometer [2]. In addition, phospholipids can cause accelerated column breakdown through enrichment in the mobile phase, which leads to changes in the separation properties and an increase in the baseline noise [3]. Therefore, the removal of the phospholipids is an important step in the sample preparation. The main methods of removing phospholipids are solid phase extraction (SPE) with strong cation exchangers, SPE and liquid-liquid extraction (LLE). However, performing these methods increases the time it takes to complete the analytical measurement process.

To reduce sample preparation time, different manufacturers have developed plates that combine the advantages of protein precipitation, removal of phospholipids and extraction, so that only a single plate or cartridge can be used [4]. In this application note we automated a sample preparation method for vitamin D determination that uses Phree phospholipid removal cartridges and solid phase extraction.

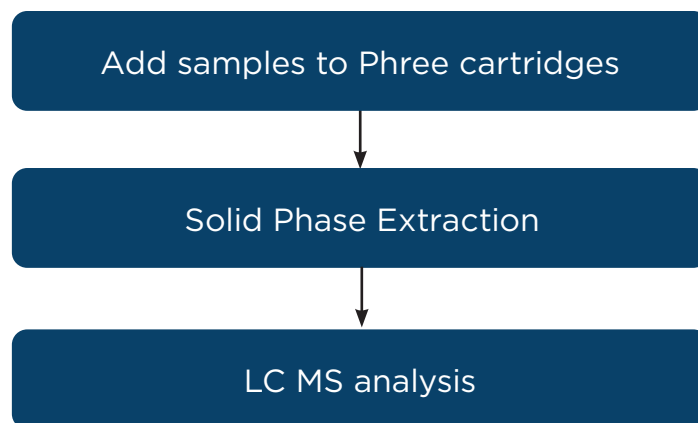


Figure 1. The sample preparation workflow. All steps except LC-MS analysis was done on a Biomek i7 hybrid workstation with integrated devices.

Materials and Methods

The automated method was based on a published vendor protocol on the usage of Phree Phospholipid Removal Cartridges (Phenomenex, Torrance, US). Pig blood serum was used for the method development and evaluation (Source: Leibniz Institute for Farm Animal Biology, Dummerstorf, DE). Materials used in the test procedures were purchased from commercial sources as listed in the tables 2-4. The Biomek i7 hybrid deck layout shown in Figure 2 is used for the automated sample processing. The deck was optimized to process 144 samples during a single run. Sample cleanups were carried out using the Positive Pressure Extractor (amplius GmbH, Rostock, DE) integrated on the right side of the Biomek i7 deck. A self-refilling reservoir (amplius GmbH, Rostock, DE) was integrated to facilitate the delivery of large volumes (e.g. 127 mL acetonitrile for protein precipitation).

The samples were placed on the deck in 1.5 mL Eppendorf safe-lock vials (Eppendorf, Hamburg, DE) using a special adapter (CELISCA, Rostock, DE) (see Figure 4) on the i7 deck and were processed according to Table 1. Serum samples with concentrations of 1 µg/mL 25(OH)D₂ and 25(OH)D₃ were used for the method validation. After extraction, the samples were analyzed by injecting 10 µL of the sample into an LC/TOF-MS system (Agilent Technologies, Santa Clara, US) with a flow rate of 0.5 mL / min. The system was calibrated according to the internal standard method in the range of 0.01 to 2 ppm 25(OH)D₂ and 25(OH)D₃.

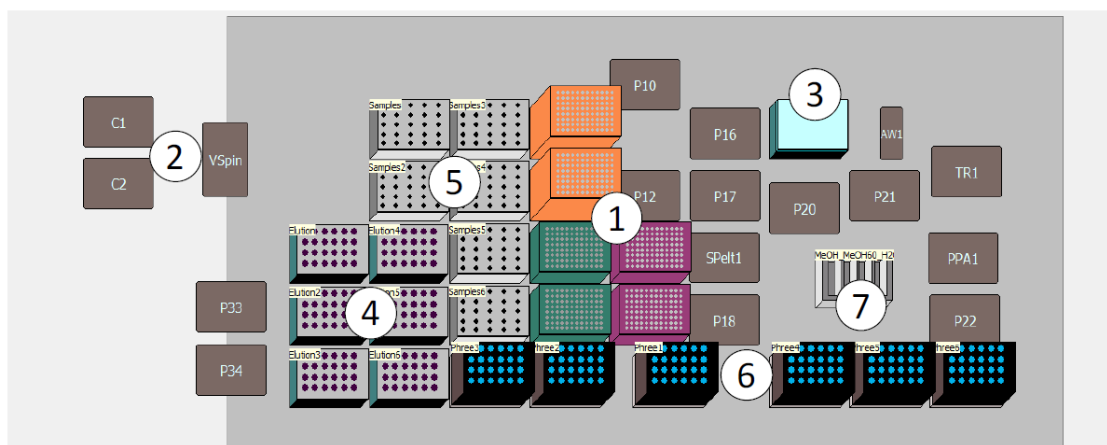


Figure 2. Deck layout for vitamin D determination using Phree phospholipid removal cartridges - (1) Tip boxes 1070 µL, 230 µL and 930 µL, (2) Centrifuge, (3) Internal Standard (ISTD), (4) Adapter made of aluminum with elution vials, (5) Samples, (6) Phree phospholipid removal cartridges, (7) Quarter self-refilling reservoir. Note: The integrated centrifuge is not used in the protocol.

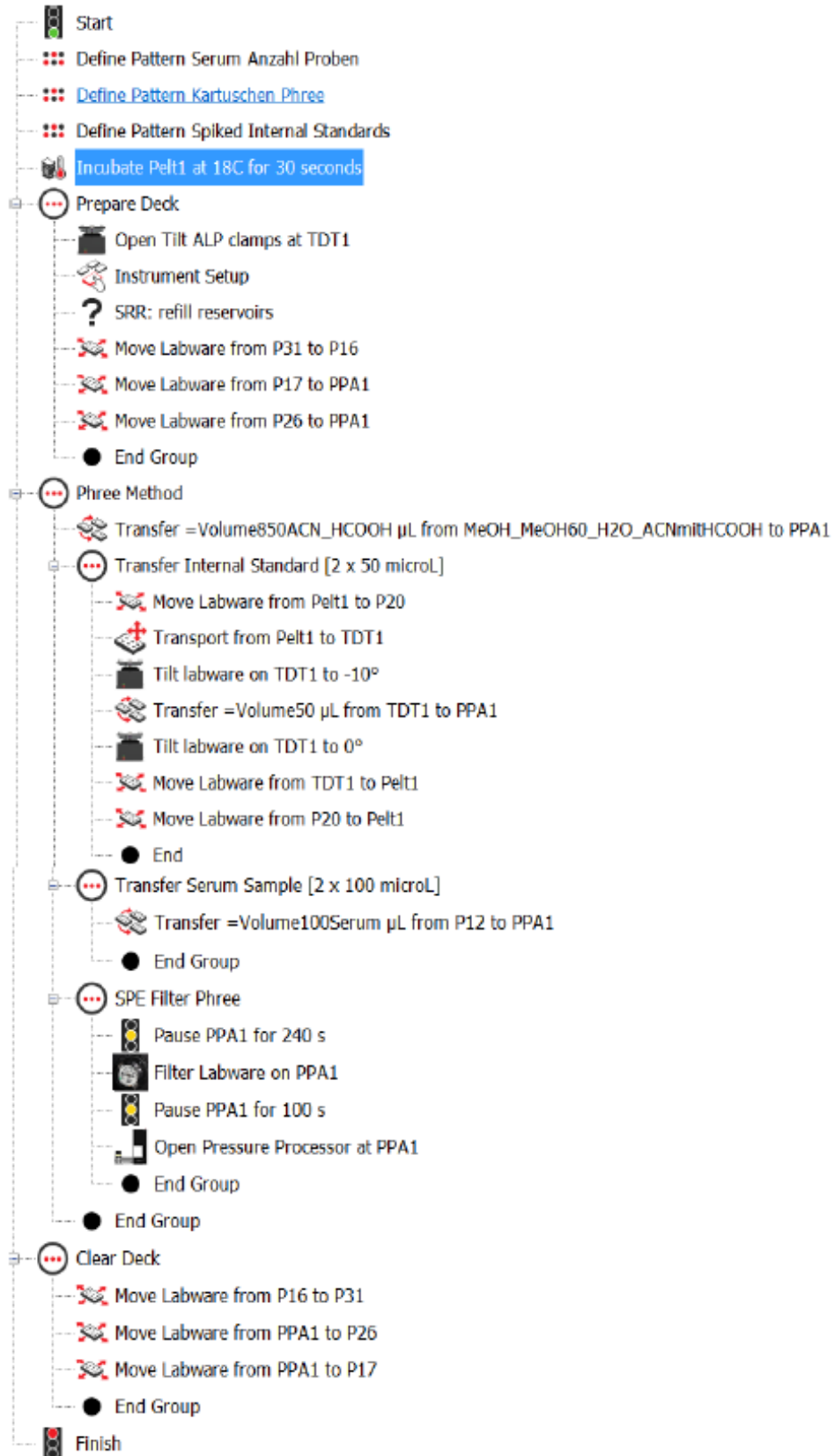


Figure 3. Biomek i7 method for vitamin D determination – Sample cleanup using Phree phospholipid removal cartridges

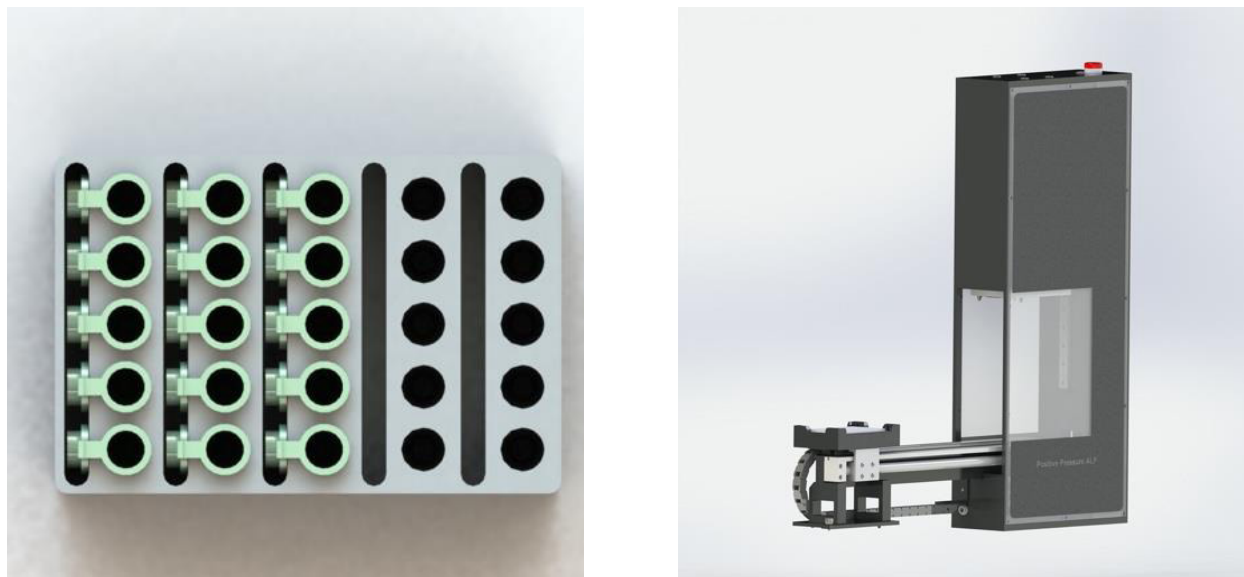


Figure 4. (a) Storage for samples vials (CELISCA, Rostock, DE), (b) Positive Pressure Extractor (amplus GmbH, Rostock, DE)

| Step Description | Manufacturer |
|------------------|--|
| Step 1 | Transfer 850 μ L acetonitrile to Phree cartridges |
| Step 2 | Transfer 50 μ L internal standard (ISTD) to Phree cartridges for all samples |
| Step 3 | Transfer 100 μ L serum sample to Phree cartridges |
| Step 4 | Mix the solution with multiple aspiration/dispensing |
| Step 5 | Incubation at room temperature for 4 min |
| Step 6 | Apply positive pressure (100 sec., 500 mbar) |
| Step 7 | Measurement using LC/MS or LC/MS/MS |

Table 1. Sample processing protocol for the vitamin D determination using Phree cartridges

Results

The measurements were calibrated between 0.01 ppm and 2 ppm for 25(OH)D₂ and 25(OH)D₃ (see Figure 5). The recovery rates were determined with values of 90.95%-106.21% for 25(OH)D₂ and 94.61%-115.50 % for 25(OH)D₃. The high recovery rates indicate optimal removal of phospholipids. The coefficient of variation (CV) of 4.36% (25(OH)D₂) and 5.1% (25(OH)D₃) and was determined with 10 samples. To determine between-experiment precision, the experiment was repeated on 5 days with 10 samples each resulting in a CV between 2.68% and 8.36% (25(OH)D₂) and 1.69-5.51% (25(OH)D₃). The limit of detection (LOD) and limit of quantification (LOQ) were determined at 5.91 ng/mL (LOD) and 12.54 ng/mL (LOQ) for 25(OH)D₂ as well as 26 ng/mL and 73.2 ng/mL for 25(OH)D₃. The detection limits can be further improved by optimizing the LC/MS/MS system. The results obtained were comparable to the results of classical methods in which protein precipitation is carried out with centrifugation.

Summary

In this application note we automated sample processing for LC-MS based vitamin D analysis on a Biomek i7 hybrid workstation. The use of Phree phospholipid removal cartridges for combined protein precipitation and sample clean-up reduced the number of processing steps and the processing time. The high recovery rate is indicative of efficient removal of phospholipids. The low CV values (2-9%) of the automated workflow indicates high repeatability.

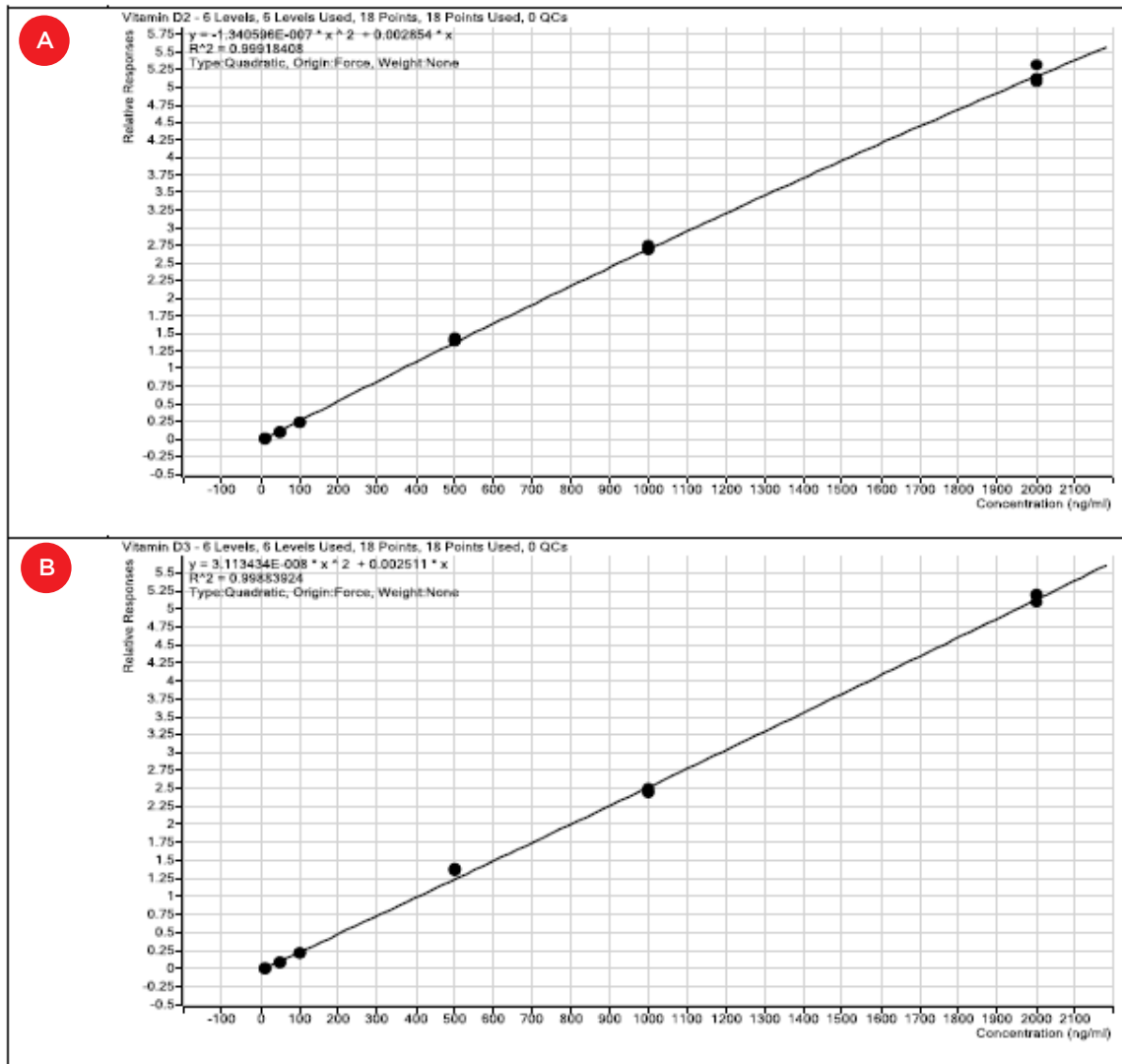


Figure 5. Calibration for (a) 25(OH)D₂ and (b) 25(OH)D₃

References

1. Ding S, Schoenmakers I, Jones K, et al. Quantitative determination of vitamin D metabolites in plasma using UHPLC-MS/MS. *Anal Bioanal Chem* 2010;378: 779-789.
2. Pucci V, Di Palma S, Alfieri A, et al. A novel strategy for reducing phospholipids-based matrix effect in LC-ESI-MS bioanalysis by means of Hybrid SPE. *J Pharm Biomed Anal* 2009;50 (5) 867-871.
3. Tulipani S, Llorach R, Urpi-Sarda M, et al. Comparative analysis of sample preparation methods to handle the complexity of the blood fluid metabolome: When less is more. *Anal Chem* 2013;85(1): 341-348.
4. Huq S, Pike Safan E, Rapid and Effective Cleanup of Vitamin D2 and D3 from Human Plasma using Phree™ Phospholipid Removal Plates and Kinetex® Core-Shell HPLC/UHPLC Columns. Application Note TN0061, Phenomenex Inc., 2013.

| Equipment | Manufacturer |
|--|--|
| Biomek i7 Workstation | Beckman Coulter Life Sciences, Indianapolis, USA |
| Shaking Peltier for Biomek® 4000, FXp, NXp | INHECO Industrial Heating & Cooling GmbH, Martinsried, Germany |
| Static Peltier Biomek® 4000, FXp, NXp | INHECO Industrial Heating & Cooling GmbH, Martinsried, Germany |
| Positive Pressure Extractor | amplus GmbH, Rostock, Germany |
| 3D Tilt ALP | amplus GmbH, Rostock, Germany |
| Self Refilling Quarter Reservoir | amplus GmbH, Rostock, Germany |

Table 2. Instruments used 3

| Reagents | Manufacturer | Part Number |
|----------------------------------|-------------------------------|-------------|
| D6-25(OH)D ₃ (50 ppm) | Sigma Aldrich, St. Louis, USA | H-074 |
| 25(OH)D ₂ (50 ppm) | Sigma Aldrich, St. Louis, USA | H-073 |
| 25(OH)D ₃ (100 ppm) | Sigma Aldrich, St. Louis, USA | H-083 |
| ACN mit HCOOH 0.1% | Sigma Aldrich, St. Louis, USA | 1590024000 |

Table 3. Reagents Method 3

| Consumables used per 144 samples | | | |
|-------------------------------------|--------|--|--------------|
| Consumables | Number | Manufacturer | Part number |
| Biomek i-series tips 90 µL | 144 | Beckman Coulter Life Sciences, Indianapolis, USA | B85881 |
| Biomek i-series tip 230 µL | 144 | Beckman Coulter Life Sciences, Indianapolis, USA | B85903 |
| Biomek i-series tips 1070 µL | 144 | Beckman Coulter Life Sciences, Indianapolis, USA | B85971 |
| Phree Phospholipid Removal 1mL Tube | 144 | Phenomenex, Torrance, USA | 8B-S133-TAK |
| GC Vials 2 mL | 144 | Agilent Technologies, Santa Clara, USA | 5182-0716 |
| Eppendorf Vials 1.5 mL (Safe Lock) | 144 | Eppendorf AG, Hamburg, Germany | EP0030121880 |
| Septum Lid | 144 | Agilent Technologies, Santa Clara, USA | 5182-0731 |

Table 4. Consumables Method 3

Biomek i-Series Automated Workstations are not intended or validated for use in the diagnosis of disease or other conditions.

©2021 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the Stylized Logo, and Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are the property of their respective owners. This protocol is for demonstration only and is not validated by Beckman Coulter. Beckman Coulter makes no warranties of any kind whatsoever express or implied, with respect to this protocol, including but not limited to warranties of fitness for a particular purpose or merchantability or that the protocol is non-infringing. All warranties are expressly disclaimed. Your use of the method is solely at your own risk, without recourse to Beckman Coulter. Not intended or validated for use in the diagnosis of disease or other conditions.



For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at beckman.com
21.01.750.AUTO