

# Quantitative Analysis of Mevalonate in Plasma Using LC-MS/MS

Flavio Giavarini<sup>1</sup>, Omar Maschi<sup>1</sup>, Samuele Scurati<sup>2</sup>, Donatella Caruso<sup>1</sup>

<sup>1</sup>Department of Pharmacological Sciences, Università degli Studi di Milano, Milano, Italy;

<sup>2</sup>Thermo Fisher Scientific, Rodano, Italy

For Research Use Only. Not for use in diagnostic procedures.

## Introduction

Cholesterol is synthesized *in vivo* through a multiple step pathway. Because mevalonate is the key intermediate of this process, its plasmatic levels are an indirect measure of *in vivo* cholesterol synthesis and, therefore, facilitate clinical research into pharmacological activity of anti-hypercholesterolemic drugs such as statins.

## Goal

To develop a reliable and fast analytical method for the quantitative determination of mevalonate in plasma using a Thermo Scientific LTQ linear ion trap mass spectrometer.

## Experimental

### Sample Preparation

The plasma sample (500  $\mu$ L) was spiked with 20 ng of Mevalonate-D<sub>7</sub>. Samples were acidified with hydrochloric acid allowing the conversion of mevalonate to mevalonolactone (Figure 1). After purification through solid phase extraction (SPE), samples were dried and dissolved in 400  $\mu$ L of 0.2% ammonium hydroxide to restore the non-lactonic form. Then 10  $\mu$ L were injected.

Quantitative analysis was performed on the basis of calibration curves, ranging from 2.5 to 250 ng/mL.

### HPLC Conditions

High performance liquid chromatography (HPLC) analysis was performed using a Thermo Scientific Surveyor autosampler and pump. The 10  $\mu$ L sample was injected directly on a Thermo Scientific BioBasic AX column (150  $\times$  2.1 mm, 5  $\mu$ m). A gradient LC method used mobile phases A (10 mM ammonium formate, pH 8) and B (acetonitrile) at a flow rate of 200  $\mu$ L/min.

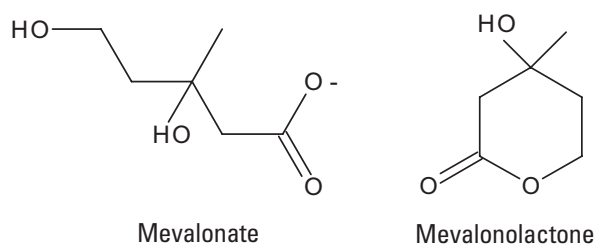


Figure 1. Structure of mevalonate and mevalonolactone

## Mass Spectrometry

MS analysis was carried out on a LTQ™ linear ion trap mass spectrometer equipped with a Thermo Scientific Ion Max source with an electrospray ionization (ESI) probe.

Ion polarity:	Negative
Spray voltage:	2 kV
Sheath/Auxiliary gas:	Nitrogen
Sheath gas pressure:	40 (arbitrary units)
Auxiliary gas pressure:	10 (arbitrary units)
Sweep gas pressure:	5 (arbitrary units)
Ion transfer tube temperature:	300 °C
Scan type:	Full Scan MS/MS
Collision gas:	Helium
Collision energy:	30%
Divert valve:	3.0-6.5 min to source
Selected ions for quantification:	$m/z$ 147 $\rightarrow$ 59 for mevalonate $m/z$ 154 $\rightarrow$ 59 for mevalonate-D <sub>7</sub>

## Results and Discussion

Figure 2 shows the ion chromatograms of a lower sample of the calibration curve. Excellent linearity ( $r^2 = 0.999$ ) fits for the calibration curve were observed over the range of 2.5 - 250 ng/mL plasma (Figures 3 and 4). The intraday CV% ( $n=3$ ) was in the range 0.5% - 4%. The limit of detection (LOD) was 2 pg, and the limit of quantification (LOQ) was 2.5 ng/mL.

Figure 5 reports an ion chromatogram of a plasma sample of a healthy volunteer (24 ng/mL plasma), extracted and analyzed as described.

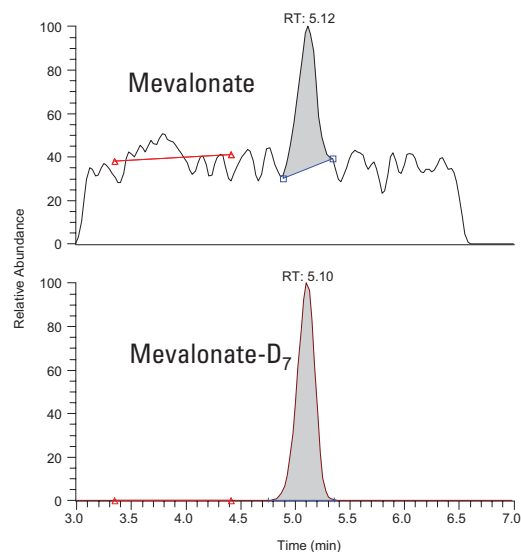


Figure 2. Ion chromatograms of 2.5 ng/mL calibration standard

## Key Words

- LTQ Ion Trap
- Clinical Research
- Cholesterol Synthesis

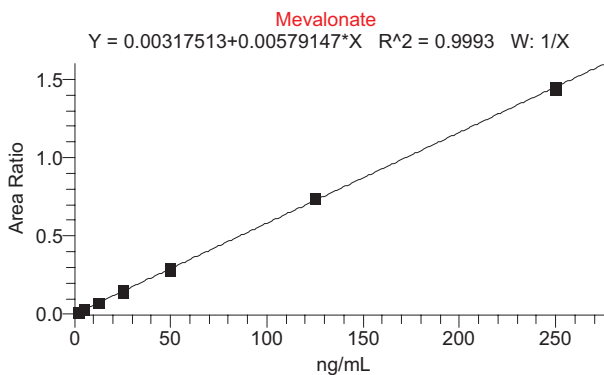


Figure 3. Calibration curve of mevalonate

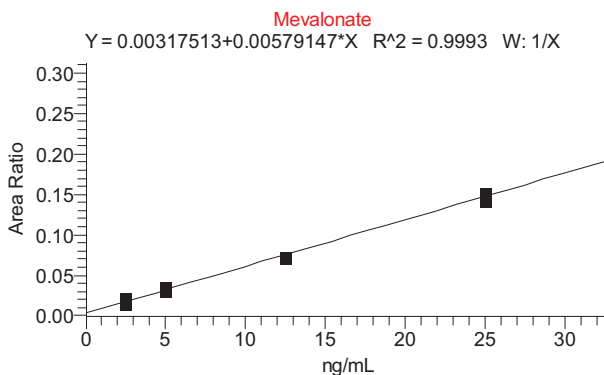


Figure 4. Zoom on low calibration points

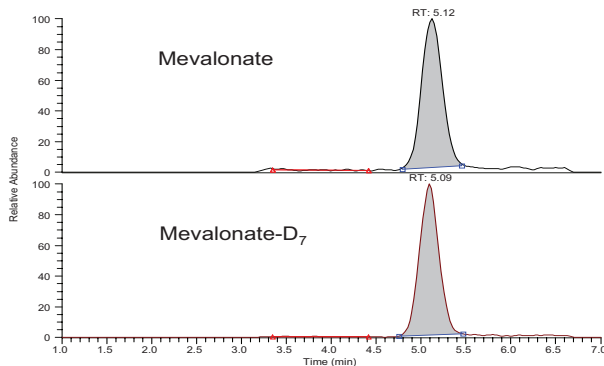


Figure 5. Ion chromatograms of plasma sample containing 24 ng/mL

### Conclusion

A robust 10-minute method for the quantification of mevalonate with a dynamic range of 2.5 - 250 ng/mL plasma has been developed for clinical research using fast SPE purification and the LTQ linear ion trap mass spectrometer.

In addition to these offices, Thermo Fisher Scientific maintains a network of representative organizations throughout the world.

- Africa-Other**  
+27 11 570 1840
- Australia**  
+61 3 9757 4300
- Austria**  
+43 1 333 50 34 0
- Belgium**  
+32 53 73 42 41
- Canada**  
+1 800 530 8447
- China**  
+86 10 8419 3588
- Denmark**  
+45 70 23 62 60
- Europe-Other**  
+43 1 333 50 34 0
- Finland/Norway/Sweden**  
+46 8 556 468 00
- France**  
+33 1 60 92 48 00
- Germany**  
+49 6103 408 1014
- India**  
+91 22 6742 9434
- Italy**  
+39 02 950 591
- Japan**  
+81 45 453 9100
- 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
- Russia/CIS**  
+43 1 333 50 34 0
- South Africa**  
+27 11 570 1840
- Spain**  
+34 914 845 965
- Switzerland**  
+41 61 716 77 00
- UK**  
+44 1442 233555
- USA**  
+1 800 532 4752

[www.thermofisher.com](http://www.thermofisher.com)

Legal Notices: ©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific Inc. 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.

AN63350\_E 06/16S