

Author:
PerkinElmer, Inc.
Waltham, MA

High Sensitivity Methotrexate Quantitation in Human Serum Utilizing PerkinElmer QSight® 220 Triple Quadrupole Mass Spectrometer

Introduction

This note outlines a rapid LC-MS/MS research method utilizing the QSight® 220 triple quadrupole mass spectrometer. The developed method provides exceptional results for the quantitation of methotrexate in serum especially in terms of LLOQs and linearity.

2. Method

Methotrexate was purchased from Toronto Research Chemicals (Toronto, ON). The stock solution 25 mg/mL of Methotrexate in methanol was prepared and stored at -15°C. Human serum was purchased from Sigma -Aldrich (Milwaukee, WI) and cleaned up by protein precipitation by mixing one volume of human serum, one volume of 0.1% formic acid and two volumes of acetonitrile. The mixture was vortexed for 1 min followed by centrifugation for 15 min. The supernatant was transferred to clean container and used as a matrix for methotrexate quantitation. A series of methotrexate standards from 0.03 to 125ng/mL were prepared by spiking the stock solution into the clean protein precipitated human serum. All the solvents used in this application are HPLC grade.

2.1. Mass Spectrometry Conditions

The QSight® 220 equipped with an ESI source was used to perform the analysis. Table 1 highlights the instrument settings used during the method.

Table 1: Settings used on the QSight® 220 during the method

ESI Voltage (V)	5000
HSID Temp (°C)	300
Nebulizer Gas Setting	400
Drying Gas Setting	120
Heating Gas	350
Source Temp. (°C)	250
Dwell Time (ms)	100
Pause Time (ms)	5

Table 2: Selected MRM operating conditions for methotrexate

MRM	CE	CCL2	CCL4	EV
455.1/308.0	28	-100	-90	30
455.1/134.0	48	-100	-90	30

Quick Facts:

- Method for high sensitivity quantitation of methotrexate using the QSight® 220 triple quadrupole mass spectrometer.
- LLOQ achieved is 0.03 ng/mL with 5µL injection
- Liquid Chromatography: Shimadzu® Prominence UFLC®
- ESI Source
- Calibration Curve $R^2 = 0.99999$

2.2. LC Conditions

The separation was performed on a Shimadzu® Prominence UFLC® system which includes binary pumps, autosampler, degasser, column oven. A sample volume of 5 µL is injected into a Imtakt C8 column (2.0 x 75 mm, 3 µm) at room temperature. The LC time program is illustrated as Table 3. The total LC cycle time is 3 min and liquid flow rate is 0.4 mL/min.

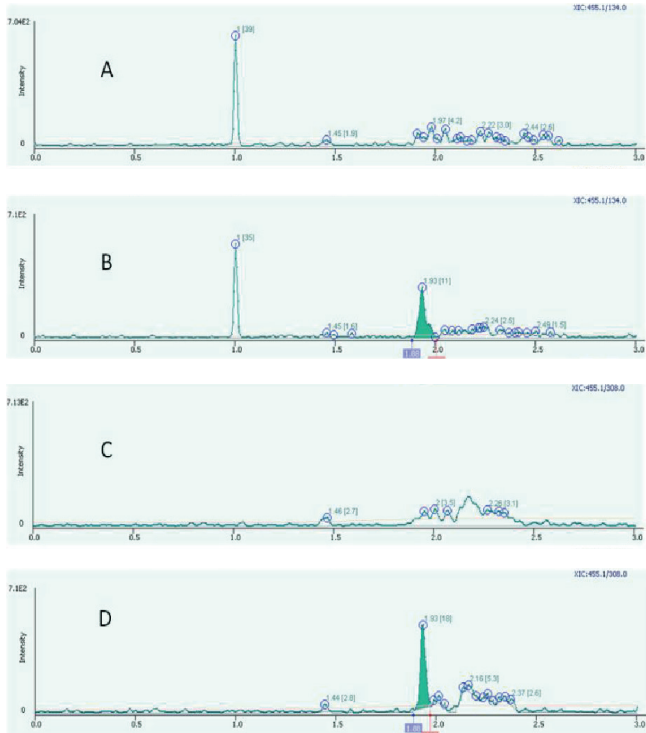
Table 3: LC Gradient Conditions; Solvent A is water with 0.1% formic acid, Solvent B is methanol with 0.1% formic acid

Time (min)	Procedure Solvent B composition (%)
0	5
0.1	5
0.6	95
1	95
1.1	5
3	5

3. Results

3.1. Extracted Ion Chromatogram (EIC)

Figure 1: (A) 455.1/134.0 Serum Blank, (B) 455.1/134.0 at a Concentration of 0.03 ng/mL of Methotrexate in Human Serum, (C) 455.1/308.1 Serum Blank, (D) 455.1/308.1 at a Concentration of 0.03 ng/mL of Methotrexate in Human Serum.



3.2. Linearity

The calibration curves generated for both MRM transitions using 1/x weighting are displayed in Figure 2 and Figure 3. Excellent linearity values, ($R^2=0.9999$) were obtained for concentrations ranging from 0.03 to 125 ng/mL for both of the monitored MRM transitions.

Table 4: Intra-day variability (%RSD)

MRM	QC1(low)	QC2(med)	QC3(high)
455.1/134.0	5.1	1.2	1.0
455.1/308.0	6.9	1.8	0.5

Table 5: Inter-day variability (%RSD)

MRM	QC1(low)	QC2(med)	QC3(high)
455.1/134.0	2.6	3.0	3.5
455.1/308.0	7.0	3.6	3.8

Figure 2: Calibration Curve for 455.1/134.0

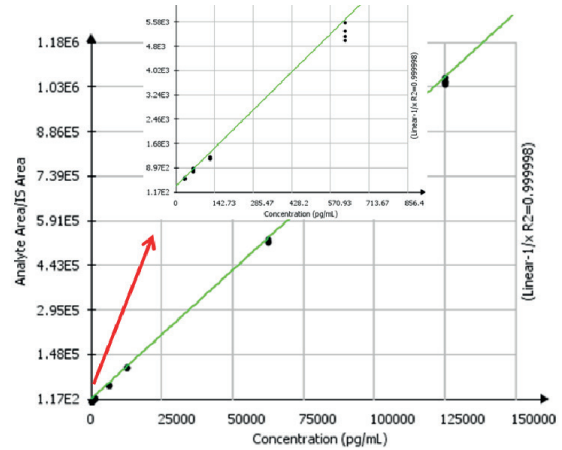
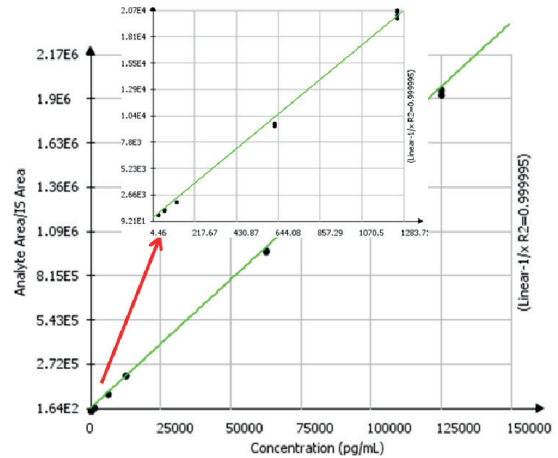


Figure 3: Calibration curve for 455.1/308.0



3.3. Quantitation Results

The inter and intra-day variability for this experiment were excellent as shown in Table 4 and 5. The intra-day variability is determined by processing 4 replicates of each QC sample and the inter-day variability is determined with 4 replicates in 3 batches.

4. Conclusion

A fast, sensitive, and accurate LC-MS/MS method was developed for methotrexate in serum using the PerkinElmer QSight® 220 triple quadrupole mass spectrometer. With an LLOQ of 0.03 ng/mL (using a 5 µL injection), the method falls well within the range necessary for methotrexate related clinical research applications. The linearity achieved by this method ($R^2=0.9999$) further illustrates the exceptional performance of the QSight® 220 for methotrexate quantitation.

5. Contact Information

To learn more about PerkinElmer Mass Spectrometry, our products or services please visit our website or contact us directly.

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
Phone: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com

PerkinElmer, Inc.
Wallac Oy
PO Box 10
20101 Turku, Finland
Phone: + 358 2 2678 111
Fax: + 358 2 2678 357

ISO 13485
ISO 9001
CMDCAS
ISO 14001
OHSAS 18001



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2018, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

1599-9832