

Reliable Quantitation on 52 L-Amino Acids in Human Plasma for Clinical Research by LC-MS/MS

¹Stephanie Samra, ²Valérie Thibert, ²Claude Netter ¹Thermo Fisher Scientific, San Jose, CA, USA; ²Thermo Fisher Scientific, Courtaboeuf, France

ABSTRACT

Purpose: An analytical method was developed that quantitatively detects 52 L-amino acids in human plasma for clinical research. An offline protein precipitation of the plasma sample without derivatization was performed prior to dilution with internal standard and followed by LC-MS/MS analysis with a triple quadrupole mass spectrometer. Separation and quantitation of compounds were performed with an 18-minute reverse phase gradient and internal or external calibration of eight standards ranging from 0.1 to 500 µM. Problematic issues such as observed carryover of the basic amino acids as well as retention time shifts were addressed by optimizing the LC method.

Methods: Plasma samples were protein precipitated and diluted with an internal standard solution prior to injection onto an LC-MS/MS triple quadrupole mass spectrometer. Separation of compounds was performed with an 18-minute aqueous/organic reverse phase gradient. Issues with carryover that originated from the mobile phases were observed in blank samples and were addressed with the use of a cation exchange trap column for their removal. Retention time shifts were kept to a minimum for excellent reproducibility by limiting pH differences to < 0.1 batch to batch.

Results: The lower limit of quantification (LLOQ) for each amino acid was set to the lowest calibration standard analyzed that yielded < 20 % inaccuracy and < 20% CV for 3 replicate injections of the same standard.

Figure 1. Thermo Scientific™ Vanquish™ Flex and TSQ Quantis™ Triple Quadrupole Mass Spectrometer.



INTRODUCTION

Amino acids are the building blocks of proteins and intermediaries to many biochemical pathways. They play an important role in biochemical regulation and any abnormalities in abundance may be indicative of greater metabolic issues or inherited disorders. Here we present a novel workflow that simplifies sample preparation by eliminating the use of a derivatization reagent to detect 52 L-amino acids for clinical research in human plasma with excellent reproducibility in 18 minutes.

Current methods most widely accepted in the routine clinical research laboratory require several steps which are highly cumbersome and time consuming. The method presented is simple, straightforward, and reproducible.

MATERIALS AND METHODS

Sample Preparation

All labeled and unlabeled standards were purchased from Cambridge Isotope Laboratories Inc., Sigma Aldrich, or Wako Pure Chemical Industries. Standard solutions were prepared at the appropriate concentrations and used for dilution of the sample prior to injection onto the LC-MS/MS system.

100 µL of calibrator, control, or plasma were mixed with 100 µL of precipitation reagent and then vortexed for 60 seconds, refrigerated for 30 minutes, and centrifuged at 12000 rpm for 5 minutes. The supernatant (50 µL) was added to 200 µL of the dilution solution containing both labeled and unlabeled standards. Samples were vortexed and then injected onto the LC-MS/MS system.

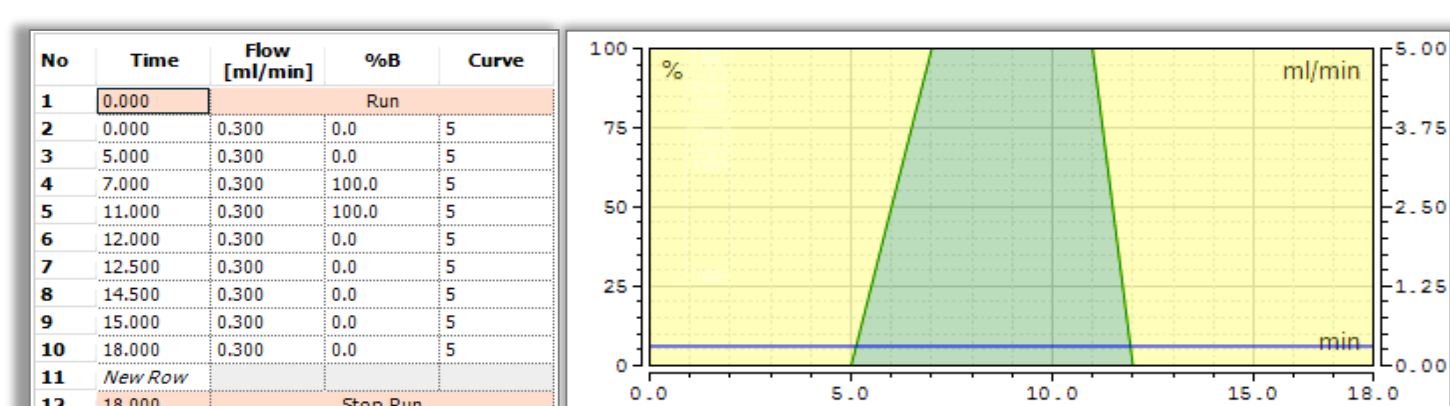
Table 1. List of 52 quantifiable L-Amino Acids. The listed amino acids are detectable in human plasma using this method.

Internal Standards		Internal Standards	
1 Alanine 13C3 15N1	27 Homocysteine	1 Alanine 13C3 15N1	
2 β-Alanine	28 Hydroxyproline	2 Arginine 13C6 15N4	
3 α-Amino adipic Acid	29 Hydroxyproline	3 Asparagine 13C4 D3 15N2	
4 α-Amino isobutyric Acid	30 Isoleucine	4 Aspartic Acid 13C4 15N1	
5 β-Amino isobutyric Acid	31 Allo-Isoleucine	5 Cysteine 13C6 15N2	
6 γ-Aminobutyric Acid	32 Lysine	6 Cystathionine 13C3 15N1	
7 α-Aminovaleric Acid	33 Lysine	7 Glutamic Acid 13C5 15N3	
8 Arginine	34 Lysine	8 Glutamine 13C5 D5 15N2	
9 Arginine	35 Methionine	9 Glycine 13C2 15N1	
10 Argininosuccinic Acid	36 1-Methylhistidine	10 Histidine 13C6 15N3	
11 Asparagine	37 3-Methylhistidine	11 Isoleucine 13C5 15N1	
12 Aspartic Acid	38 Ornithine	12 Leucine 13C6 15N1	
13 Carnosine	39 Phenylalanine	13 Lysine 13C6 15N2	
14 Citrulline	40 Phosphoethanolamine	14 Methionine 13C5 15N1	
15 Cystathionine	41 Phosphoserine	15 Ornithine D6	
16 Cysteine	42 Picopalic Acid	16 Phenylalanine 13C9 15N1	
17 Cysteine-Homocysteine Disulfide	43 Proline	17 Picopalic Acid 13C6 15N1	
18 Cysteine	44 Saccharopine	18 Proline 13C5 15N1	
19 Ethanolamine	45 Sarcosine	19 Serine 13C3 15N1	
20 Glutamic Acid	46 Serine	20 Taurine 13C2 15N1	
21 Glutamine	47 Sulfocysteine	21 Threonine 13C4 15N1	
22 Glutathione Reduced	48 Taurine	22 Tryptophan 13C11 15N2	
23 Glycine	49 Threonine	23 Tyrosine 13C3 15N1	
24 Histidine	50 Tyrosine	24 Valine 13C5 15N1	
25 Homocitrulline	51 Tyrosine	25 γ-Aminobutyric Acid 13C4	
26 Homocysteine	52 Valine		

Liquid Chromatography Conditions

- Mobile phases consisted of an organic/aqueous composition.
- Columns used for this method consisted of a trap column to remove basic amino acids seen consistently in the background, a guard column to increase column lifetime, and an analytical column for amino acid separation.
- Column Temperature: 30 ° C
- Injection volume: 4 µL

Figure 2. Liquid chromatography gradient used for analysis.



Conditions for TSQ Quantis Mass Spectrometer

Table 2. HESI Source Parameters.

HESI Source Parameters			
Polarity	Positive	Sheath Gas	45
Spray Voltage	3.5 kV	Auxiliary Gas	15
Capillary Temp	270 °C	Sweep Gas	0
Vaporizer Temp	370 °C	Source Fragmentation	15.0 V

Table 3. Scan Properties used for SRM experiment type.

SRM Scan Properties			
Use Cycle Time	True	Collision Gas Pressure	1.5 mTorr
Cycle Time	0.4 sec	Use Calibrated RF	False
Chrom Filter	3.0 sec	Q1 Resolution (FWHM)	0.7
Data Mode	centroid	Q3 Resolution (FWHM)	0.7

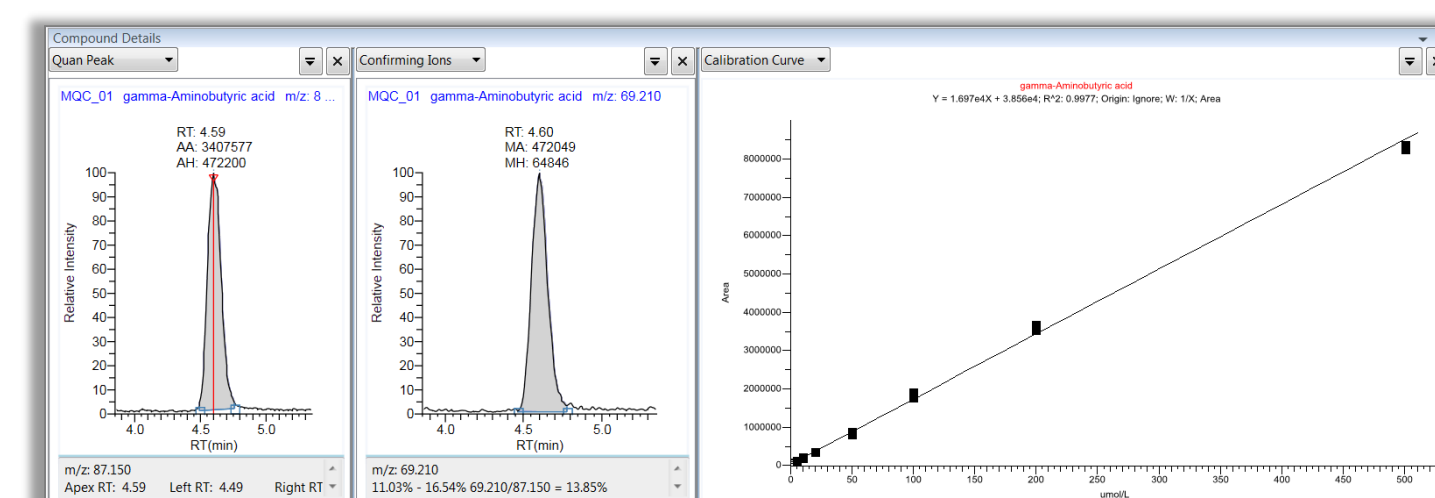
RESULTS

Table 4. SRM Table and Calibration Curves.

Amino Acid	Precursor (m/z)	Product (m/z)	RT	R ²	Range µmol/L	Type	Weighting
1-methylhistidine	170.12	124.11	10.41	0.9915	2-200	Linear	1/X
3-methylhistidine	170.12	126.1	10.27	0.9945	2-100	Linear	1/X
alanine	90.274	44.44	2.03	0.9969	20-500	Linear	1/X
α-amino adipic	162.152	98.22	2.35	0.9949	5-250	Linear	1/X
α-aminobutyric	104.213	58.33	2.43	0.9819	10-250	Linear	1/X
anserine	241	109	10.93	0.9966	5-200	Linear	1/X
arginine	175.12	70.26	12.12	0.9912	5-200	Linear	1/X
asparagine	133.2	87.15	1.9	0.9977	20-500	Linear	1/X
aspartic acid	134.11	74.22	2.03	0.9949	10-500	Linear	1/X
β-aminoisobutyric acid	104.252	86.15	4.22	0.9942	5-500	Linear	1/X
carnosine	227.183	110.17	11.21	0.9953	5-200	Linear	1/X
citrulline	176.213	159.04	2.87	0.9962	5-200	Linear	1/X
cystathionine	223.152	134.11	8.94	0.9985	2.5-250	Linear	1/X
cystine	241.1	152	8.71	0.9973	5-500	Linear	1/X
ethanolamine	62.365	62.37	3.75	0.9938	10-200	Linear	1/X
γ-aminobutyric acid	104.252	87.15	4.69	0.9968	5-200	Linear	1/X
glutamic acid	148.1	84.15	2.11	0.9968	5-500	Linear	1/X
glutamine	147.12	130.04	1.98	0.9926	5-500	Linear	1/X
glycine	76.3	76.3	2.16	0.9989	50-500	Linear	1/X
histidine	156.13	110.1	10.46	0.9594	5-100	Linear	1/X ²
hydroxylysine	163	128	10.15	0.9909	2-100	Linear	1/X ²
isoleucine	132.18	69.26	3.52	0.9974	5-500	Linear	1/X
leucine	132.18	86.22	3.74	0.9917	20-200	Linear	1/X
lysine	147.17	84.22	10.49	0.9906	5-500	Linear	1/X
methionine	150.098	133.07	3.08	0.9981	5-500	Linear	1/X
ornithine	133.243	70.26	10.22	0.9933	20-500	Quadratic	1/X
phenylalanine	166.12	103.17	6.85	0.9885	2-500	Linear	1/X ²
phosphoethanolamine	142.07	44.44	1.56	0.9946	2.5-100	Linear	1/X
phosphoserine	186	88	3.72	0.9957	2.5-250	Linear	1/X
pipecolic acid	130.16	84.22	2.98	0.9987	10-500	Quadratic	1/X
proline	116.243	70.26	2.13	0.9942	5-500	Quadratic	1/X
serine	106.191	60.33	1.88	0.9982	20-500	Linear	1/X
sulfocysteine	202.091	120.06	6.44	0.9989	2-500	Linear	1/X
taurine	126.152	126.15	1.54	0.9917	10-100	Linear	1/X
threonine	120.16	74.28	1.89	0.9965	10-500	Linear	1/X
tryptophan	205.078	187.99	9.64	0.998	5-500	Linear	1/X
tyrosine	182.078	136.04	5.95	0.9995	5-500	Linear	1/X
valine	118.17	72.26	2.62	0.9942	20-500	Linear	1/X

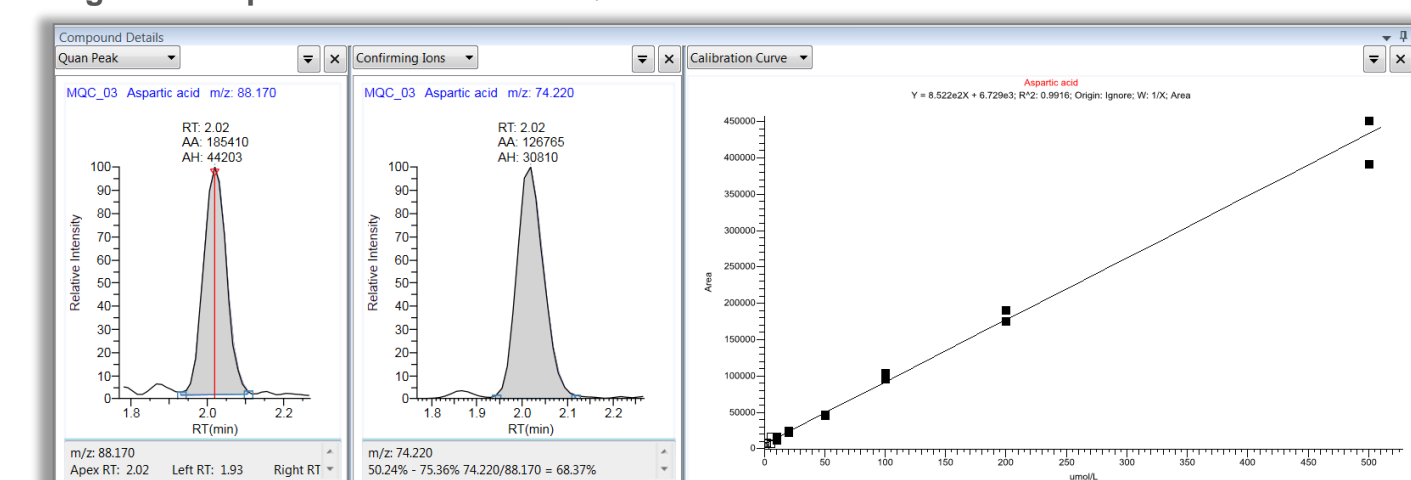
Amino acids quantitated with retention times, calibration R² values, calibration ranges, calibration type, and calibration weighting. Limit of Quantification (LOQ) was determined as the lowest value in the calibration curve giving <20%Diff and <20% CV for 3 injections of calibrators in neat solution.

Figure 3. Aspartic Acid Mid Level QC with calibration curve.



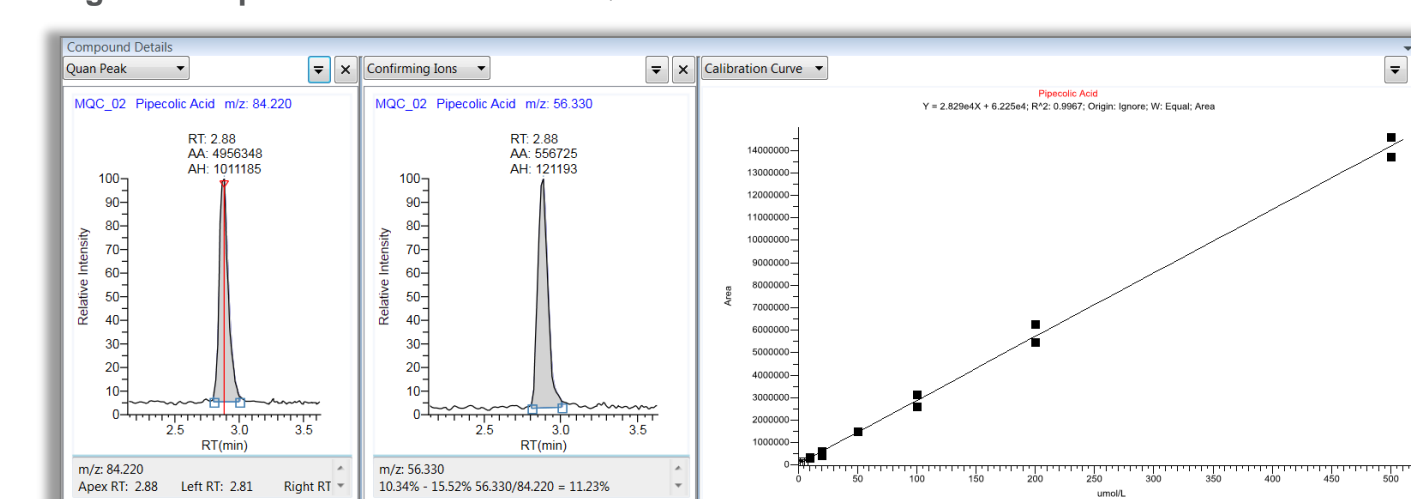
Example Chromatograms and Calibration Curves

Figure 3. Aspartic Acid Mid Level QC with calibration curve.



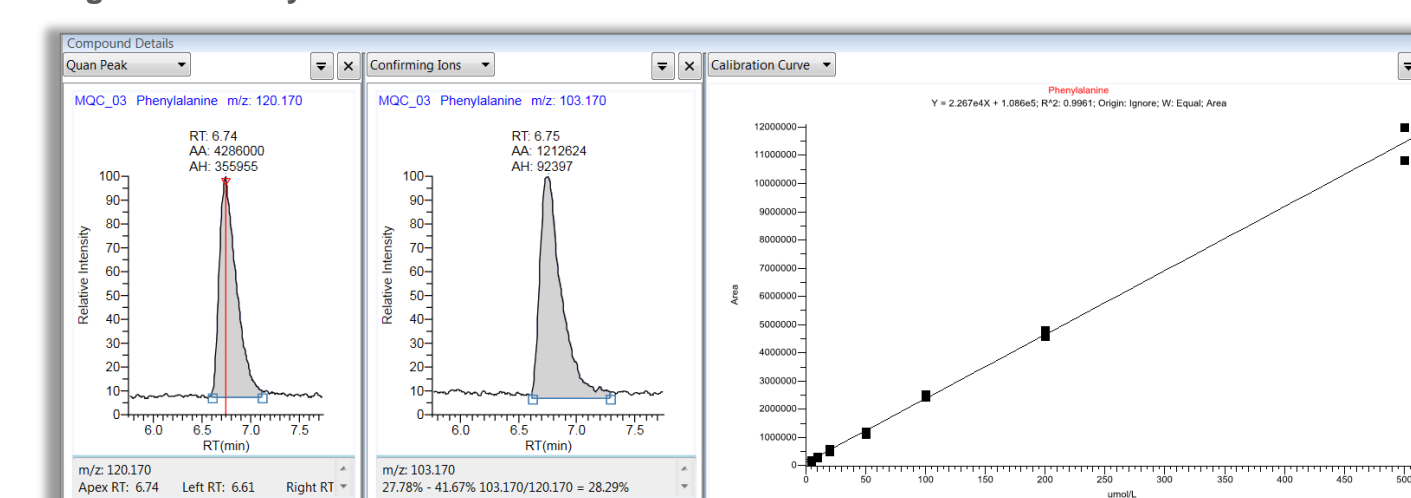
Chromatogram of quantitative and confirming ion peaks. Mid Level QC (200 µmol/L) was extracted using the protein precipitation method.

Figure 4. Pipecolic Acid Mid Level QC with calibration curve.



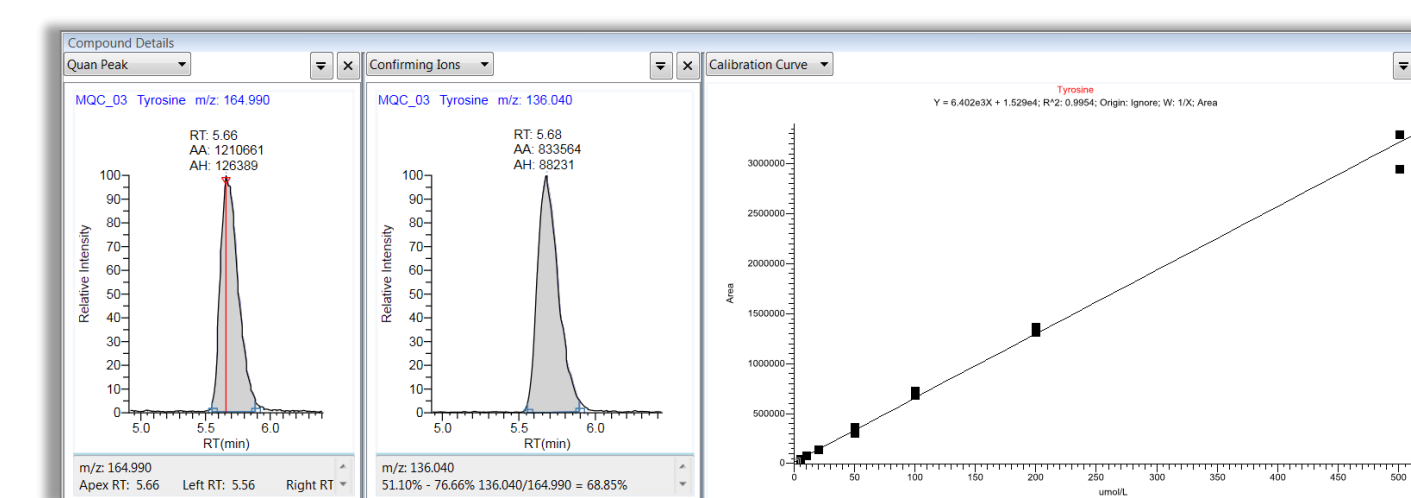
Chromatogram of quantitative and confirming ion peaks. Mid Level QC (200 µmol/L) was extracted using the protein precipitation method.

Figure 5. Phenylalanine Mid Level QC with calibration curve.



Chromatogram of quantitative and confirming ion peaks. Mid Level QC (200 µmol/L) was extracted using the protein precipitation method.

Figure 6. Tyrosine Mid Level QC with calibration curve.



Chromatogram of quantitative and confirming ion peaks. Mid Level QC (200 µmol/L) was extracted using the protein precipitation method.

Repeatability

Table 5. Repeatability of Mid Level QC (Control L1), High Level QC (Control L2), and Plasma. Controls, quality controls, and plasma samples were prepared using the protein precipitation method. Calibration curves were determined as described in Table 4. Controls and Plasma samples were analyzed using five replicate injections with % RSD and % Difference reported here for ten of the amino acids.

Amino Acid	Control L1				Control L2				Plasma	
	Theoretical µmol/L	AVG µmol/L	%RSD	%Diff	Theoretical µmol/L	AVG µmol/L	%RSD	%Diff	AVG µmol/L	% RSD
Aspartic Acid	200	194	4.23	-3.06	500	474	2.67	-5.17	44.4	17.7
Pipecolic Acid	200	178	3.98	-11.0	500	493	3.35	-1.46	N/F	-
Phenylalanine	200	180	3.63	-9.86	500	460	3.39	-8.10	37.1	5.78
Tyrosine	200	176	4.77	-12.0	500	443	2.37	-11.4	44.0	3.39
γ-aminobutyric acid	200	188	4.28	-6.08	500	474	2.67	-5.17	N/F	-
alanine	200	180	6.13	-10.2	500	467	2.71	-6.63	30.3	12.3
α-Amino adipic acid	100	86.7	5.79	-13.3	250	223	8.37	-10.6	5.20	8.23
α-Aminobutyric acid	100	88.9	9.33	-11.1	250	236	4.55	-5.52	N/F	-
β-Aminoisobutyric acid	200	175	2.18	-12.3	500	478	3.91	-4.31	N/F	-
citrulline	200	186	4.53	-7.03	500	442	3.97	-11.5	21.2	9.12

CONCLUSIONS

The elimination of derivatization simplifies the sample preparation and increases overall throughput with shorter preparation and analysis time, and also minimizes the need for costly reagents.

The LC-MS/MS method was able to simultaneously quantify 52 amino acids within 18 minutes, representing a significant improvement over conventional methods.

The method has been evaluated for 37 amino acids in neat standard and in human plasma, and will be further assessed for more extensive reproducibility and repeatability for all 52 amino acids.

TRADEMARKS/LICENSING

© 2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.