

A Comparison of Amino Acid Analysis (AAA) using UHPLC-ToF and CE-ToF

Bob Giuffre, Agilent Technologies, 550 Clark Dr, Budd Lake NJ, 07825
Dawn Stickle, Agilent Technologies, 40 Shattuck Rd, Andover , MA 01810
Dat Phan, Agilent Technologies, 770 Stanford Dr, Columbia MD, 20580

Purpose of Work

- Detection and quantification of amino acids plays an important role in protein analysis, food analysis, and physiological fluids.
- Sample preparation is an important issue in amino acid analysis and the sample matrix.
- Samples are often complex and the separation can prove challenging.
- Amino acids are often difficult to detect so amino acid analysis (AAA) typically involves chemical derivatization to improve detection by UV or fluorescence.
- Time of Flight (ToF) Mass Spectrometry coupled with either Ultra-High Pressure Liquid Chromatography (UHPLC) or Capillary Electrophoresis (CE) provides a sensitive and discriminating detection technique which will preclude the need for derivatization and avoid some of the sample matrix interferences.

- **Goal:** To compare these two sample introduction systems using a time-of-flight mass spectrometer as the detection system.

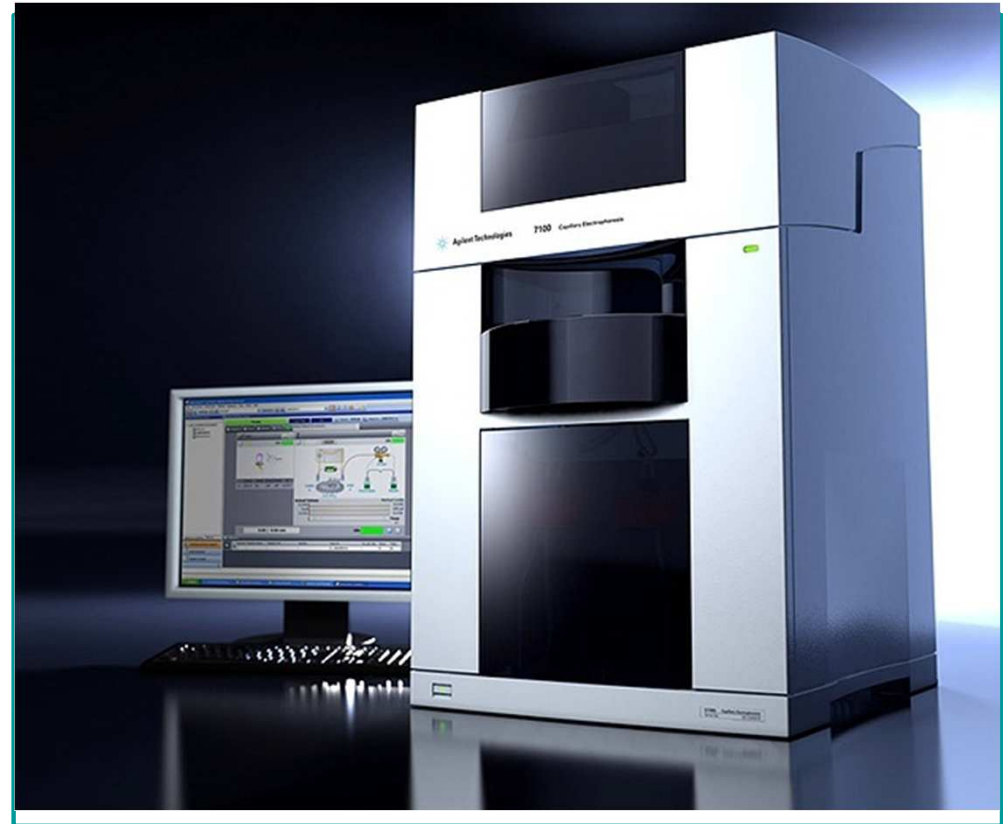
Agilent 1290 Infinity LC



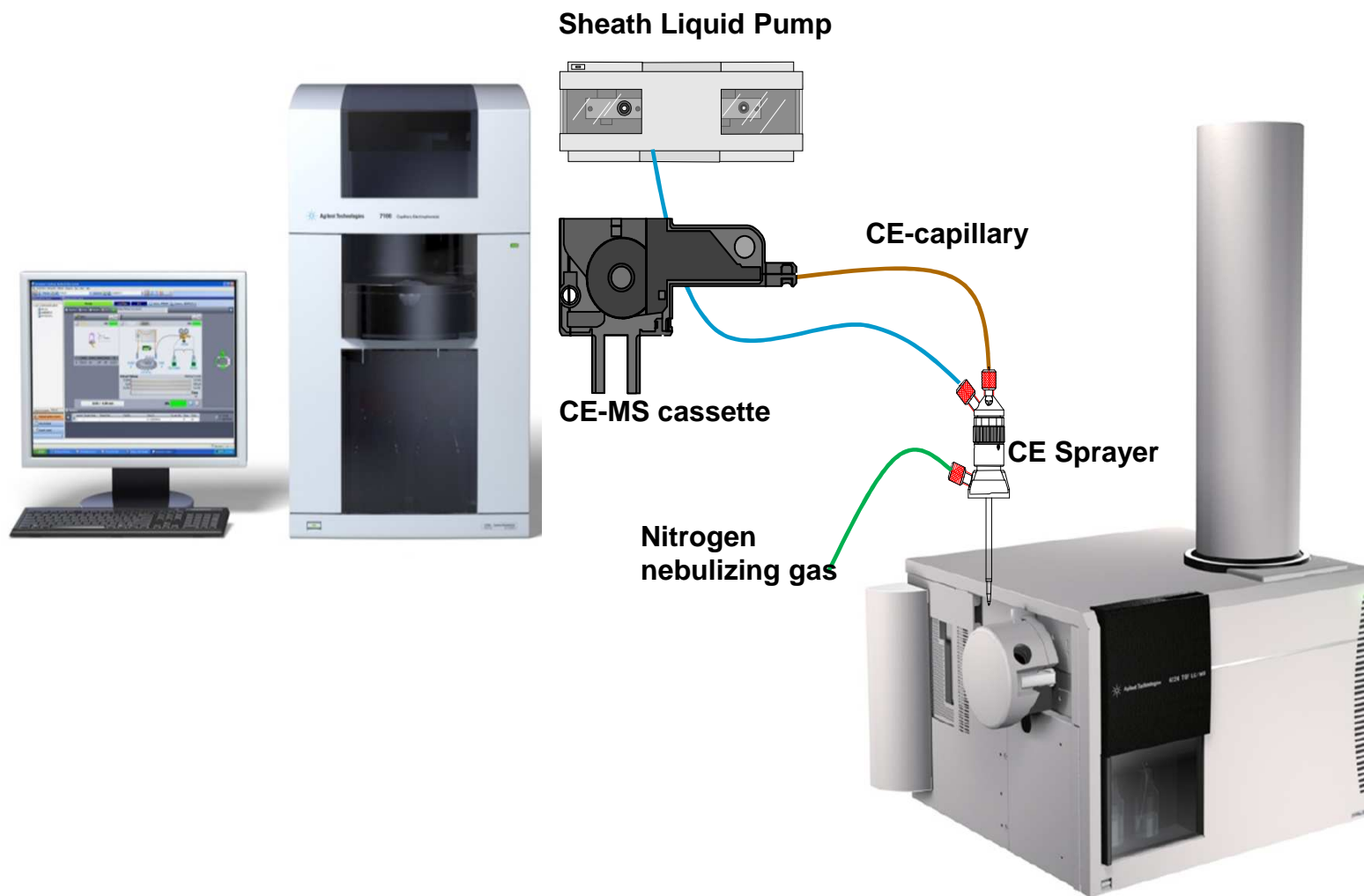
- Power range 1200 bar (2 ml/min);
800 bar 5 ml/min)
- Delay volume (Pump (w/o) mixer: 10 µl)
- Pump: e.g RT stability < 0.02 % (1.5 min runs)
- ALS precision for small volumes
- Injection principles: flow-through
- TCC: column switching, easy valve exchange
- DAD: lowest detector noise < +/- 3 µAU/cm
Max-Light High Sensitivity cell

Agilent 7100 Capillary Electrophoresis System Summary

- Updated UV detector with 40Hz acquisition rate
- Smaller footprint
- Flexible detection options
 - UV, MS, LIF, CCD
- Plug & play CE-MS
 - Grounded source allows easy interfacing



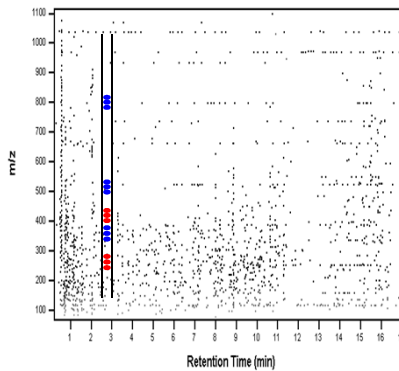
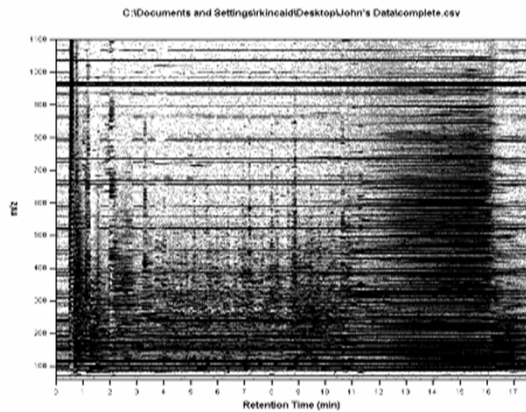
Agilent CE-ESI-MS Setup



Molecular Feature Extractor (MFE)

- How the MFE works

- Map signals in the 3-dim. space in time and mass at the MS level.
IMPORTANT: High resolution in mass and time!
- Remove areas which only contain noise, and NO signals.
- Identify all mass signals with a common RT (narrow time window)
- Combine mass signals with common RT and chemical relation (isotope, adduct, dimer, higher charge state) => **Molecular feature** or **Compound**.
- Create **Extracted Compound Chromatograms (ECC)** and **Compound Mass Spectra** from ions associated to a molecular feature (compound)
- Feature consists of an accurate neutral mass, a retention time, and a peak volume



RT	m/z	Abund
...
...
2.11	195.1745	21000
2.11	257.2566	33550
2.34	224.2134	11784
...
...



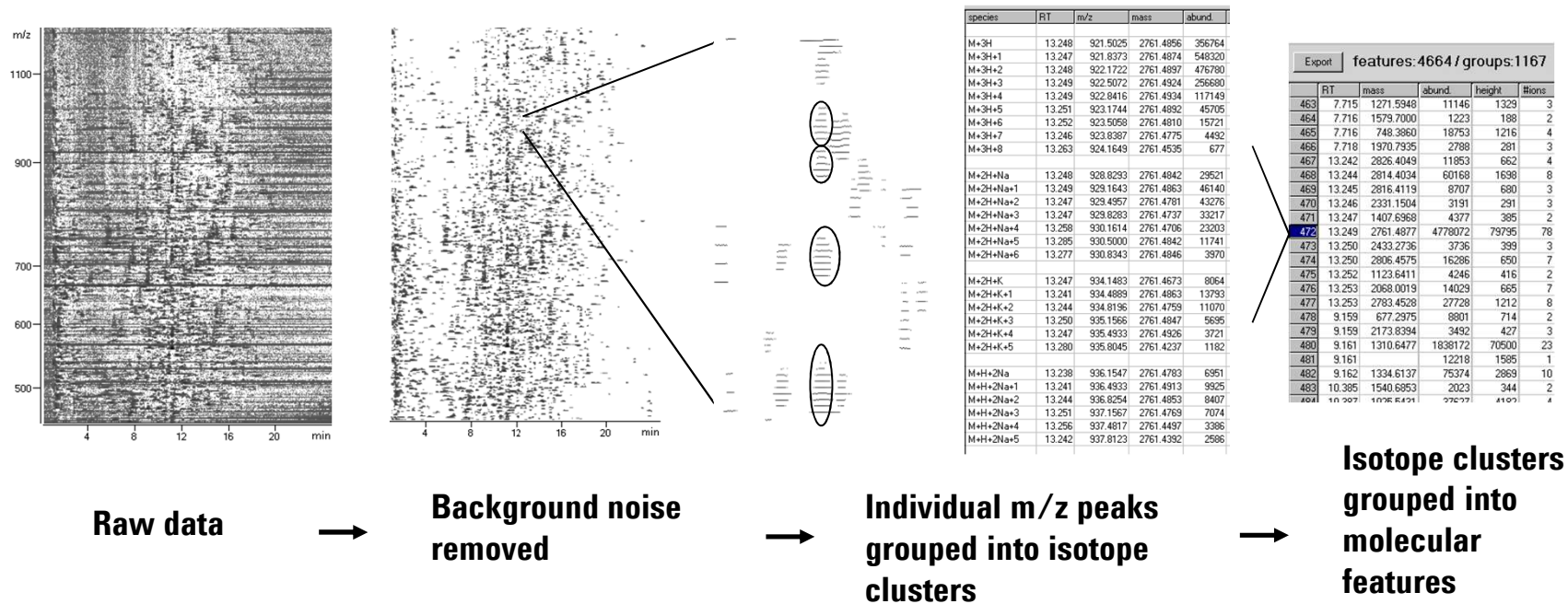
Export features:4664 / groups:1167

RT	mass	abund.	height	#ions
463	7.715	1271.5348	11146	1329
464	7.716	1579.7000	1223	188
465	7.716	748.3860	18753	1216
466	7.718	1970.7935	2788	281
467	13.242	2626.4049	11853	662
468	13.244	2814.4034	60168	1638
469	13.245	2816.4119	8707	680
470	13.246	2331.1504	3191	291
471	13.247	1407.6368	4377	385
472	13.249	2761.4877	4778072	79795
473	13.250	2433.2736	3736	399
474	13.250	2896.4575	16286	650
475	13.252	1123.6411	4246	416
476	13.253	2068.0019	14029	665
477	13.253	2783.4528	27728	1212
478	9.159	677.2975	8901	714
479	9.159	2173.8384	3492	427
480	9.161	1310.6477	1838172	70500
481	9.161		12218	1595
482	9.162	1334.6137	75374	2869
483	10.385	1540.6853	2023	344
484	10.397	1076.8491	97277	4103

Compound list or Molecular feature list

Isotope clusters grouped into molecular features

Data Analysis: Molecular Feature Extractor



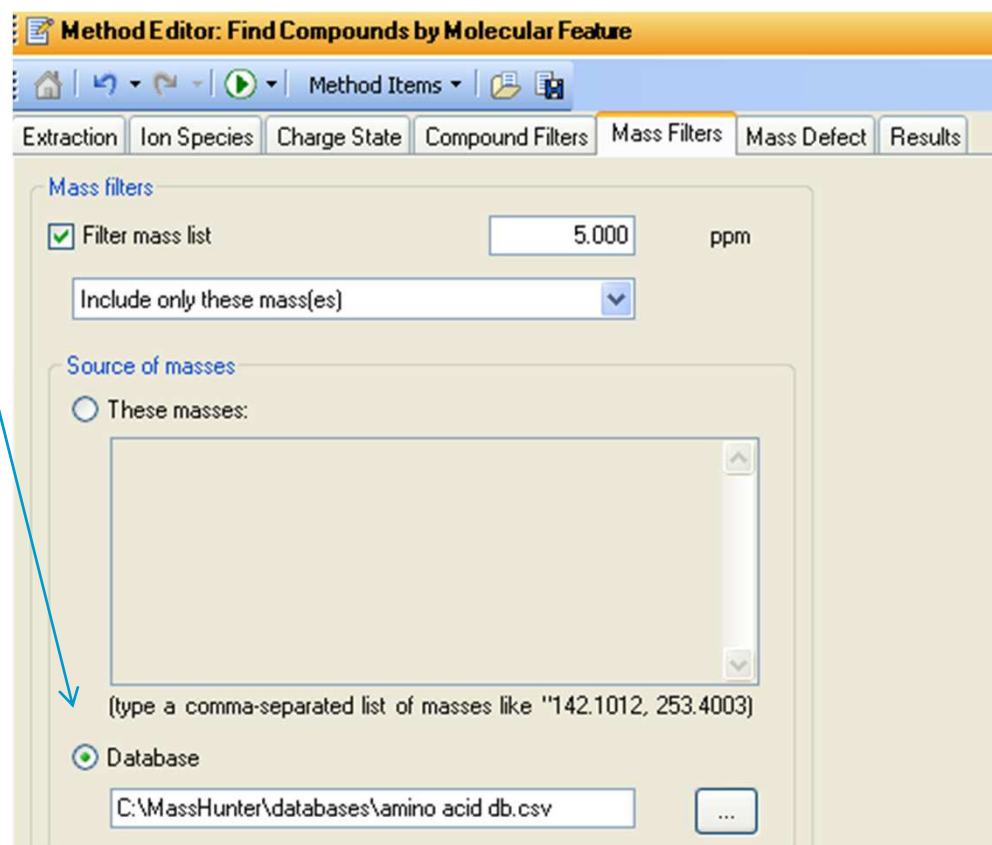
Identification, Quantification, Differential Analysis are performed on *chemically qualified compound data*

- Looks at Mass spectral data first → Groups co-eluting isotopes & converts to neutral mass
- Checks that there is a chromatographic response
- Groups all charge states, adducts, etc. associated with a given feature into a “peak volume”
- Feature consists of an accurate neutral mass, a retention time, and a peak volume

Data Analysis: Filter Molecular Feature Extractor Results by Amino Acid Database

Contents of Database (.csv file)

Formula	Mass	Compound name
C3H7NO2	89.0477	L-Alanine
C6H14N4O2	174.1117	L-Arginine
C4H7NO4	133.0375	L-Aspartic Acid
C6H12N2O4S2	240.0238	L-Cystine
C5H9NO4	147.0532	L-Glutamic Acid
C2H5NO2	75.032	Glycine
C6H9N3O2	155.0695	L-Histidine
C6H13NO2	131.0946	L-Isoleucine
C6H13NO2	131.0946	L-Leucine
C6H14N2O2	146.1055	L-Lysine
C5H11NO2S	149.051	L-Methionine
C9H11NO2	165.079	L-Phenylalanine
C5H9NO2	115.0633	L-Proline
C3H7NO3	105.0426	L-Serine
C4H9NO3	119.0582	L-Threonine
C9H11NO3	181.0739	L-Tyrosine
C5H11NO2	117.079	L-Valine



*Database contents only contain amino acids found in the standard mix

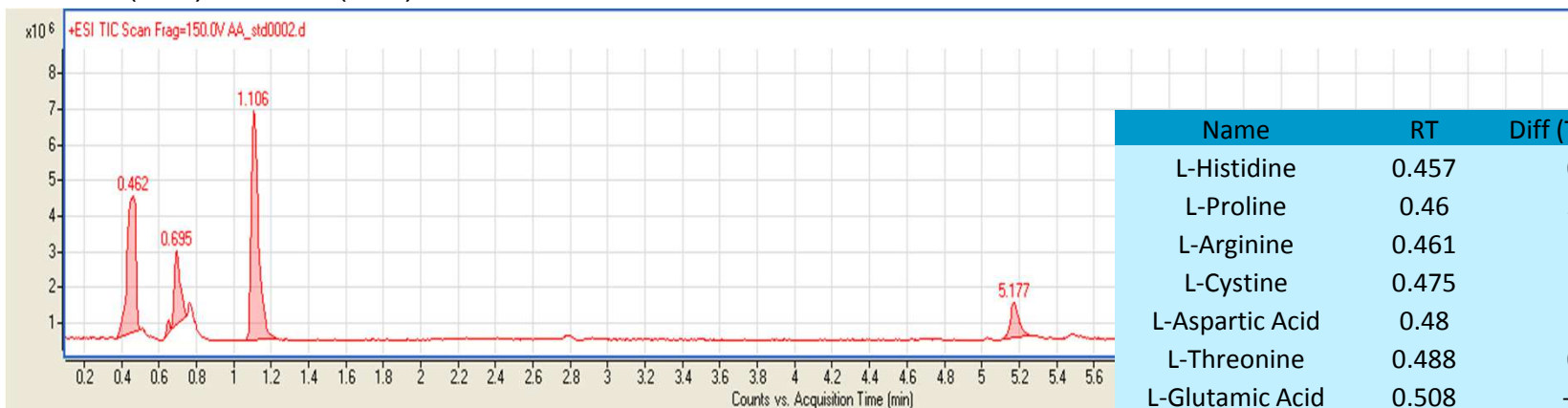
TOF Instrument Parameters for LC method:

Parameter	Setting
Source Type	Dual spray ESI in Positive mode
Gas temperature	325 °C
Drying Gas Flow rate	12 L/min
Nebulizer Pressure	45 psi
Capillary Voltage	4000 V
Fragmentor Voltage	150 V
Acquisition rate	3 scans/sec
Reference correction	1 point @ 922.009798
Acquisition Mode	2 GHz, extended dynamic range 50 to 1000 M/Z

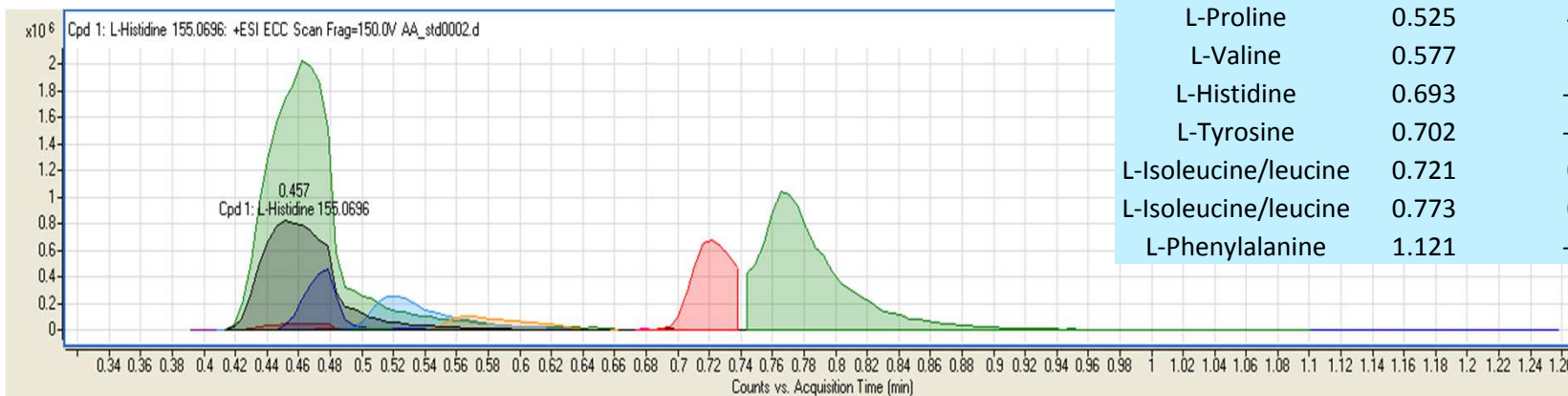
Initial LC Method Development (Amino Acid Standard)

Traditional chromatography using typical MS compatible mobile phases

- Zorbax HD SB C18 (1.8µm, 2.1x150mm) with 0.1% formic acid in water and 0.1% formic acid in ACN
- 95/5 (A/B) to 5/95 (A/B) over 10 minutes



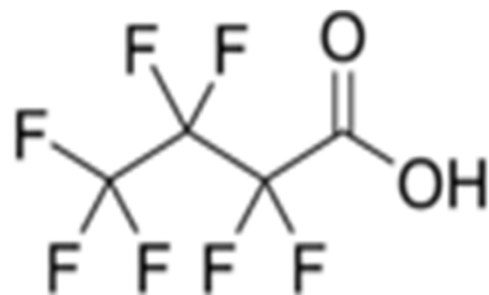
Name	RT	Diff (Tgt, ppm)	Vol %
L-Histidine	0.457	0.99	16.95
L-Proline	0.46	3.08	0.88
L-Arginine	0.461	1.14	39.12
L-Cystine	0.475	1.63	4.65
L-Aspartic Acid	0.48	2.82	0.17
L-Threonine	0.488	0.87	0.29
L-Glutamic Acid	0.508	-1.09	0.29
L-Proline	0.525	4.71	5.43
L-Valine	0.577	2.56	2.52
L-Histidine	0.693	-0.79	0.23
L-Tyrosine	0.702	-0.19	0.26
L-Isoleucine/leucine	0.721	0.19	7.94
L-Isoleucine/leucine	0.773	0.25	21.12
L-Phenylalanine	1.121	-2.28	0.17



- But underivatized amino acids are more polar than derivatized . . .
 - minimal retention and separation
 - Ion suppression (from co-elution) results in lack of sensitivity
 - Possible co-elution with sample matrix (eluting near void volume)

Change Modifier from Formic Acid to Heptafluorobutyric acid

HFBA has a variety of niche applications in analytical and synthetic chemistry. It is an ion pair reagent for reverse phase HPLC. It is used in the sequencing, synthesis, and solubilizing of proteins and peptides.



Final 1290 Infinity UHPLC Parameters

HFBA offers better retention/peak shape and is MS compatible

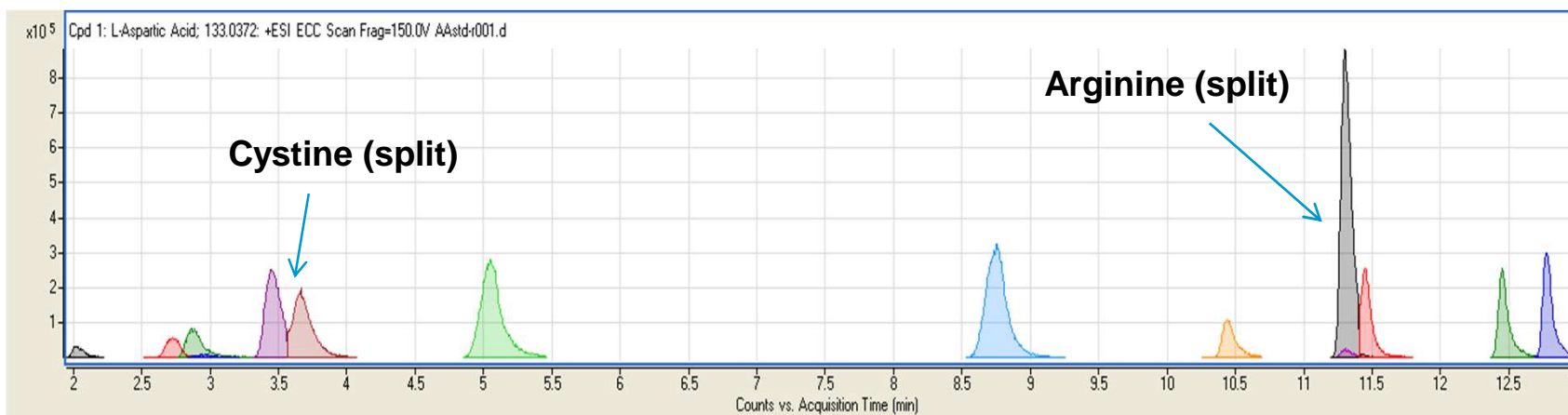
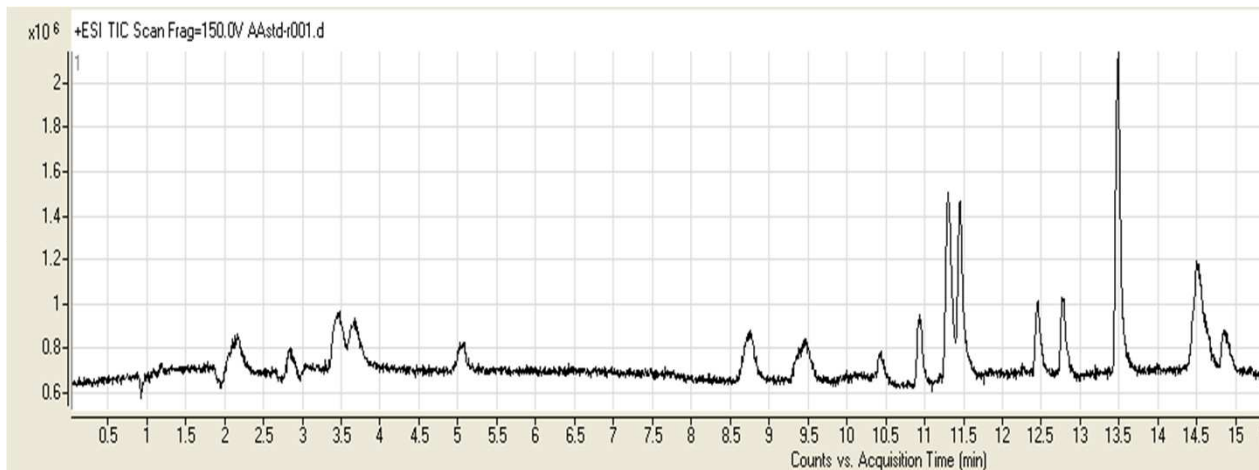
Parameter	Setting
Mixer	35 μ L jet weaver
Column	Zorbax HD SB C18 (1.8 μ m, 2.1x150mm)
Column Temperature	45 $^{\circ}$ C
Flow rate	0.3 mL/min
Mobile phase A	Water, 0.1 % HFBA
Mobile phase B	Acetonitrile, 0.1% HFBA
Injection Volume	3 μ L

	Time (min)	%A	%B
Gradient Program	0.0	100	0
	5.0	100	0
	15.0	70	30
	15.1	40	60
	18.0	40	60
	18.1	100	0
	21.0	100	0



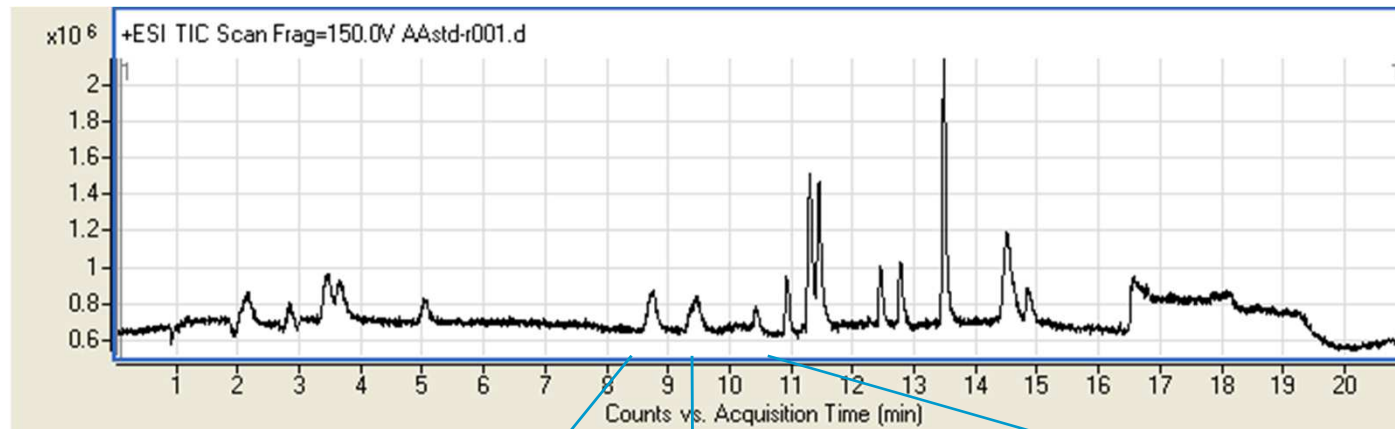
Final LC Conditions (Amino Acid Standard)

Name	RT	Diff (Tgt, ppm)	Vol %
L-Aspartic Acid	2.032	-2.48	0.86
L-Threonine	2.729	-3.13	1.92
L-Glutamic Acid	2.879	-1.97	3.26
L-Alanine	2.971	1.91	0.32
L-Cystine	3.456	0.83	9.14
L-Cystine	3.671	0.93	8.42
L-Proline	5.05	0.68	13.54
L-Histidine	8.747	-0.62	15.67
L-Valine	10.444	-1.15	3.42
L-Proline	11.306	-0.03	0.57
L-Arginine	11.307	-0.31	23.35
L-Arginine	11.455	0.4	6.27
L-Isoleucine/Leucine	12.455	-1.34	6.05
L-Isoleucine/Leucine	12.779	-1.28	7.21

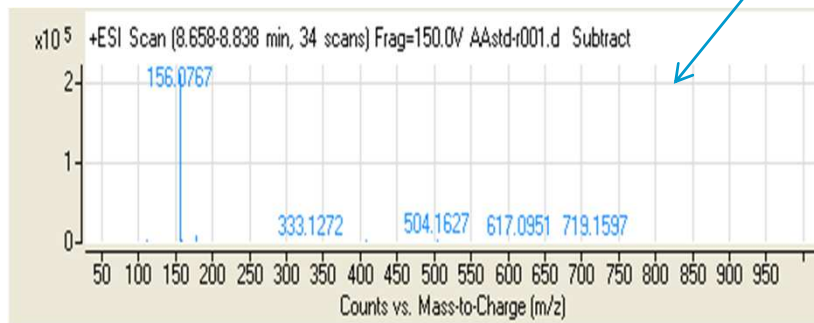


- Sufficient retention and separation
- Ion suppression from co-elution should be minimal
- Possible co-elution with sample matrix (near void volume) less likely

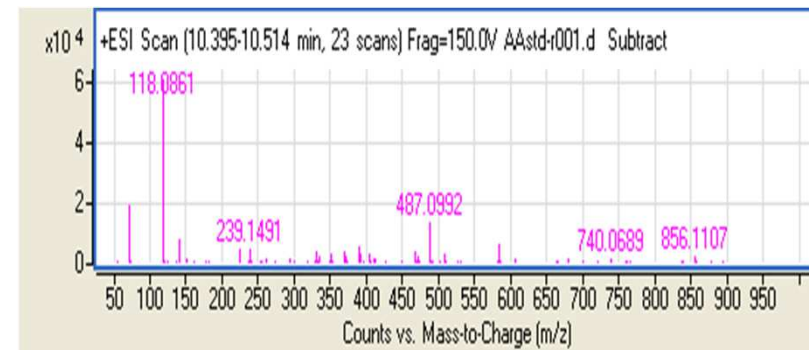
LC Method Example Spectra (Amino Acid Standard)



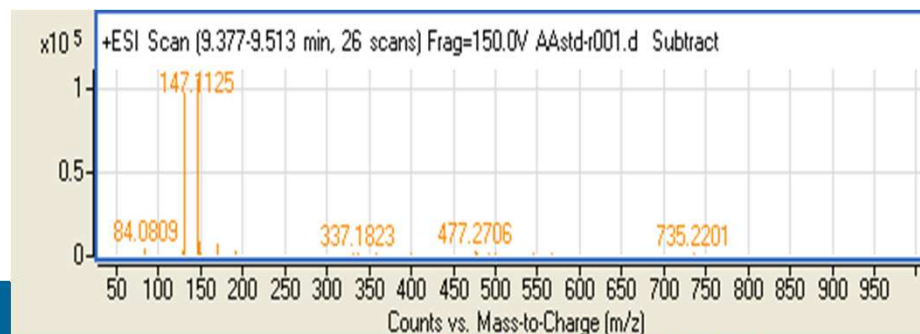
Histidine



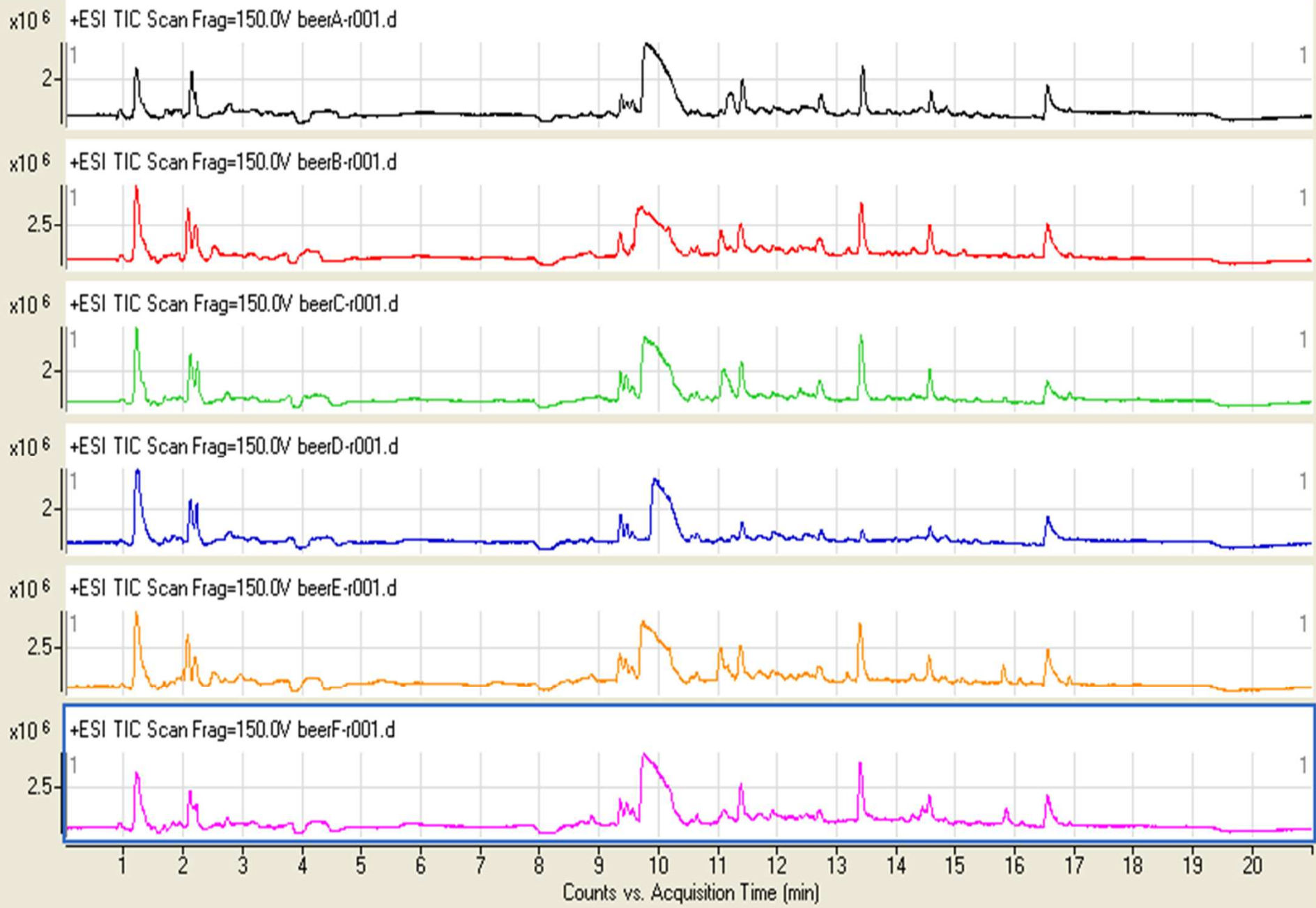
Valine



Lysine



TICs of Different Beer Brands by LC

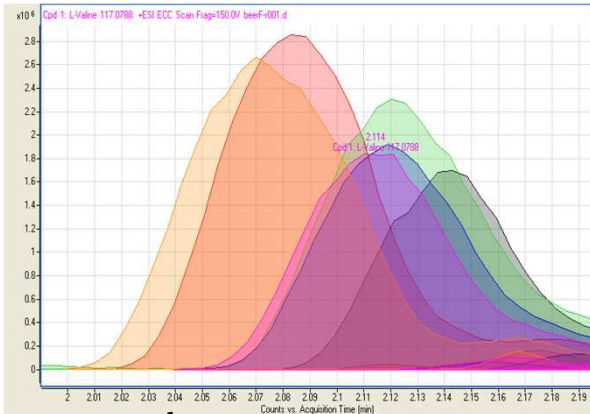


Overlay of Amino Acids In Different Beer Brands By LC

Present in all 6
Misidentified as Valine!
Valine is at 10.3min

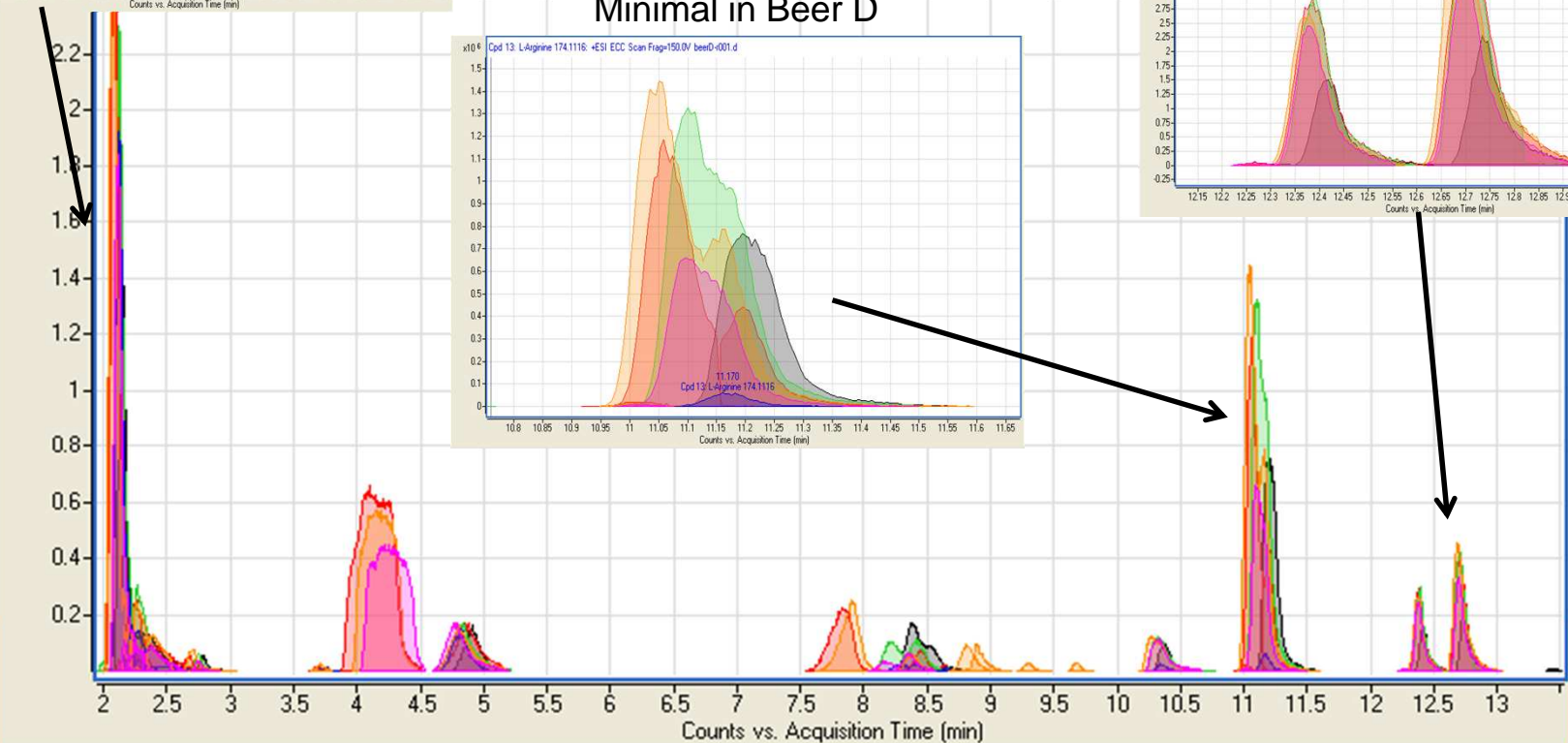
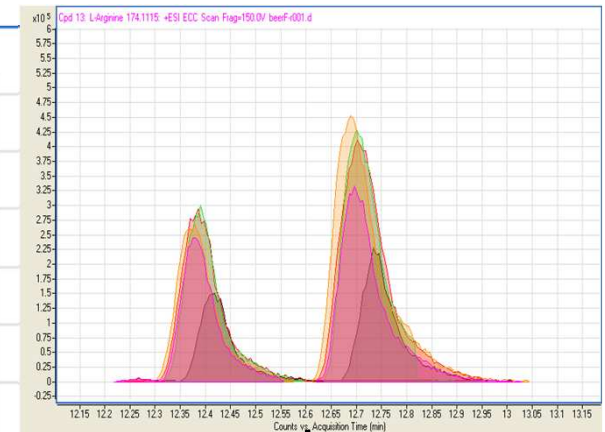
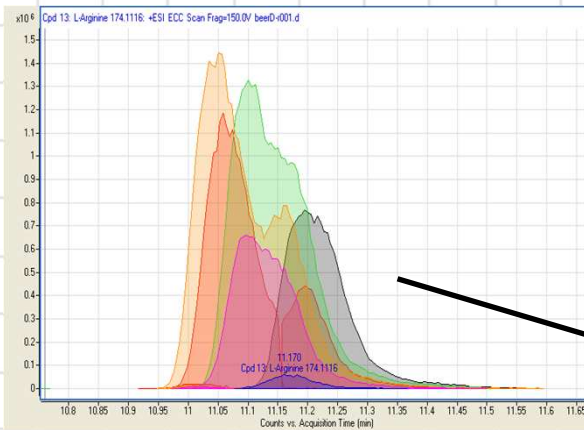
- Arginine has a bad peak shape
- Leucine and Isoleucine is baseline separated
- Overall peaks appeared tailed

Leucine and Isoleucine
present in 5 or 6 beers
Missing in Beer D



Scan Frag=150.0V beerF-r001.d

Arginine Present in all 6
Minimal in Beer D



7100 CE Parameters

Parameter	Setting
Capillary Type	50 μ m uncoated fused silica, 105 cm long
Cassette Temperature	20 $^{\circ}$ C
Voltage	20 kV
Buffer	1N Formic Acid
Injection	50 mBar for 10 sec
Sheath Buffer	50:50 water:Methanol 0.1% Formic Acid
Sheath Flow rate	5 μ L/min

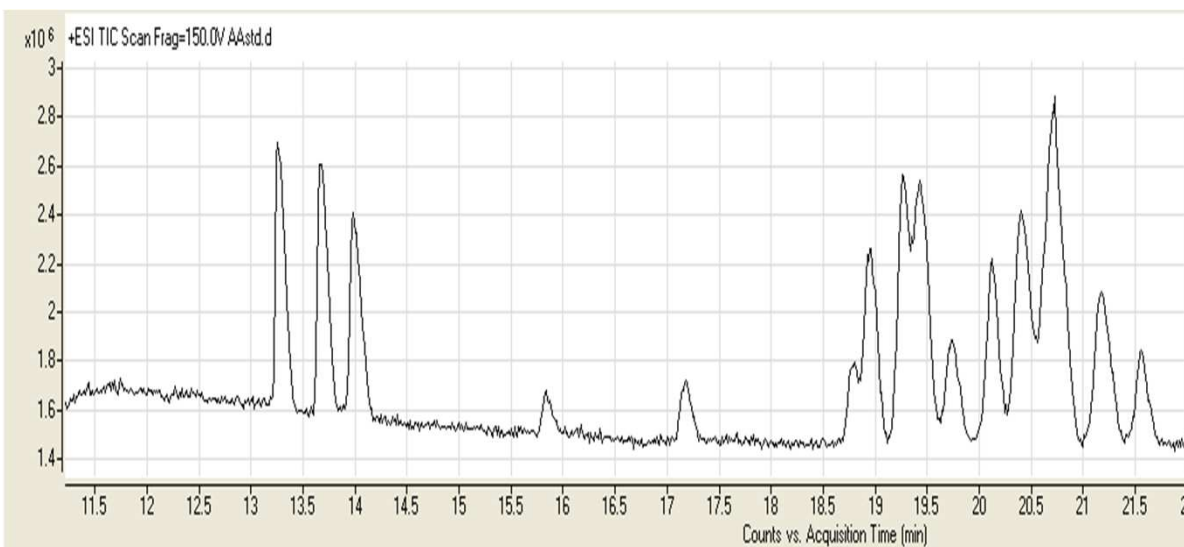


TOF Instrument Parameters for CE Method:

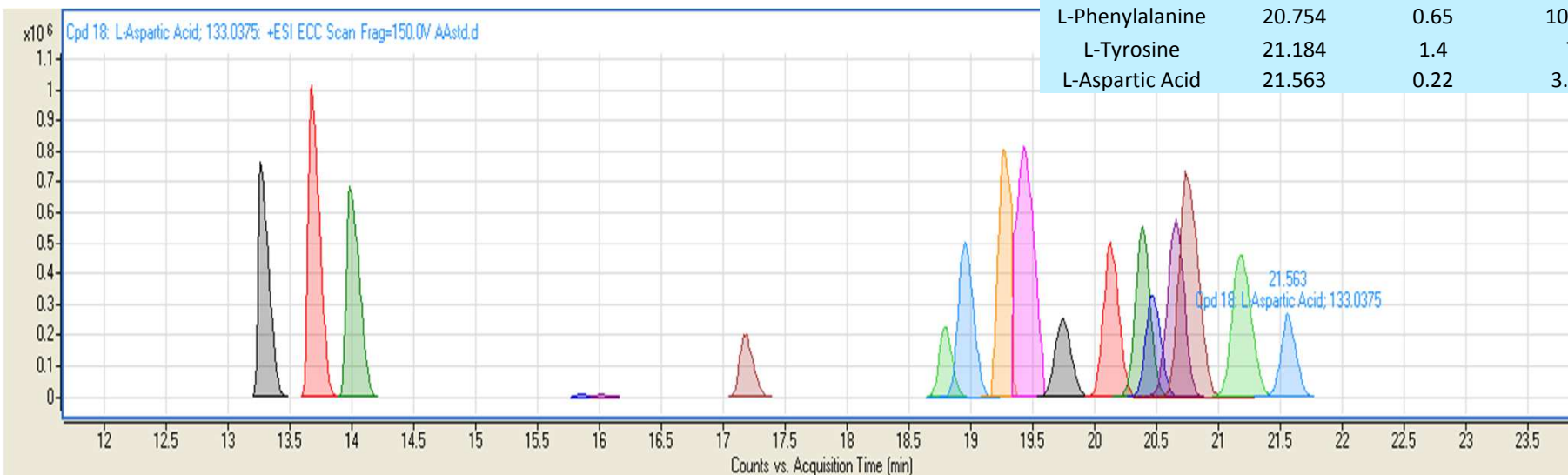
Parameter	Setting
Source Type	ESI in Positive mode
Gas temperature*	250 °C
Drying Gas Flow rate*	5 L/min
Nebulizer Pressure*	5 psi
Capillary Voltage	4000 V
Fragmentor Voltage	150 V
Acquisition rate	3 scans/sec
Reference correction	1 point @ 922.009798
Acquisition Mode	2 GHz, extended dynamic range 50 to 1000 M/Z

***These parameters need to be adjusted lower compared to the source conditions used for the LC method due to the lower flow rate (5 µL/min from the sheath pump vs 0.3 mL/min for the LC method)**

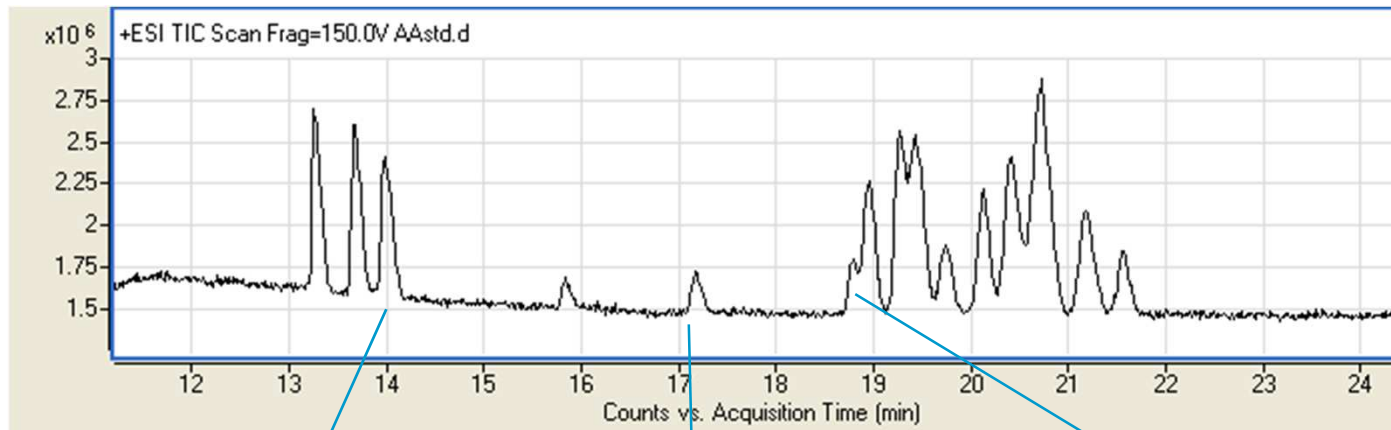
CE data: Amino Acid Standard



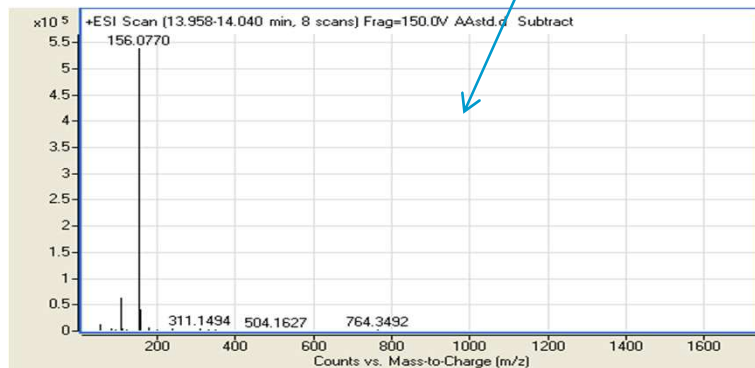
Name	RT	Diff (Tgt, ppm)	Vol %
L-Lysine	13.276	1.64	6.23
L-Arginine	13.684	2.08	8.88
L-Histidine	14	1.64	6.97
L-Alanine	17.18	4.85	2.11
L-Serine	18.792	1.32	2.28
L-Valine	18.955	1.4	5.92
L-Isoleucine/leucine	19.269	1.76	7.98
L-Isoleucine/leucine	19.436	1.82	11.33
L-Threonine	19.742	-2.87	3.21
L-Methionine	20.13	1.21	5.95
L-Proline	20.387	1.77	6.38
L-Glutamic Acid	20.47	0.86	3.98
L-Cystine	20.654	1.59	7.59
L-Phenylalanine	20.754	0.65	10.89
L-Tyrosine	21.184	1.4	7
L-Aspartic Acid	21.563	0.22	3.16



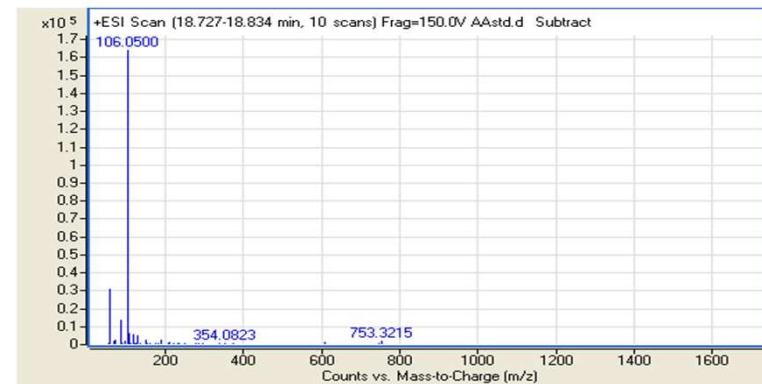
CE Method Example Spectra (Amino Acid Standard)



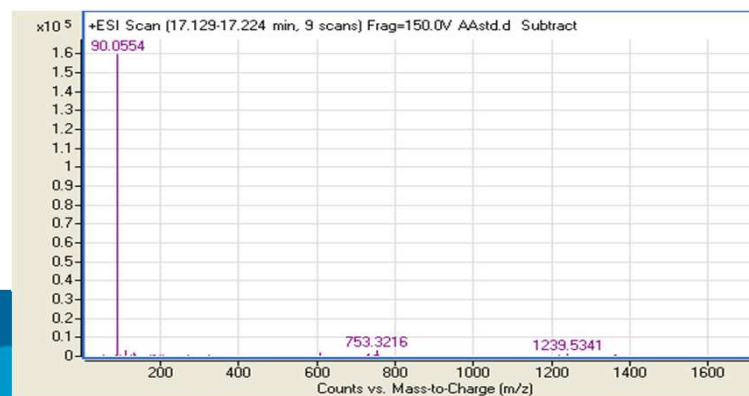
Histidine



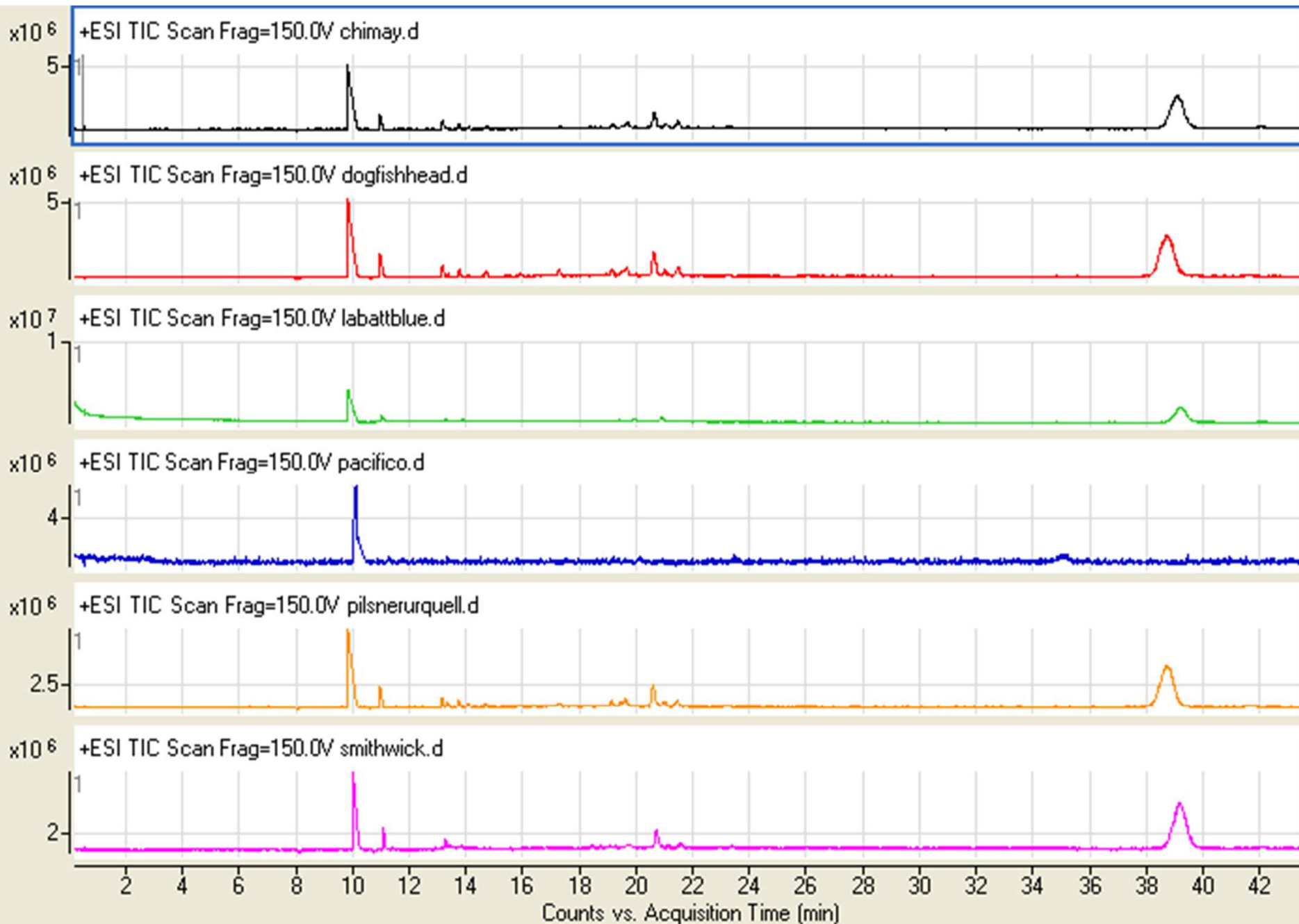
Serine



Alanine



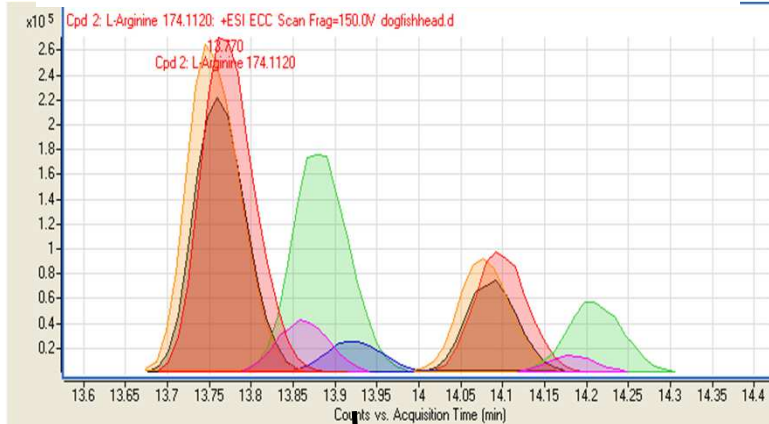
TIC of Different Beer Brands by CE



Overlay of Amino Acids In Different Beer Brands by CE

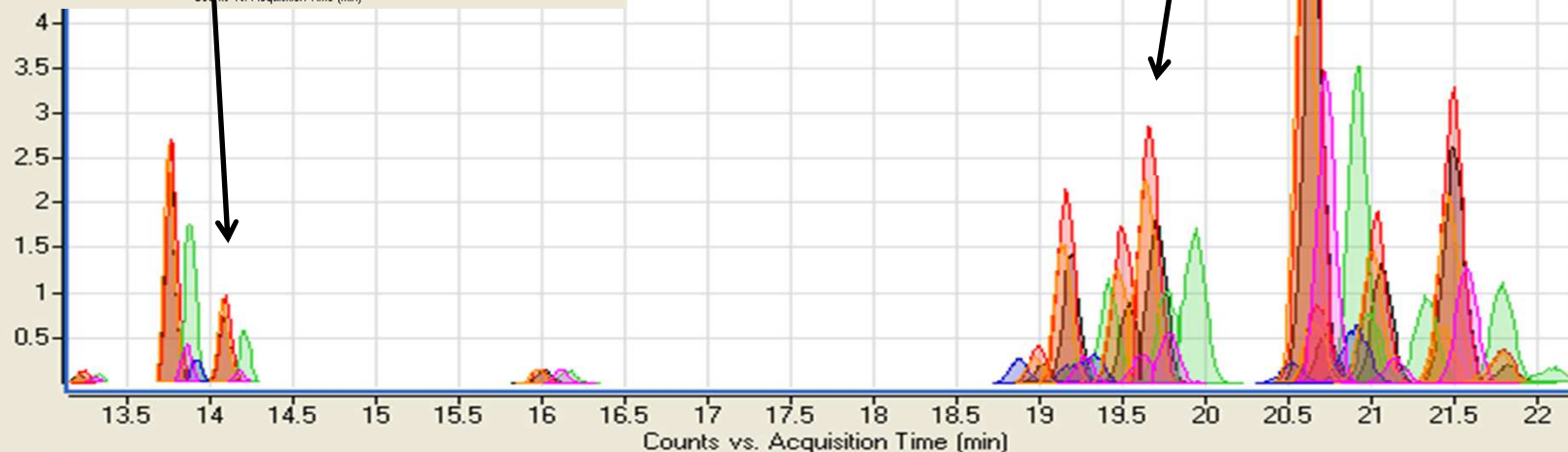
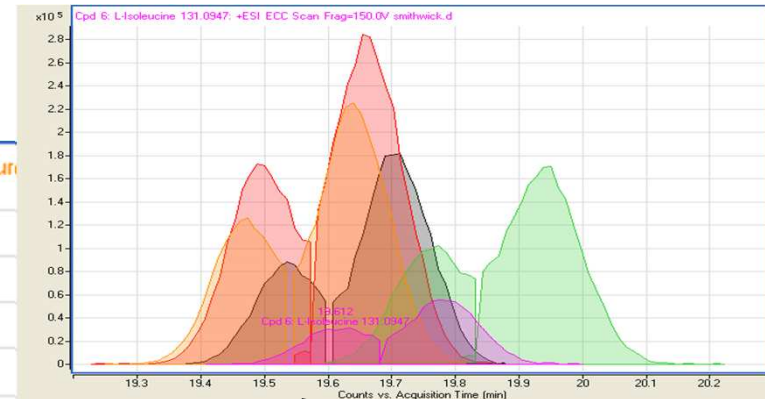
- Peak shapes are more symmetrical
- Leucine and Isoleucine not fully baseline separated
- Retention times have a higher %RSD

Arginine (present in all 6, minimal in D)
Histidine (present in 5 of 6 beers,
missing in Beer D)



Scan Frag=150.0V pilsnerun

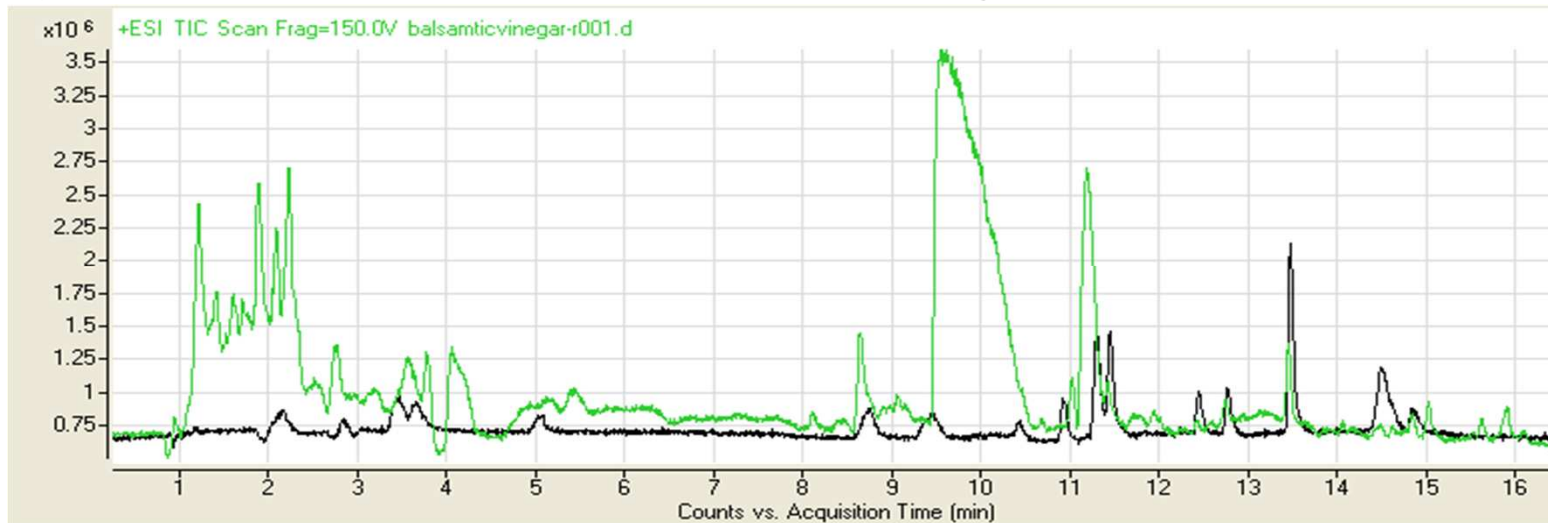
Leucine and Isoleucine
present in 5 or 6 beers
Missing in Beer D



Comparison of TICs for Balsamic Vinegar

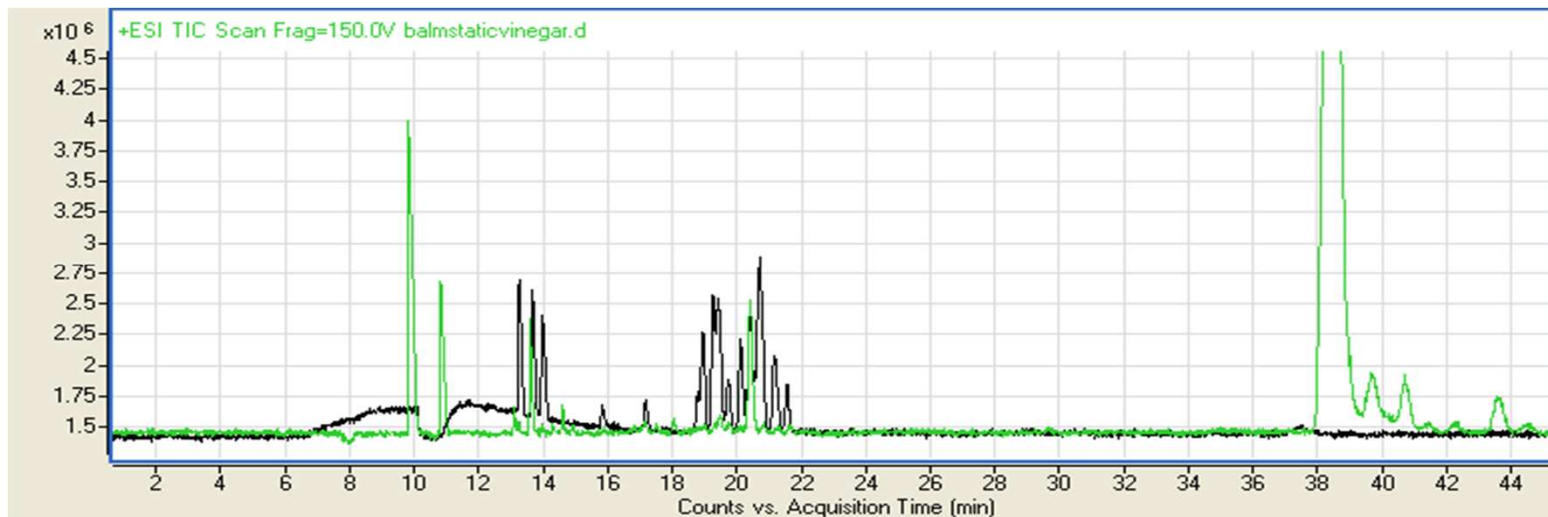
Overlay of Amino Acid Standard (Black) with Balsamic Vinegar (Green)

LC



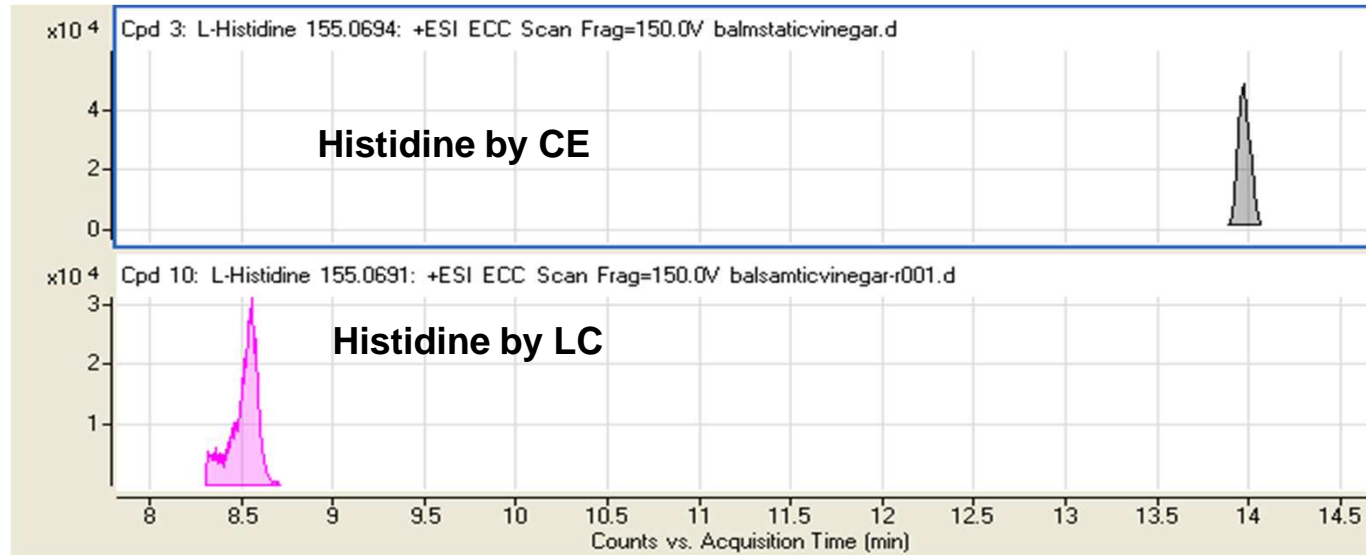
For LC the interferences are spread out across the run where amino acids are eluting
For CE there are less interferences from the sample matrix (comes early and late in the run)

CE

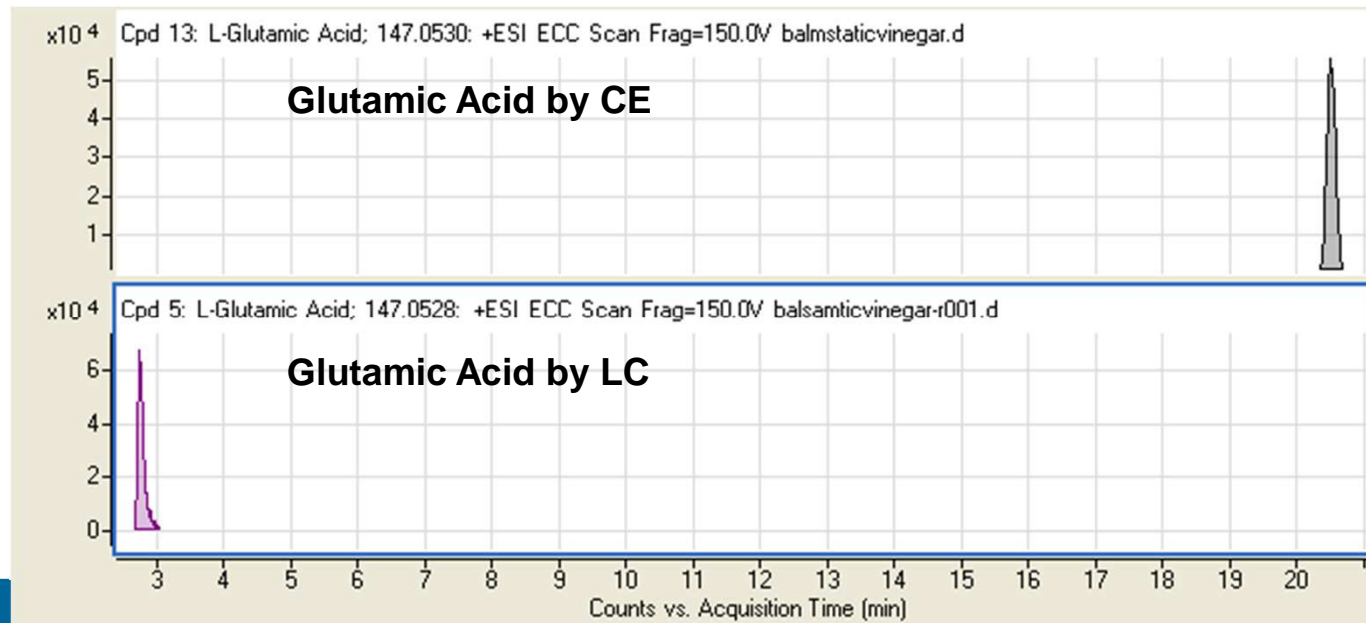


Comparison of Peak Shapes for Balsamic Vinegar

Example 1



Example 2



Comparison of Leucine/Isoleucine Separation

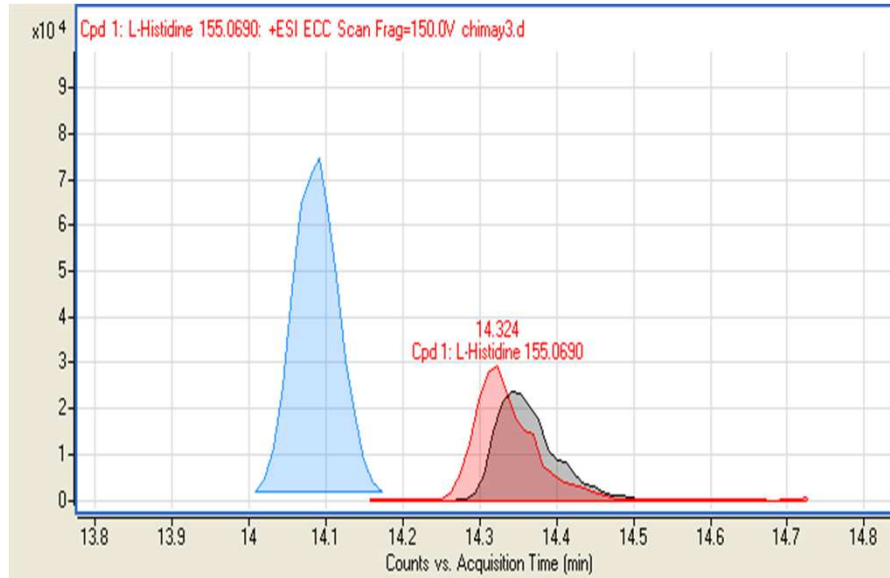


LC baseline
separated

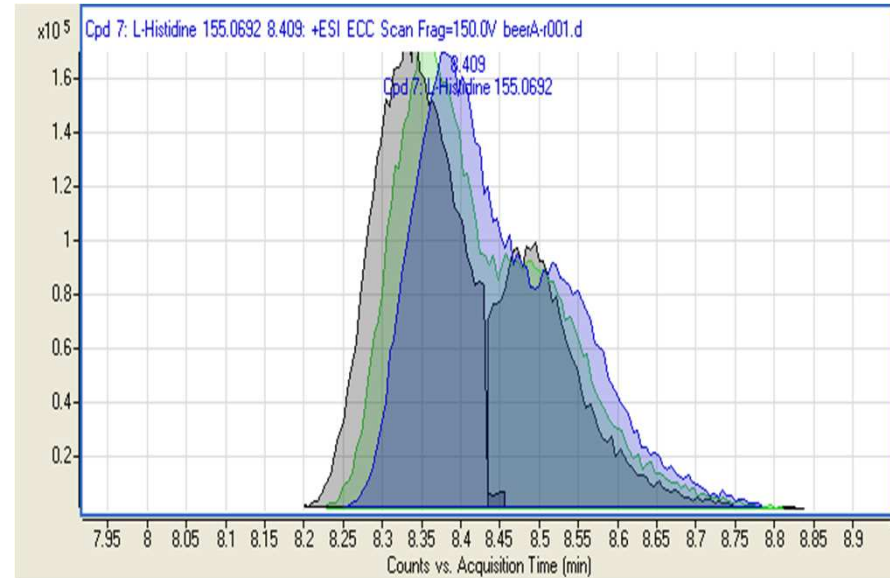
CE partially
resolved

Comparison of Migration/Retention Time Precision

CE: Histidine for 3 replicate injections



LC: Histidine for 3 replicate injections



	CE RT (min)	LC RT (min)
Run 1	14.324	8.409
Run 2	14.321	8.382
Run 3	14.087	8.336
Average	14.244	8.375667
Std Dev	0.135974	0.03691
% RSD	0.954607	0.440679

Conclusion

LC method Advantages

- Easier to setup and implement
- User's have more familiarity with LC
- Faster run and cycle time
- Better retention time precision
- Leucine and Isoleucine are fully baseline separated

CE method Advantages

- Amino acids are better separated from matrix effects
- Less false positive identifications
- Sample cleanup/preparation is less involved
- Peak shapes are more symmetrical