



Tips & Tricks to Achieve Fast, Sensitive, and Reproducible Separation of Amino Acids

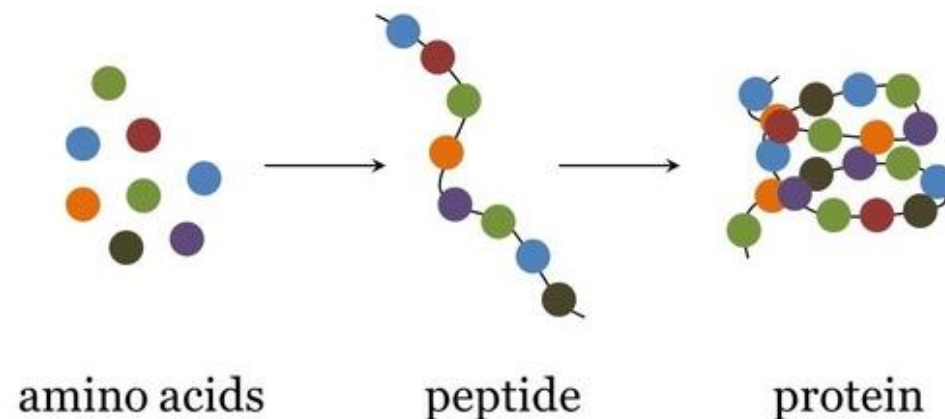
Anne Blackwell, Ph.D.

Biocolumns Product Support Scientist

September 12, 2017

Some Basics

Amino Acids are the building blocks of proteins



Amino Acids require derivatization to be detected by UV or FL

- OPA/FMOC, Ninhydrin, Dansyl chloride, and PITC are common reagents used

Derivatization can be done pre-column or post column

- OPA/FMOC, Dansyl chloride, and PITC are common reagents used for pre-column
- Ninhydrin is common for post column methods

Analysis of AA can be done by several methods:

- GC, CE, HPAE-PAD
- LC/UV/FL, LC/MS

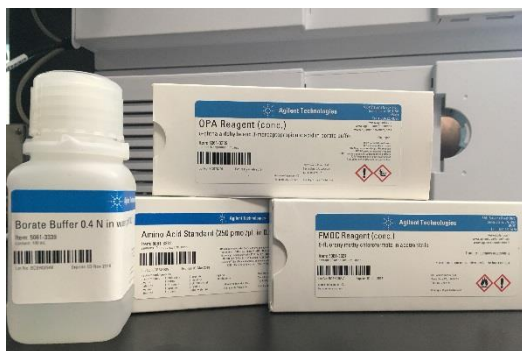


Why is Amino acid analysis important?

- Important for protein and peptide identification and quantitation
- Part of reverse-phase characterization in biopharma
- Required by the FDA
- Important for monitoring cell culture media
- Used for the analysis of metabolic intermediates - “Bound vs. Free”



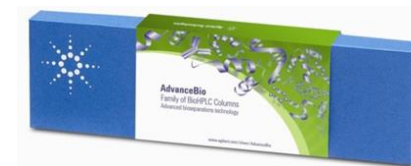
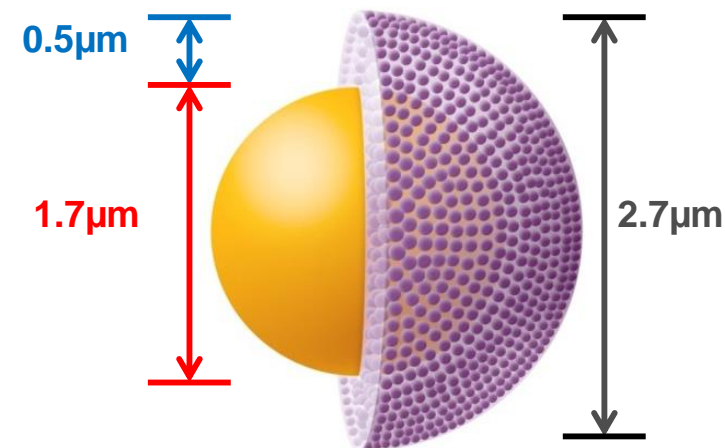
The Agilent Amino Acid Analysis solution



Ready to use
AdvanceBio AAA kit
(Standards and
Reagents)



All Agilent LC
systems including
Infinity II systems



AdvanceBio AAA
Columns
Fast and rugged



Agilent Technologies

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

Agilent AdvanceBio AAA

Previous Agilent AAA Method

Agilent has a well established solution for Amino Acid Analysis

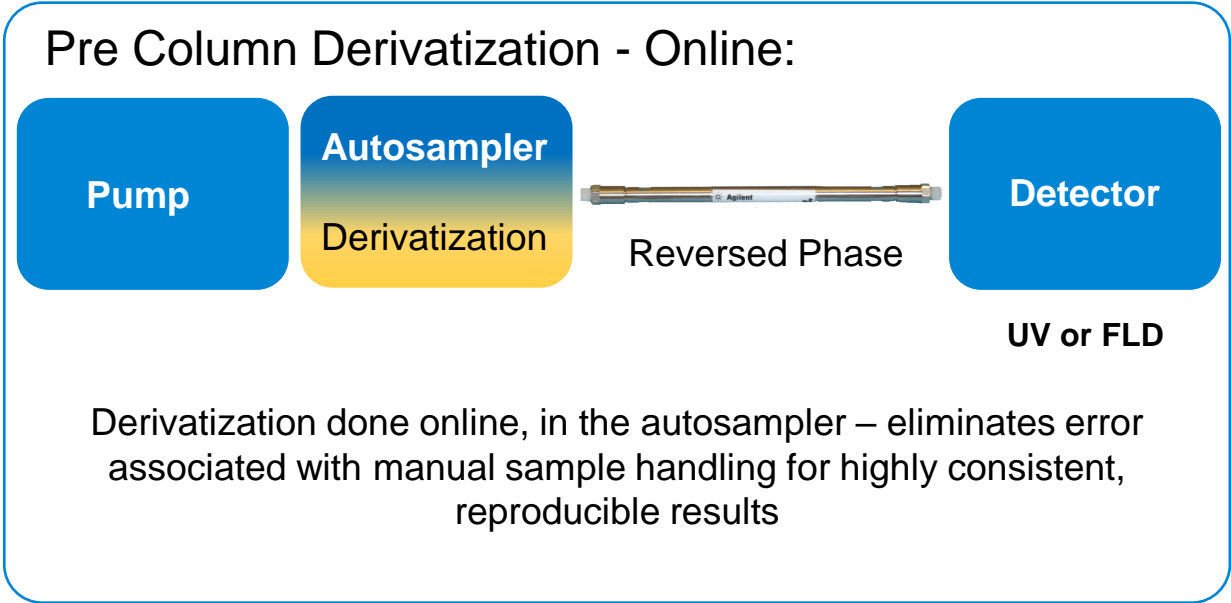
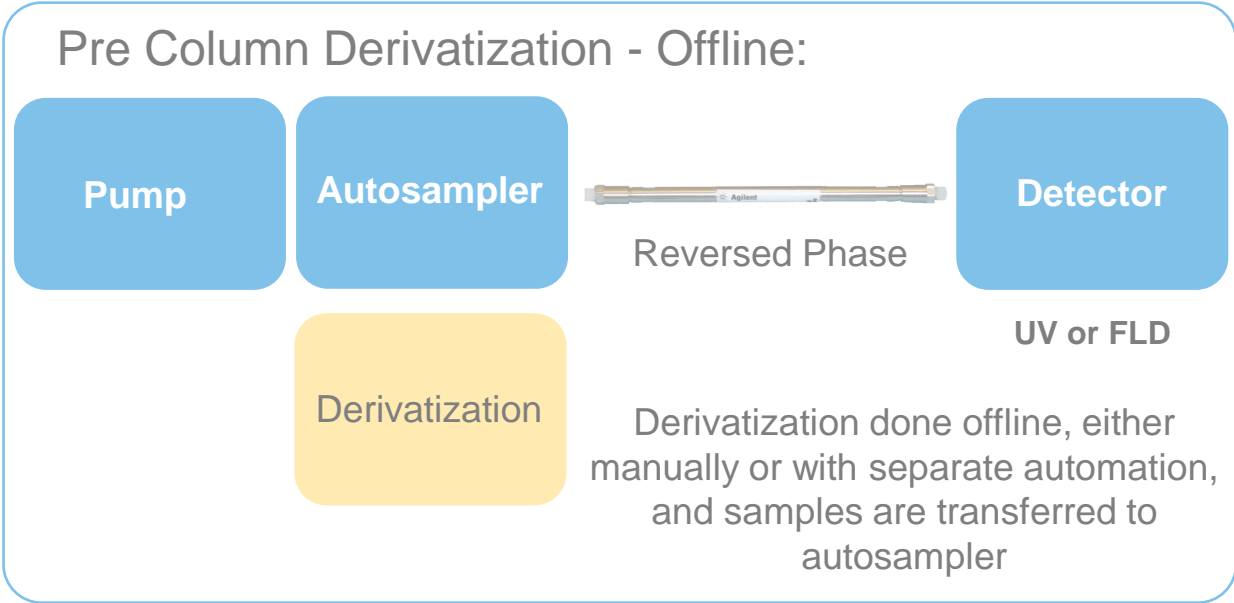
- based on automated pre-column derivatization capabilities of Agilent Autosamplers
- Uses ZORBAX Eclipse AAA column
- Well established method using reagents and standards from Agilent

What's New?

- All reagents conveniently kitted together under a single part number
- Introduced an HpH chemistry on a Poroshell particle for improved column lifetime
 - Traditional silica columns dissolve above neutral pH, but HpH chemistry stabilizes column
 - AA derivatization and separation are most efficient at higher pH
 - Poroshell column with 2 μm frits is less susceptible to clogging

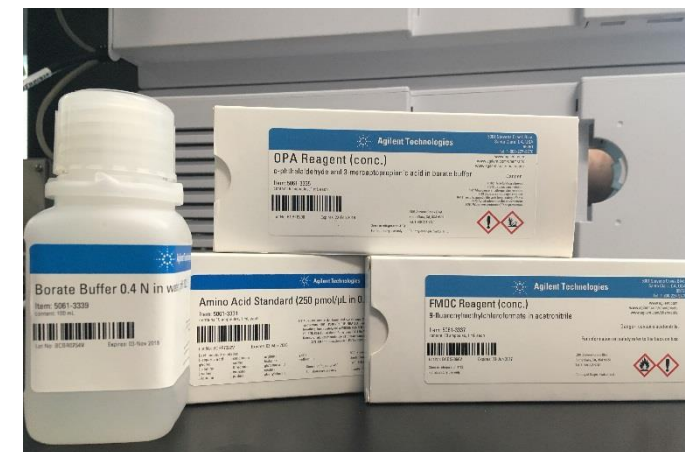


Pre- vs Post-Column Derivatization



AdvanceBio AAA Reagent Kit

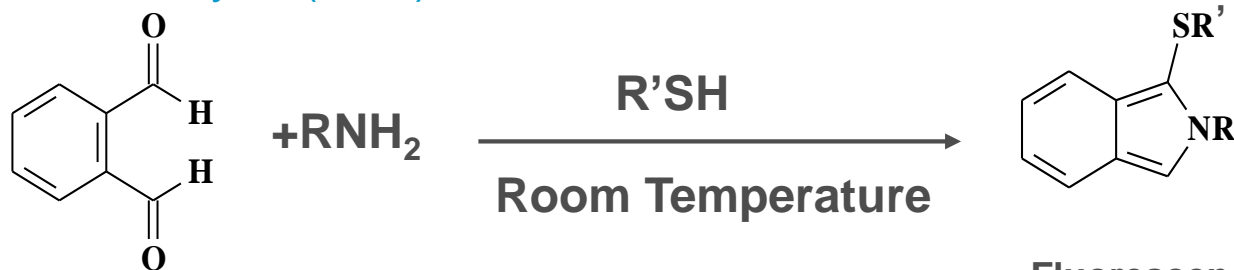
Part Number	Material
5061-3339	100mL Borate Buffer
5061-3337	FMOC reagent - 10 ampoules, 1 mL each
5061-3335	OPA reagent, 10 mg/mL, 6 ampoules
5062-2479	Dithiodipropionic acid (DTDPA)
5061-3330	AA, standard 1nmol 10/PK
5061-3331	AA standards, 250 pmol 10/PK
5061-3332	AA standards, 100 pmol 10/PK
5061-3333	AA standards, 25 pmol 10/PK
5061-3334	AA standards, 10 pmol 10/PK
5062-2478	AA supplements, 1g each



Order components individually, or together as part of a kit with a single part number (5190-9426)

Automated Derivatization in the Autosampler

Ortho Phthalaldehyde (OPA)



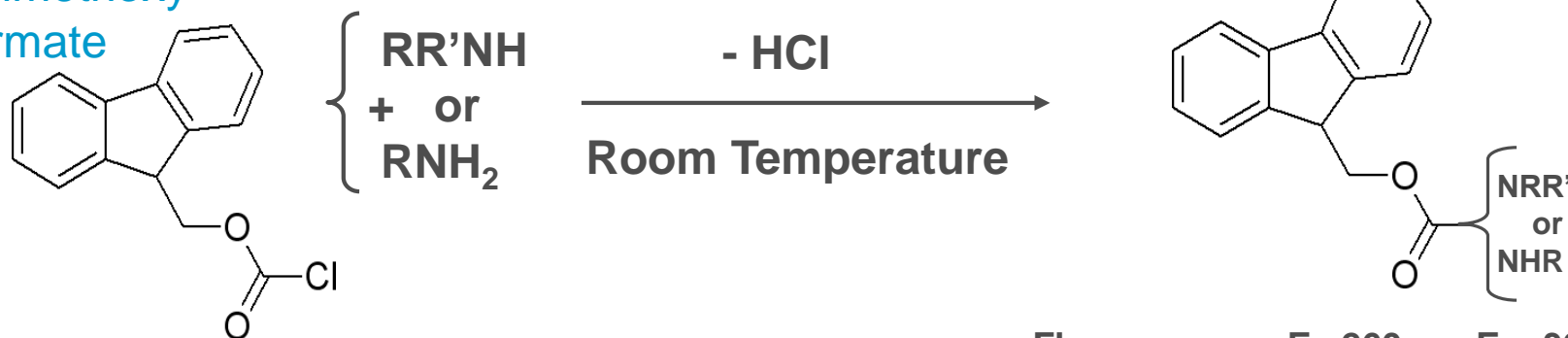
Non-fluorescent
Does not absorb at 338nm

Fluorescence: Ex 340nm, Em 450nm
DAD: 338, 10nm; Ref. 390, 20nm

1. Allows visualization by UV or FL
2. Helps retain very polar compounds



Fluorenylmethoxy chloroformate (FMOC)



Fluorescent
Absorbs at 262nm and
Fluorescences at 324nm

Fluorescence: Ex 260nm, Em 325nm
DAD: 262, 16nm; Ref. 324, 8nm

Optimal pH for reaction with AA: ~10.0



Agilent Technologies

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

Online derivatization/Injection program

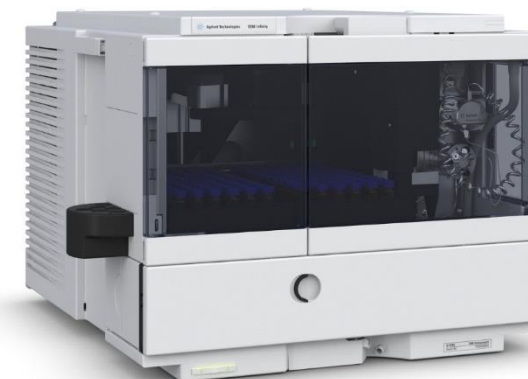
- Draw 2.5 μL from borate vial (Agilent p/n 5061-3339)
- Draw 1.0 μL from sample vial
- Mix 3.5 μL in wash port 5 times
- Wait 0.2 min
- Draw 0.5 μL from OPA vial (Agilent p/n 5061-3335)
- Mix 4.0 μL in wash port 10 times default speed
- Draw 0.4 μL from FMOC vial (Agilent p/n 5061-3337)
- Mix 4.4 μL in wash port 10 times default speed
- Draw 32 μL from injection diluent vial
- Mix 20 μL in wash port 8 times
- Inject
- Wait 0.1 min
- Valve bypass

1290 Infinity II Multisampler



Method can be programmed into ANY Agilent autosampler –

- Eliminates manual labor and variability
- Enables highly precise data

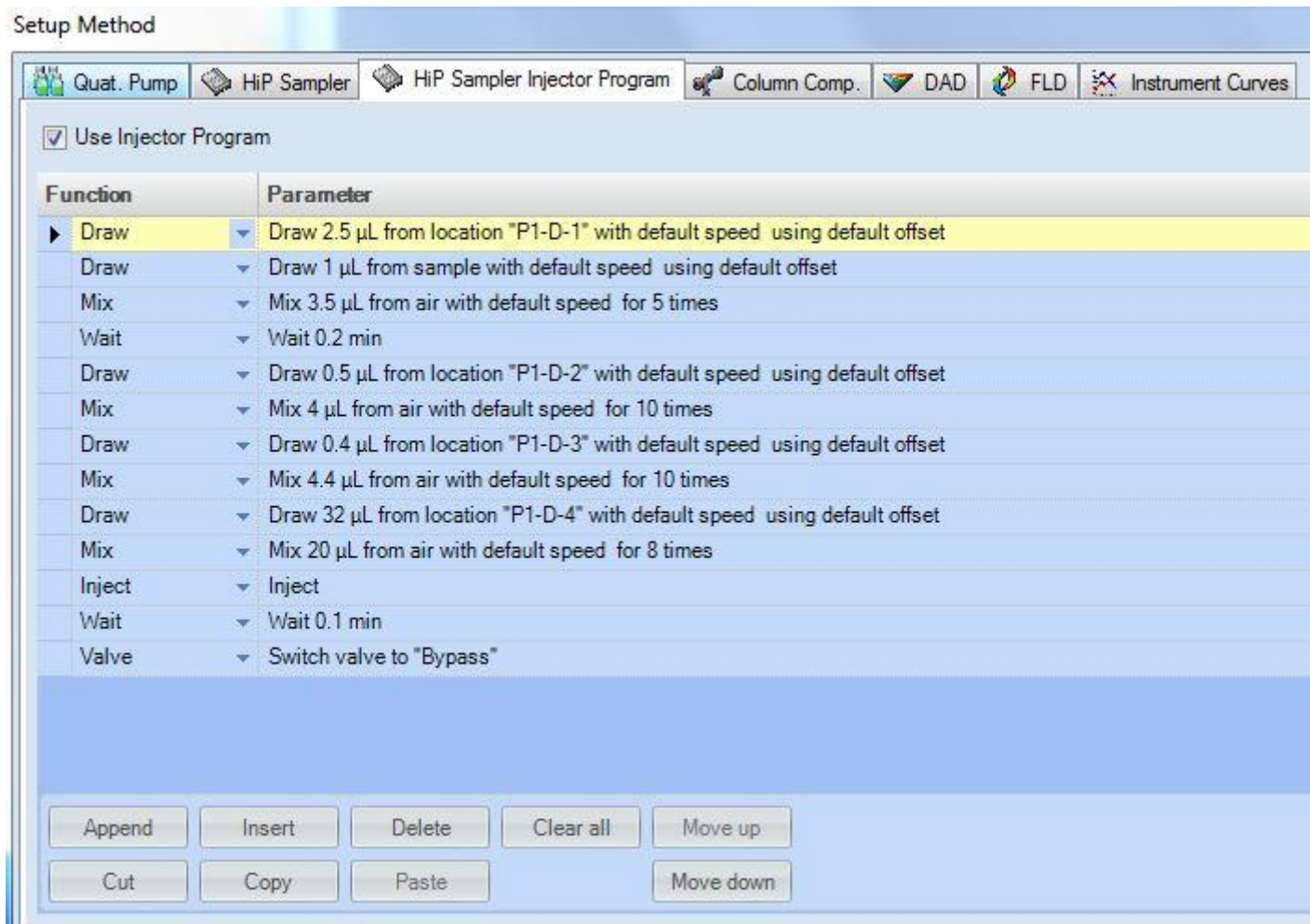


1260 Infinity II Vialsampler



Online derivatization/Injection program

OpenLab ChemStation C.01

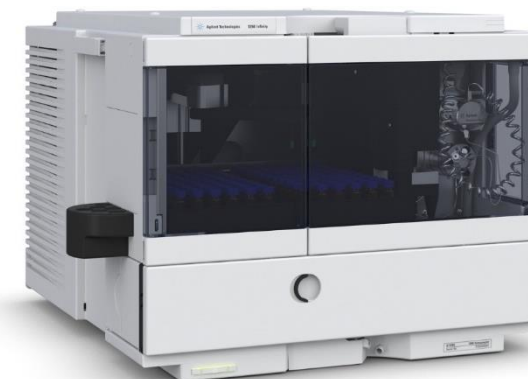


1290 Infinity II Multisampler



Method can be programmed into ANY Agilent autosampler –

- Eliminates manual labor and variability
- Enables highly precise data



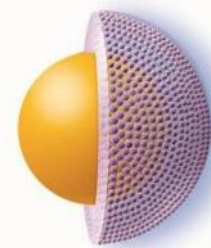
1260 Infinity II Vialsampler



Agilent Technologies

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

Robust Columns for AAA



A robust, high efficiency Fast LC column with resistance to elevated pH and temperature offering users performance comparable to that of sub-2 μm alternatives but with up to 50% less back pressure.

Utilizes a proprietary technology for particle synthesis



Core



P120 Particle



Treated P120

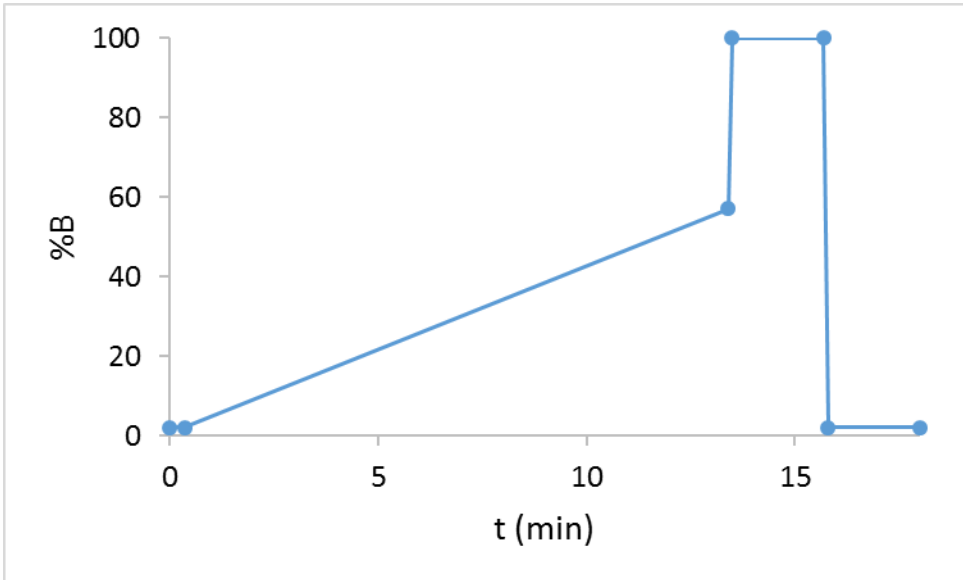
- 2.7 μm particles, 110 Å pore size
- Two dimensions available: 3.0 x 100 mm, 4.6 x 100 mm
 - Guard columns also available in each i.d.
- Each individual column is tested for efficiency
- Each batch is tested with amino acid standards to ensure performance



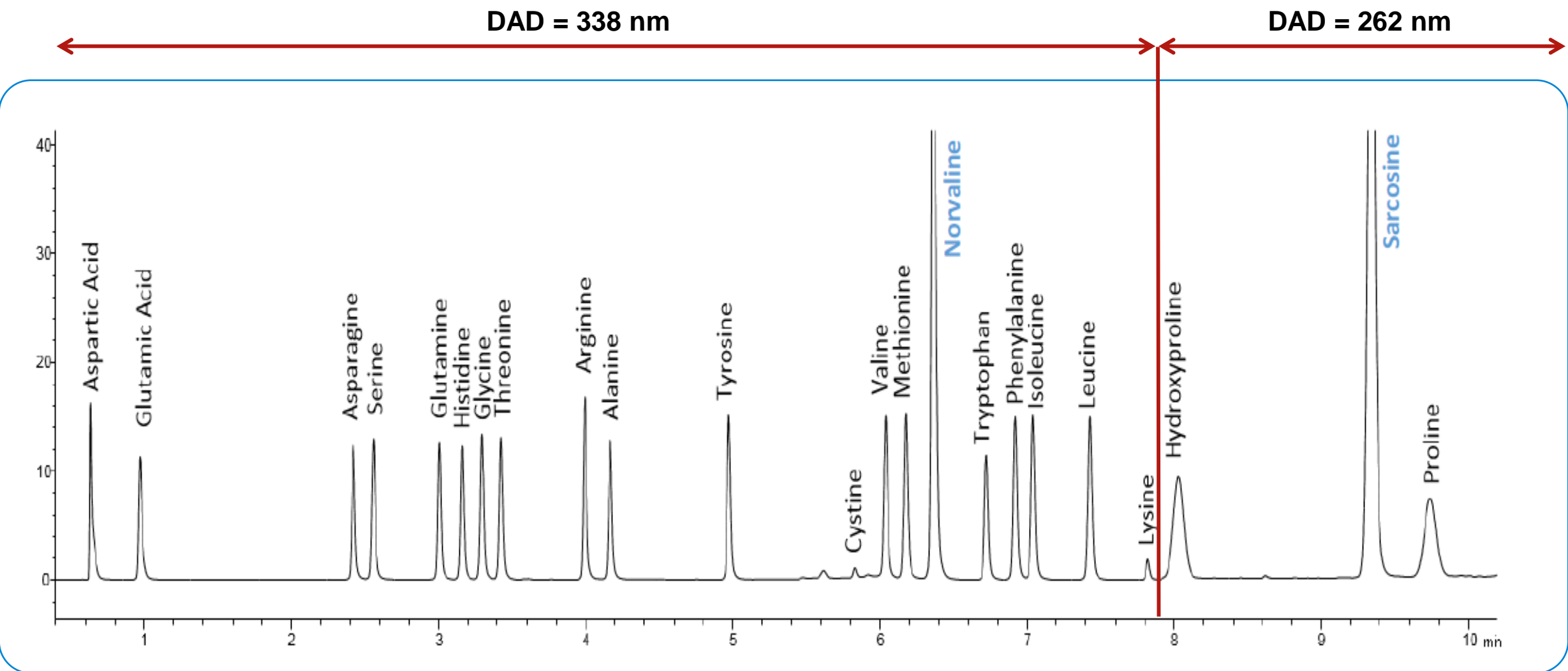
Chromatographic Method

- Flow rate – 1.5 mL/min for 4.6 mm and 0.62 mL/min for 3 mm i.d.
- Injection volume – 1µL with needle wash at the wash port for 7s
- Column temperature – 40 °C
- Detection wavelength – 338 and 262nm
- Samples- Agilent AAA standards, media samples and protein hydrolysate standards

Time (min)	%B
0	2
0.35	2
13.4	57
13.5	100
15.7	100
15.8	2
18	stop



Fast and Rugged Amino Acids Separation

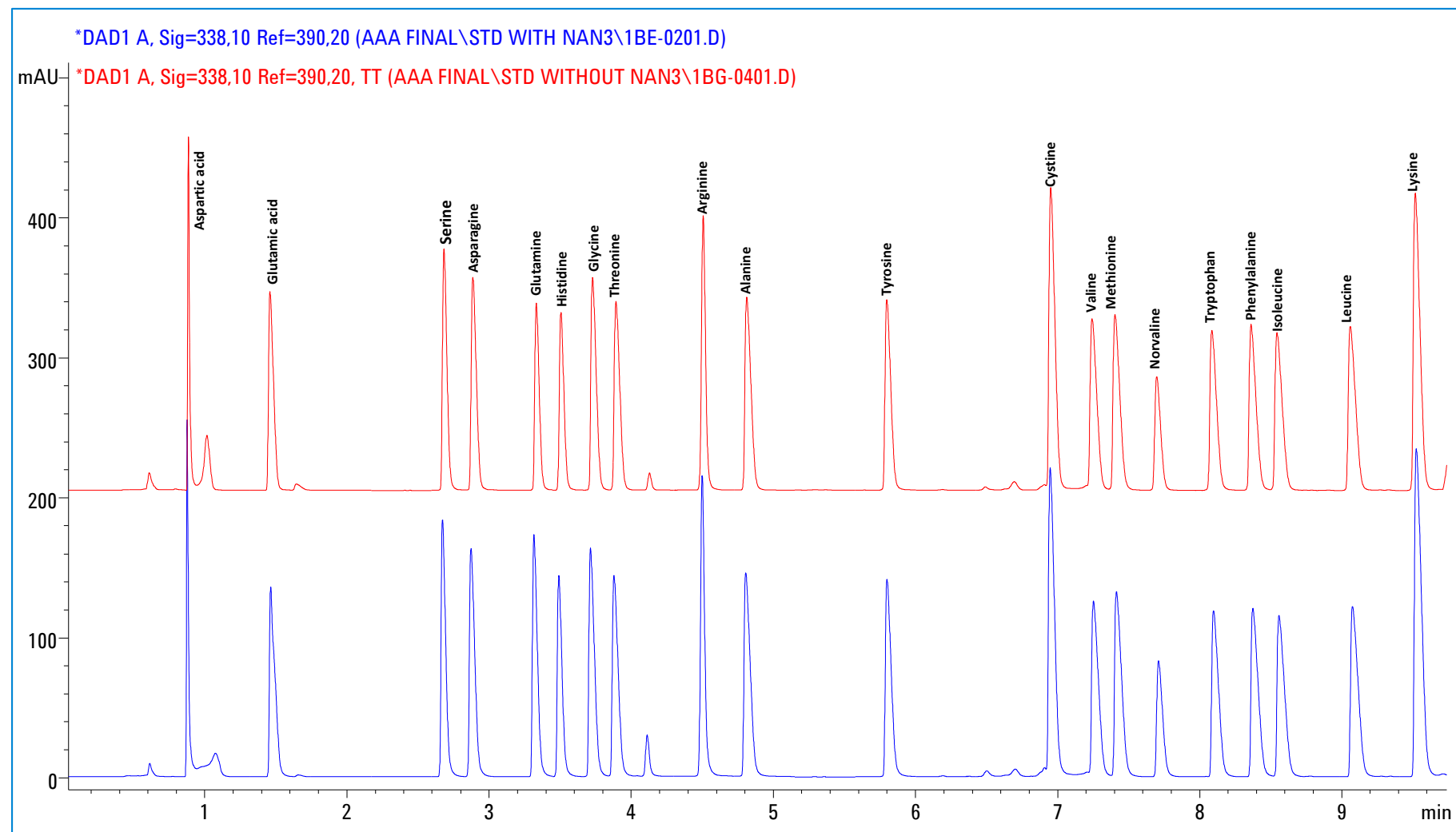


Order of Elution for OPA and FMOC derivatives

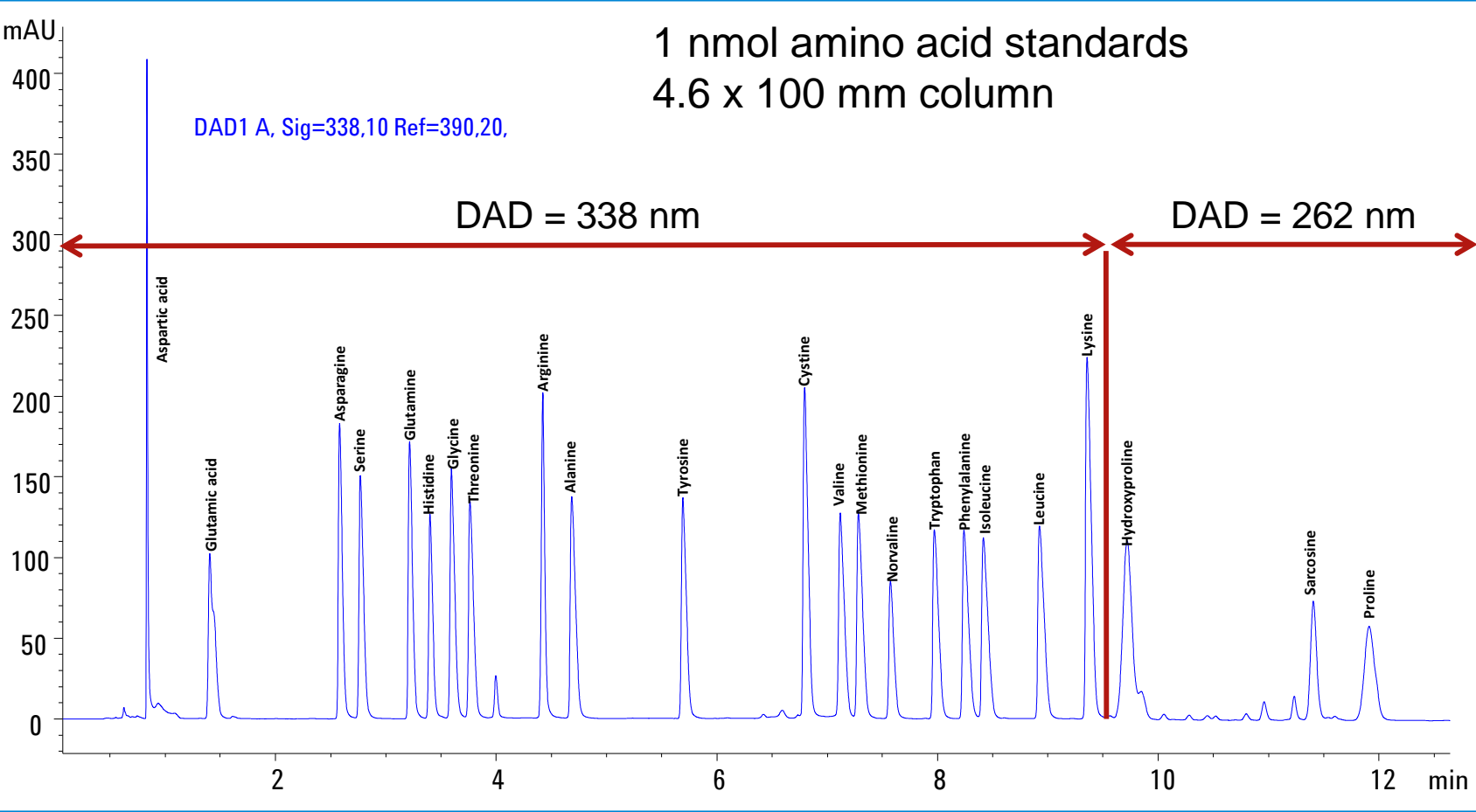
Peak #	AA Name	AA Abbreviation	Derivative Type	
1	Asparic Acid	ASP	OPA	Primary AA
2	Glutamic Acid	GLU	OPA	
3	Aspraragine	ASN	OPA	
4	Serine	SER	OPA	
5	Glutamine	GLN	OPA	
6	Histidine	HIS	OPA	
7	Glycine	GLY	OPA	
8	Threonine	THR	OPA	
9	Arginine	ARG	OPA	
10	Alanine	ALA	OPA	
11	Tyrosine	TYR	OPA	
12	Cysteine	CYS-CYS	OPA	
13	Valine	VAL	OPA	
14	Methionine	MET	OPA	
15	Norvaline*	NVA	OPA	
16	Tryptophan	TRP	OPA	
17	Phenylalanine	PHE	OPA	
18	Isoleucine	ILE	OPA	
19	Leucine	LEU	OPA	
20	Lysine	LYS	OPA	
21	Hydroxyproline	HYP	FMOC	Secondary AA
22	Sacrosine (IS)	SAR	FMOC	
23	Proline	PRO	FMOC	

Elution Profile with and without Sodium Azide

- Historically NaN_3 has been added to aqueous mobile phase to reduce bacterial growth.
- NaN_3 is highly toxic.
- No effect on the separation.
- Highly recommend filtering mobile phases (0.45 or 0.2 μm) to reduce bacterial growth.



Reproducible Separations



- Retention time %RSD mostly under 1%
- Peak area %RSD mostly under 3%

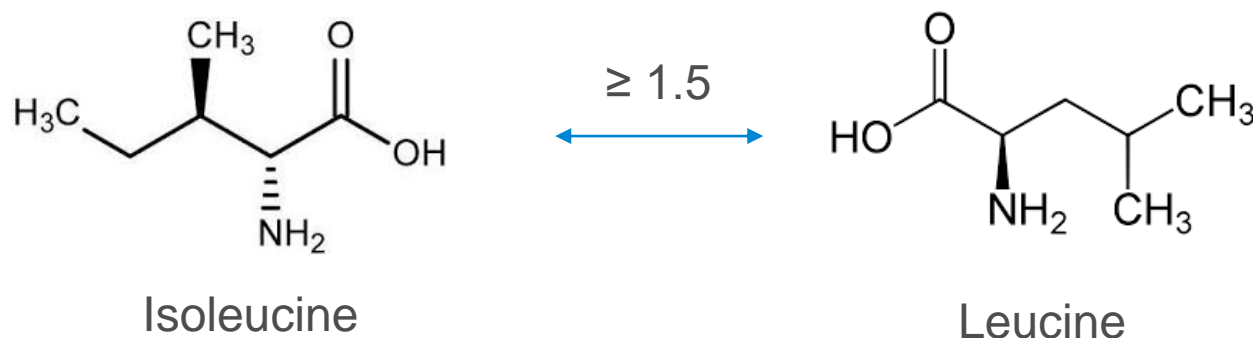
Amino Acids	RT RSD (%)	Area RSD (%)
1. Aspartic acid	1.270	1.066
2. Glutamic acid	0.973	1.85
3. Asparagine	0.605	1.79
4. Serine	0.629	1.82
5. Glutamine	0.470	1.56
6. Histidine	0.430	1.22
7. Glycine	0.477	1.92
8. Threonine	0.440	1.95
9. Arginine	0.251	2.15
10. Alanine	0.280	3.06
11. Tyrosine	0.128	1.65
12. Cystine	0.067	1.9
13. Valine	0.084	2.47
14. Methionine	0.073	1.82
15. Norvaline	0.073	1.72
16. Tryptophan	0.054	1.57
17. Phenylalanine	0.051	1.66
18. Isoleucine	0.047	1.72
19. Leucine	0.03	1.7
20. Lysine	0.028	1.66
21. Hydroxyproline	0.021	4.13
22. Sarcosine	0.026	1.15
23. Proline	0.021	4.36



System suitability as per European Pharmacopoeia (Ph.Eur)

The European Pharmacopoeia (Ph. Eur.) defines requirements for the qualitative and quantitative composition of amino acids and mixtures of amino acids. The requirements for allowed impurities are also defined. Manufacturers of amino acids are legally bound to prove that their amino acids meet these specifications before they can distribute their products in Europe.

Leucine (Leu) is a branched-chain α -amino acid and is produced by the fermentation process. During this process, isoleucine can be produced as a by-product. The European Pharmacopoeia states that leucine and isoleucine should have a resolution of not less than 1.5



Ref: Ph.Eur.9.0 (2.2.56) Amino Acid Analysis

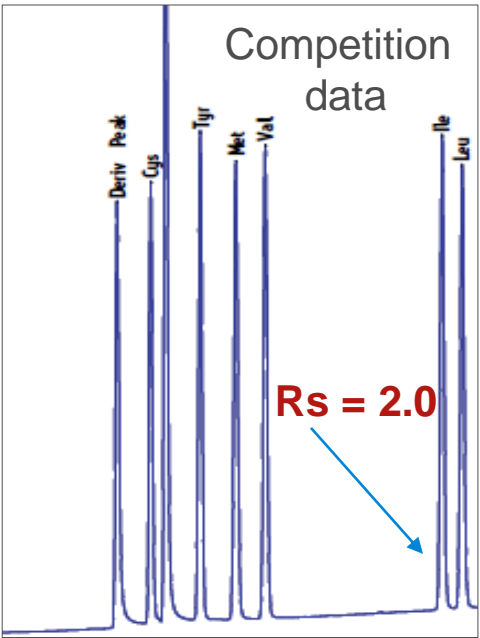
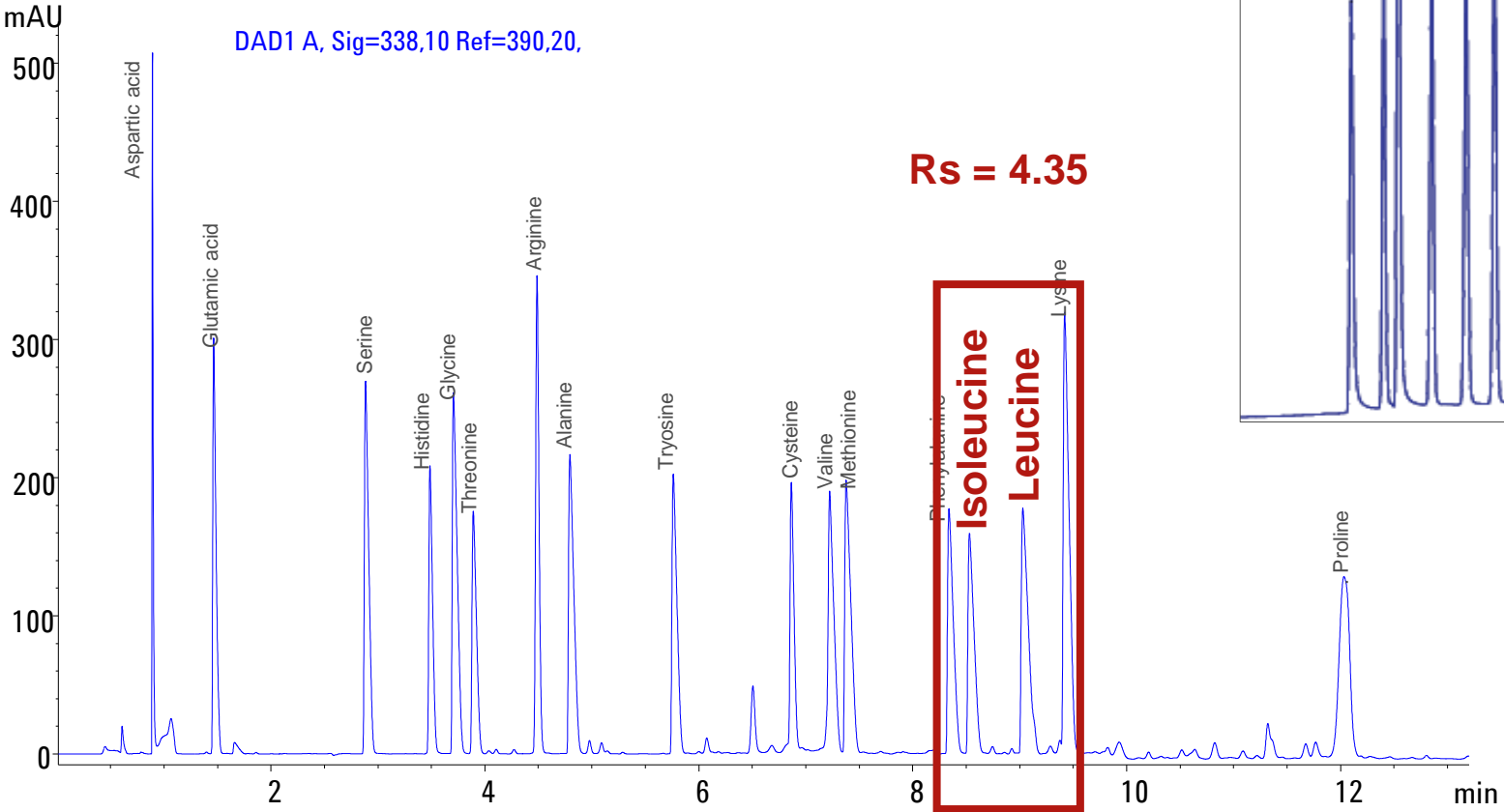


Agilent Technologies

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

Ample Resolution of Leucine & Isoleucine

Baseline resolution of isoleucine and leucine ($R_s = 4.35$) meeting the regulatory requirements for these components.

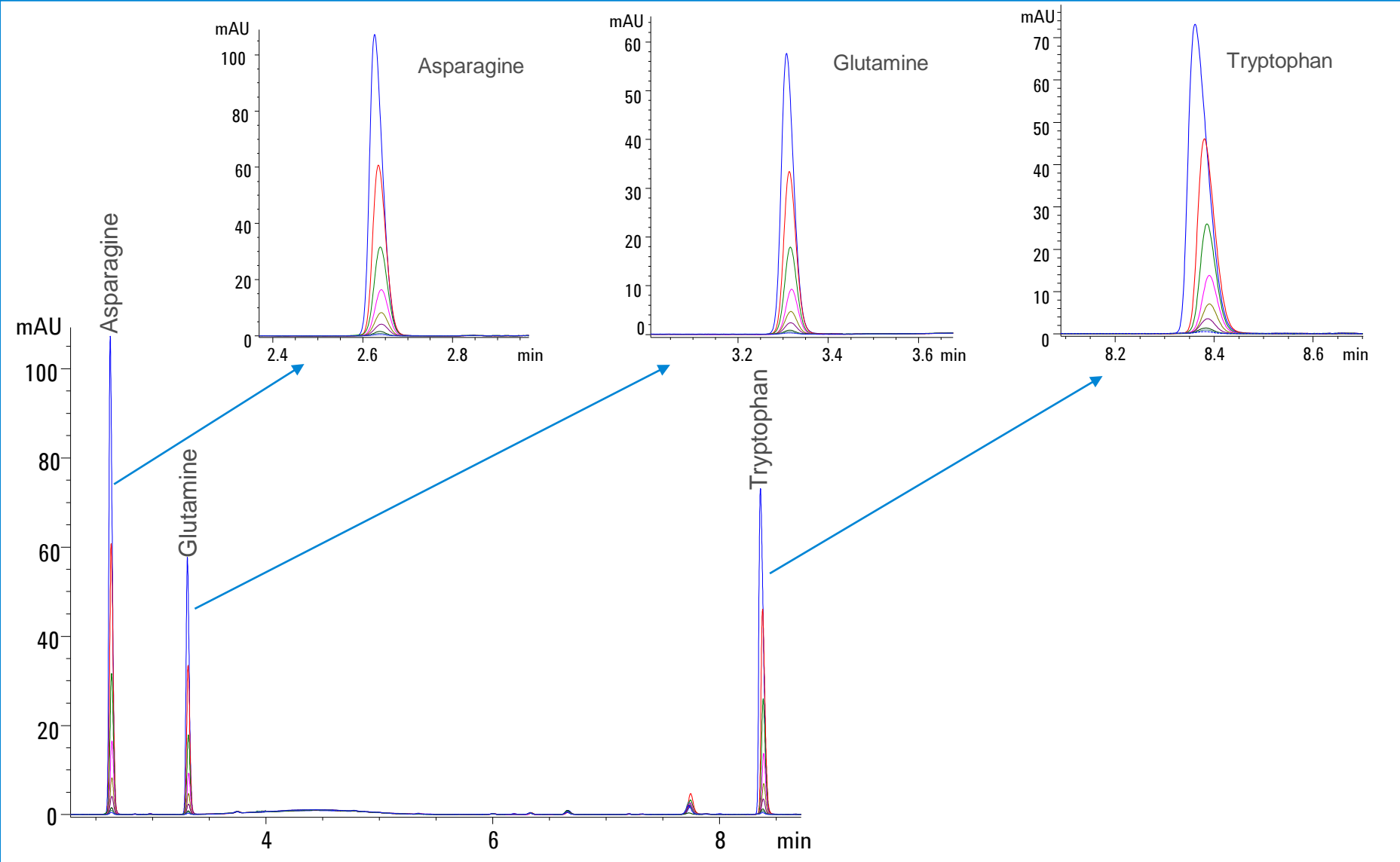


Protein hydrolysate sample

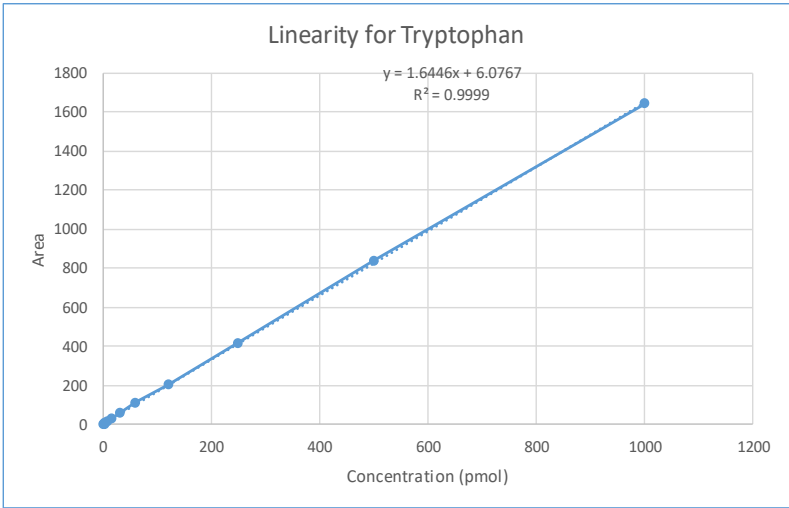
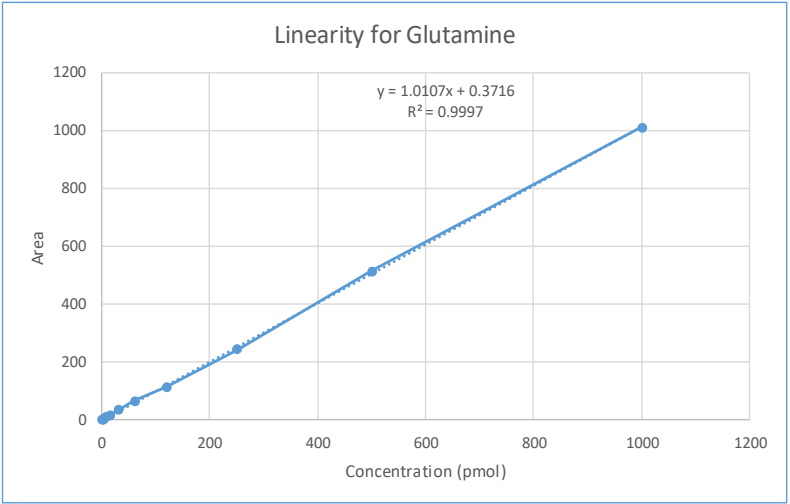
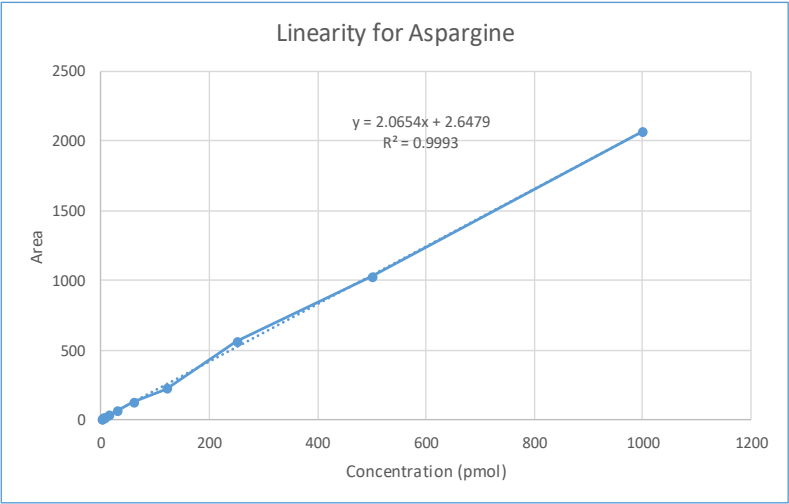
Column	Leu/Ile R_s (≥ 1.5)
Agilent AdvanceBio AAA 4.6 × 100 mm	4.5
Agilent AdvanceBio AAA 3 × 100 mm	4.6



Linearity and Limits of Detection & Quantitation



Linearity and Limits of Detection & Quantitation

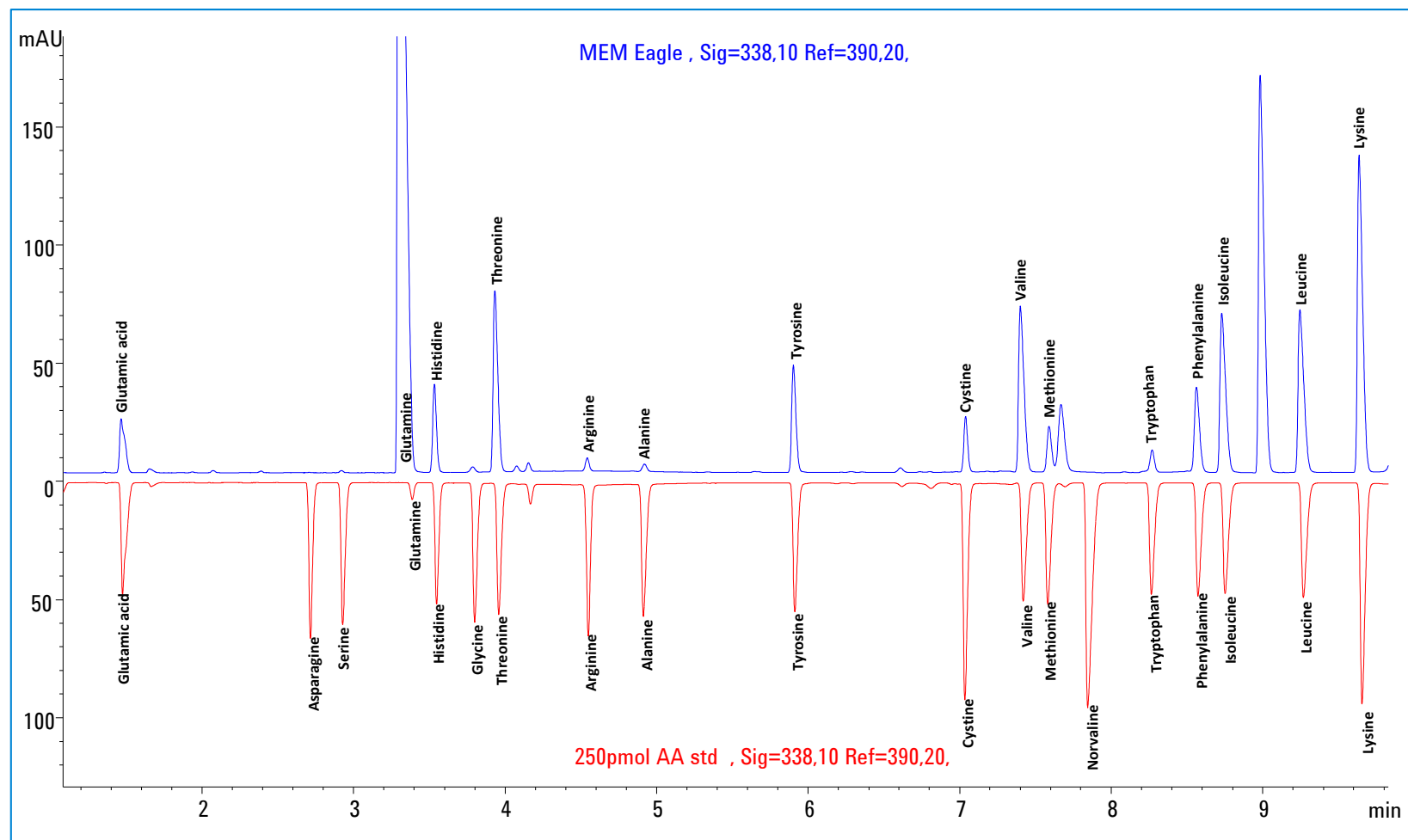


Concentration (pmol)	S/N ratio
Asparagine	
0.9 (LOD)	5.3
1.9 (LOQ)	10.8
Glutamine	
0.9 (LOD)	3.0
3.8 (LOQ)	13.8
Tryptophan	
0.9 (LOD)	4.5
3.8 (LOQ)	20.5

S/N >3.0 = LOD
S/N > 10 = LOQ

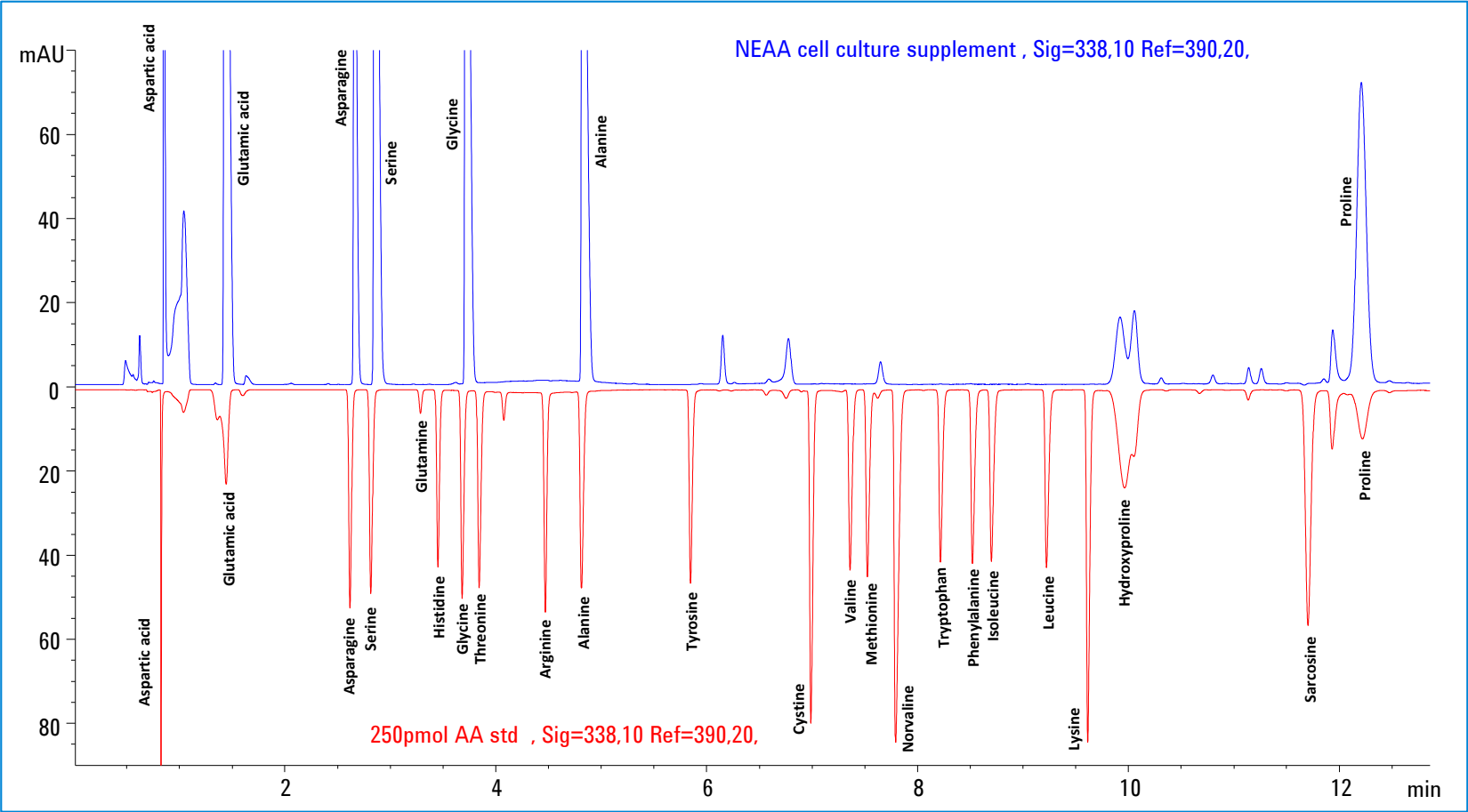
AAA of Cell Culture Media – MEM

L-Arginine, L-Cystine, L-Glutamine, L-Histidine, L-Isoleucine, L- Leucine, L-Lysine, L-Methionine, L- Phenylalanine, L- Threonine, L-Tryptophan, L- Tyrosine and L-Valine, L-Glutamic acid



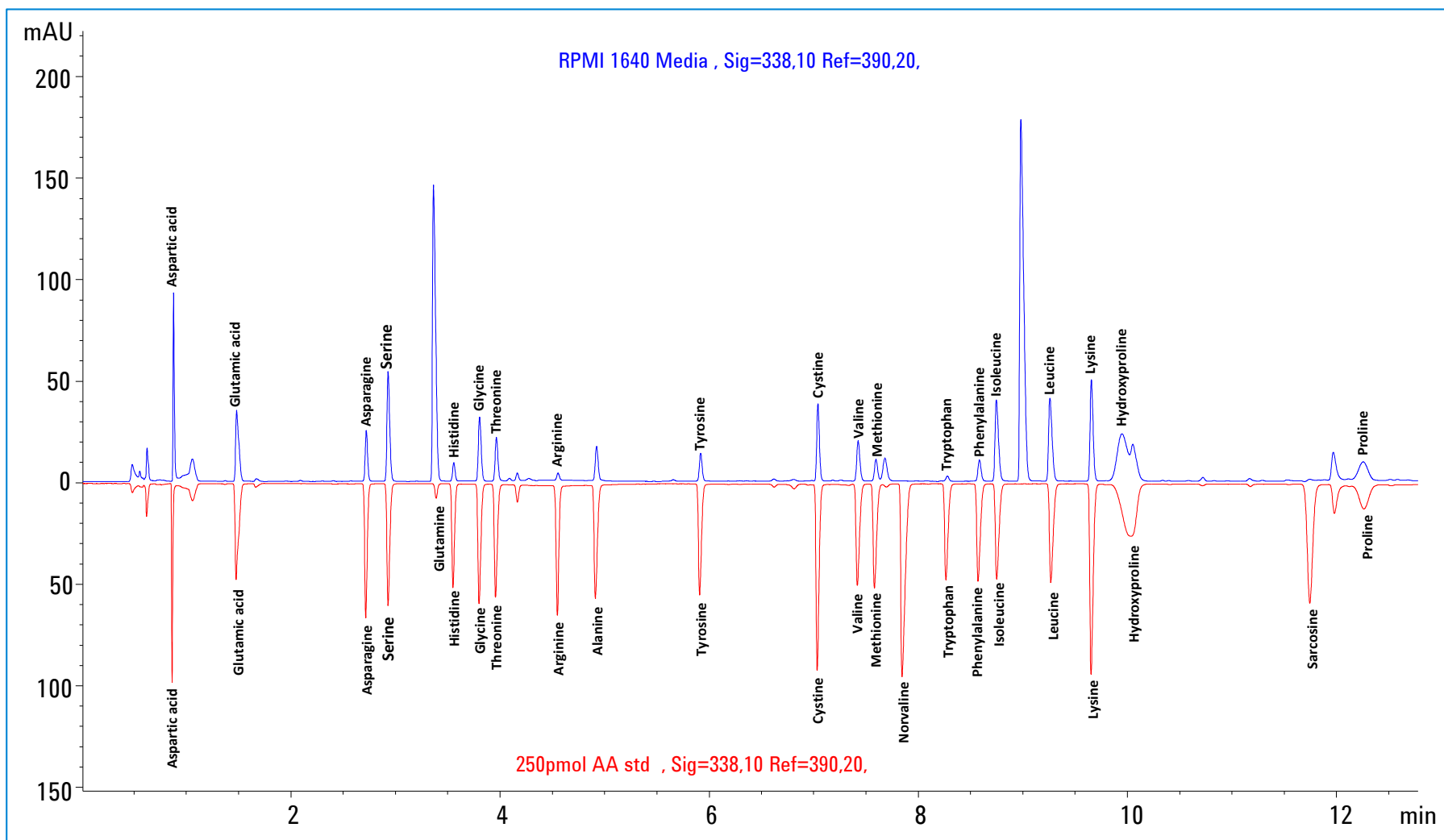
AAA of Cell Culture Media – NEAA cell culture supplement

L-Alanine, L-Asparagine, L-Aspartic acid, L-Glutamic acid, Glycine, L-Proline and L-Serine



AAA of Cell Culture Media – RPMI 1640

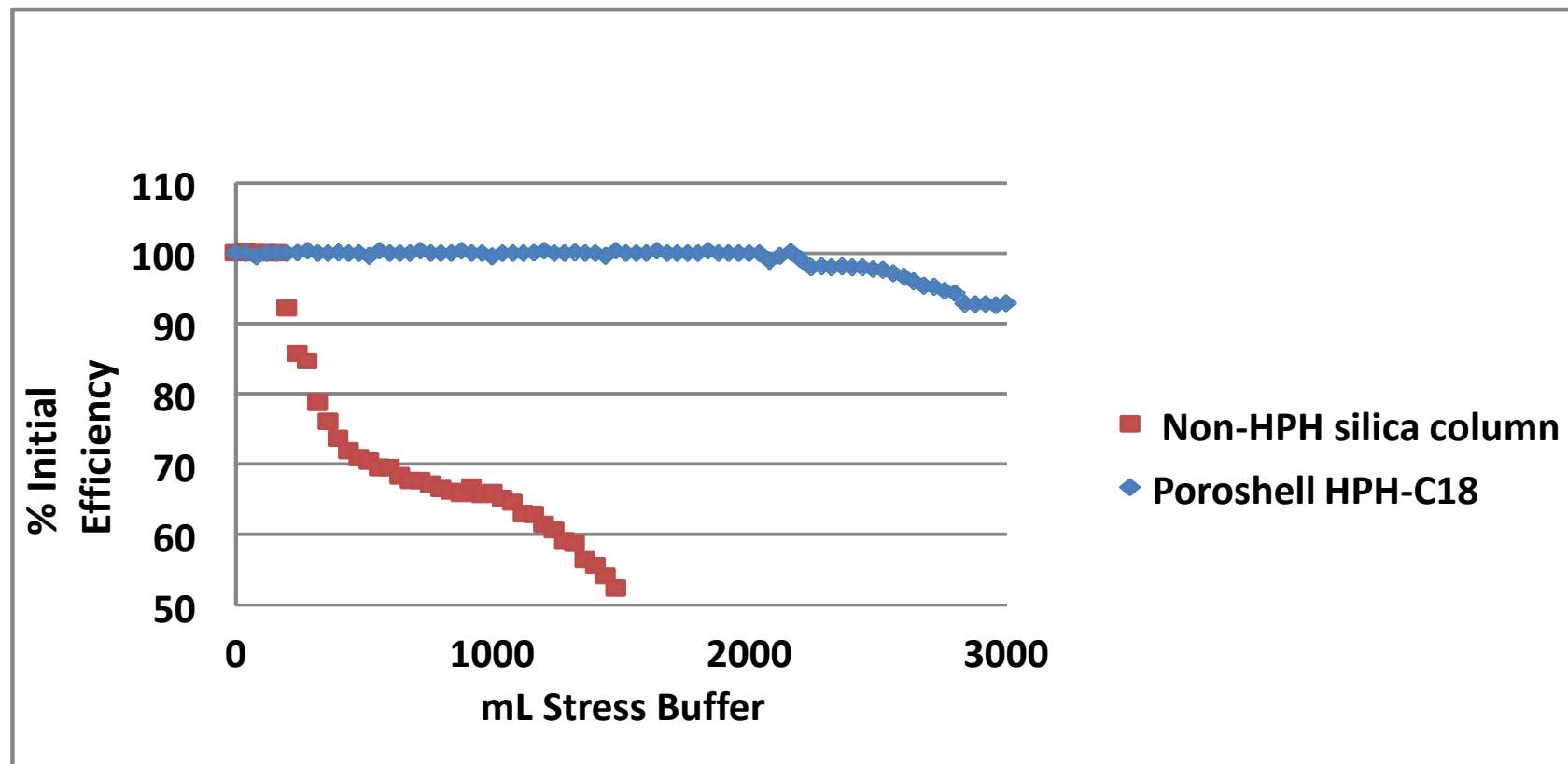
L-Arginine, L-Glutamic acid, L-Asparagine, L-Cystine, Glycine, L-Histidine, Hydroxy-L-Proline, L-Isoleucine, L-Leucine, L-Lysine, L-Methionine, L-Phenylalanine, L-Proline, L-Serine, L-Threonine, L-Tryptophan, L-Tyrosine and L-Valine



Agilent Technologies

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

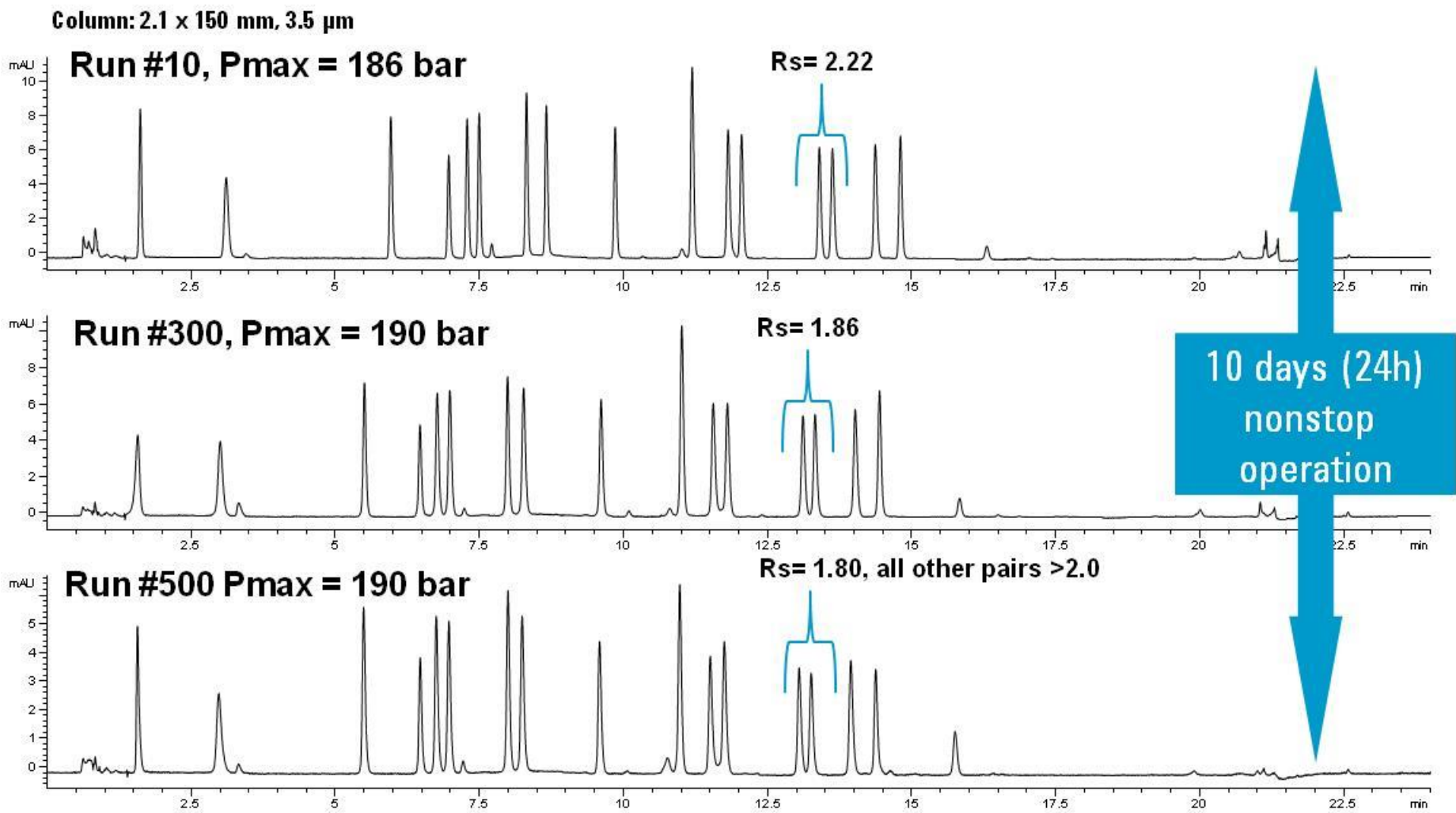
Poroshell HPH C18 last longer in phosphate buffer



Lifetime of SPP columns in phosphate buffer, **pH 8**, at elevated temperature. Mobile phase: Premixed 60% 30 mM sodium phosphate buffer at pH 8 and 40% acetonitrile; Flow rate 0.4 mL/min; UV absorbance 254 nm; 65 °C; Columns: 2.1 x 50 mm, 2.7 µm; Analyte: Naphthalene.

Traditional Lifetime Testing

Using conventional column testing Eclipse Plus C18 column test looks great and will perform well for most customers. However it is not a realistic test for most lab workflow.



Improvements in Lifetime Testing for AA Analysis

Past lifetime testing consisted of 500 injections without stopping

- This test, while it could differentiate long life columns, was not indicative of a customer workflow

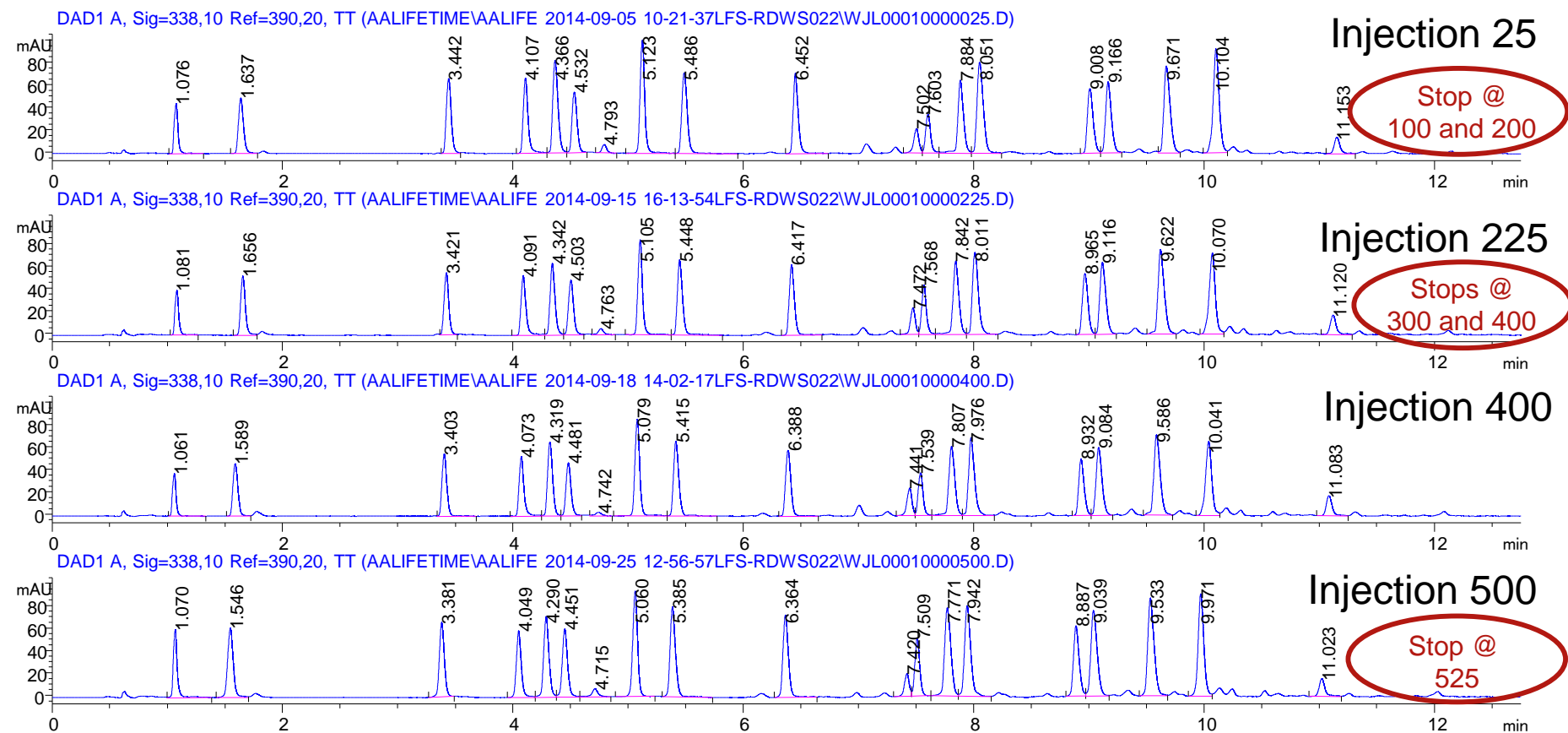
Customers will typically run ~20 to 50 or 100 samples, then stop, and resume a day or more later

- This is much more stressful on the column, and will cause columns to fail sooner versus running 500 injections straight through.

The lifetime testing for the Poroshell HPH-C18 was adapted to more customer focused workflow, with breaks between series of injections to give a true indicator of the lifetime in the customers' hands

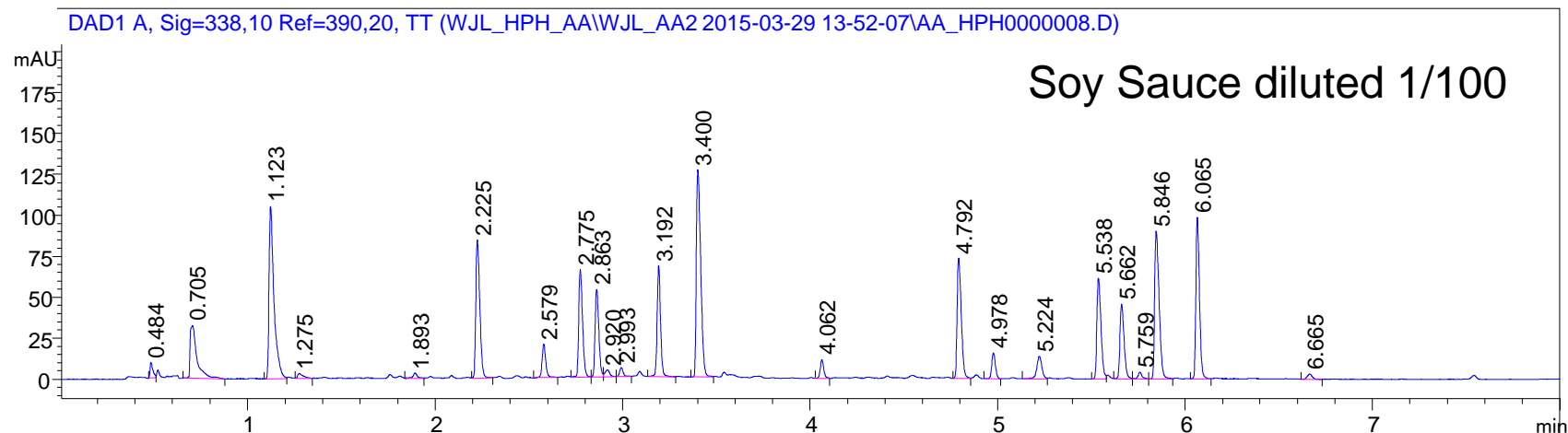
Test like a customer runs: 100 injections (3 days), Store (4 days)

Most labs run a batch of samples (25-100) and then shut off for a few days so we did too



Looks great after 500 injections run 100 injections shut down 4 days repeat

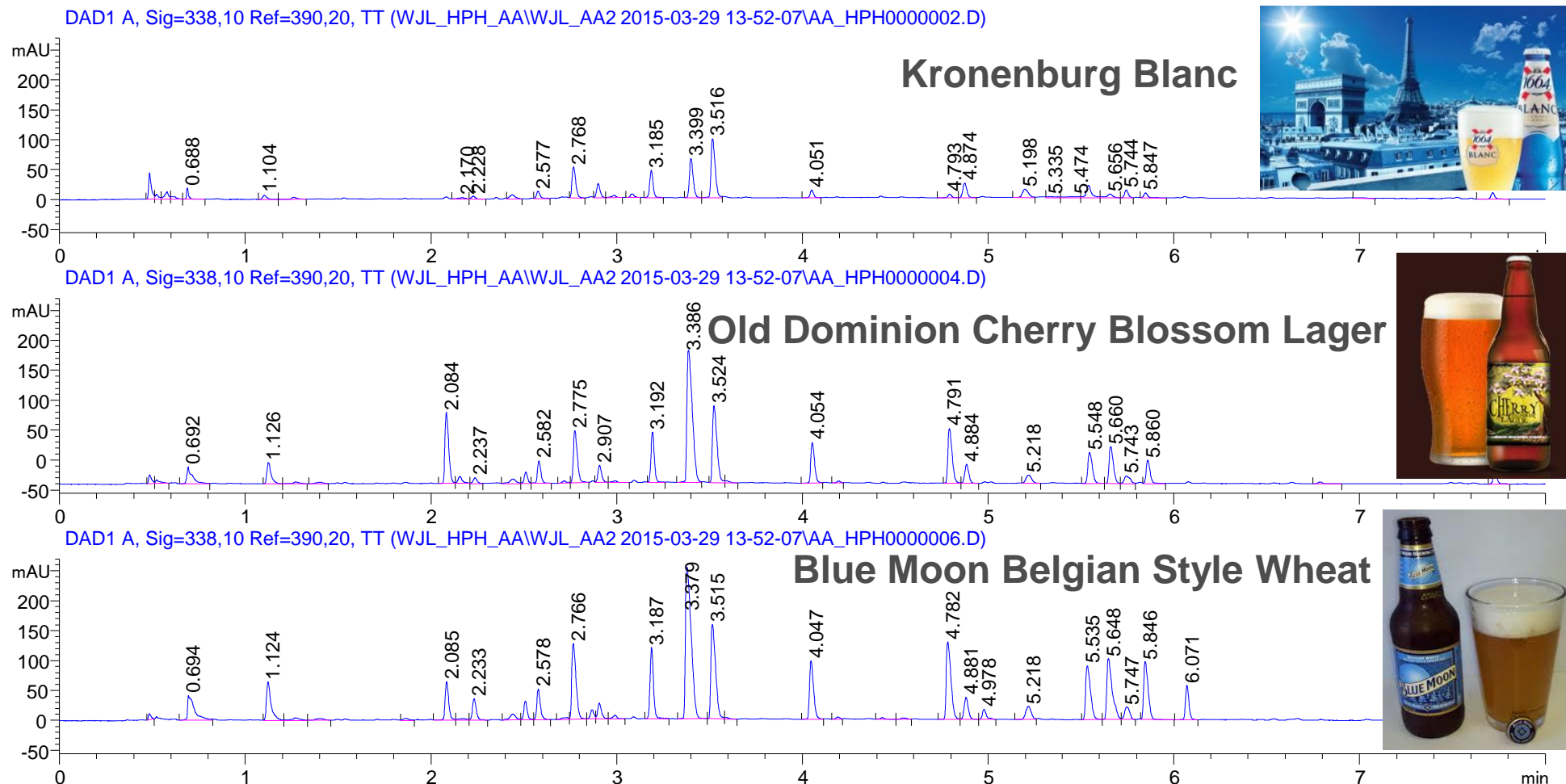
Amino Acid Analysis in Fermentation Applications



Many other foods (such as soy sauce) and pharmaceuticals that are produced using fermentation processes are monitored by AAA



Amino Acids Analysis for Batch Comparison



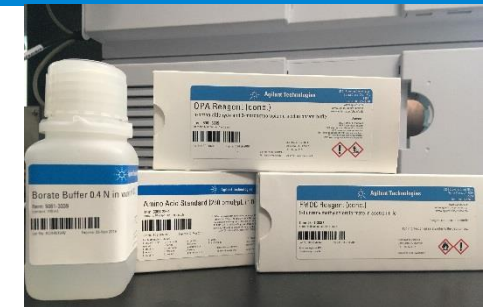
Quantity and diversity of amino acids is evident, can be used to monitor reactions and compare batches



Agilent Technologies

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

Tips & Tricks - Maintenance



- Replace derivatization reagent, borate buffer, amino acid standard daily
- Recalibrate for retention times and response factors daily
- Check column and guard column performance by following specs (Rs for 2 pairs of AA)
- Replace mobile phase A and B with fresh ones every other day
- Exchange guard column if high back pressure develops
- Avoid using MAX mixing speed during sample derivatization
 - The max speed on newer LCs is much faster than older LCs (1100s, 1200s), and can cause excessive wear on the autosampler.



Tips & Tricks - Troubleshooting

Poor chromatographic resolution?

- Cell culture media does not require any sample preparation, however appropriate dilutions have to be made to suit detector response
- In all cases, use the low-volume heat exchanger with short red tubing to minimize extra column volume
- Ensure proper connections
- Damaged guard or analytical column

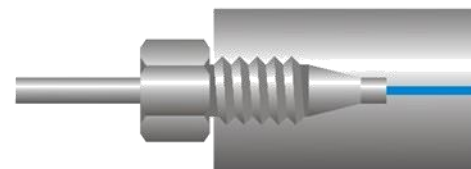
Low intensity chromatogram?

- OPA/FMOC reagent deteriorated
- Air bubble in vial insert

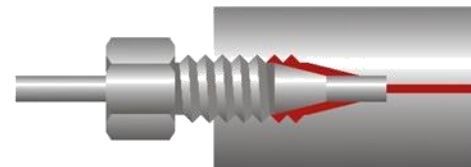


Column storage?

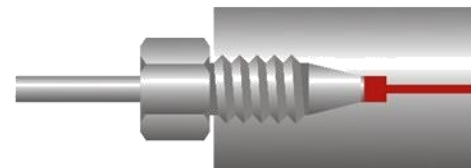
- Never leave the column in mobile phase A even if it's just overnight
- For short term always store the column in mobile phase B
- For long term, store column in 50/50 acetonitrile/H₂O



Good Connection



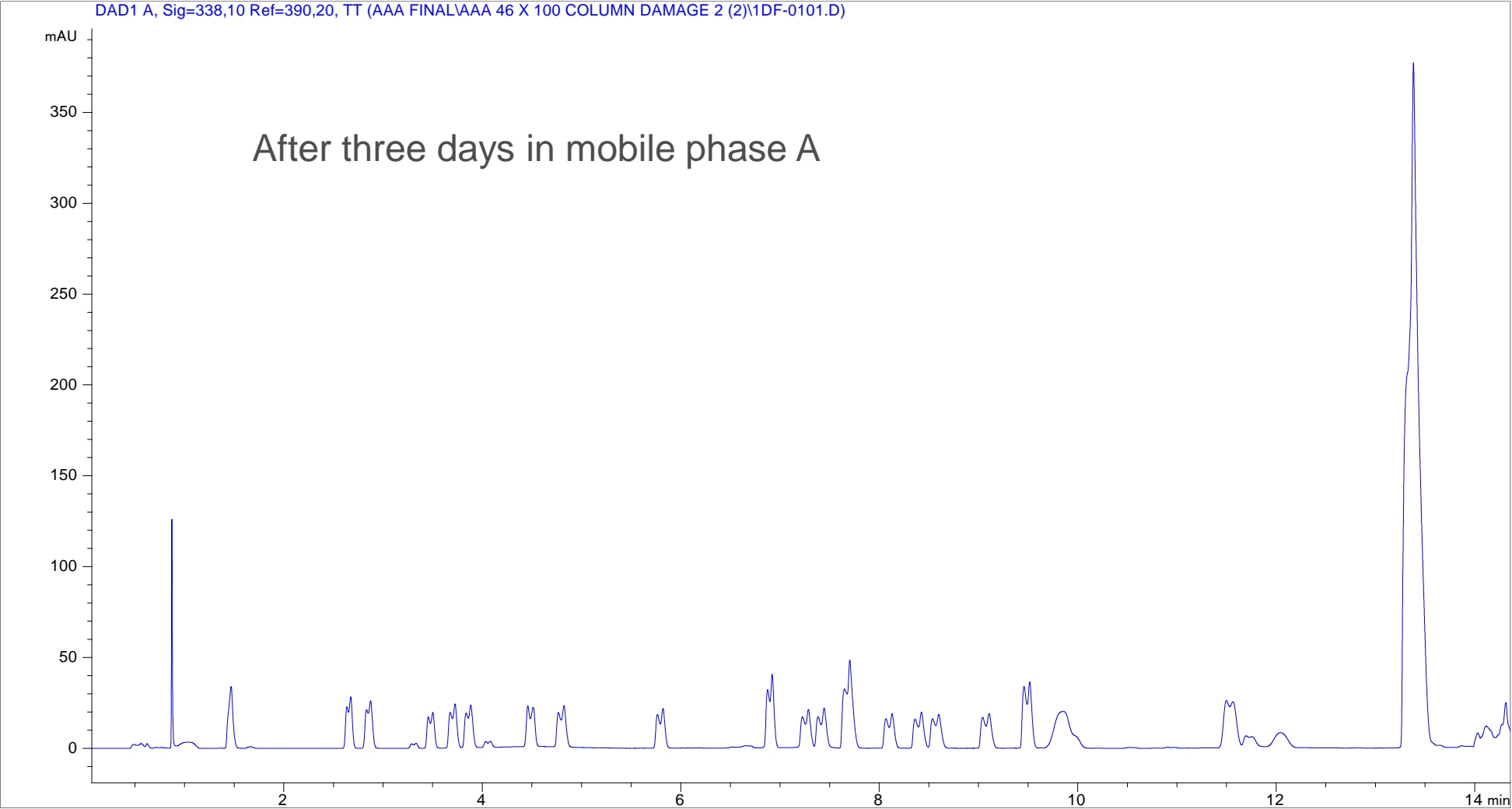
Bad Connection – capillary too long so ferrule not seated; leaking will occur



Bad Connection – capillary too short causing extra dead volume and potential mixing chamber

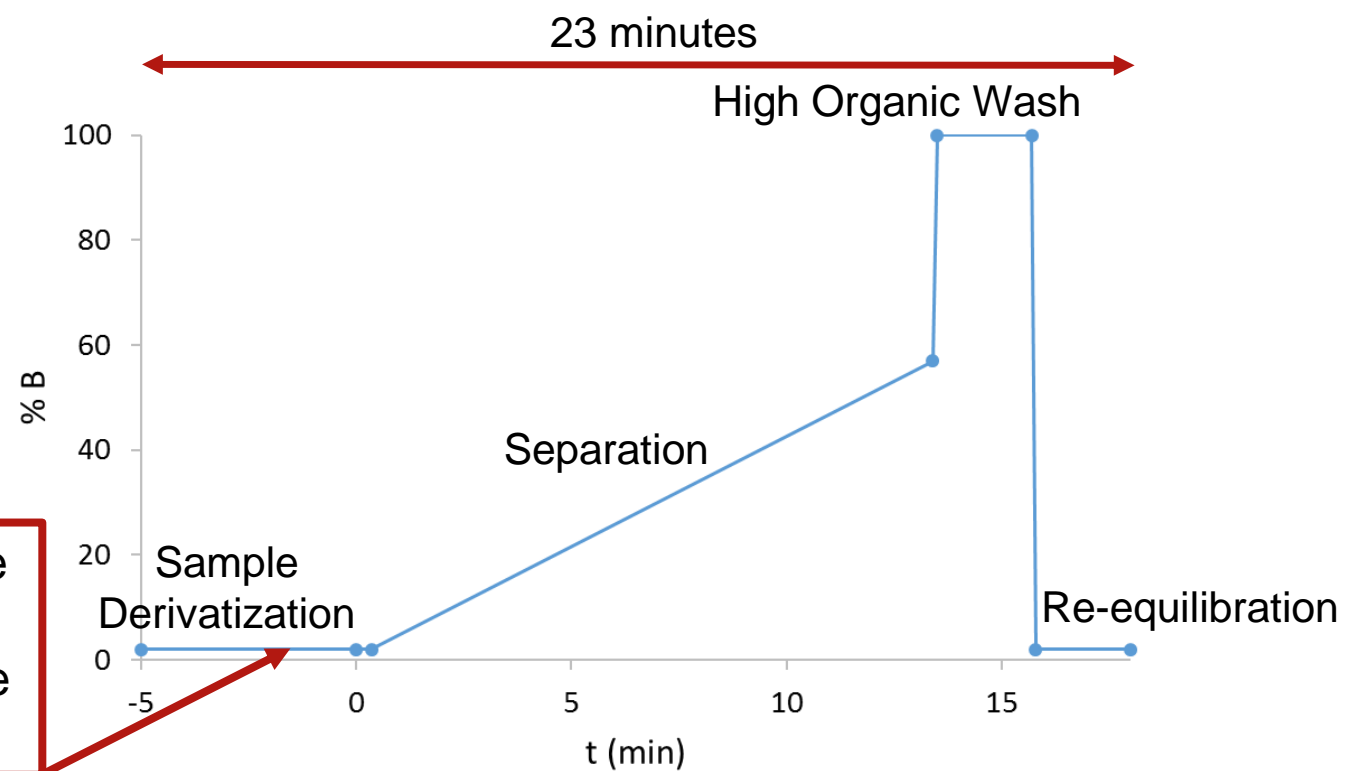


Damaged Column



Tips & Tricks – Saving Time

- After Injection 1, during the sample derivatization for Injection 2, the initial mobile phase condition is flowing through the LC and the column
- Save time by shortening the length of the re-equilibration time at the end of the method by the amount of time consumed by sample derivatization

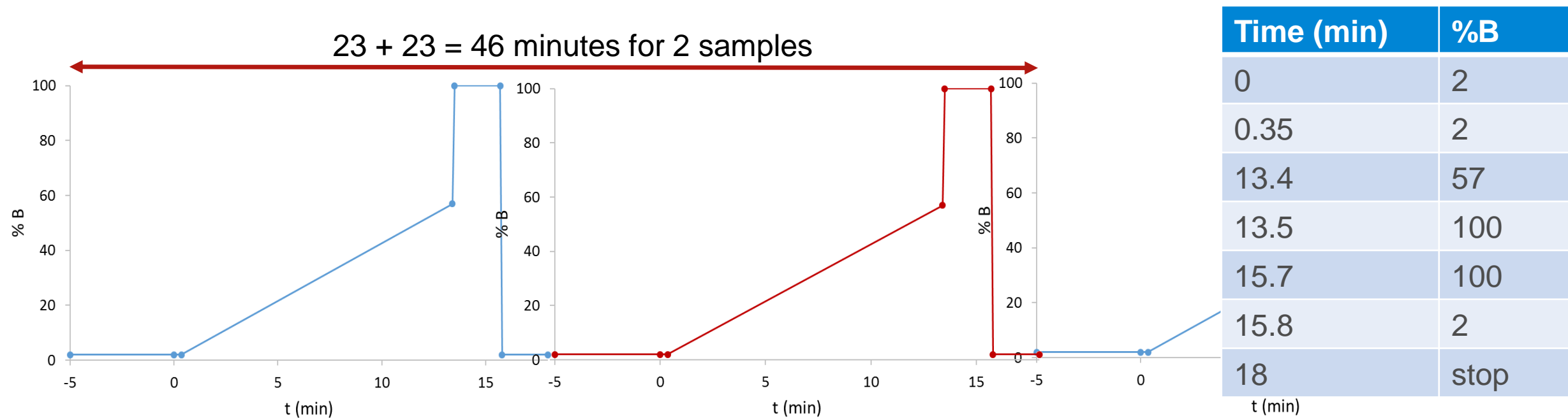


Initial mobile phase conditions are running through the column!

Time (min)	%B
0	2
0.35	2
13.4	57
13.5	100
15.7	100
15.8	2
18	stop

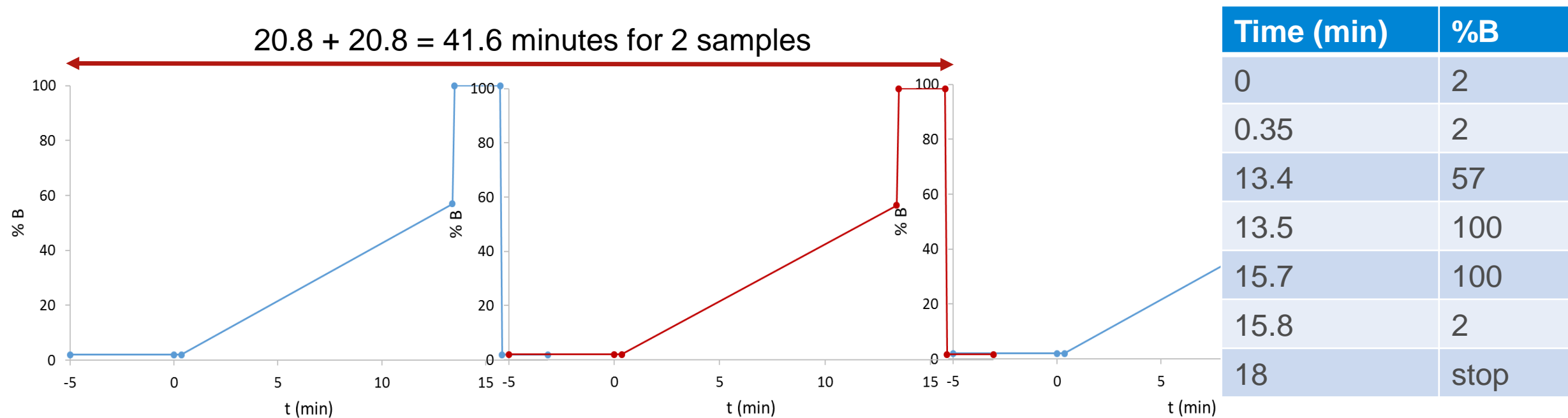
Tips & Tricks – Saving Time

- After Injection 1, during the sample derivatization for Injection 2, the initial mobile phase condition is flowing through the LC and the column
- Save time by shortening the length of the re-equilibration time at the end of the method by the amount of time consumed by sample derivatization

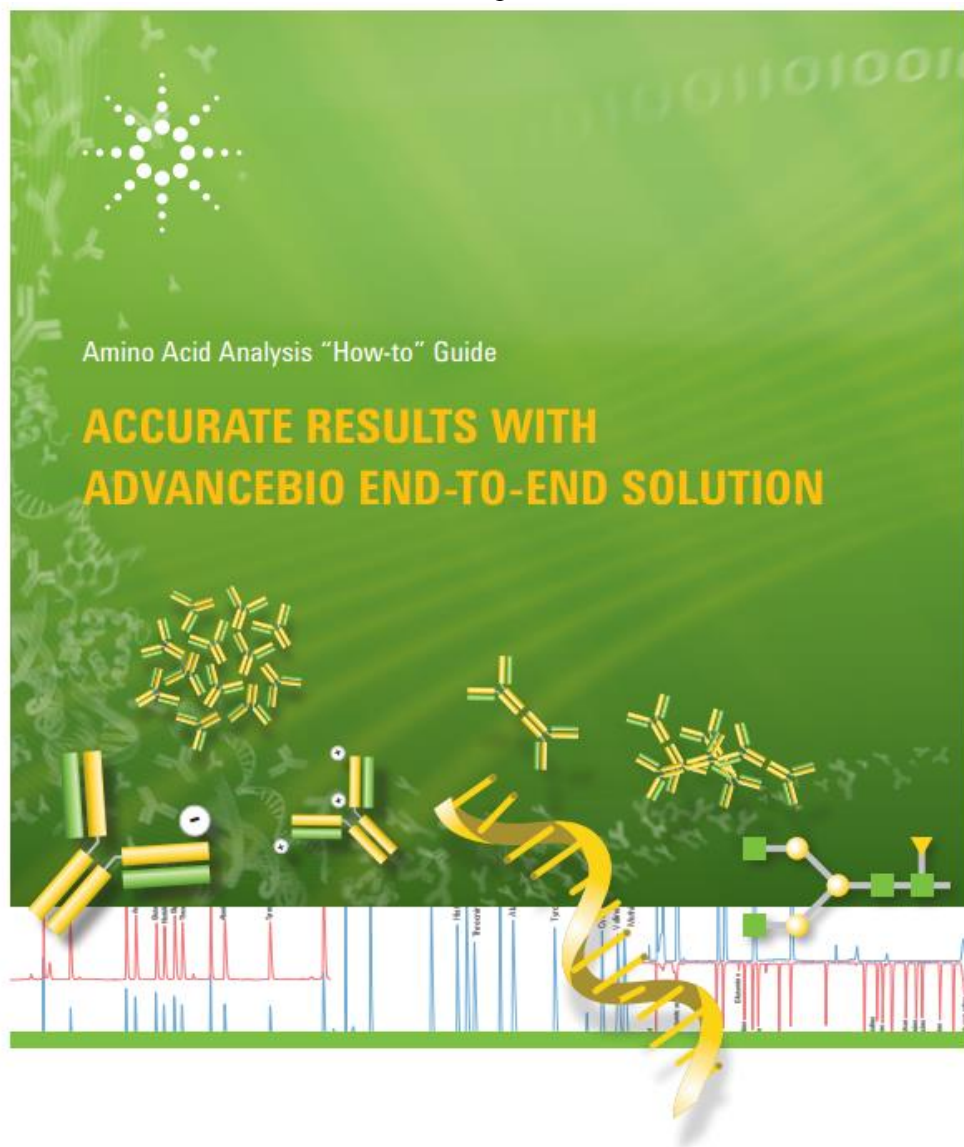


Tips & Tricks – Saving Time

- After Injection 1, during the sample derivatization for Injection 2, the initial mobile phase condition is flowing through the LC and the column
- Save time by shortening the length of the re-equilibration time at the end of the method by the amount of time consumed by sample derivatization



How-To Guide – Step by Step Instructions & Method Details



All the information you need

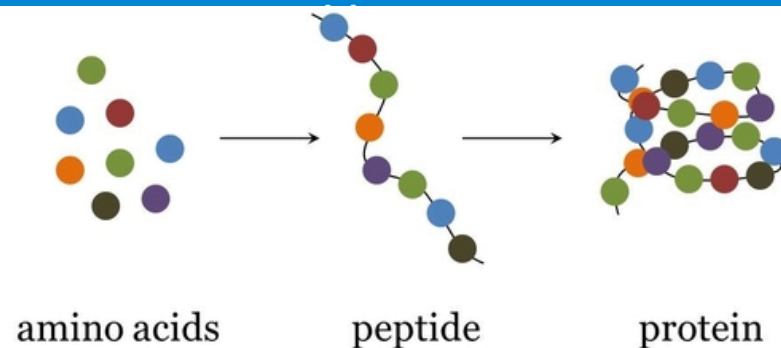
Document number 5991-7694EN



Agilent Technologies

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

Summary



- We used the Agilent AdvanceBio AAA solution for the automated online derivatization and separation of amino acids.
- Area and RT precision of the method were excellent, and Leu/Ile resolution met the system suitability requirement.
- Linearity curves with ten standard concentrations of three amino acids, ranging from 0.9pmol to 1nmol, had excellent coefficient of linearity values, indicating that the method was quantitative and accurate.
- The LOD and LOQ for the amino acids were 0.9pmol and 3.8 pmol respectively, indicating that the method was sensitive.
- This method was able to separate and detect, amino acids from a variety of samples, including cell culture media, protein hydrolysate, and fermentation reactions.

