



ThermoFisher
S C I E N T I F I C

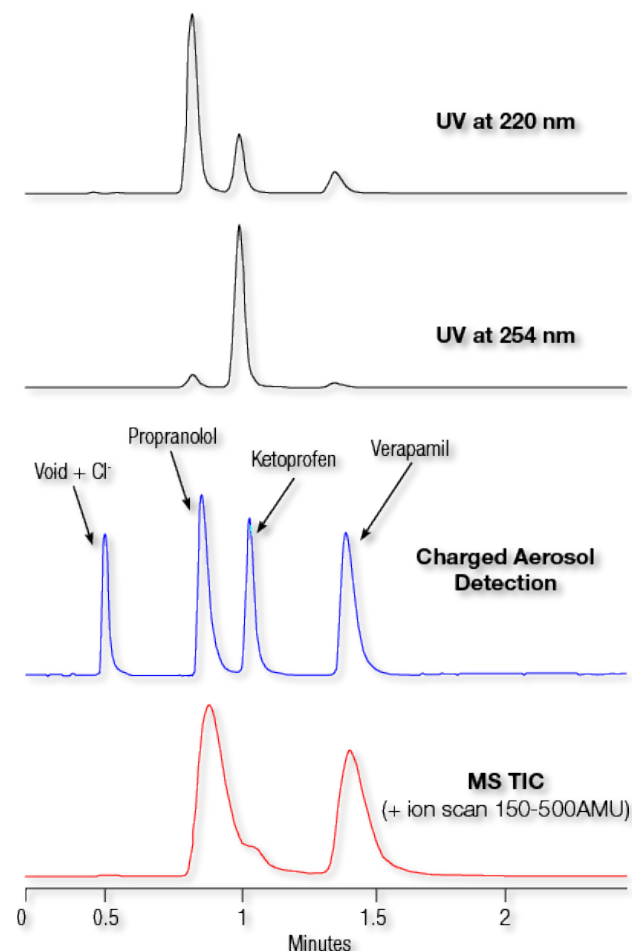
Charged Aerosol Detection 101

Dr. Alexander Schwahn
European Sales Support Expert for Biopharma Industry
Thermo Fisher Scientific, Reinach, Switzerland

- Introduction to charged aerosol detection (CAD)
- How charged aerosol technology works
- Comparison of CAD to ELSD
- CAD product evolution
- Example applications
- Summary

Introduction to Charged Aerosol Detection (CAD)

- Used to quantitate any non-volatile and many semi-volatile analytes with LC
- Provides consistent analyte response independent of chemical structure and molecule size
- Neither a chromophore, nor the ability to ionize, is required for detection.
- Dynamic range up to four orders of magnitude from a single injection (*sub-ng to μg quantities on column*)
- Mass sensitive detection – CAD provides relative quantification without the need for reference standards
- Compatible with gradient conditions for HPLC, UHPLC, and micro LC

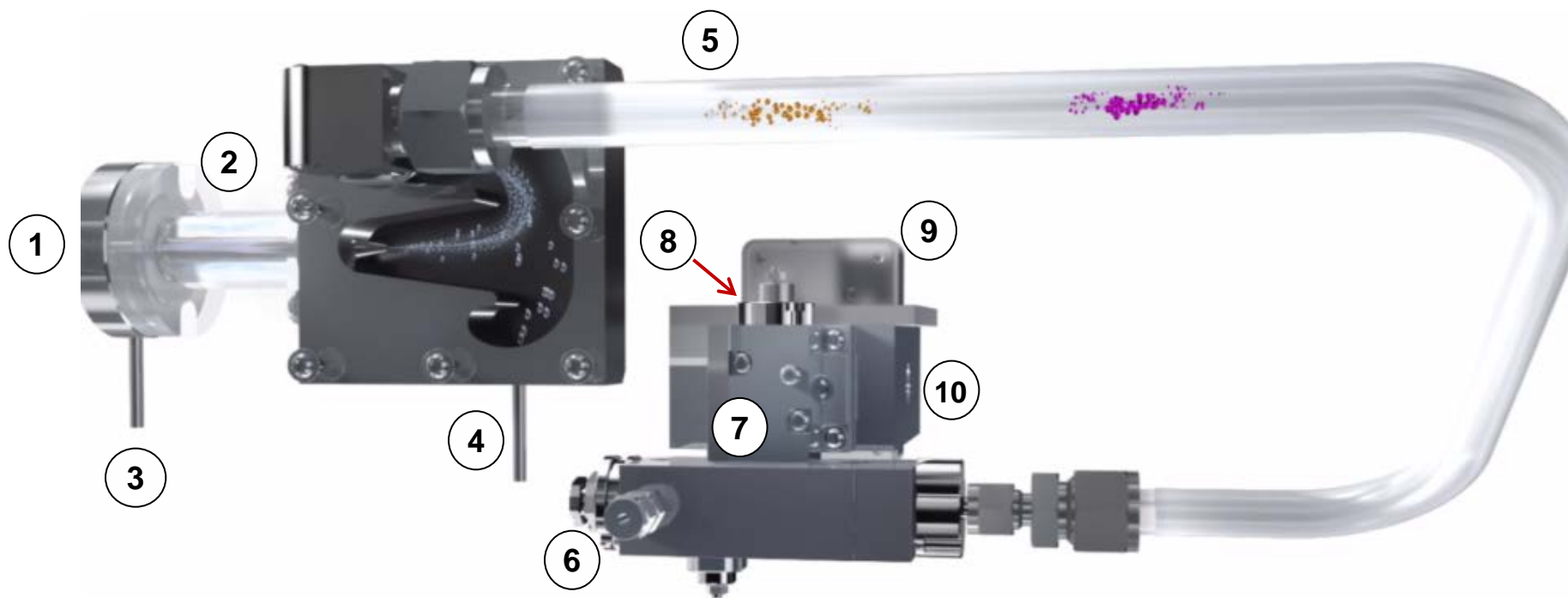


Comparison of charged aerosol detection to UV and MS

Charged Aerosol Detection – How It Works

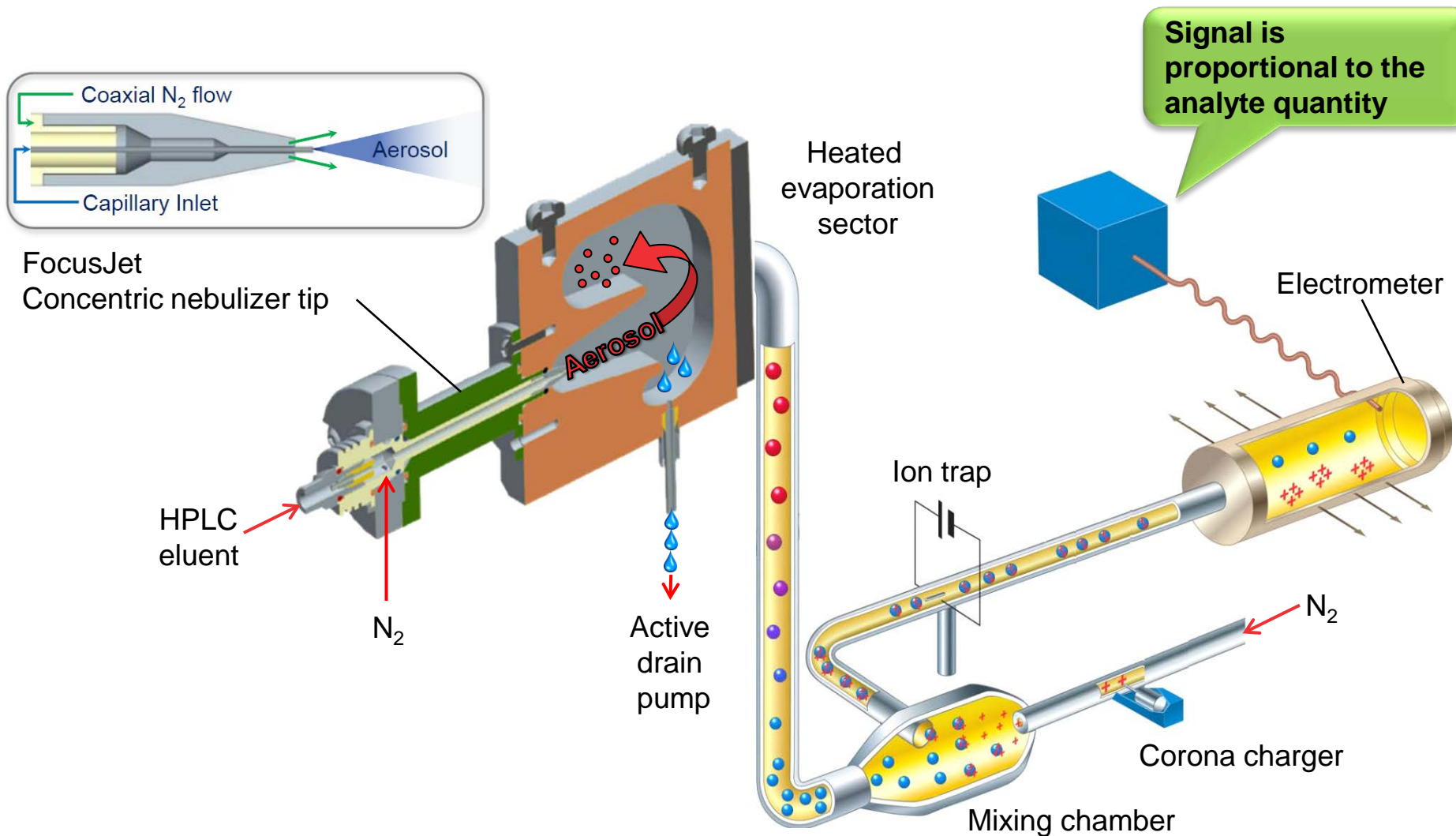
Flow path through a Thermo Scientific™ Dionex™ Corona™ Veo™ charged aerosol detector

- | | | | |
|---|--------------------------------|----|----------------|
| 1 | Inlet from column | 6 | Corona charger |
| 2 | FocusJet™ concentric nebulizer | 7 | Mixing chamber |
| 3 | Gas inlet | 8 | Ion trap |
| 4 | Micro drain pump | 9 | Electrometer |
| 5 | Evaporation tube | 10 | Gas exhaust |



Charged Aerosol Detection – How It Works

Flow path through a Thermo Scientific™ Dionex™ Corona™ Veo™ charged aerosol detector



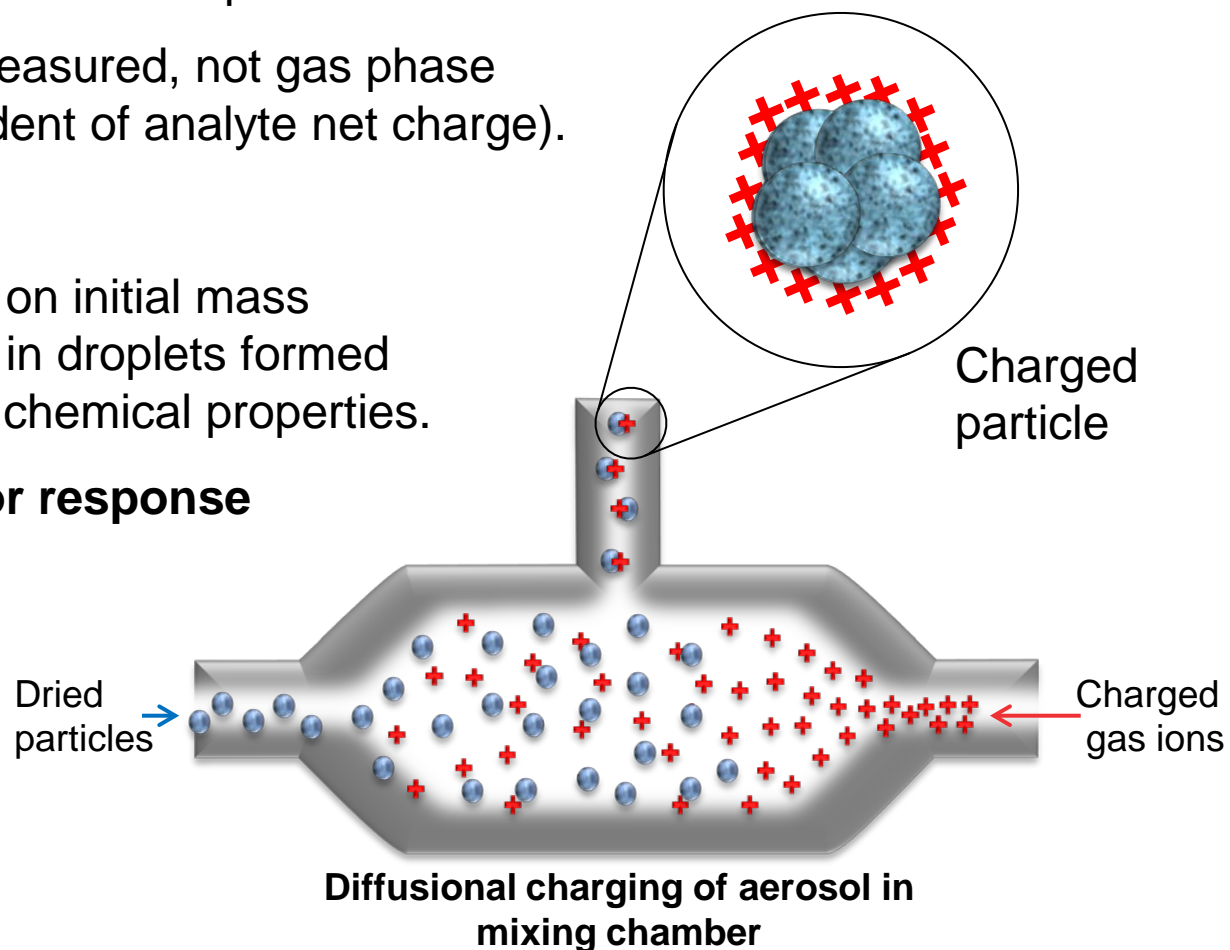
Particle Charging for Charged Aerosol Detection

- Particle size proportional to mass of analyte
- Charge per particle proportional to particle size
- Charged particles are measured, not gas phase ions as in MS (Independent of analyte net charge).

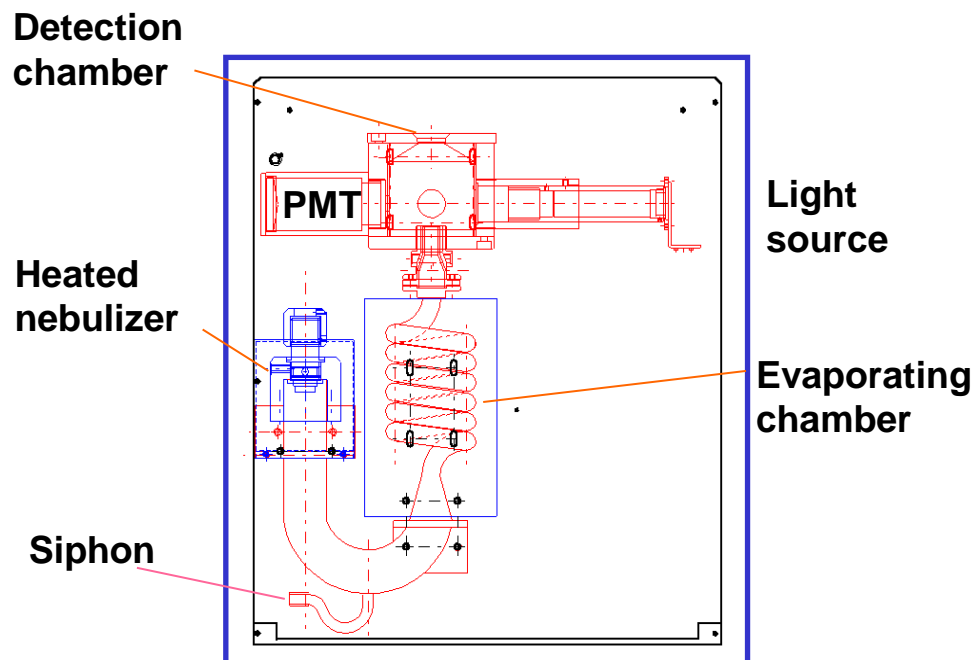
➔ CAD response depends on initial mass concentration of analyte in droplets formed but is independent of its chemical properties.

➔ **Nearly uniform detector response**

However, sample needs to be non-volatile or at least only semi-volatile.

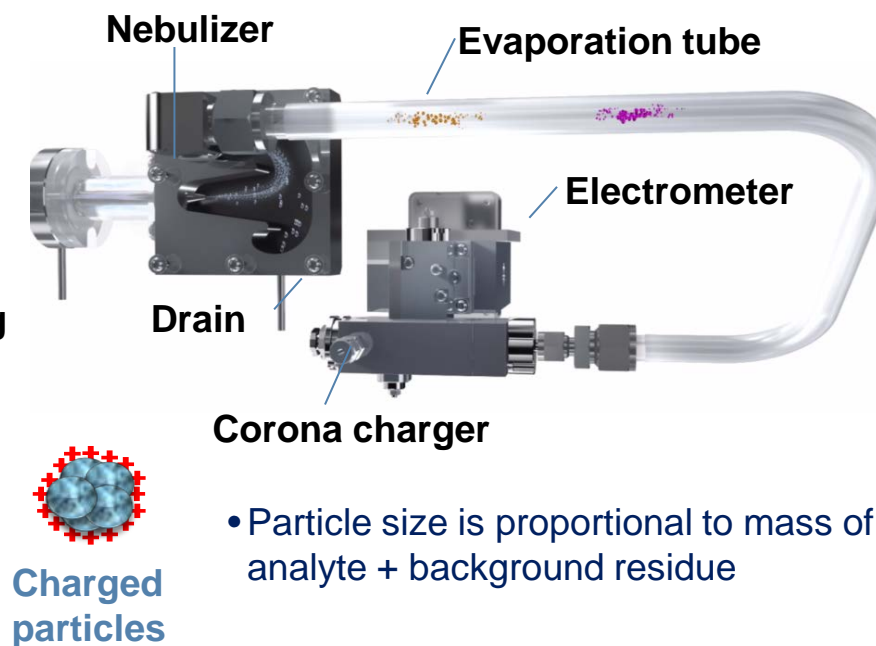


Evaporative light scattering detector (ELSD)



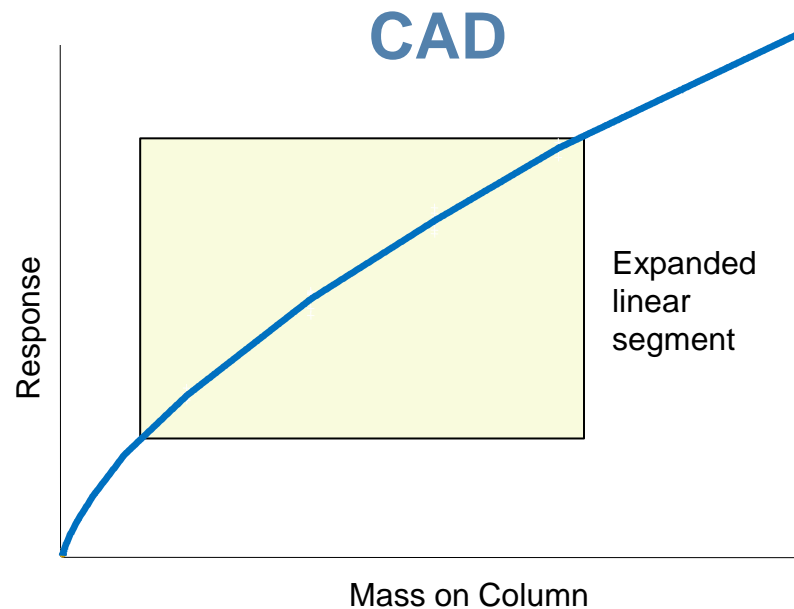
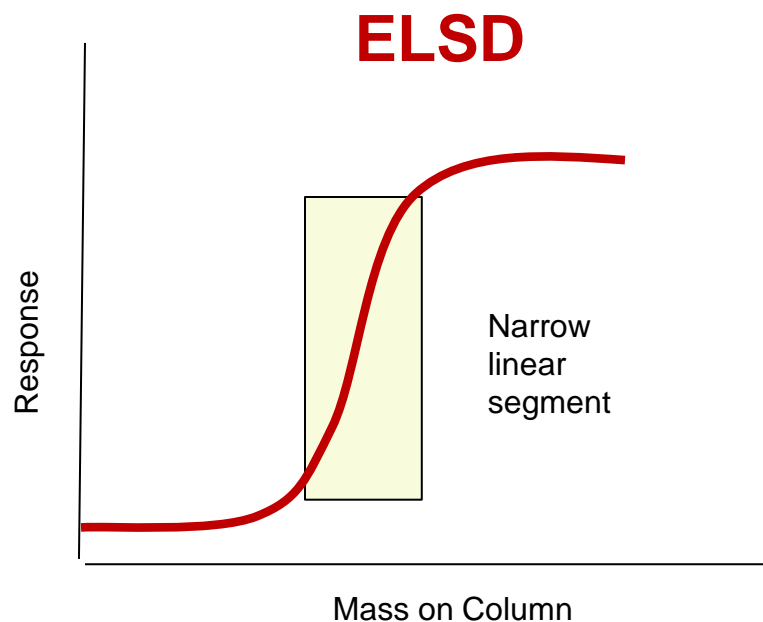
Measures the **optical reflection** of solute particles after the sample has been passed through a nebulizer

Charged aerosol detector



Measures **charged particles** by an electrometer generating a signal that is proportional to particle size (Mass of analyte) after nebulization

Detector Response Characteristics



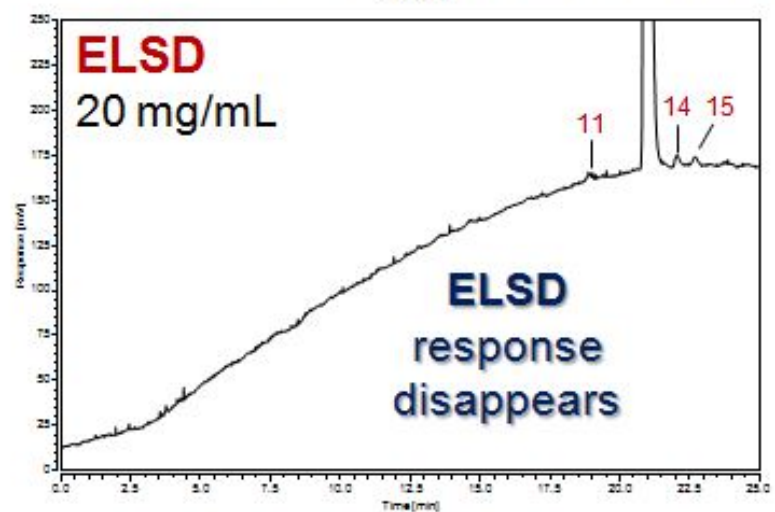
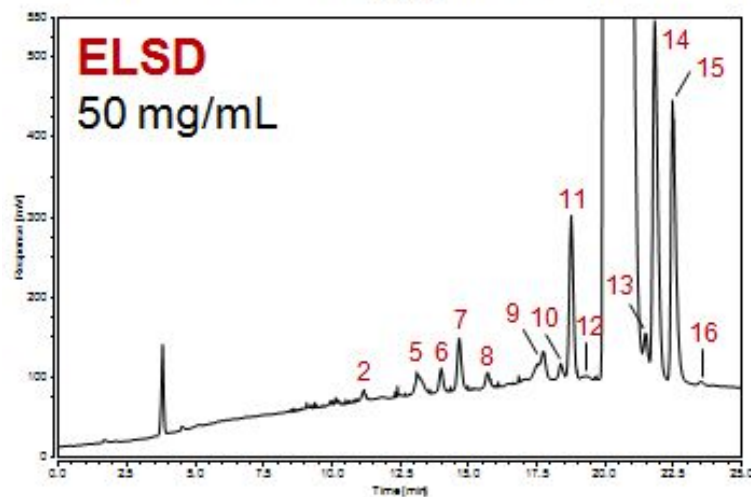
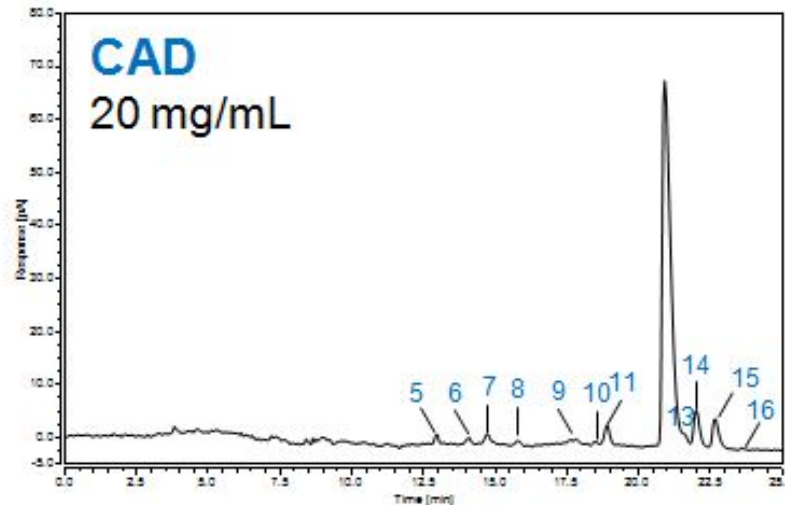
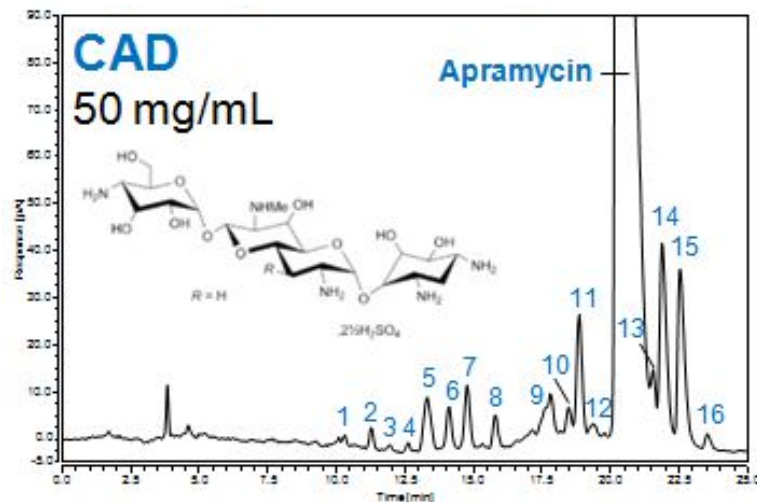
- For Rayleigh scattering: $b = 2$
- For Mie scattering: $b = 1\frac{1}{3}$
- For Refraction and reflection scattering: $b = \frac{2}{3}$

- Nonvolatiles - Decreasing slope with increasing mass ($b \sim \frac{2}{3}$)

ELSD exhibits a narrower linear calibration range than CAD.

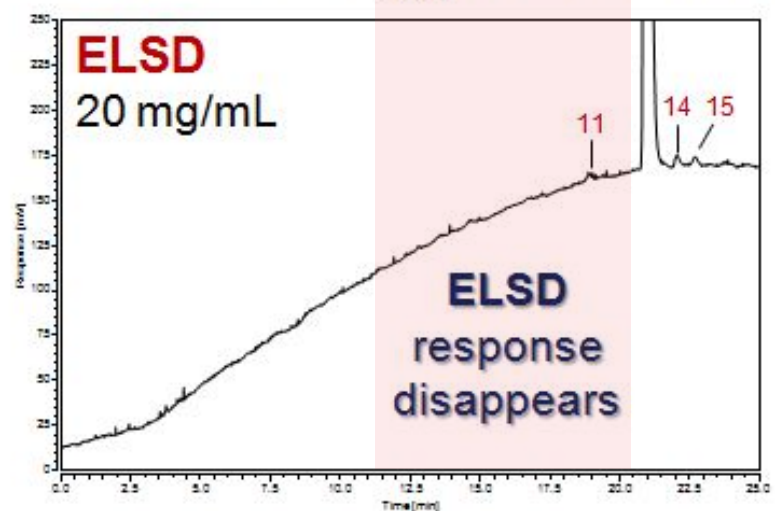
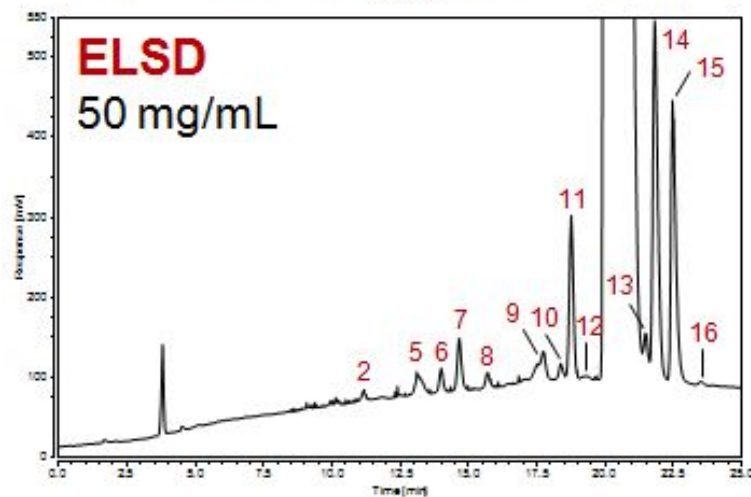
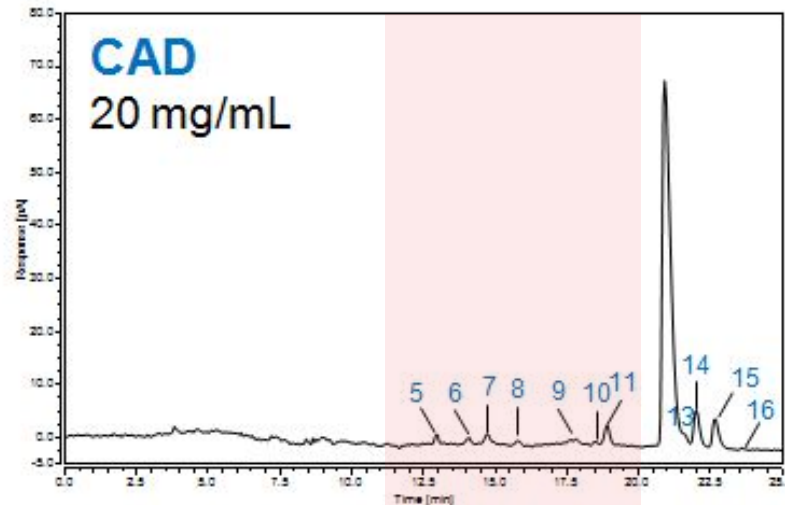
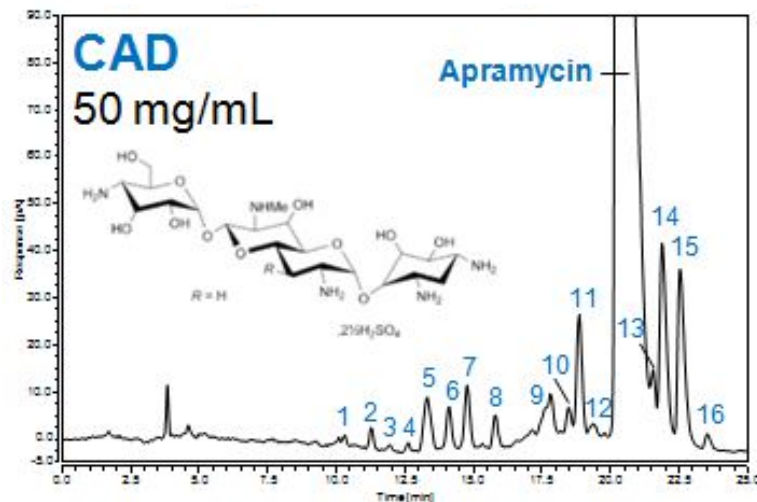
Detector Response Characteristics

Apramycin and impurities



Detector Response Characteristics

Apramycin and impurities



Calibration Consideration with Using Universal Detectors

- Over short ranges, both ELSD and charged aerosol detector can offer linear response.
- All aerosol-based detectors exhibit a non-linear response over large concentration ranges.
- Several calibration curve approaches are available:
 - Selections –

ELSD	CAD
Sigmoidal response behavior	Parabolic response behavior
Log-log	Log-log
Point to point	Quadratic
	Power function

The most appropriate approach depends upon the data.

Comparison Review

Feature	Evaporative light scattering	Charged aerosol
Response	Sigmoidal	Curvilinear
Dynamic range	2–3 orders	>4 orders
LoQ and LoD	LoQ and LoD often higher (Worse) than estimated by SNR	LoQ and LoD often lower (Better) than estimated by SNR
Sensitivity (LoD)	>10 ng	<1 ng
Semi-volatility range	Similar	Similar
Analyte response	Variable - Dependent on compound	Independent of structure
Flow rate range (0.2 – 2 mL/min)	Possibly several nebulizers	One nebulizer
Ease of operation	Can be complex	Simple

Evolution of Charged Aerosol Detectors

2015 Thermo Scientific™ Vanquish™ charged aerosol detector Full integration with Thermo Scientific™ Vanquish™ UHPLC platform, slide-in module design, reduced flow path for optimum operation



2013 Corona Veo RS CAD Extended micro flow rate range; total redesign with concentric nebulization and optimized spray chamber for enhanced sensitivity, heated evaporation and electronic gas regulation



Dionex Corp. acquired by Thermo Fisher Scientific Inc.

2011 Corona *ultra* RS CAD Unified with Dionex™ UltiMate™ 3000 UHPLC+ system, added on-board diagnostics / monitoring, automated flow diversion capability and selection of linearization parameters



ESA Biosciences, Inc. acquired by Dionex Corp.

2009 Corona *ultra* CAD UHPLC compatible, stackable design, enhanced sensitivity, touch-screen user interface with real-time chromatogram display, incorporated precision internal gas regulation system



2006 Corona *Plus* CAD Expanded solvent compatibility with heated nebulization, software drivers for popular CDS systems and external gas conditioning module for improved precision.



2005 Corona CAD Introduction of the first commercial charged aerosol detector for HPLC with full control via front panel interface. Designed for near-universal detection on any HPLC system using isocratic or gradient separations



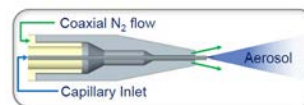
Corona Veo and Vanquish Charged Aerosol Detectors

- Concentric nebulization system improves sensitivity and precision.
- Thermally controlled evaporation scheme widens the scope of applications.
- Corona Veo and Vanquish RS model includes low flow capabilities for micro LC, as well as UHPLC.
- Usability and serviceability have been enhanced.
- CDS drivers available for use with all Thermo Scientific and many other vendor systems.



	Corona Veo	VeoRS
Nebulization	FocusJet	FocusJet
Flow rates	0.2 – 2.0 mL/min	0.01 – 2.0 mL/min
Data collection	100 Hz	200 Hz
Evaporation Temp.	35°C or 50°C	RT+5°C – 100°C
Gas pressure control	manual	electronic

Vanquish CAD



FocusJet



Corona Veo RS

Vanquish CAD model	Flow rate range (ml/min)	Data rate (Hz)	Evap temp (°C)	Positioning
Horizon	0.01 – 2.0 (Microflow)	2 – 200	Settable from ambient +5–100	Ideal for R&D and methods development Labs
Flex	0.2 – 2.0	2 – 100	Selectable 35, 50 or 70	Suitable for routine analysis in QC/QA Labs

Pharma and Biopharma Application Areas

- Drug composition
 - Impurity testing
- Formulation
 - Counterions
 - Surfactants / Excipients
- Degradation / Stability testing
- Characterization
 - Glycan analysis
 - Adjuvant analysis
- Excipient raw material analysis and lot-to-lot variability

Pharma and Biopharma Application Areas

- Drug composition
 - Impurity testing
- Formulation
 - Counterions
 - Surfactants / Excipients
- Degradation / Stability testing
- Characterization
 - Glycan analysis
 - Adjuvant analysis
- Excipient raw material analysis and lot-to-lot variability
- Cleaning validation
- Mass balance
- Extractables / Leachables
- PEGylation and antibody-drug conjugates
- siRNA lipid delivery vehicles
- QbD
- MIST (Metabolites in safety testing)

Visit the charged aerosol detection website and the free Thermo Scientific AppsLab library of analytical applications, to see more examples of HPLC-CAD solutions:

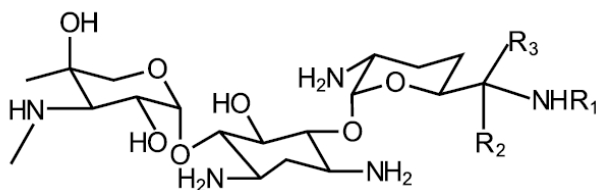
www.thermofisher.com/cad

www.thermofisher.com/appslab

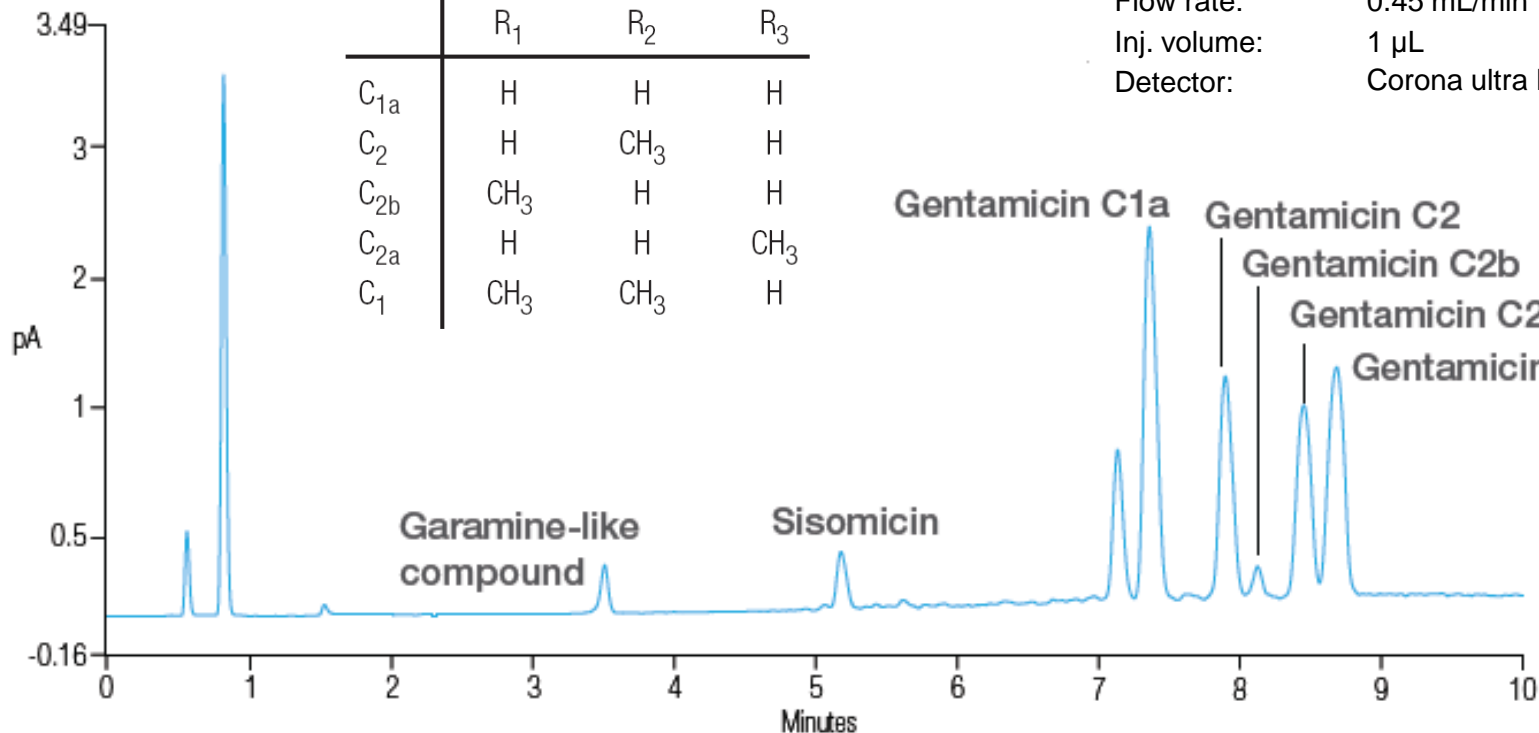
Drug Composition – Example: Aminoglycoside Antibiotic

Analysis of Gentamicin standard

(200 µg/mL)



	R ₁	R ₂	R ₃
C _{1a}	H	H	H
C ₂	H	CH ₃	H
C _{2b}	CH ₃	H	H
C _{2a}	H	H	CH ₃
C ₁	CH ₃	CH ₃	H



Column: Thermo Scientific™ Acclaim™ RSLC PolarAdvantage II, 2.2 µm, 2.1 × 100 mm

Mobile phase A: 0.025:95:5 HFBA:water:acetonitrile

Mobile phase B: 0.3:95:5 TFA:water:acetonitrile

Gradient: 0 to 1.5min, 1 to 10%B
1.5 to 7min, 10 to 100% B
7 to 10min, 100% B
4 min. pre-injection equilibration

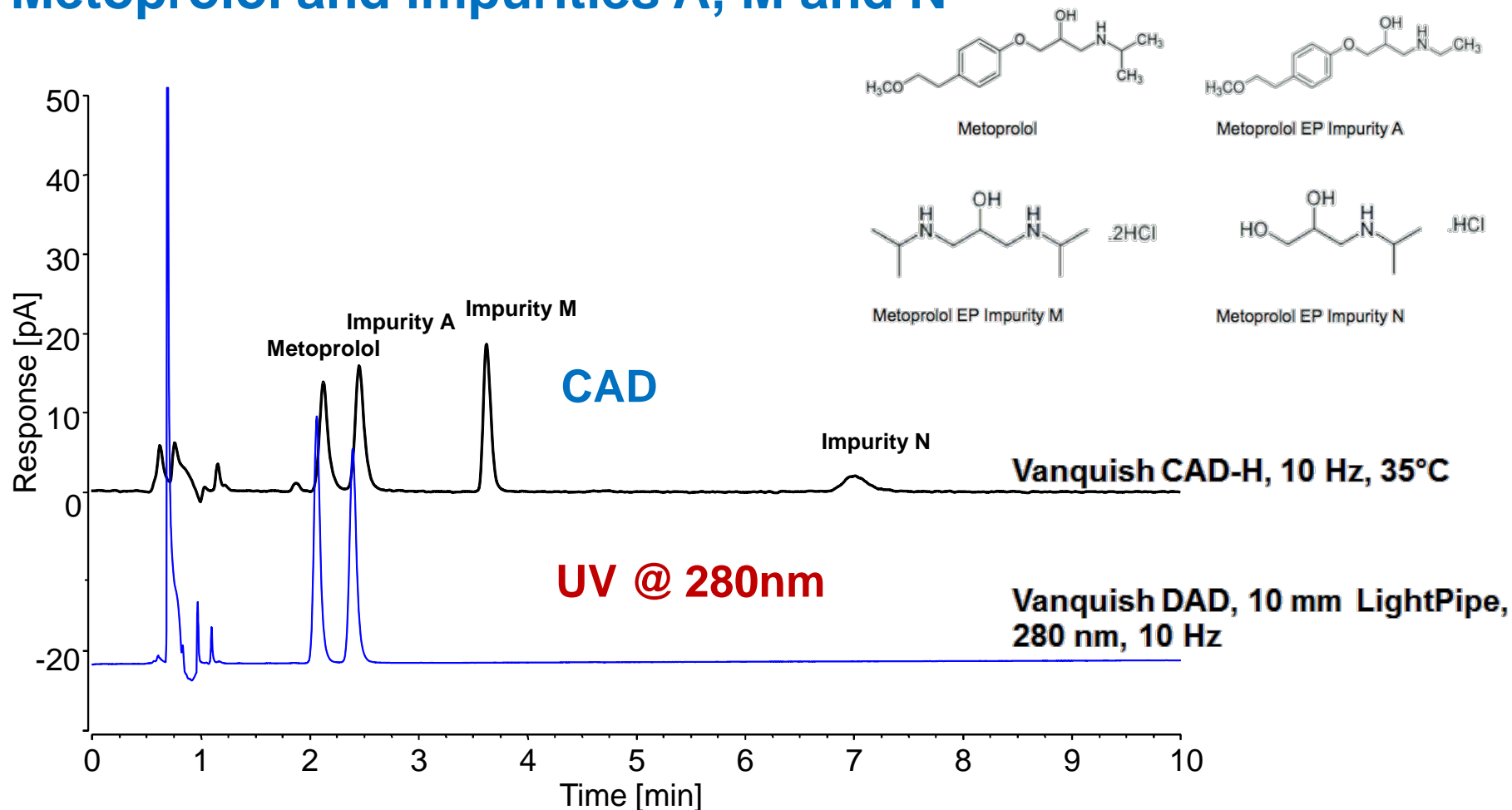
Flow rate: 0.45 mL/min

Inj. volume: 1 µL

Detector: Corona ultra RS, 15 °C, 60 Hz

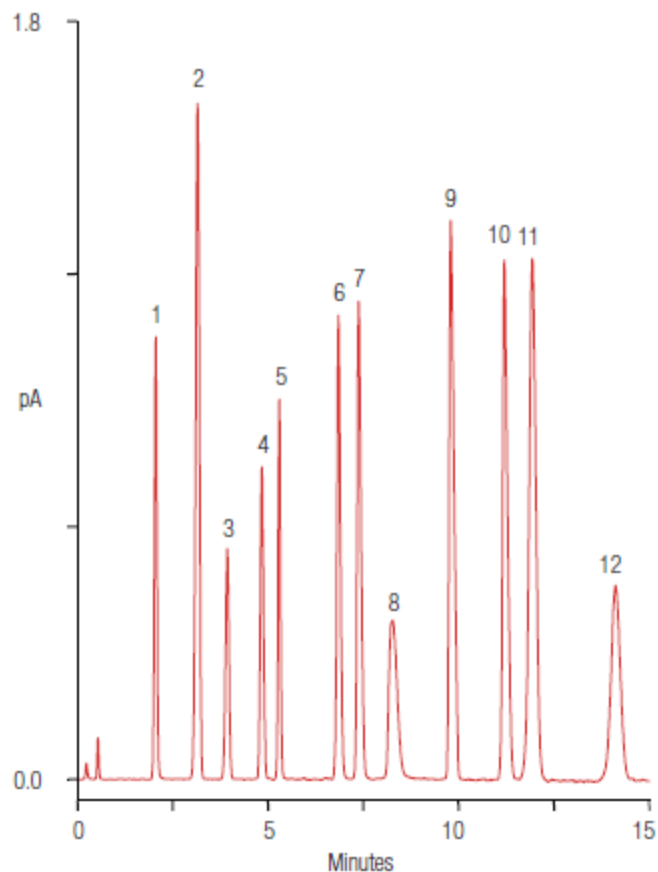
Orthogonal and Complimentary Detection with DAD and CAD

Metoprolol and impurities A, M and N



Isocratic HILIC chromatographic method using both UV and charged aerosol detection

Drug counterions



Instrumentation: Thermo Scientific™ Dionex™ UltiMate™ 3000 RSLC system
Column: Acclaim Trinity P2, 3 μm, 3 × 50 mm
Col. temp: 30 °C
Flow rate: 0.6 mL/min
Inj. volume: 2 μL
Mobile phase A: Water
Mobile phase B: 100 mM ammonium formate, pH 3.65
Gradient:

Time (min)	-8.0	0.0	1.0	11	15
%A	90	90	90	0	0
%B	10	10	10	100	100

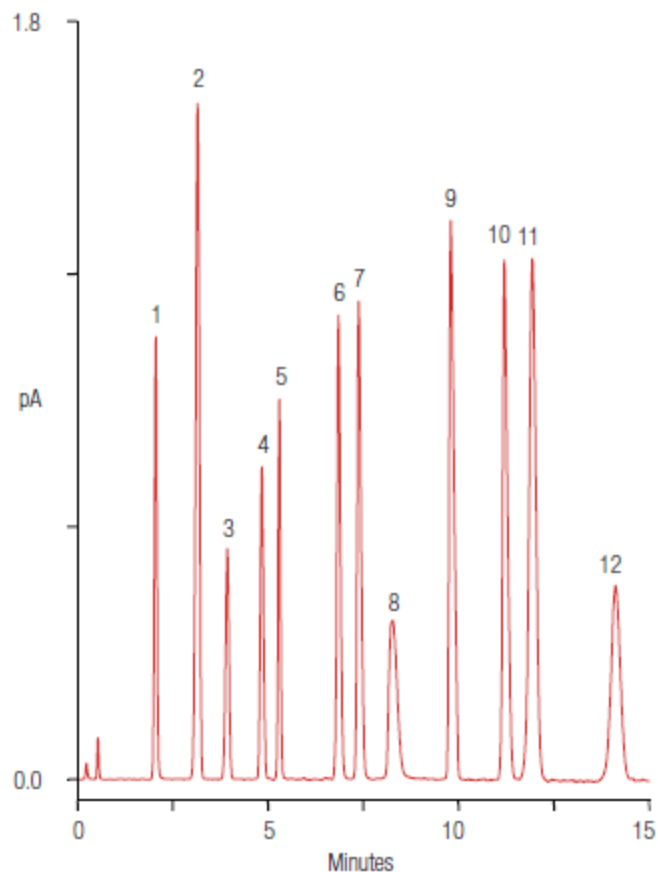
Charged aerosol: Corona Veo
RS; 55 °C, 5 Hz, 2 s, PF 1.5
Sample: 20 to 100 ng/μL each in deionized water

Peaks

- | | |
|--------------|---------------|
| 1. Phosphate | 7. Nitrate |
| 2. Sodium | 8. Citrate |
| 3. Potassium | 9. Fumarate |
| 4. Chloride | 10. Sulfate |
| 5. Malate | 11. Magnesium |
| 6. Bromide | 12. Calcium |

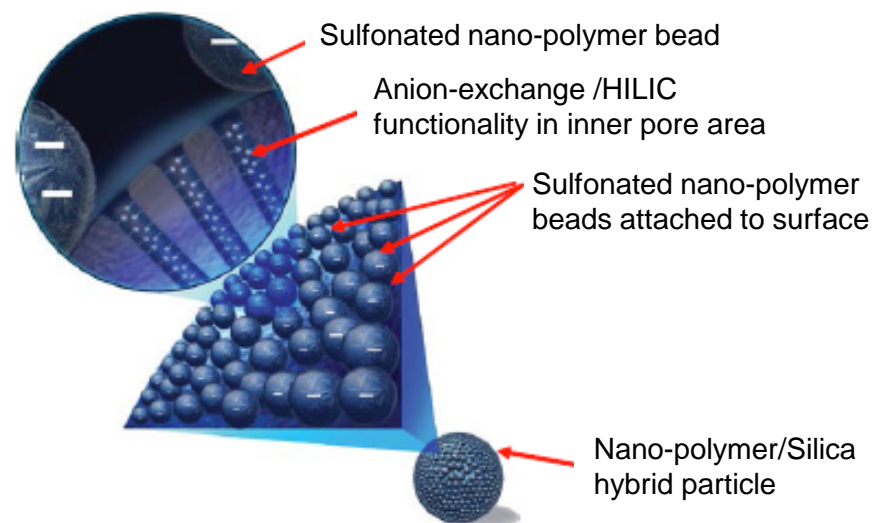
Anions, cations, organic and inorganic ions simultaneously

Drug counterions



Instrumentation: UltiMate 3000 RSLC system

Column: **Acclaim Trinity P2, 3 μ m, 3 \times 50 mm**



Peaks

- 1. Phosphate
- 2. Sodium
- 3. Potassium
- 4. Chloride
- 5. Malate
- 6. Bromide

- 7. Nitrate
- 8. Citrate
- 9. Fumarate
- 10. Sulfate
- 11. Magnesium
- 12. Calcium

Anions, cations, organic and inorganic ions simultaneously

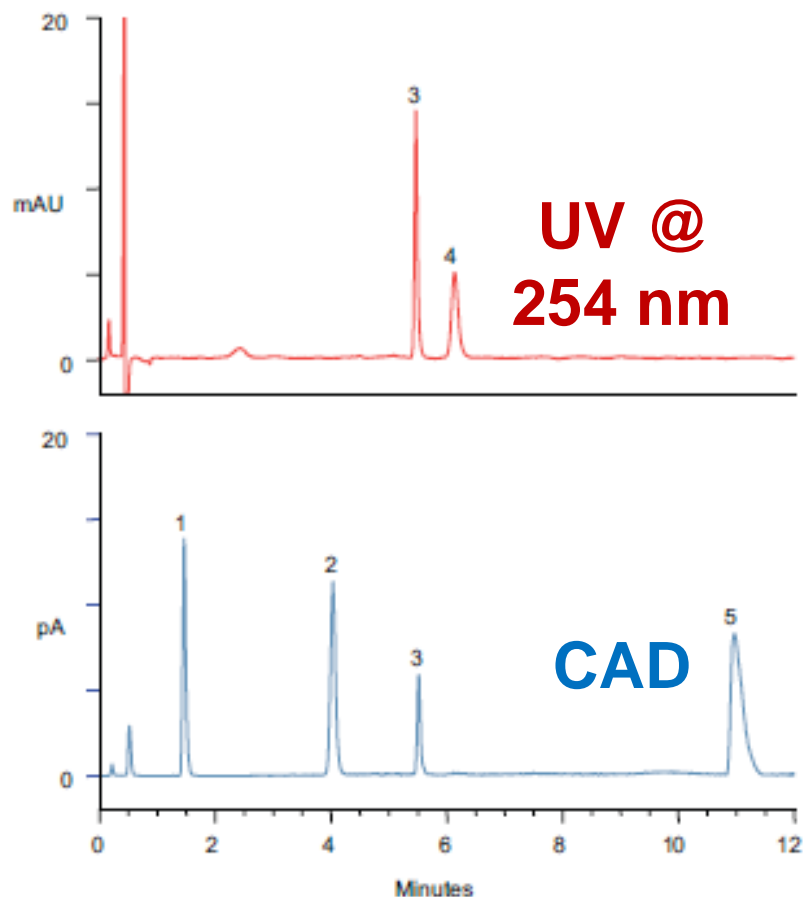
Adderall® (Shire Pharmaceuticals) and counterions

Instrumentation: UltiMate 3000 RSLC system
 Column: Acclaim Trinity P2, 3 µm, 3 × 50 mm
 Col. temp: 30 °C
 Flow rate: 0.6 mL/min
 Inj. volume: 5 µL
 Mobile phases: A: Acetonitrile
 B: Water
 C: 100 mM ammonium formate, pH 3.65

Gradient:

Time (min)	A	B	C
-8.0	35	59	6
0.0	35	59	6
0.5	35	59	6
5.0	35	0	65
10	20	0	80
12	20	0	80

UV detector: UV diode array; 254 nm, 5 Hz, 0.5 s
 Charged aerosol: Corona Veo RS; 55 °C, 5 Hz, 2 s, PF 1.5
 Peaks: 1 aspartate 24 µg/mL
 2 sodium
 3 saccharin 24 µg/mL
 4 amphetamine 122 µg/mL
 5 sulfate 26 µg/mL
 Ref: AN20870

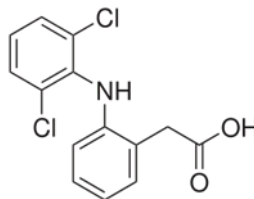
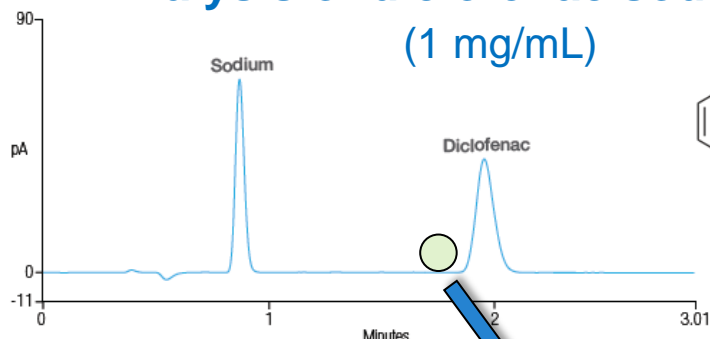


Complimentary detection by CAD and UV/Vis

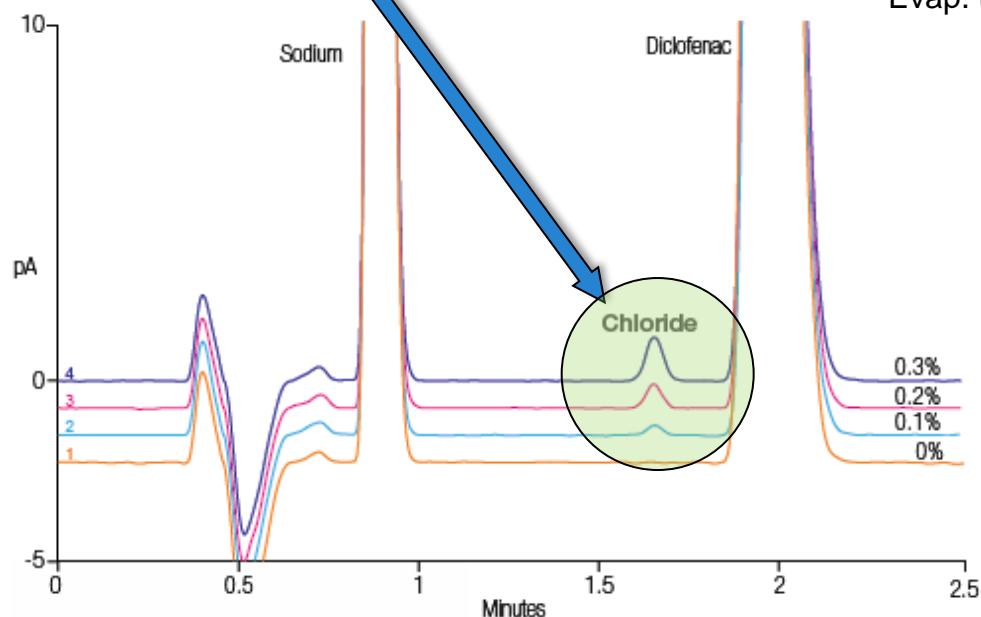
Formulation – API, Counter-Ions and Impurities

Analysis of diclofenac sodium salt

(1 mg/mL)

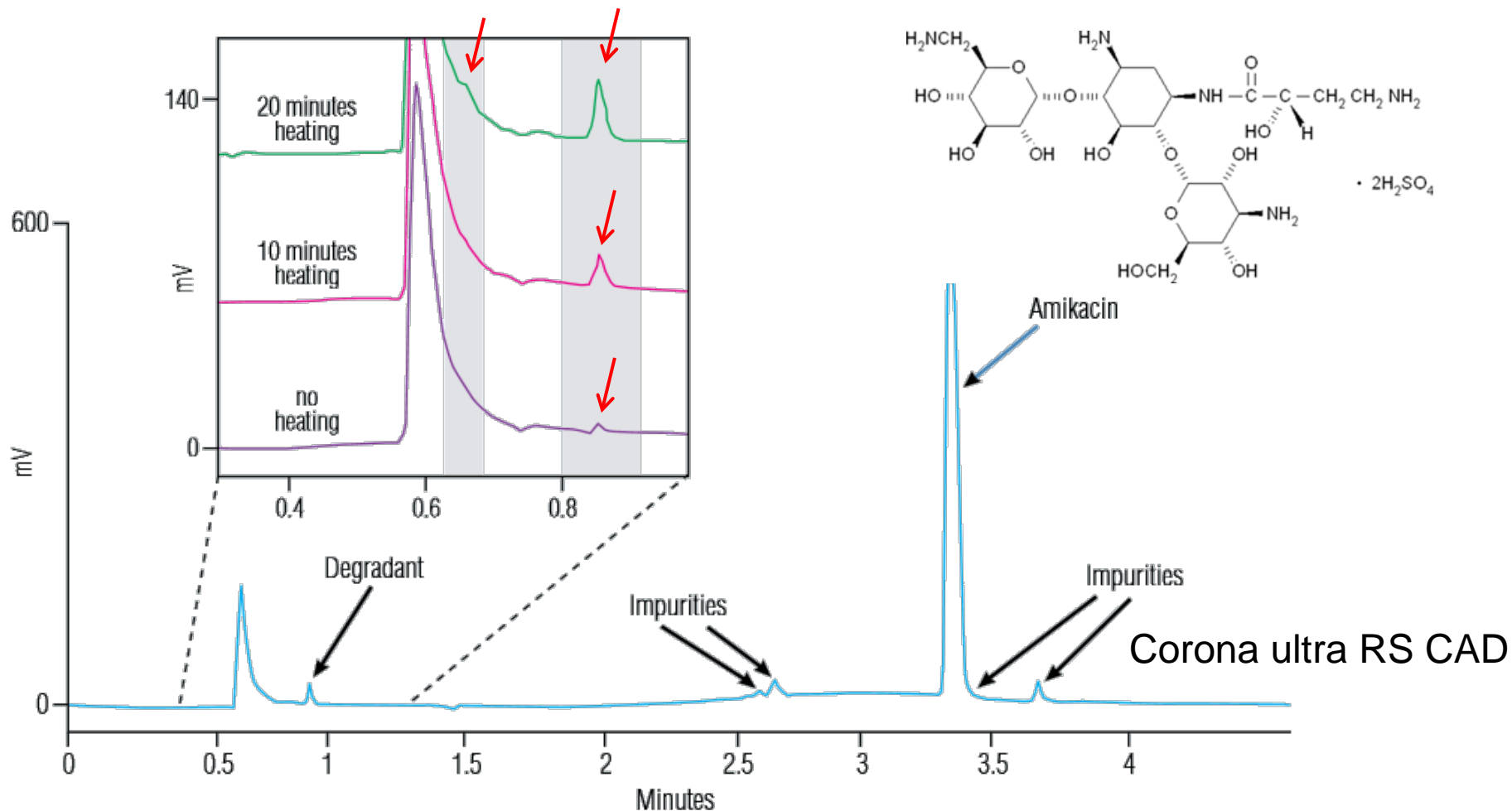


Column: Acclaim Trinity P1, 3 μ m, 3.0 \times 50 mm
Mobile phase A: 75% Acetonitrile
Mobile phase B: 25% 200 mM Ammonium acetate pH 4
Flow rate: 0.8 mL/min
Inj. volume: 5 μ L
Col. temp: 30 $^{\circ}$ C
Detector: Corona Veo RS CAD
Evap. temp: 60 $^{\circ}$ C



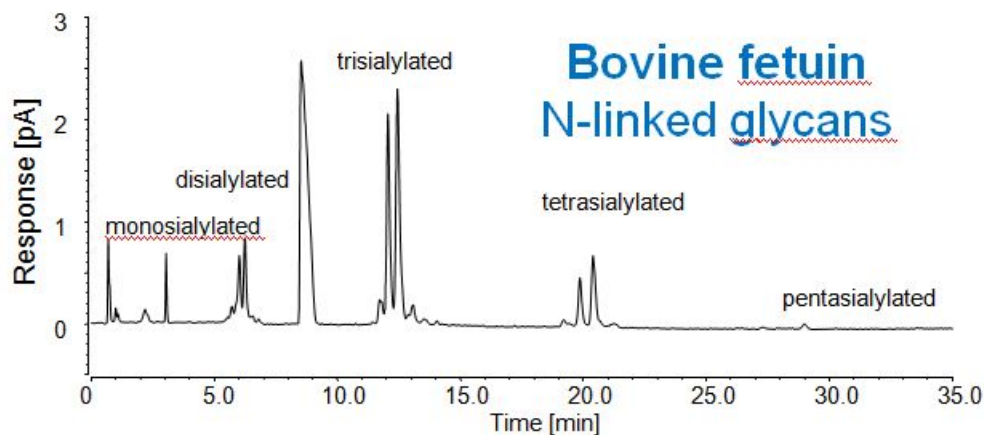
Charged aerosol even detects chloride impurity.

Stability – Forced Degradation



Follow forced degradation of Amikacin sulfate

Glycoprotein Characterization – Released Glycan Analysis



System: Thermo Scientific™ Vanquish™ UHPLC system

Column: Thermo Scientific™ GlycanPac™ AXR-1, 1.9 μ m, 2.1 \times 150 mm

Mobile phase A: Deionized water

Mobile phase B: 1.00 mM Ammonium formate, pH 4.4

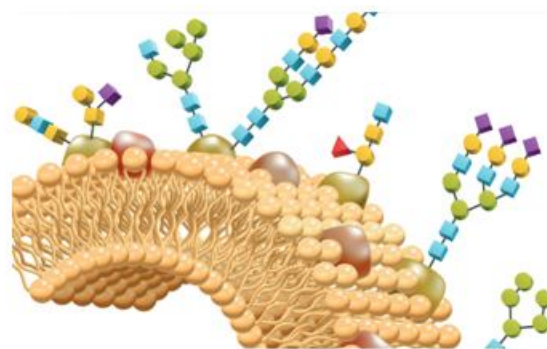
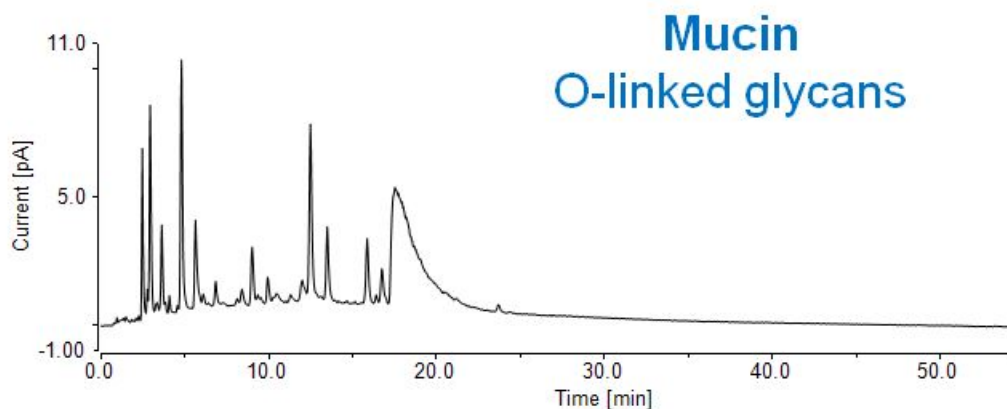
Gradient: 4 % B to 39% B in 35 min

Flow rate: 0.4 mL/min

Inj. volume: 2 μ L

Col. temp: 30° C

Detector: Vanquish Charged Aerosol Detector H
50° C, PF 1.0, 10 Hz, 5s



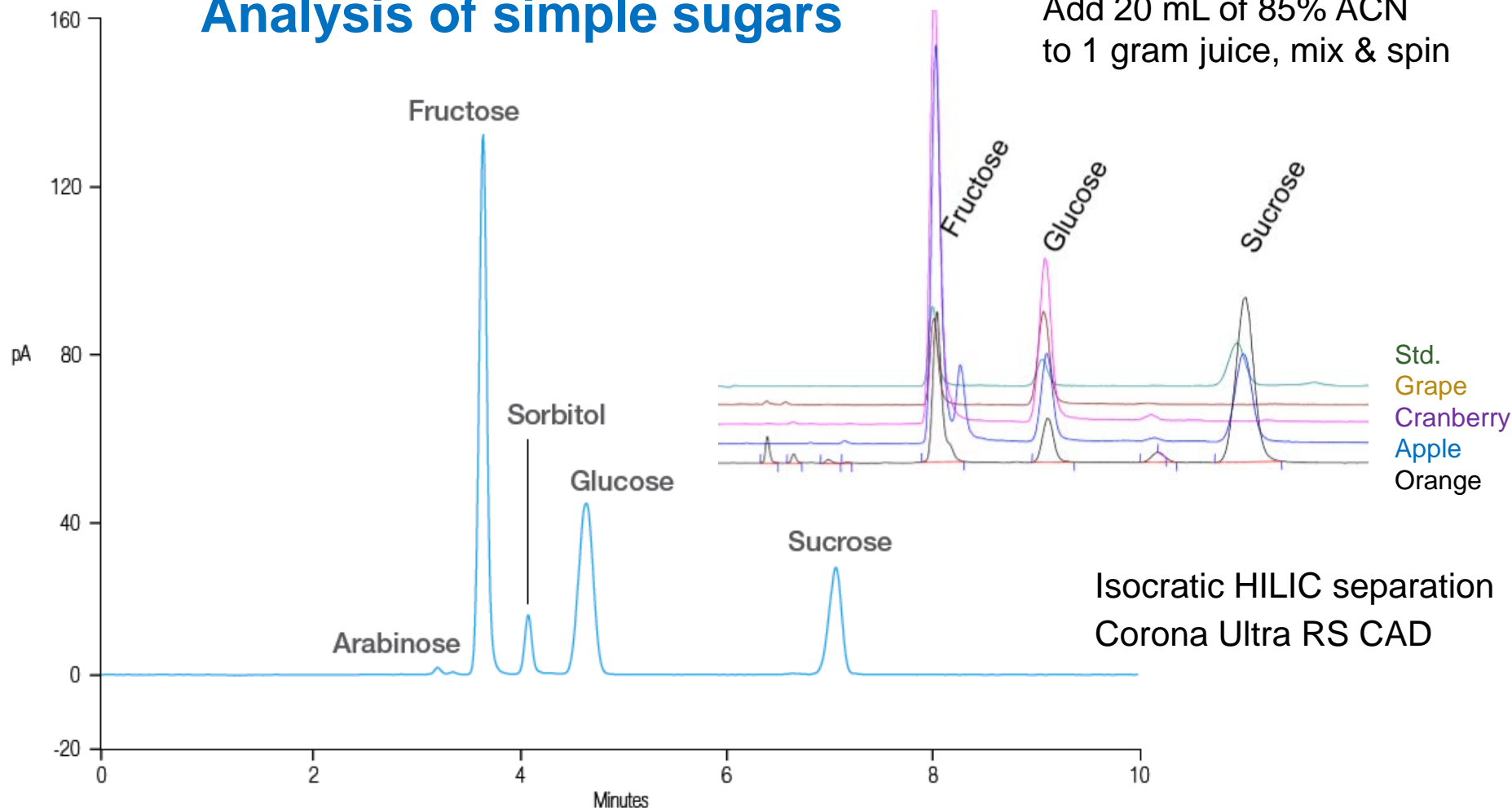
Detection of released glycans – No labeling required

Food and Beverage Application Areas

- Simple carbohydrates
- Lipids
 - Profiling methods
 - Targeted methods
- Artificial sweeteners

Analysis of simple sugars

Add 20 mL of 85% ACN
to 1 gram juice, mix & spin



Simplified sample preparation “Dilute-and-shoot” method

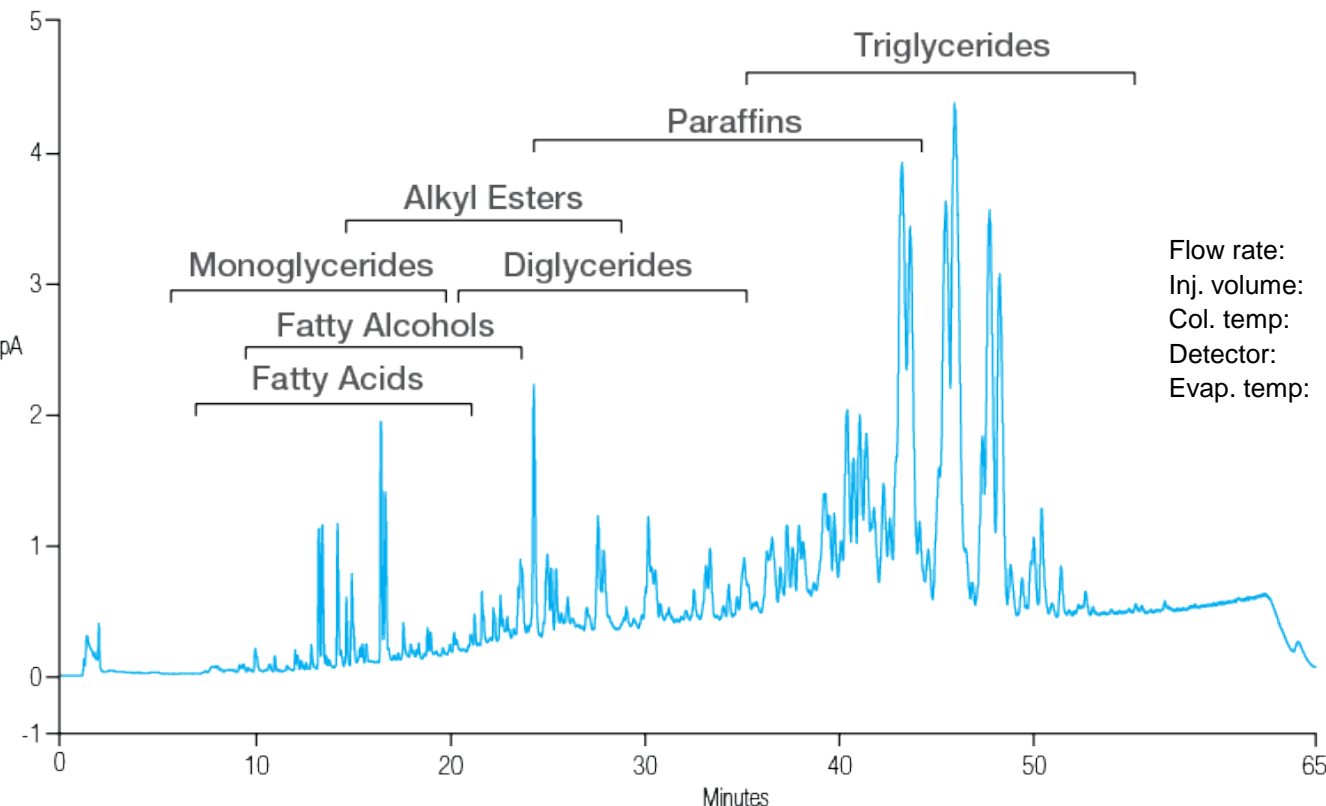
Global lipids – Algal oils

Column: Thermo Scientific™ Accucore™ C18, 2.6 μm, 3.0x150 mm
 Mobile phases: A: Methanol:water:acetic acid (600:400:4)
 B: Tetrahydrofuran:acetonitrile (50:950)
 C: Acetone:acetonitrile (900:100)

Gradient:

Time (min)	FlowRate (mL/min)	%A	%B	%C
-10.0	1.00	90	10	0
-0.1	1.00	90	10	0
0.0	0.25	90	10	0
20.0	0.50	15	85	0
35.0	0.50	2	78	20
60.0	0.50	2	3	95
65.0	0.50	90	10	0

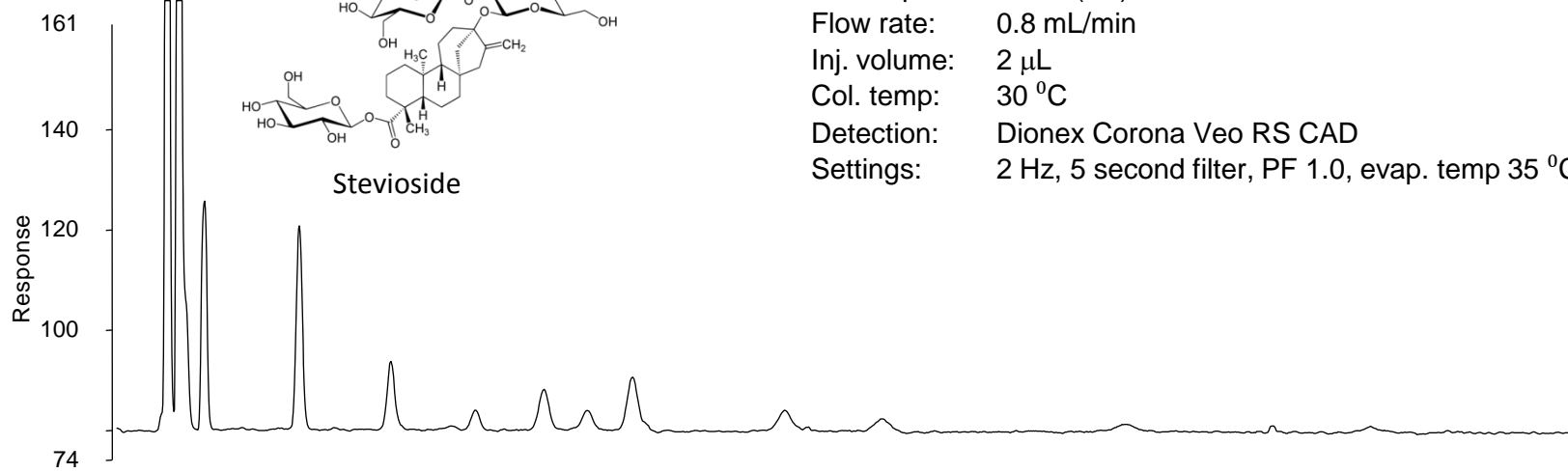
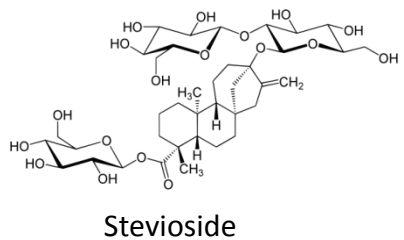
Flow rate: 1.0 mL/min
 Inj. volume: 2 μL
 Col. temp: 40 °C
 Detector: Corona Veo RS CAD
 Evap. temp: 40 °C



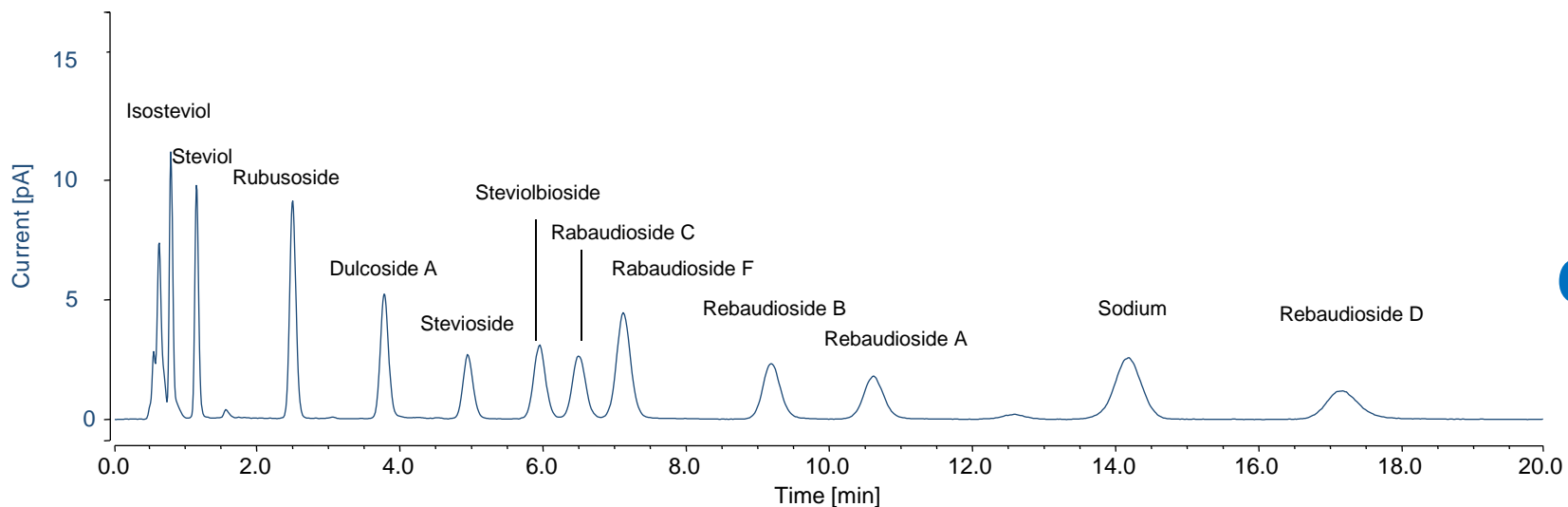
Complex sample – Minimal sample prep

Food Additives: Steviol Glycosides (Sweetener)

Column: Acclaim Trinity P1, 3 μm , 2.1 \times 150 mm
Mobile phase: 88:12 (v/v) Acetonitrile:10 mM ammonium formate, pH 3.1
Flow rate: 0.8 mL/min
Inj. volume: 2 μL
Col. temp: 30 $^{\circ}\text{C}$
Detection: Dionex Corona Veo RS CAD
Settings: 2 Hz, 5 second filter, PF 1.0, evap. temp 35 $^{\circ}\text{C}$



ELSD

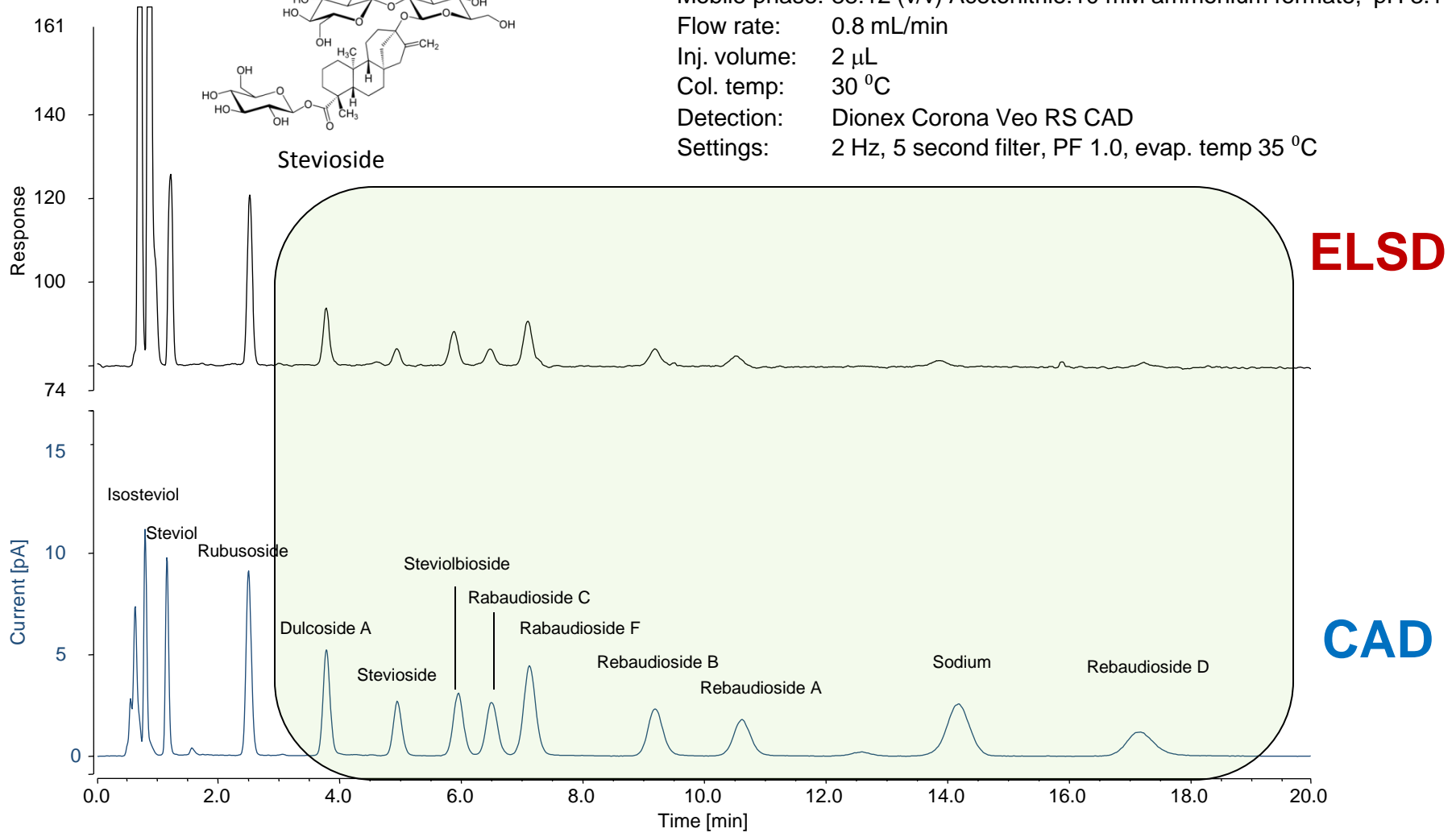
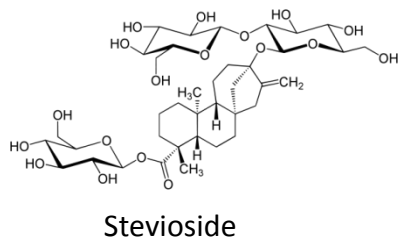


CAD

CAD exhibits more uniform response.

Artificial Sweeteners: Steviol Glycosides

Column: Acclaim Trinity P1, 3 μm , 2.1 \times 150 mm
Mobile phase: 88:12 (v/v) Acetonitrile:10 mM ammonium formate, pH 3.1
Flow rate: 0.8 mL/min
Inj. volume: 2 μL
Col. temp: 30 $^{\circ}\text{C}$
Detection: Dionex Corona Veo RS CAD
Settings: 2 Hz, 5 second filter, PF 1.0, evap. temp 35 $^{\circ}\text{C}$



CAD exhibits more uniform response.

Summary

- Charged aerosol detection delivers accurate and precise quantification of lipids, carbohydrates, surfactants, amines and counterions that UV/Vis absorbance cannot detect.
- For analytes with chromophores, charged aerosol detection provides uniform response independent of the extinction coefficient.
- Charged aerosol detection provides a good estimate of the amount of unknown impurities and degradation products.
- Charged aerosol detection is superior to ELSD in terms of sensitivity, dynamic range, response uniformity, precision and ease of use.
- More information on charged aerosol detection can be found at www.thermofisher.com/cad
- Bibliography of charged aerosol detector applications can be downloaded from <http://analyteguru.com/resources/charged-aerosol-detection-list-of-published-articles/>

Thank You Very Much for Your Attention!



Questions?

**Do you have additional questions
or do you want to talk to an expert
from Thermo Fisher Scientific?**

Please send an E-Mail to
analyze.eu@thermofisher.com
and we will get back to you.