

Pesticide Analysis FAQs

Julie Kowalski
Owner/ Technical Consultant
JA Kowalski Science Support
julie@kowalskiscience.com
www.kowalskiscience.com

DE.3001736111



goal is to provide practice information



focus on

GC and LC tandem mass spectrometry

- QQQ
- MRM or transitions

Matrix-matched calibration

important references

Document N° SANTE/12682/2019

ANALYTICAL QUALITY CONTROL AND METHOD VALIDATION PROCEDURES
FOR PESTICIDE RESIDUES ANALYSIS IN FOOD AND FEED

https://www.eurl-pesticides.eu/userfiles/file/EurlALL/AqcGuidance_SANTE_2019_12682.pdf

**Guidelines for the Validation of Chemical Methods
for the FDA FVM Program, 3rd Ed.**

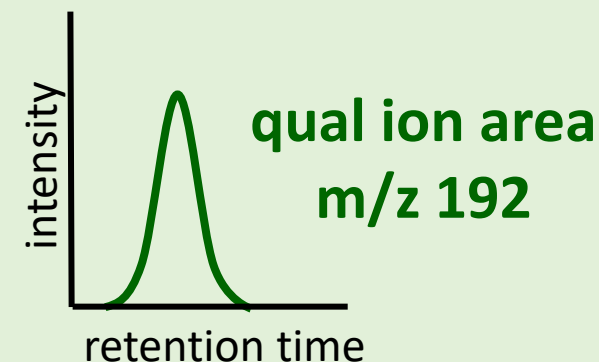
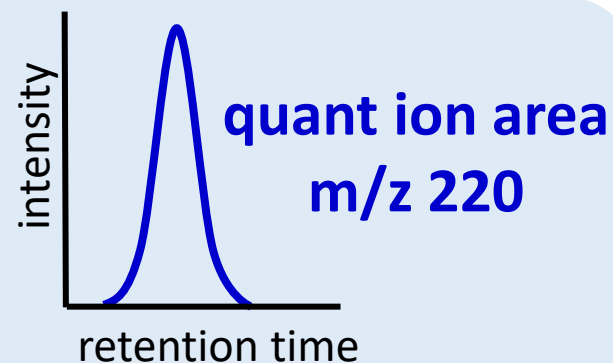
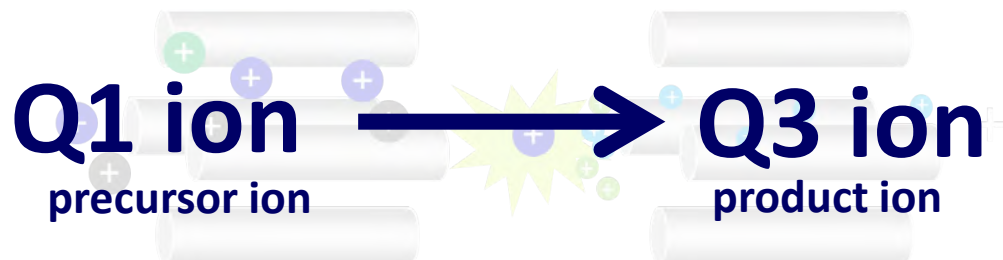
**US Food & Drug Administration
Office of Foods and Veterinary Medicine**

**Guidelines for the Validation of Chemical Methods
for the FDA FVM Program, 3rd Edition**

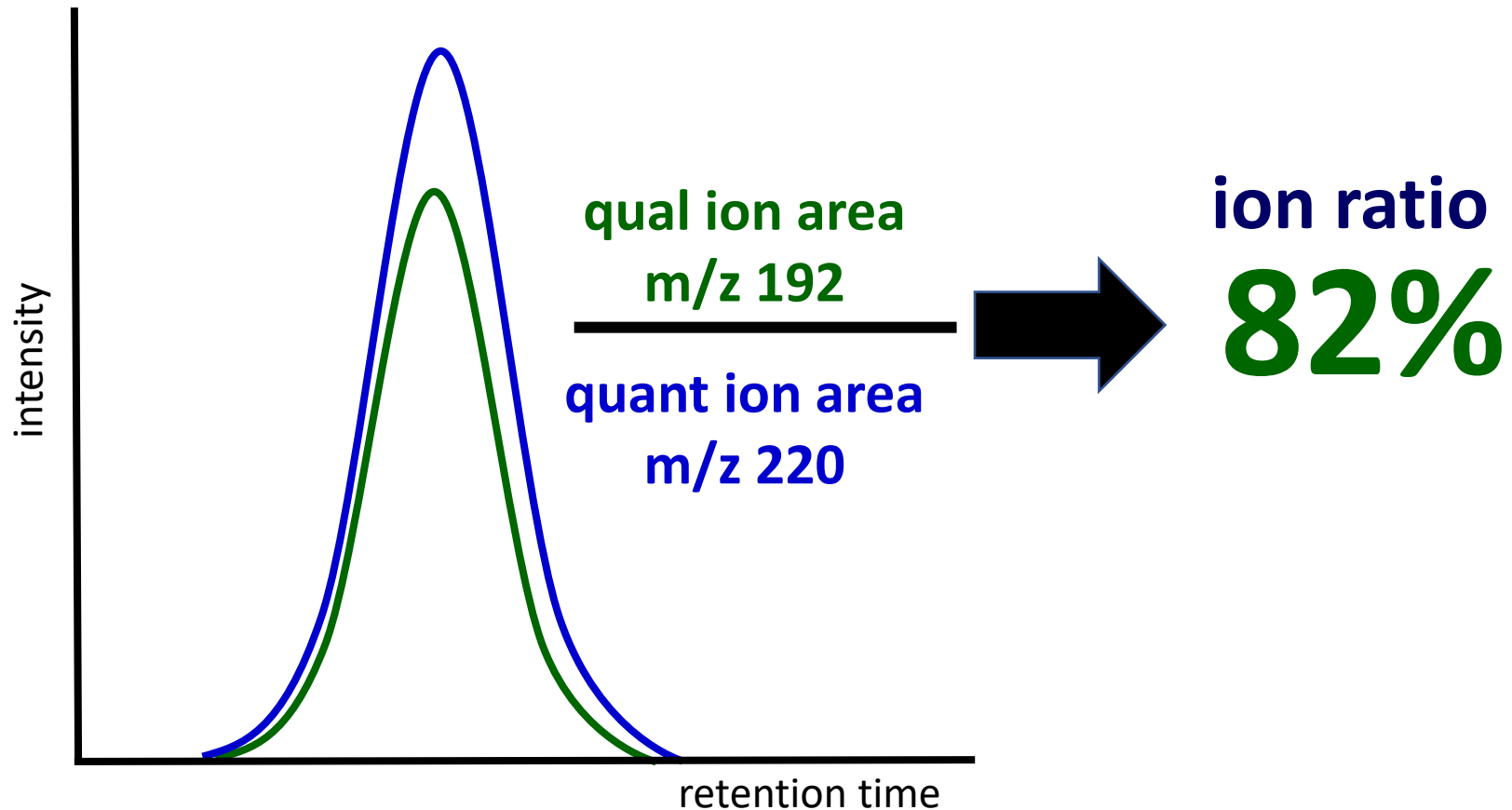
<https://www.fda.gov/media/81810/download>

DE.3001736111

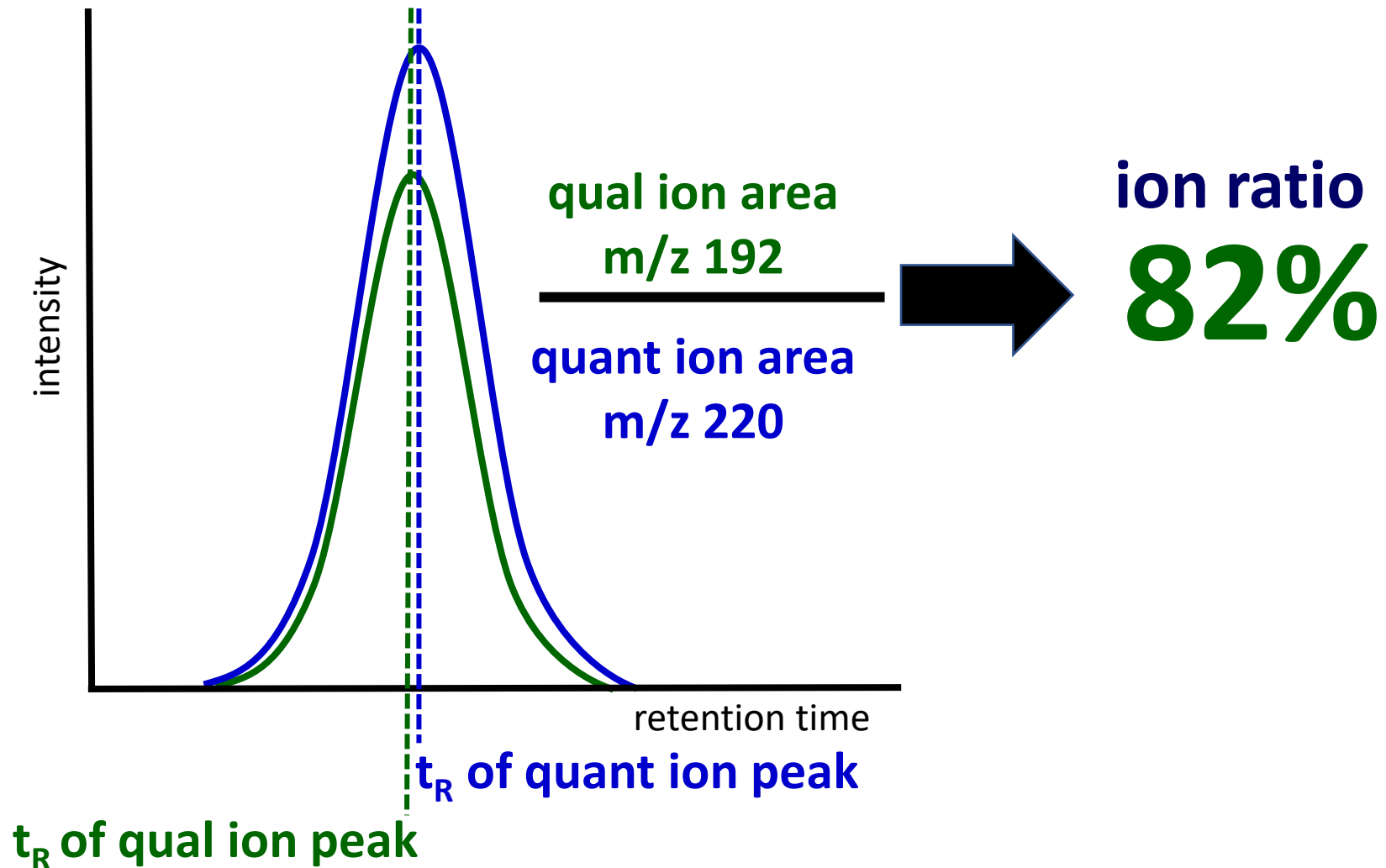
product ion peaks



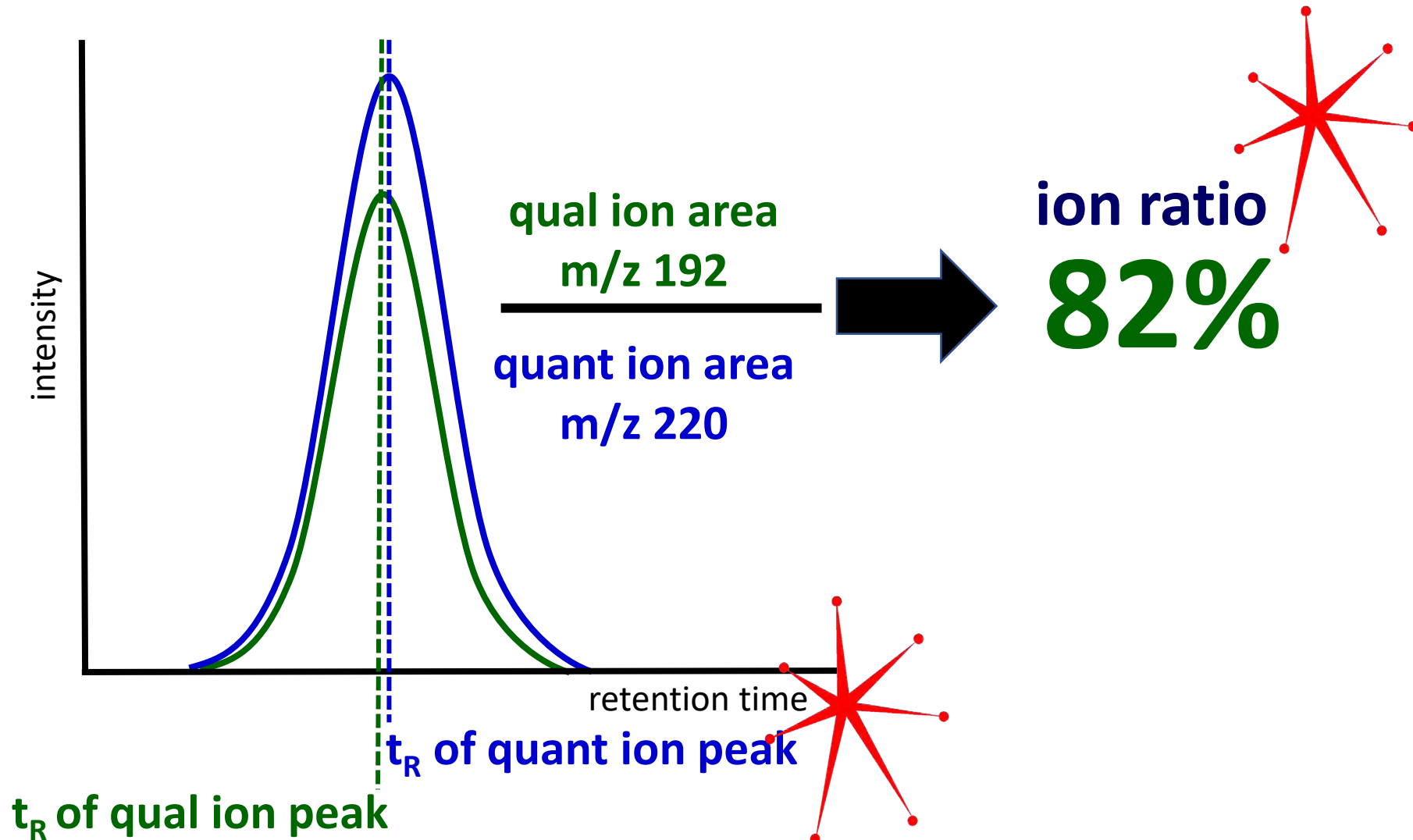
identification data



identification data

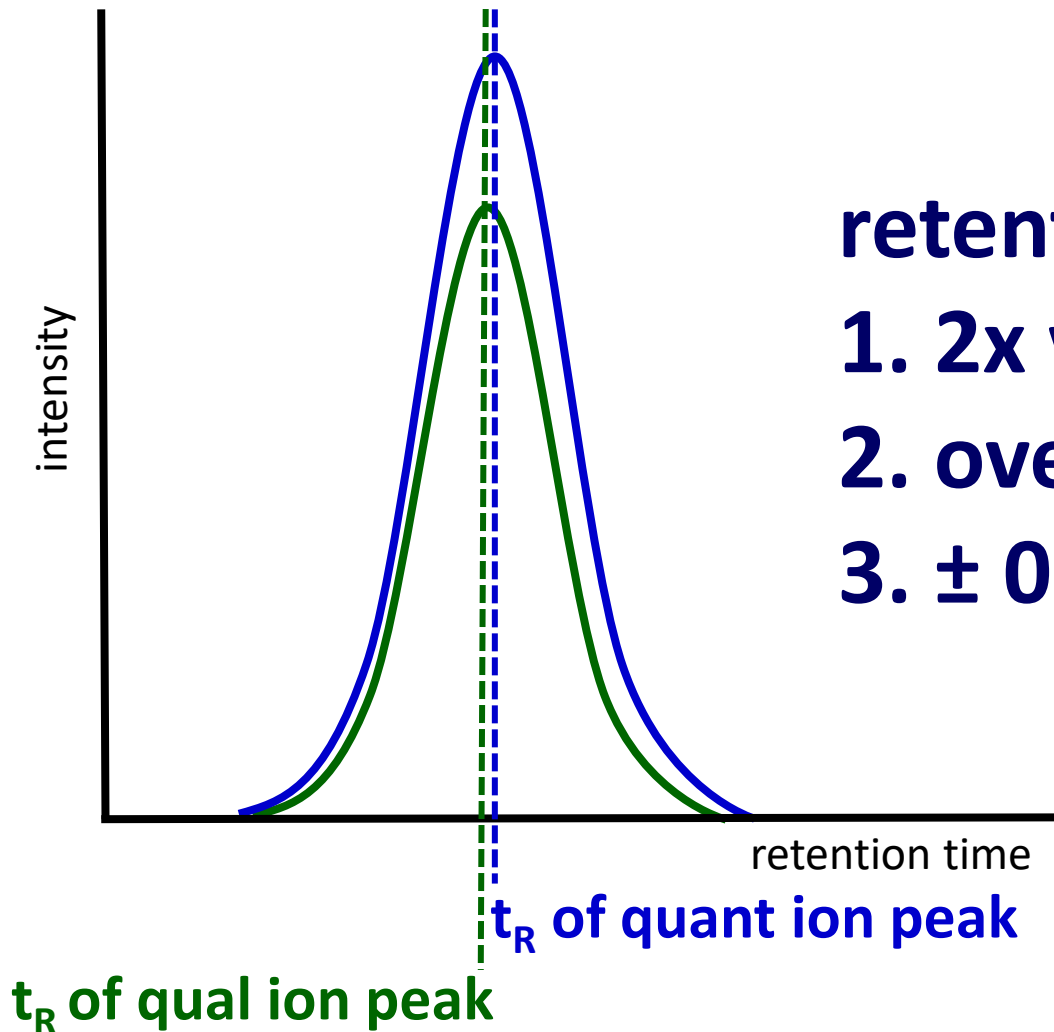


identification data



identification data –retention time

quant ion
qual ion



retention time

1. 2x void

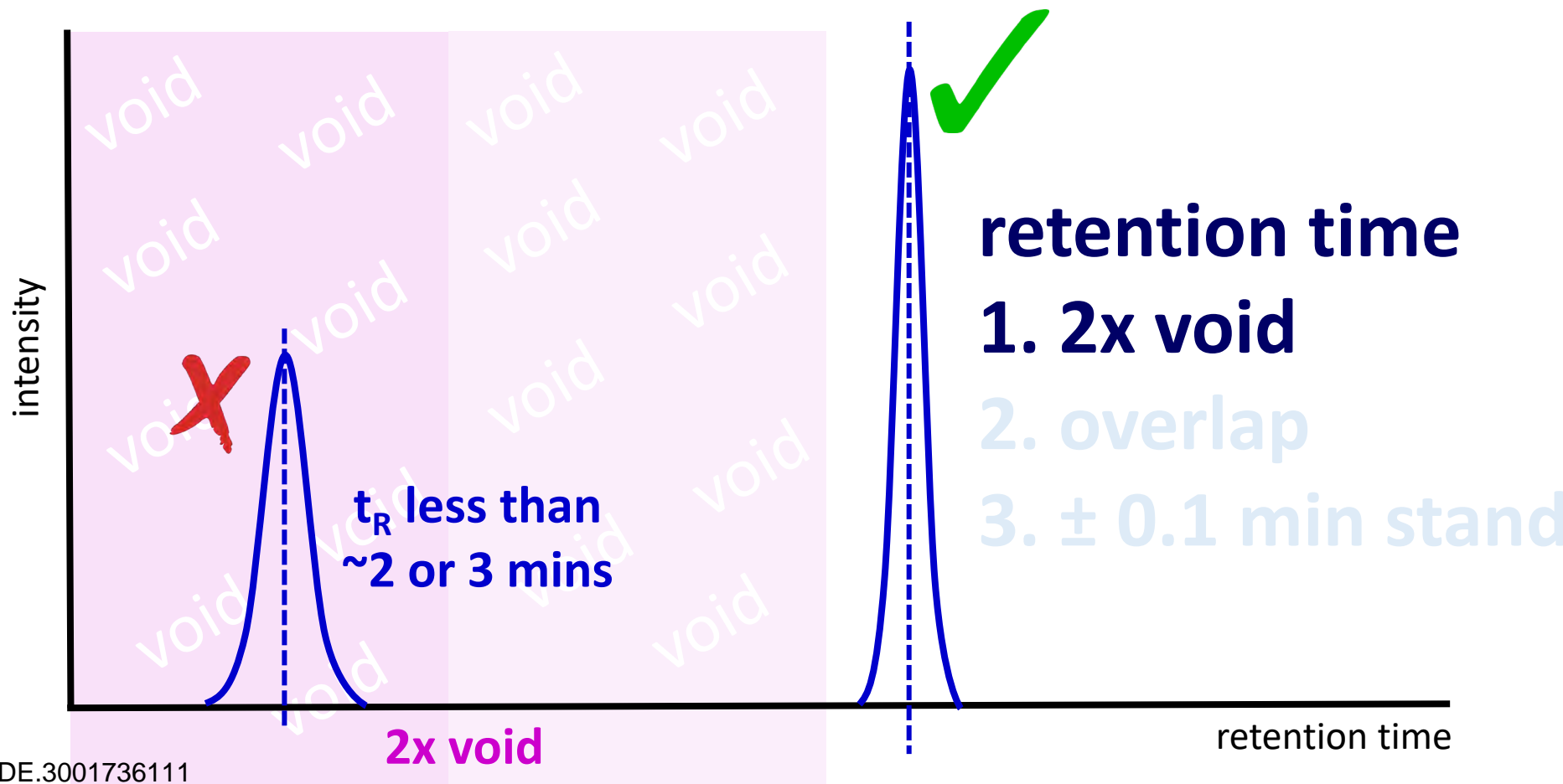
2. overlap

3. ± 0.1 min standard

identification data –retention time

“real” retention for polar pesticides

daminozide, chlormequat chloride, methamidophos, acephate



identification data –retention time

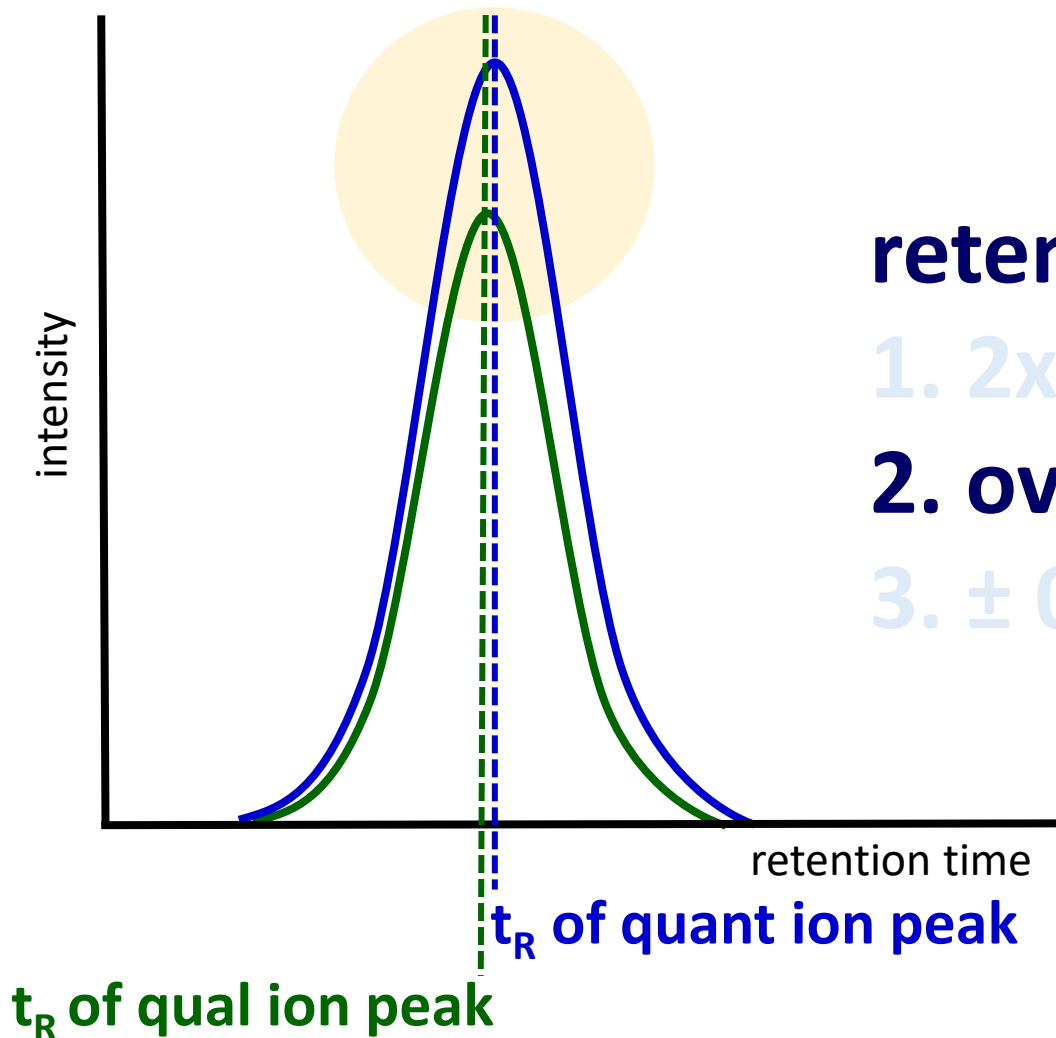
quant ion
qual ion

retention time

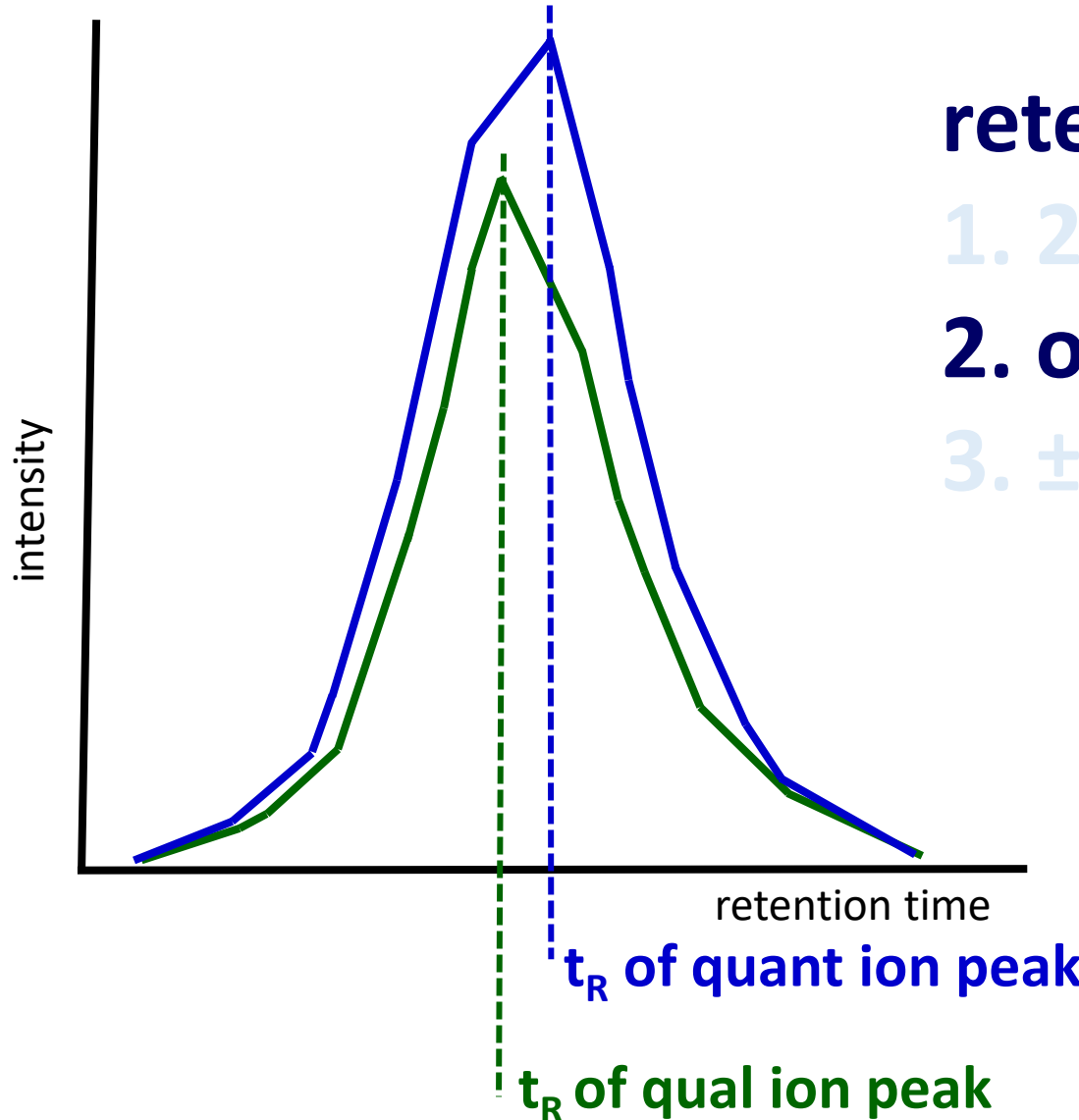
1. 2x void

2. overlap

3. ± 0.1 min standard



identification data –retention time



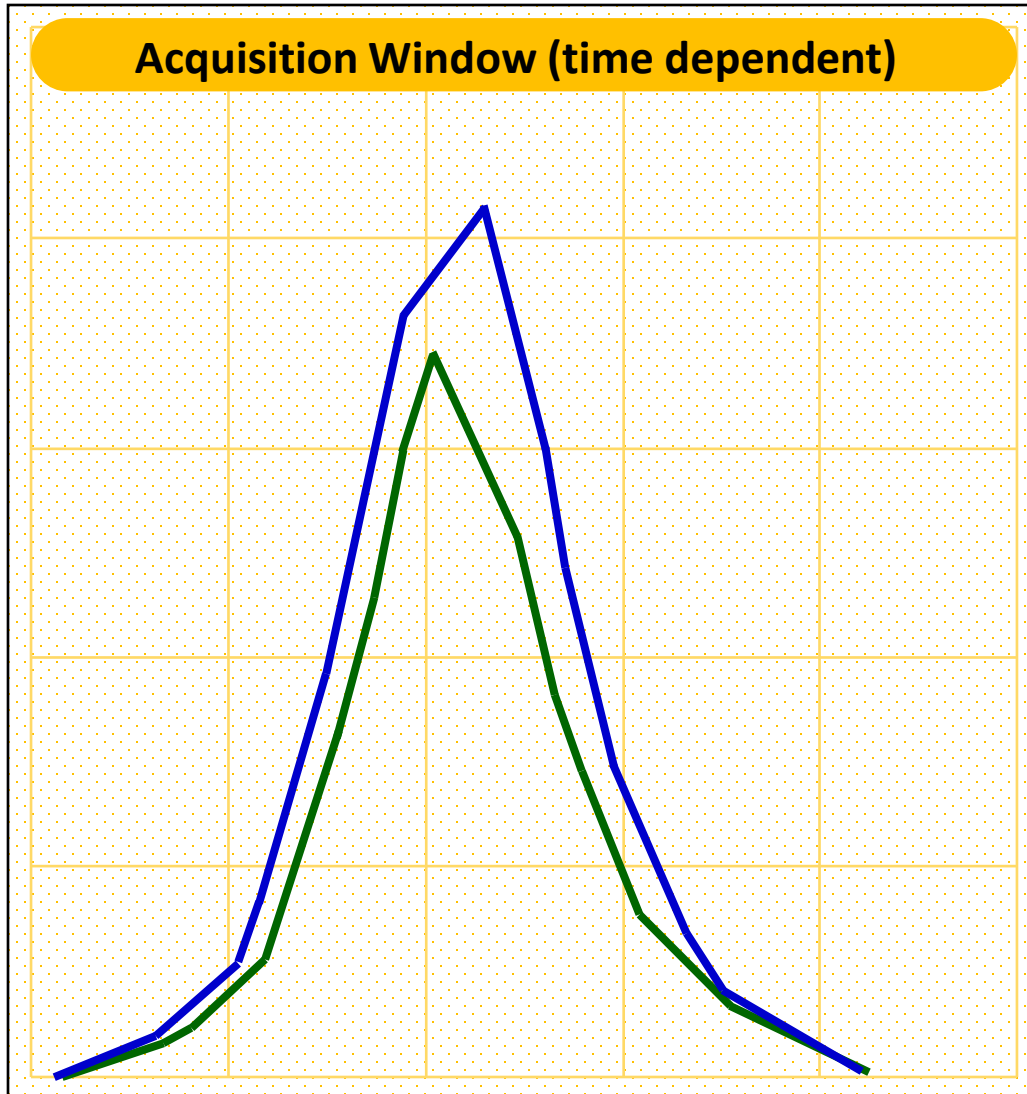
retention time

1. 2x void

2. overlap

3. ± 0.1 min standard

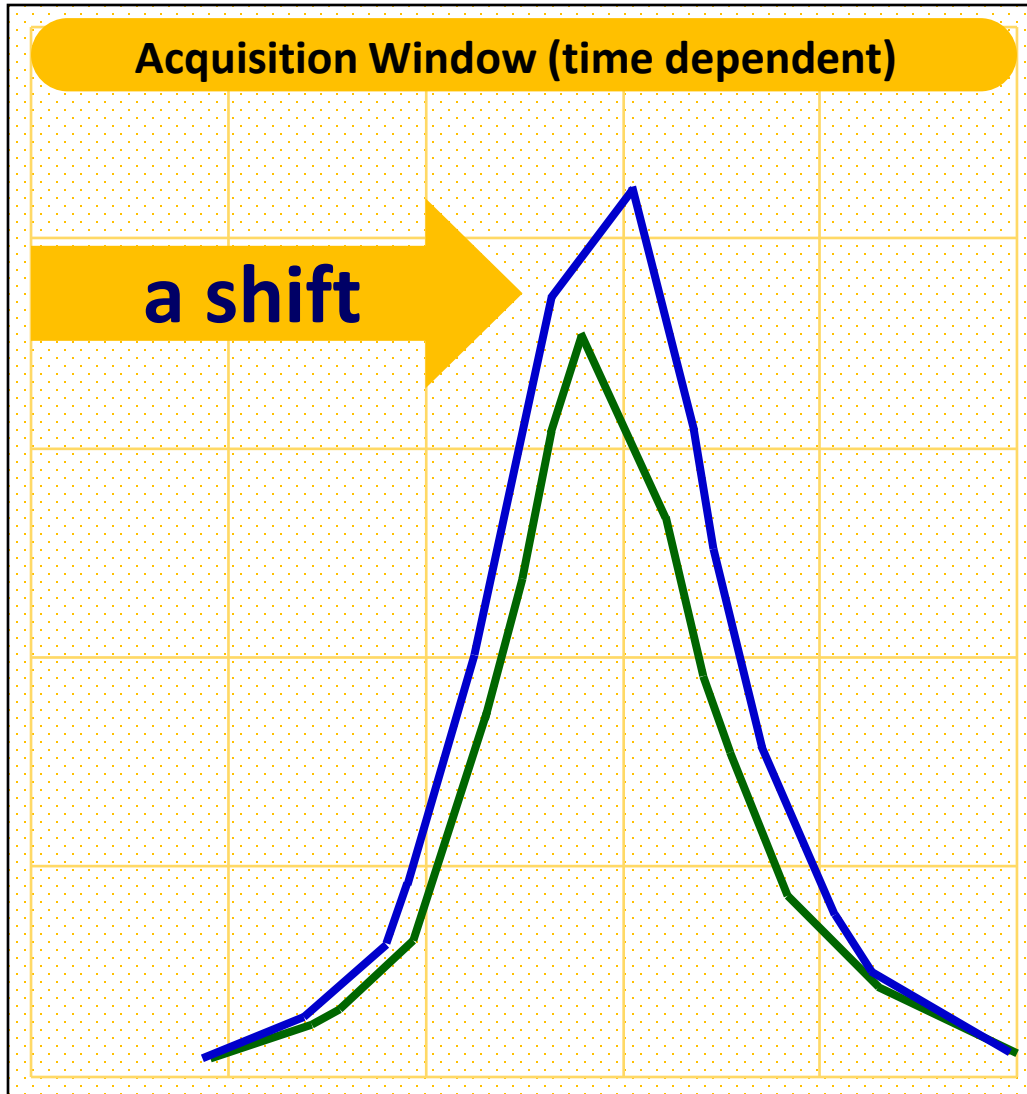
identification data –retention time



**sensitivity
&
robustness**

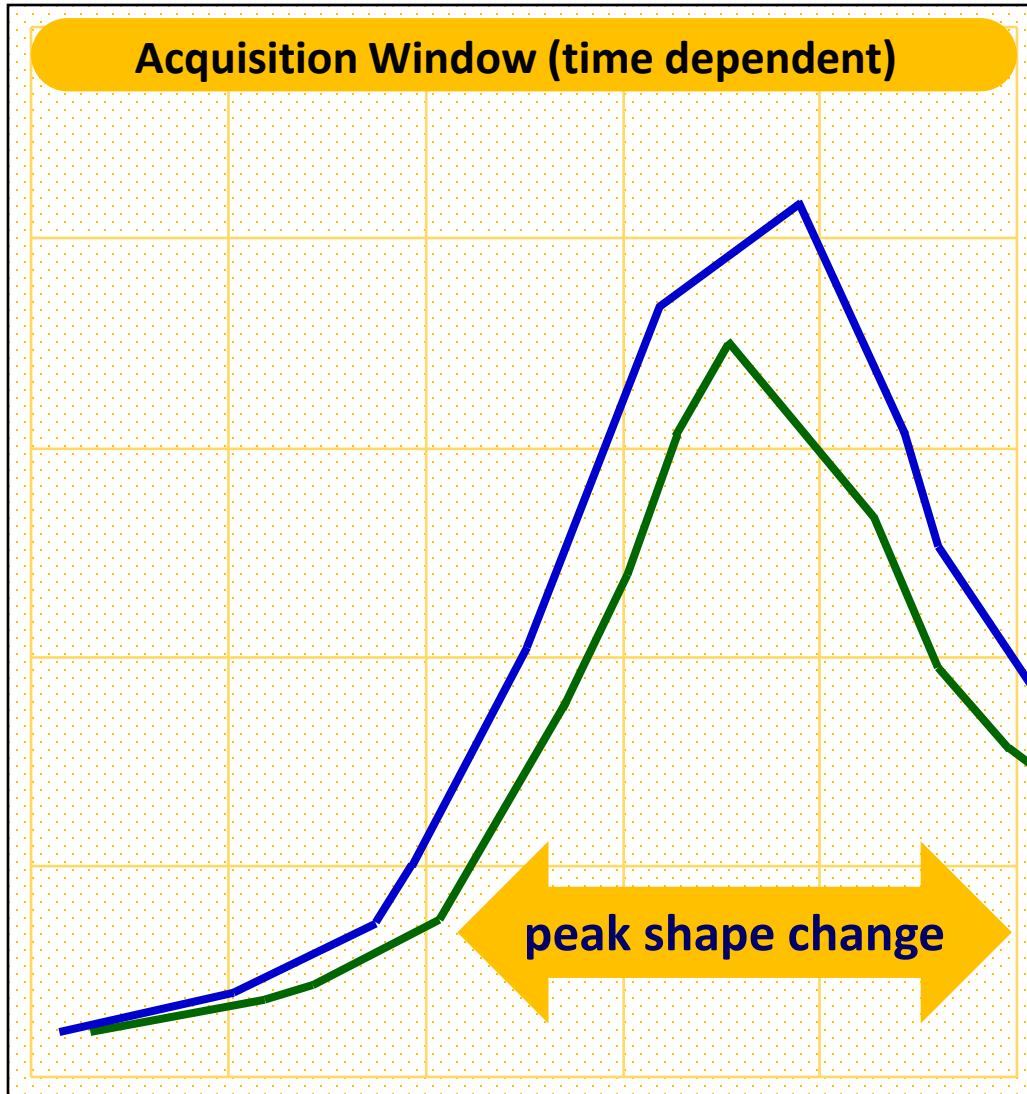
**recommend a
conservative
approach**

identification data –retention time



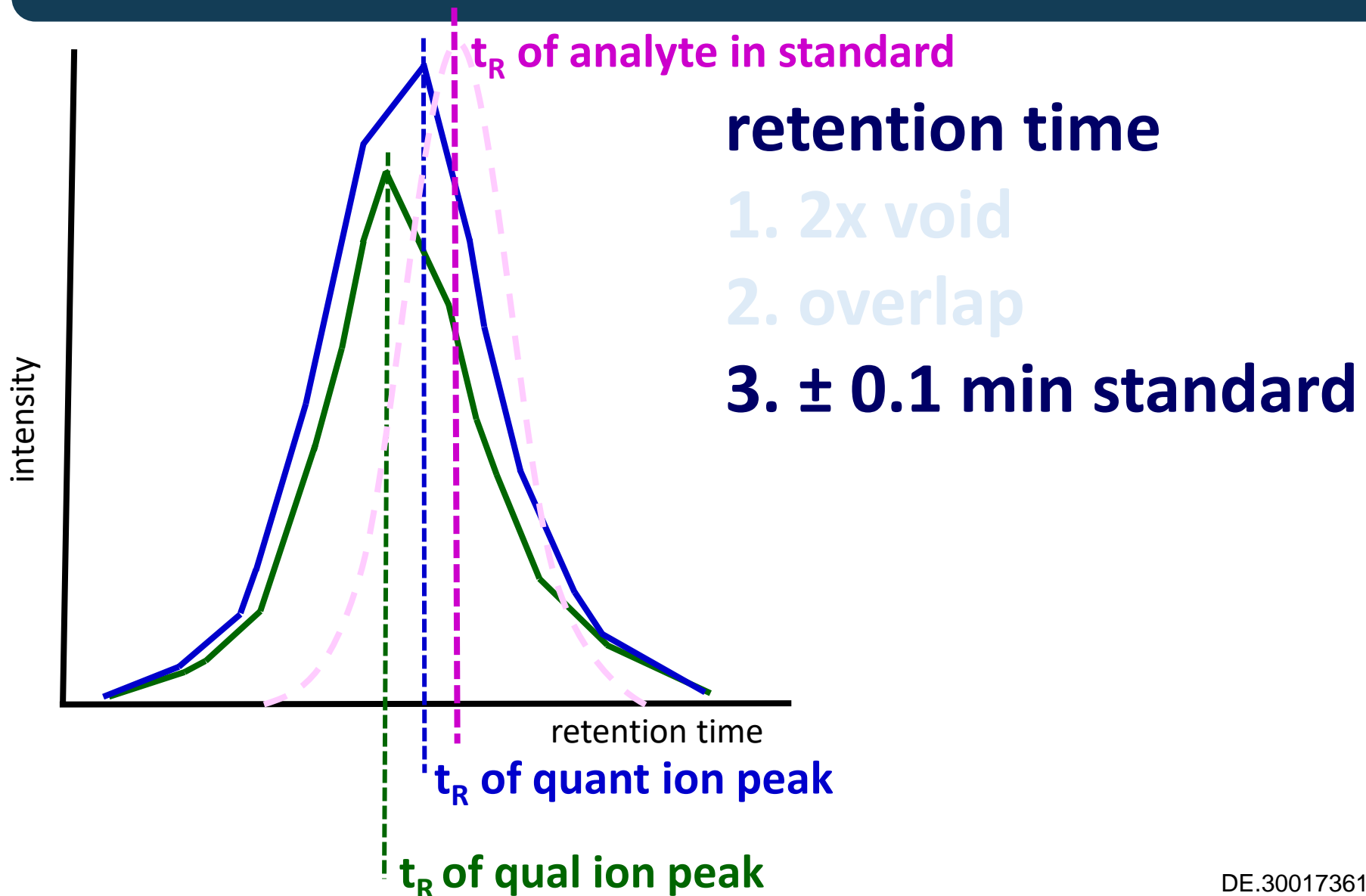
after a
few batches...

identification data –retention time

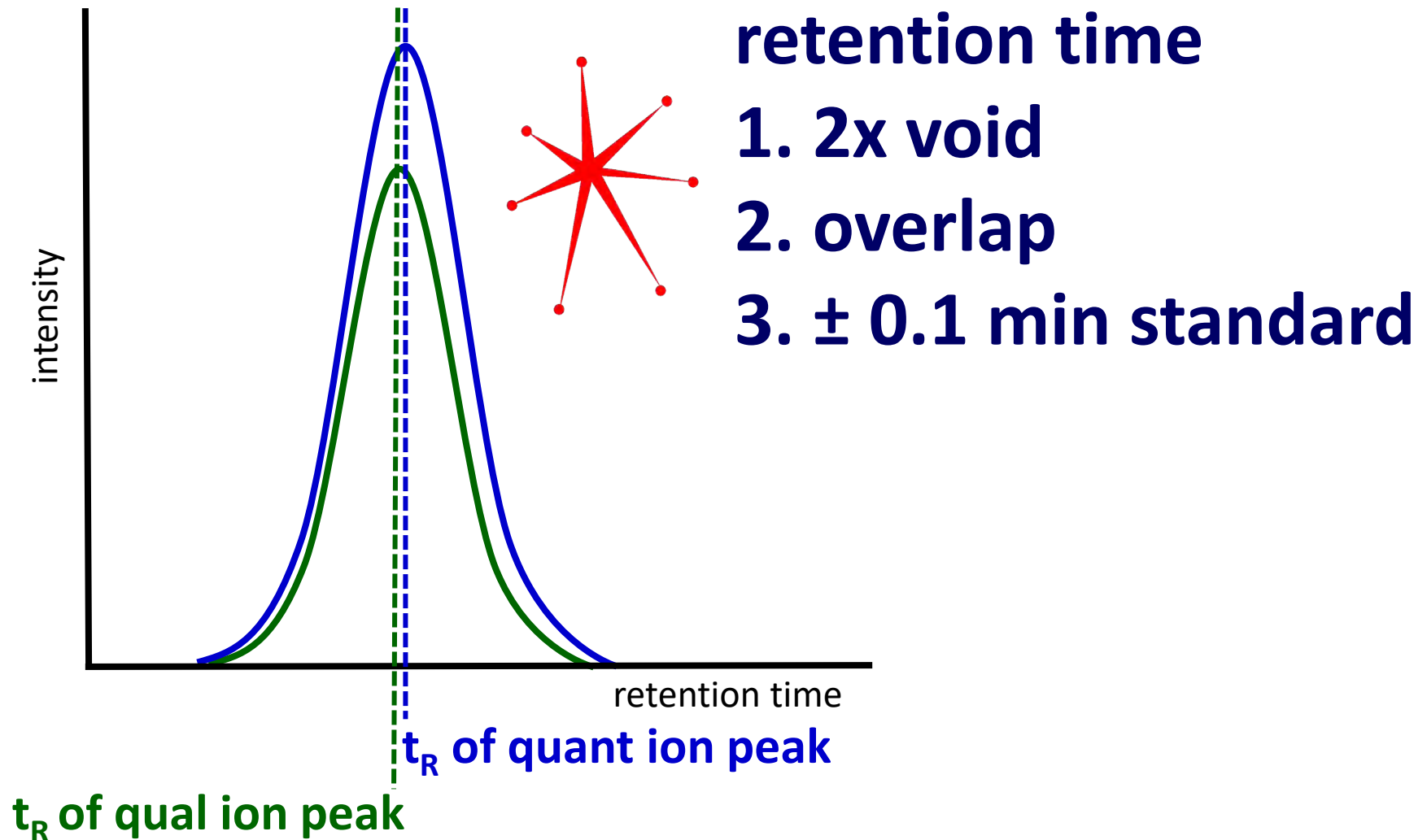


after a
few more batches...

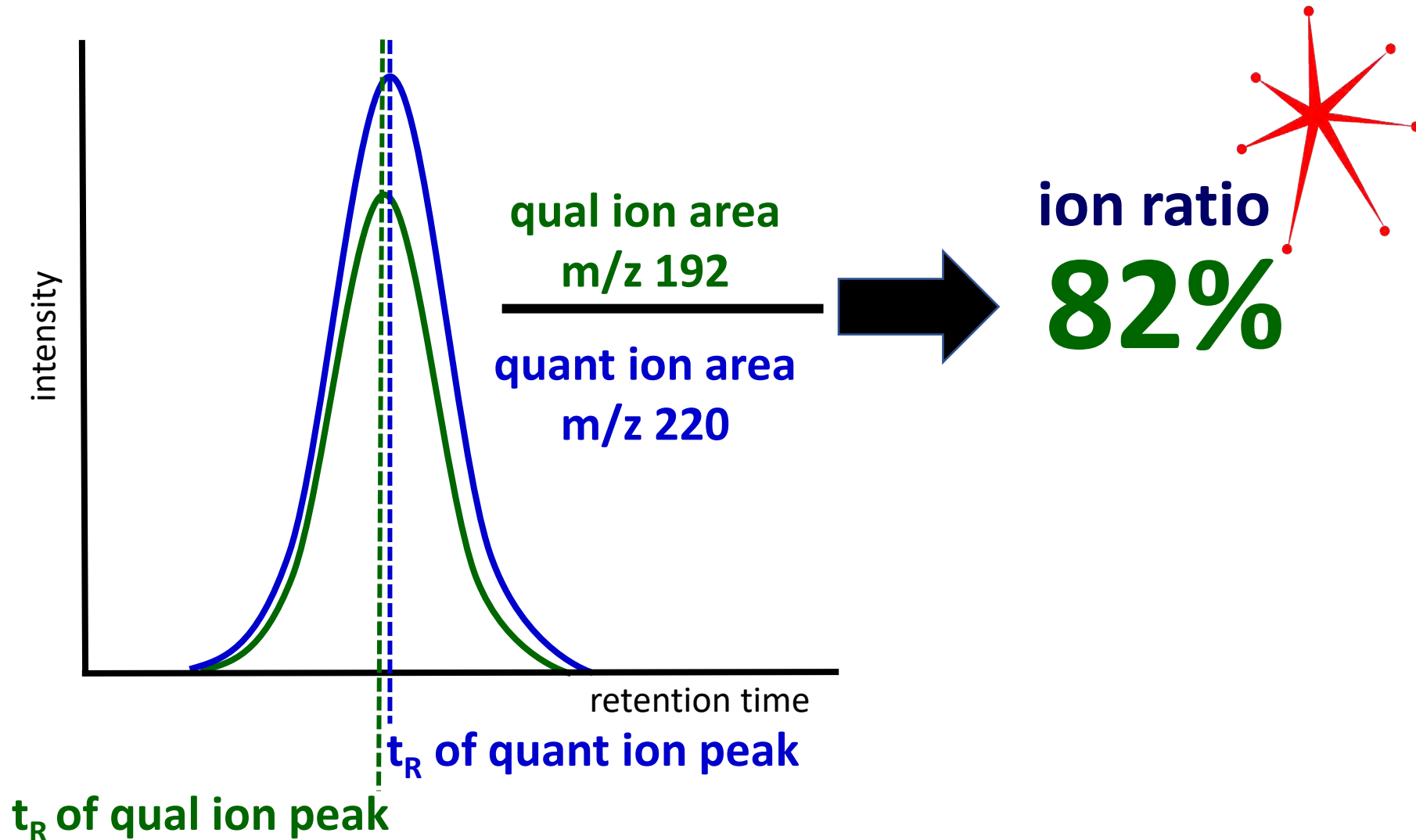
identification data –retention time



identification data- retention time



identification data- ion ratio



identification data- ion ratio

Table 3. Identification requirements for different MS techniques.¹

MS detector/Characteristics		Acquisition	Requirements for identification	
Resolution	Typical systems (examples)		minimum number of ions	other
Unit mass resolution	Single MS quadrupole, ion trap, TOF	full scan, limited m/z range, SIM	3 ions	S/N $\geq 3^d$ Analyte peaks from both product ions in the extracted ion chromatograms must fully overlap.
	MS/MS triple quadrupole, ion trap, Q-trap, Q-TOF, Q-Orbitrap	selected or multiple reaction monitoring (SRM, MRM), mass resolution for precursor-ion isolation equal to or better than unit mass resolution	2 product ions	Ion ratio from sample extracts should be within $\pm 30\%$ (relative) of average of calibration standards from same sequence
Accurate mass measurement	High resolution MS: (Q-)TOF (Q-)Orbitrap FT-ICR-MS sector MS	full scan, limited m/z range, SIM, fragmentation with or without precursor-ion selection, or combinations thereof	2 ions with mass accuracy ≤ 5 ppm ^{a, b, c)}	S/N $\geq 3^d$ Analyte peaks from precursor and/or product ion(s) in the extracted ion chromatograms must fully overlap. Ion ratio: see D12

^{a)} preferably including the molecular ion, (de)protonated molecule or adduct ion

^{b)} including at least one fragment ion

^{c)} < 1 mDa for m/z < 200

^{d)} in case noise is absent, a signal should be present in at least 5 subsequent scans

identification data- ion ratio

target ion ratio

(average from calibration for batch)

80% \pm 30%

identification data- ion ratio

Let's put the
breaks on

identification data- ion ratio

Setting up method or adding compounds

- Ion ratios determined for individual mixes
(without coeluting compounds)
- Check ion ratio for combined mix

identification data- ion ratio

Metalaxyl

Q1/Q3

quant: 234/146

qual: 132/117

ion ratio: 70%



identification data- ion ratio

Metalaxyl

Q1/Q3

quant: 234/146

qual: 132/117

ion ratio: 70%

added analytes

**ion ratio:
300%**

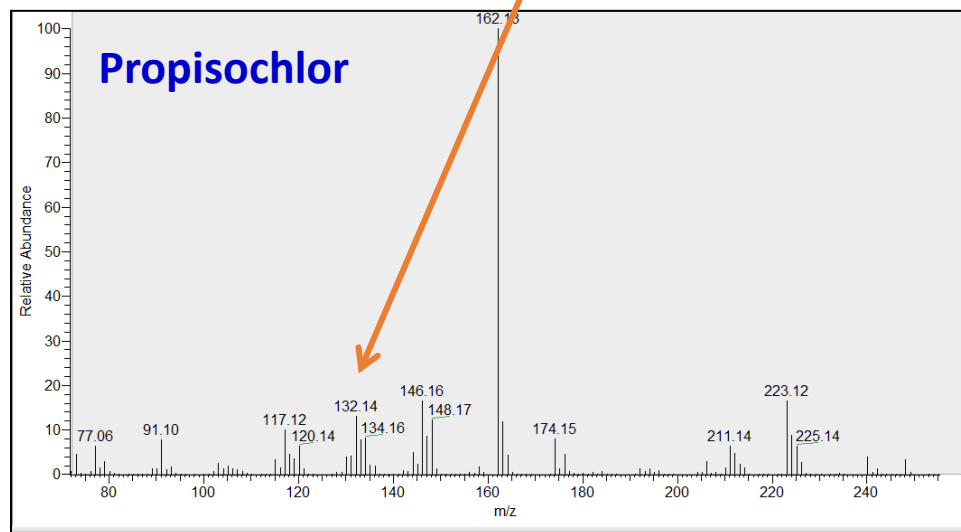


identification data- ion ratio

Metalaxyl
Q1/Q3
quant: 234/146
qual: 132/117
ion ratio: 70%



Propisochlor
0.02 min different t_R
Corruption
of 132/117 MRM



identification data- ion ratio

target ion ratio

(average from calibration for batch)

80% \pm 30%

absolute

Simply add/subtract 30

$$80\% - 30\% = 50\%$$

$$80\% + 30\% = 110\%$$

50 to 110% = “pass”

identification data- ion ratio

target ion ratio

(average from calibration for batch)

80% \pm 30%

absolute

Simply add/subtract 30

$$80\% - 30\% = 50\%$$

$$80\% + 30\% = 110\%$$

50 to 110% = “pass”

relative

$$80\% (0.3) = 24\%$$

$$80\% - 24\% = 56\%$$

$$80\% + 24\% = 104\%$$

56 to 104% = “pass”

identification data- ion ratio

target ion ratio

(average from calibration for batch)

80% + 30%

abs

Simply a

80% -

80% +



%

5%

4%

50 to 110% = “pass”

56 to 104% = “pass”

identification data- ion ratio

target ion ratio

(average from calibration for batch)

20% ± 30%

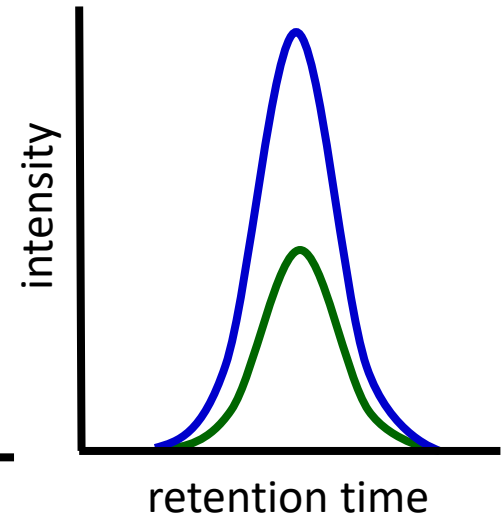
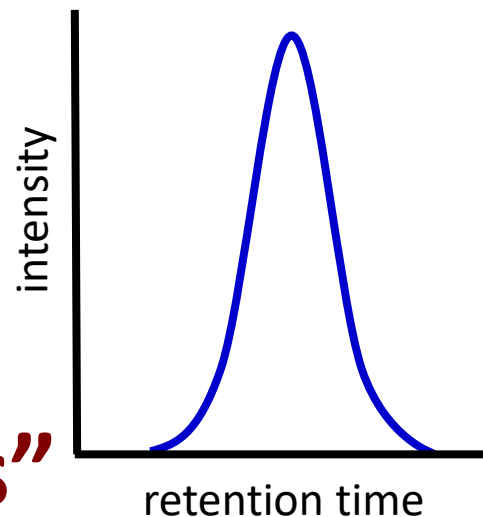
absolute

Simply add/subtract 30

$$20\% - 30\% = -10\%$$

$$20\% + 30\% = 50\%$$

“0” to 50% = “pass”



identification data- ion ratio

target ion ratio

(average from calibration for batch)

20% \pm 30%

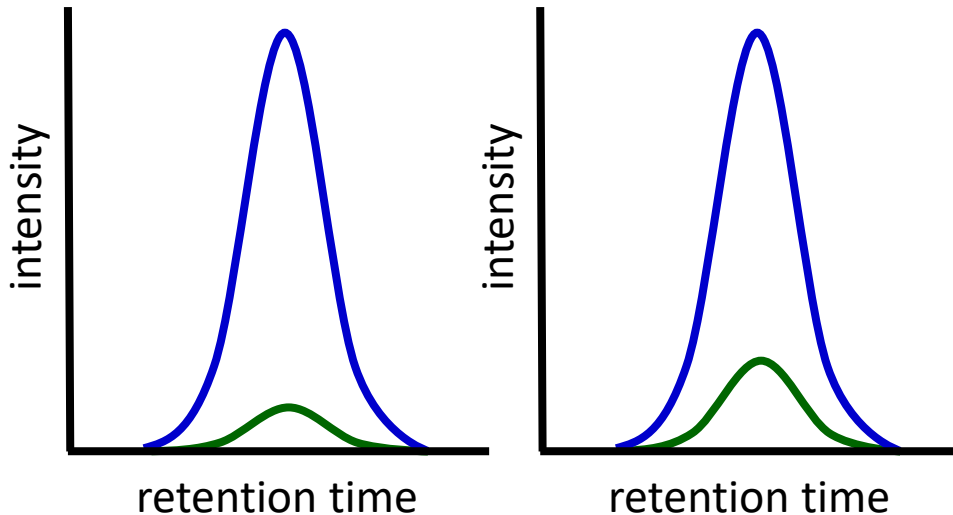
relative

$$20\% (0.3) = 6\%$$

$$20\% - 6\% = 14\%$$

$$20\% + 6\% = 26\%$$

14 to 26% = “pass”



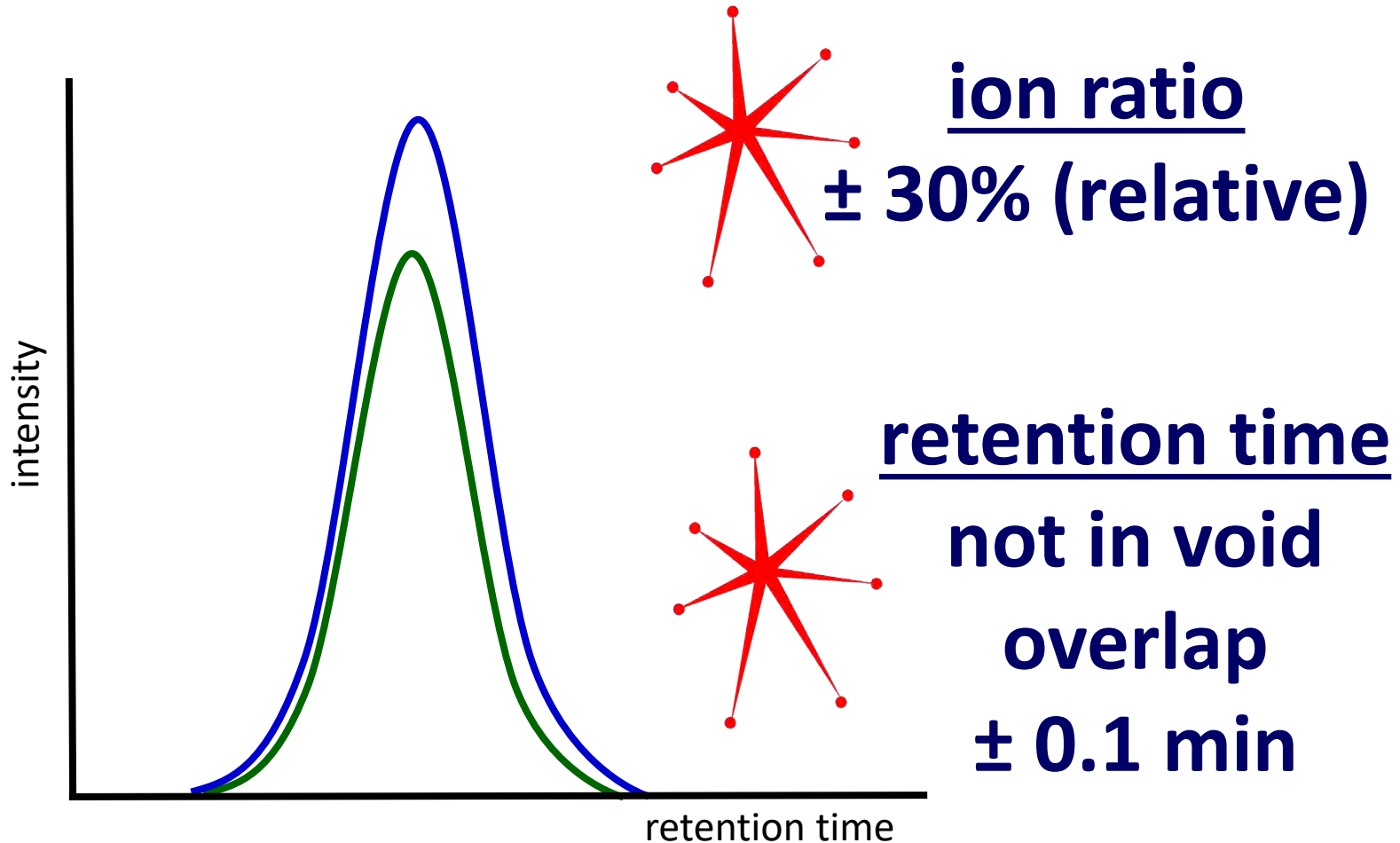
identification data- ion ratio

- ion ratio range = false – and false + control
- too wide (absolute) = false positives
too narrow = false negative


Why?

appear more sensitivity (lower LODs)
easier to say meeting regulatory limits

identification criteria



limit of detection

different ways
disagreement
 fit-for-purpose

Several:

Instrument – solvent

Matrix – matrix without prep

Method – all method steps

limit of detection

DE.3001736111

S/N popular but...

S/N is not really appropriate for estimating detection limits

Modern MS/MS systems have very low noise...

A statistical approach will yield more robust calculations of detection limits



Signal, Noise, and Detection Limits in Mass Spectrometry

Technical Note

Chemical Analysis Group

Authors

Greg Wells, Harry Prest, and
Charles William Russ IV,
Agilent Technologies, Inc.,
2850 Centerville Road
Wilmington, DE 19809-1610
USA

Abstract

In the past, the signal-to-noise of a chromatographic peak determined from a single measurement has served as a convenient figure of merit used to compare the performance of two different MS systems. Design evolution of mass spectrometry instrumentation has resulted in very low noise systems that have made the comparison of performance based upon signal-to-noise increasingly difficult, and in some modes of operation impossible. This is especially true when using ultra-low noise modes such as high resolution mass spectrometry or tandem MS; where there are often no ion



Why use Signal-To-Noise as a Measure of MS Performance When it is Often Meaningless?

Technical Overview

Authors

Greg Wells, Harry Prest, and
Charles William Russ IV,
Agilent Technologies, Inc.,
2850 Centerville Road
Wilmington, DE 19809-1610
USA

Abstract

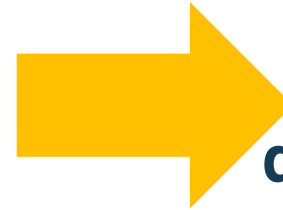
The signal-to-noise of a chromatographic peak determined from a single measurement has served as a convenient figure of merit used to compare the performance of two different MS systems. The evolution in the design of mass spectrometry instrumentation has resulted in very low noise systems that have made the comparison of performance based upon signal-to-noise increasingly difficult, and in some modes of operation impossible. This is especially true when using ultra-low noise modes such as high

limit of detection

$$DL = t_{(n-1, 1-\alpha=0.99)} (S)$$

students' t value

standard deviation
of area counts



calculated
detection limit

1. replicate determinations (n=8)
2. calculate standard deviation of area counts (S)
3. determine t value from table (for n-1=7, t=2.998)
4. convert DL area counts to DL pg

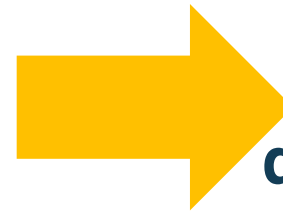
DL as pg = (DL)*(pg on column/average area counts)

limit of detection

$$DL = t_{(n-1, 1-\alpha=0.99)} (S)$$

students' t value

standard deviation
of area counts



calculated
detection limit

Verify 

1. replicate determinations (n=8)
2. calculate standard deviation of area counts (S)
3. determine t value from table (for n-1=7, t=2.998)
4. convert DL area counts to DL pg

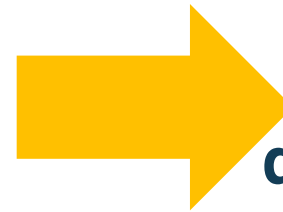
DL as pg = (DL)*(pg on column/average area counts)

limit of detection

$$DL = t_{(n-1, 1-\alpha=0.99)} (S)$$

students' t value

standard deviation
of area counts



calculated
detection limit

1. replicate determinations (n=8)

2. calculate standard deviation of area counts (S)

Must be near DL

concentration

Identification criteria

must be met

Different concentrations
for different analytes?

Quant ion or qual ion?



**What do I need
to know?**

limit of detection

Pesticides: different signal



different levels needed

Lowest level meeting ion ratio criteria

Metalaxyl

	0.1 ppb	1 ppb	10 ppb	100 ppb	1000 ppb
Injection 1	✗	✓	✓	✓	✓
Injection 2	✓	✗	✓	✓	✓
Injection 3	✗	✗	✓	✓	✓
Injection 4	✗	✓	✓	✓	✓
Injection 5	✓	✓	✓	✓	✓
Injection 6	✗	✓	✓	✓	✓
Injection 7	✓	✓	✓	✓	✓
Injection 8	✗	✓	✓	✓	✓

limit of detection

Pesticides: different signal



different levels needed

Lowest level meeting ion ratio criteria

Metalaxyl

	0.1 ppb	1 ppb	10 ppb	100 ppb	1000 ppb
Injection 1	✗	✓	✓	✓	✓
Injection 2	✓	✗	✓	✓	✓
Injection 3	✗	✗	✓	✓	✓
Injection 4	✗	✓	✓	✓	✓
Injection 5	✓	✓	✓	✓	✓
Injection 6	✗	✓	✓	✓	✓
Injection 7	✓	✓	✓	✓	✓
Injection 8	✗	✓	✓	✓	✓

limit of detection

Pesticides: different signal



different levels needed

Lowest level meeting ion ratio criteria

Metalaxyl

	0.1 ppb	1 ppb	10 ppb	100 ppb	1000 ppb
standard deviation of area counts from 10 ppb level					
Injection 1			✓		✓
Injection 2			✓		✓
Injection 3	✗	✗	✓	✓	✓
Injection 4	✗	✓	✓	✓	✓
Injection 5		✓	✓	✓	✓
Injection 6	✗	✓	✓	✓	✓
Injection 7		✓	✓	✓	✓
Injection 8	✗	✓	✓	✓	✓

DL = $t_{(n-1, 1-\alpha=0.99)} (S)$

Option 1 (arrow pointing to 0.1 ppb and 1 ppb columns)

Option 2 (arrow pointing to 100 ppb and 1000 ppb columns)

Verify (green checkmark icon)

Group and (text next to Verify)

- Perform on “large” number of concentration levels
- Set this level to lower reporting limit

recovery



Known 200 ppb



loss of analyte during
all steps of the
method procedure

(calculated)

$$\text{Recovery} = \frac{146}{200} \times 100 = 73\%$$

(known)



Calculated 146 ppb

recovery



80 – 120%

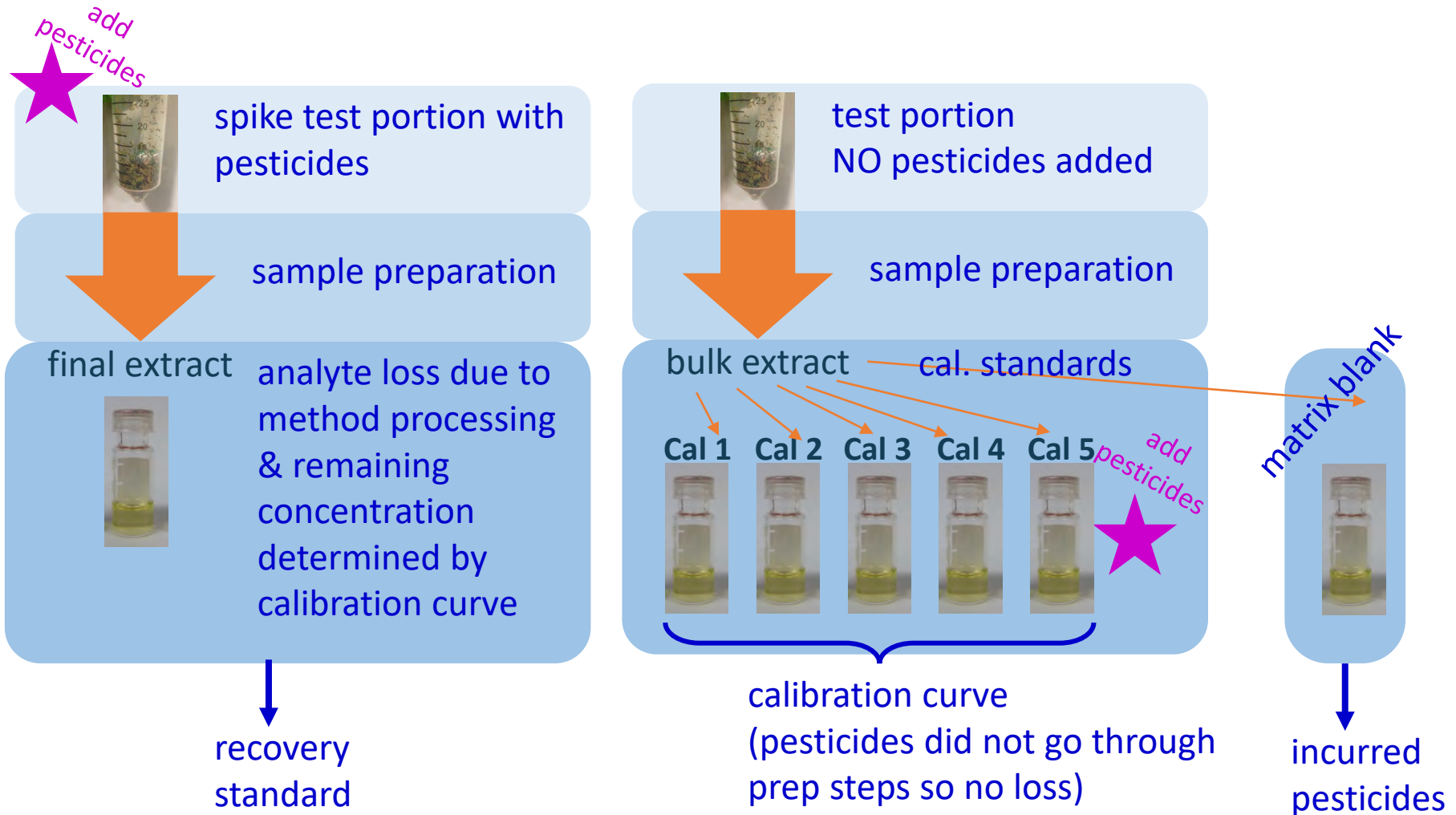
low levels

difficult pesticides

difficult matrices

See SANTE for recommendations on recovery correction
See FDA for concentration dependent ranges

recovery- matrix matching



recovery \neq procedural calibration

Typical Matrix Matched Calibration

pesticides added to aliquoted final extract

- recovery not compensated
- matrix effects are compensated
- can use for recovery calculation

1. test portion



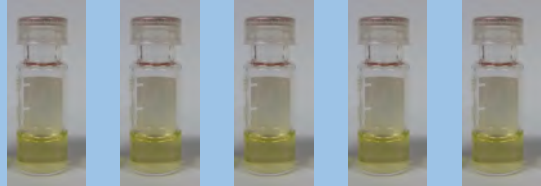
2. preparation



bulk extract

3. cal. standards

Cal 1 Cal 2 Cal 3 Cal 4 Cal 5



add pesticides

4. calibration curve

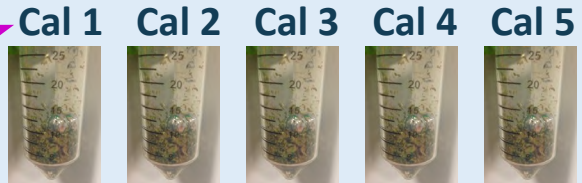
Procedural Calibration

pesticides added to individual test portions

- recovery compensated
- matrix effects are compensated

1. test portions

add pesticides

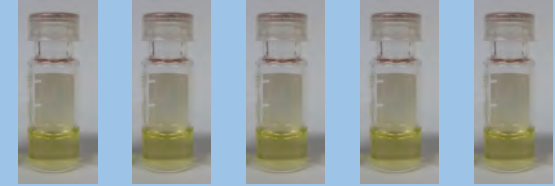


2. preparation



3. cal. standards

Cal 1 Cal 2 Cal 3 Cal 4 Cal 5



4. calibration curve

DE.3001736111

conclusions

Identification

Retention time

- 2x void
- overlap
- ± 0.1 min of standard

Ion Ratio

- $\pm 30\%$ Relative
- no coelutions (setup)

Detection Limits

S/N~~X~~

Lowest Level ID confirmed

- repeat determinations
- statistical determination
- combine for several DL
- verify

Recovery

Recovery definition

- loss due to prep

Calibration

- matrix-matched
- procedural



**FDA and
EU SANTE references**

DE.3001736111



Thank you!

Julie Kowalski

julie@kowalskiscience.com

www.kowalskiscience.com



DE.3001736111