

How to Use TOF and Q-TOF Mass Spectrometers

October 2011

What do TOF and Q-TOF offer?

TOF

- Fast scanning of full spectrum
- High resolution full scan spectra
- Accurate mass measurements

Q-TOF

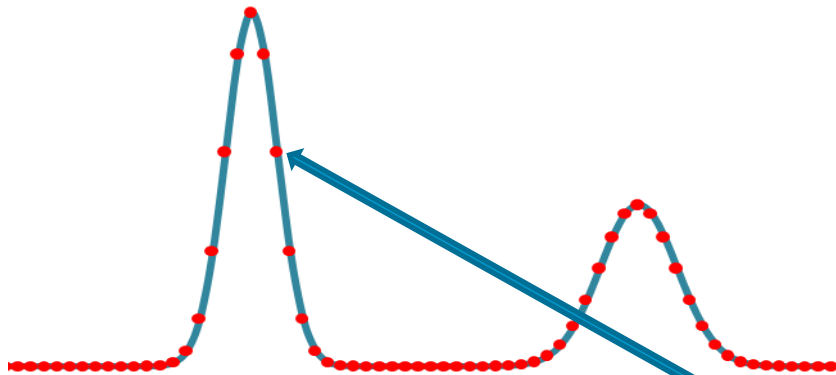
- Fast scanning of full spectrum
- High resolution full scan spectra
- Accurate mass measurements
- MS/MS product ion spectra
- Full scan product ion spectra with high resolution and accurate mass



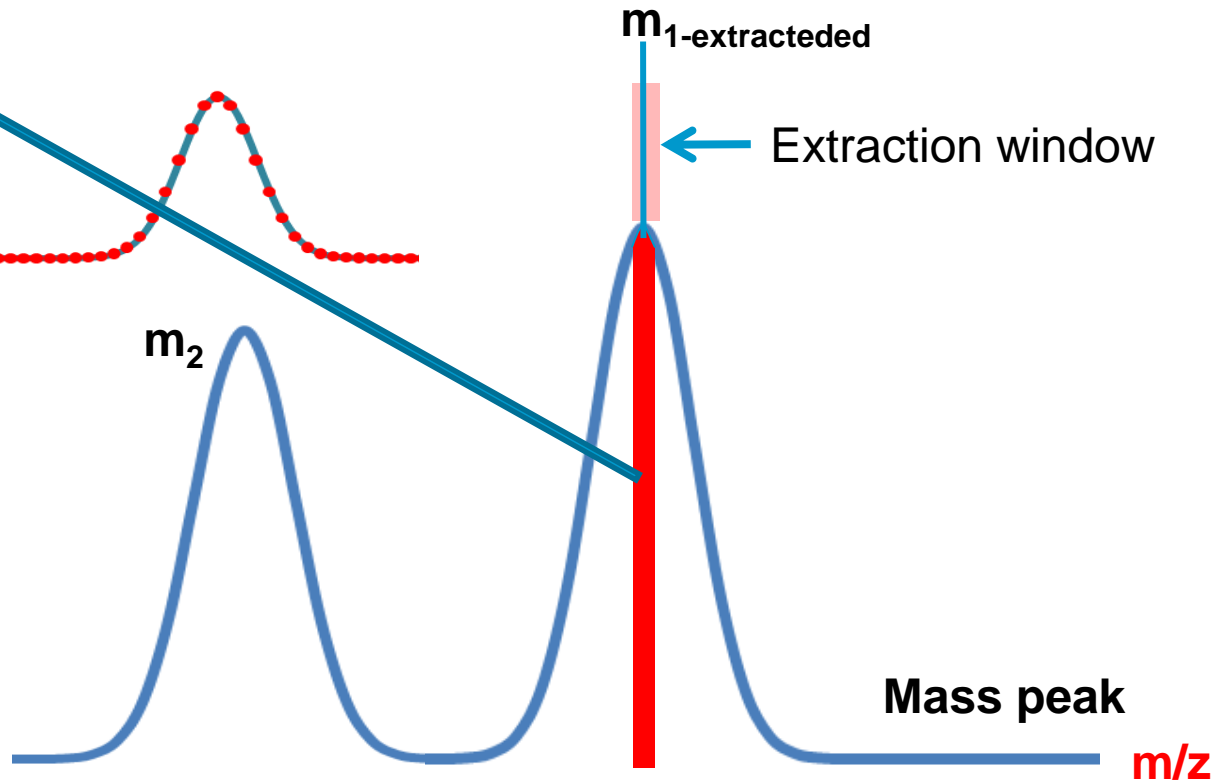
How is accurate mass used?

Extracted ion chromatograms - EIC

Chromatogram – EIC m_1

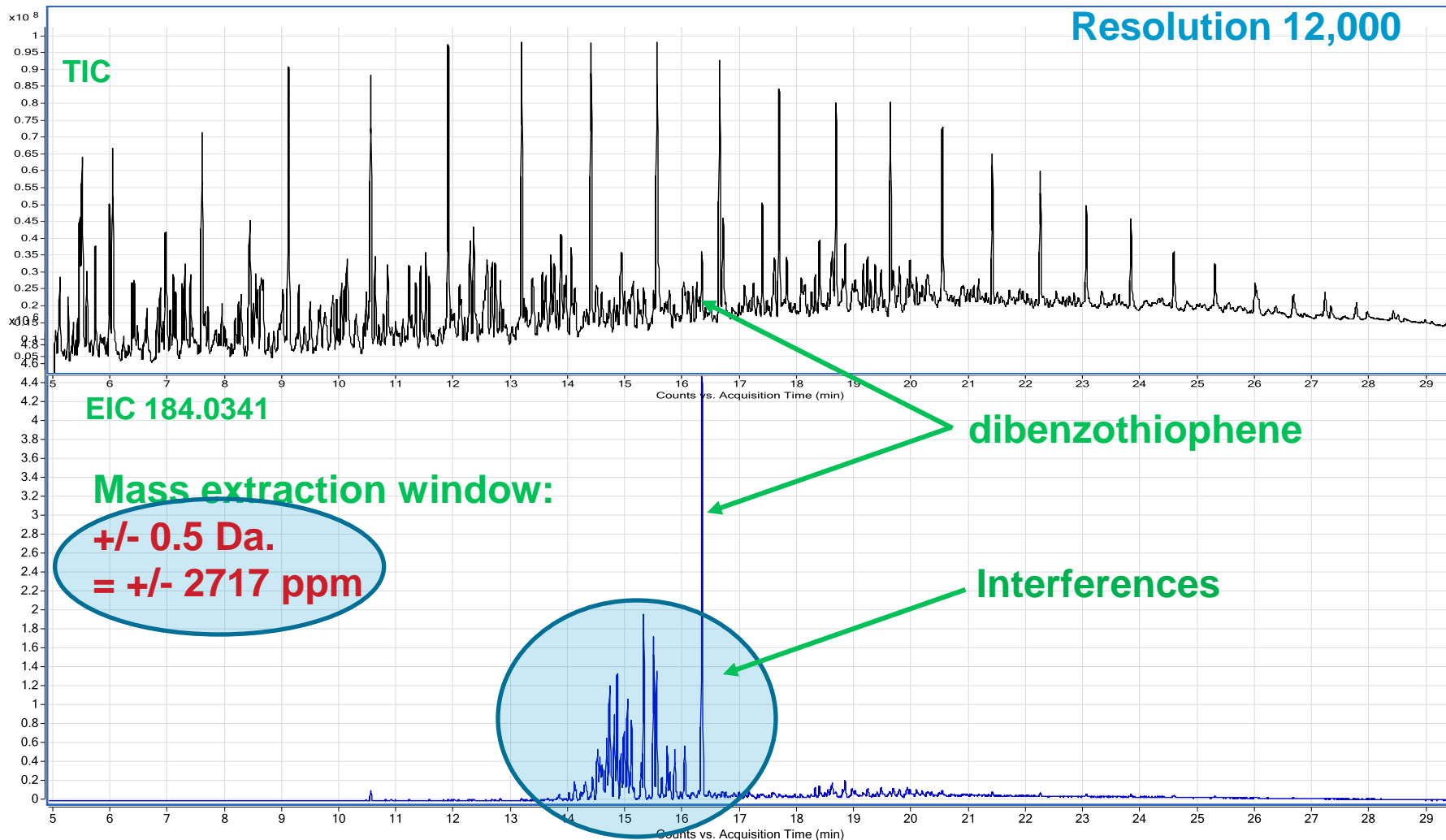


Each point of EIC represents the sum of ions in mass peak within a specified mass range



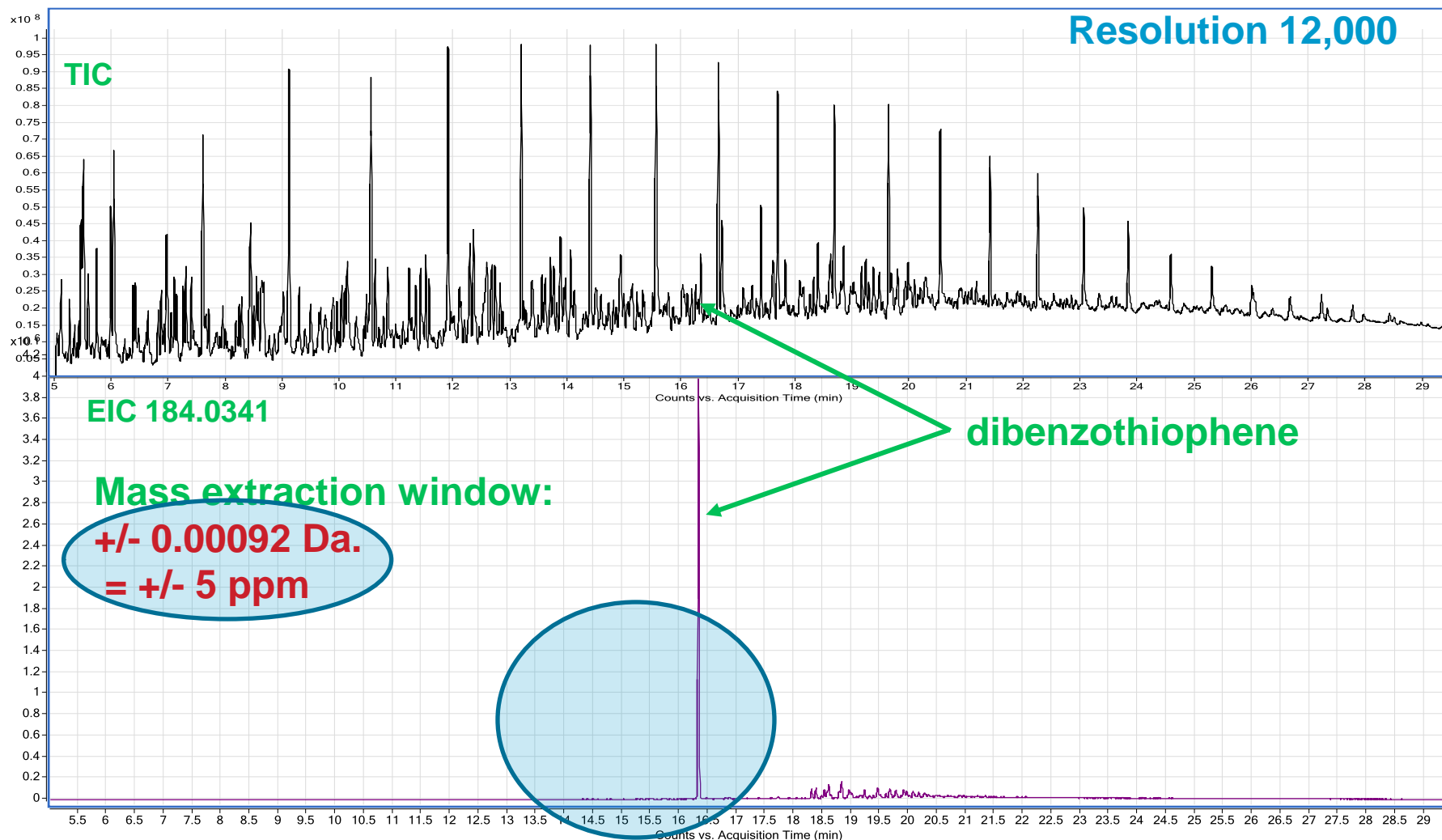
Crude oil target analysis – target analysis

Problem: identify and quantify dibenzothiophene



Crude oil target analysis – target analysis

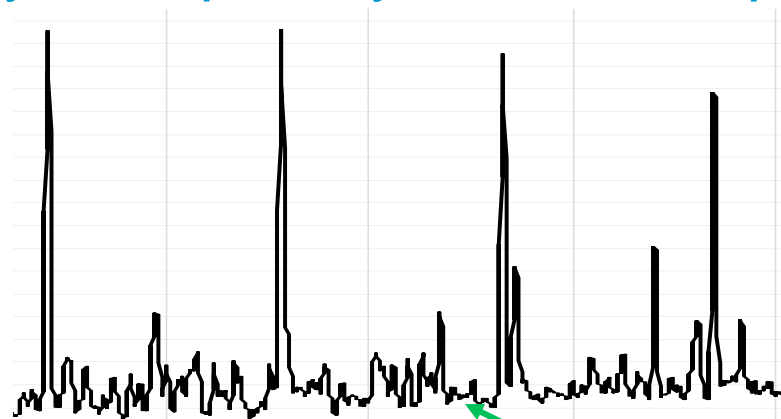
Problem: identify and quantify dibenzothiophene



Crude oil target analysis – target analysis

Problem: identify and quantify dibenzothiophene

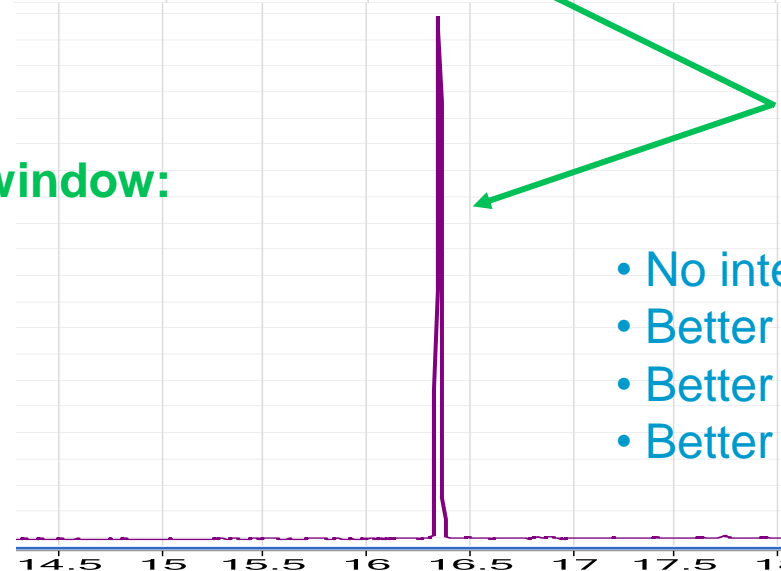
TIC



Resolution 12,000

EIC 184.0341

Mass extraction window:
+/- 5 ppm



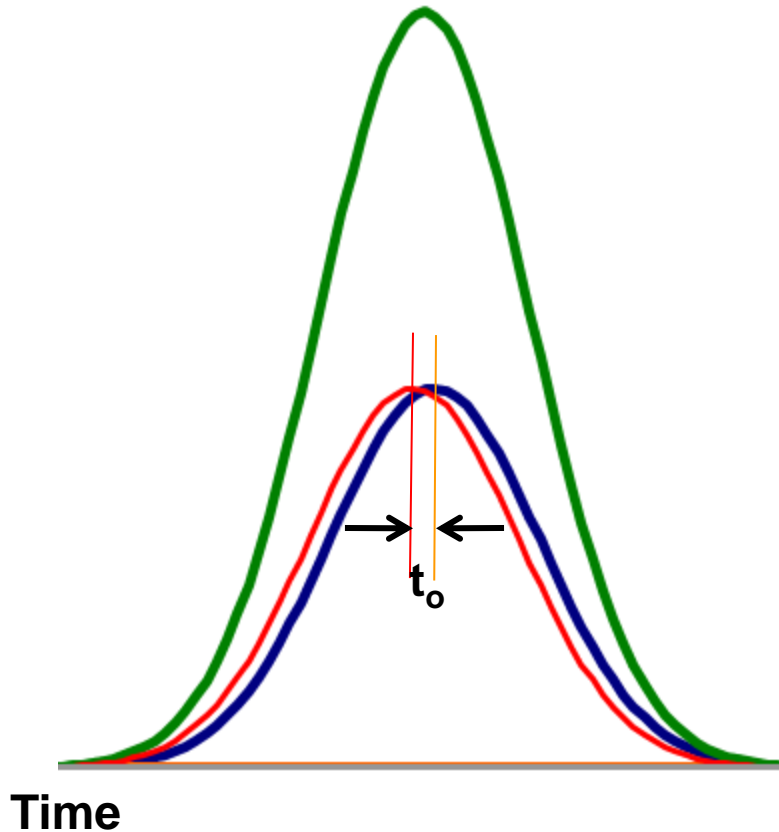
dibenzothiophene

- No interferences
- Better low level peak integration
- Better quantitation
- Better confirmation



Fast scanning of full spectrum

“Speed” enhances deconvolution



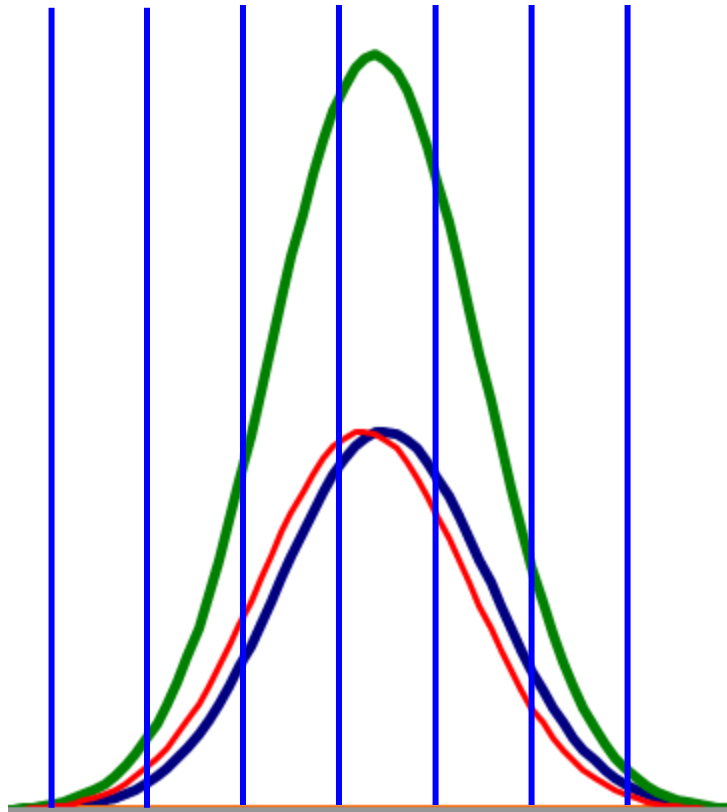
Deconvolution Requires:

- Retention time difference $t_0 > (1/10) W_{1/2}$
- At least one mass that is different for each component
- Peak sampled $< t_0/2$

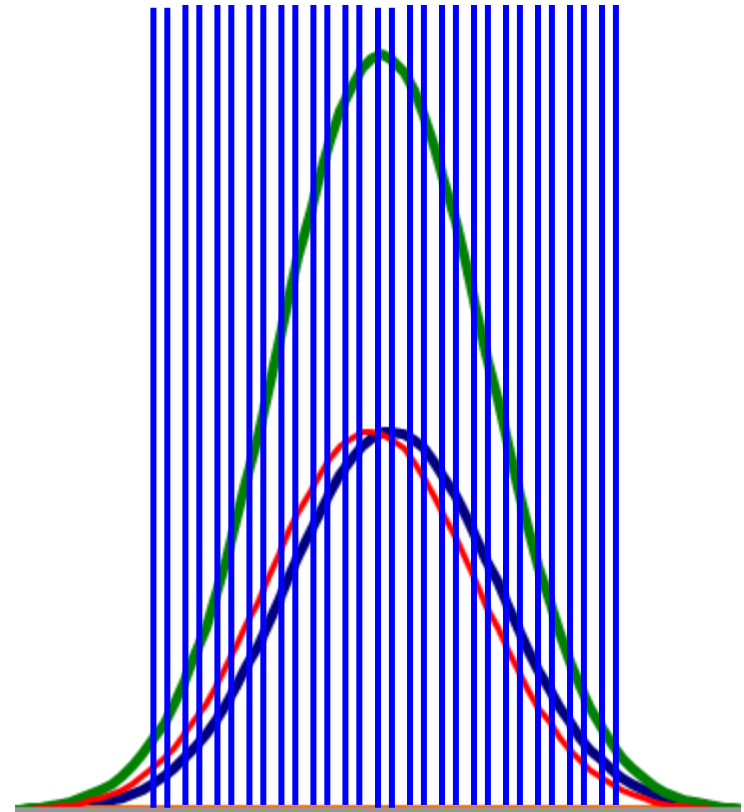
Example:

- $W_{1/2} = 1$ second
- $t_0 > 0.1$ second
- Peak sampled < 0.05 second
= 20 spectra/second

High data rate = better deconvolution



Time Slow data rate will not pick each peak apex



Time Fast data rate will allow deconvolution of closely eluting peaks

High resolution, mass accuracy and scan speed

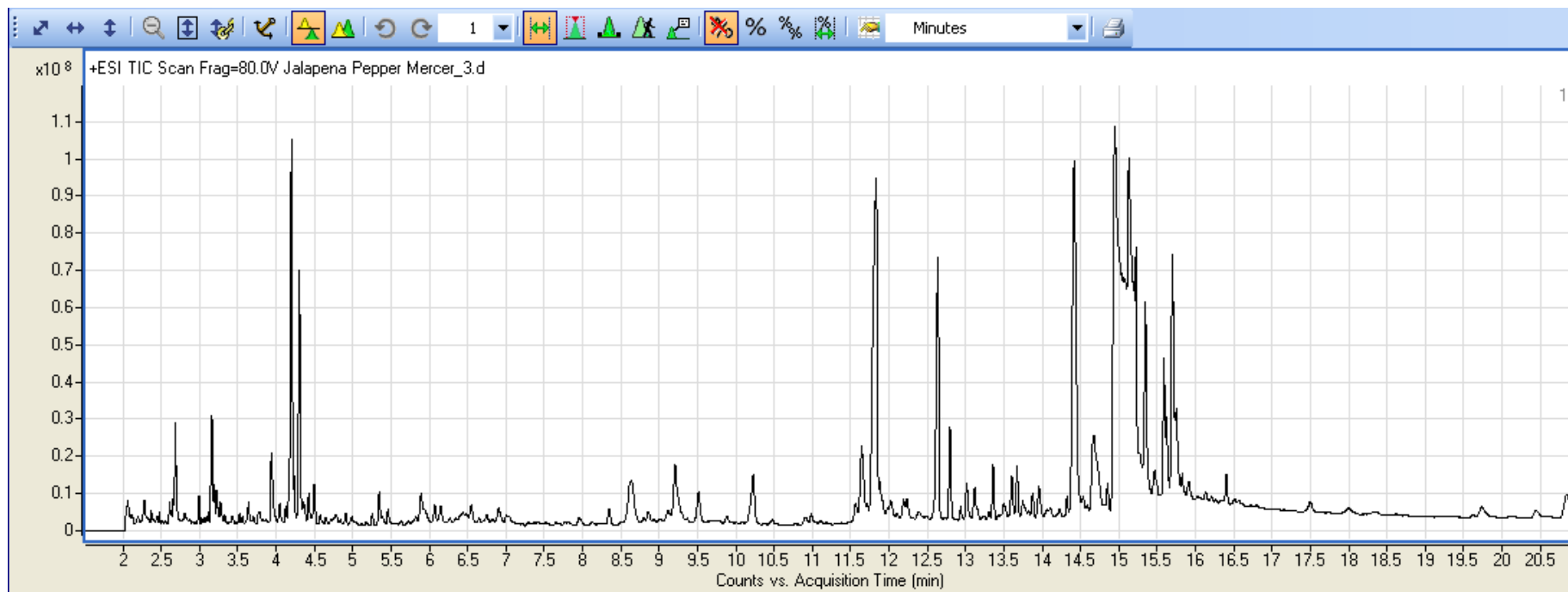
Using fractional masses for deconvolution

TOF and Q-TOF

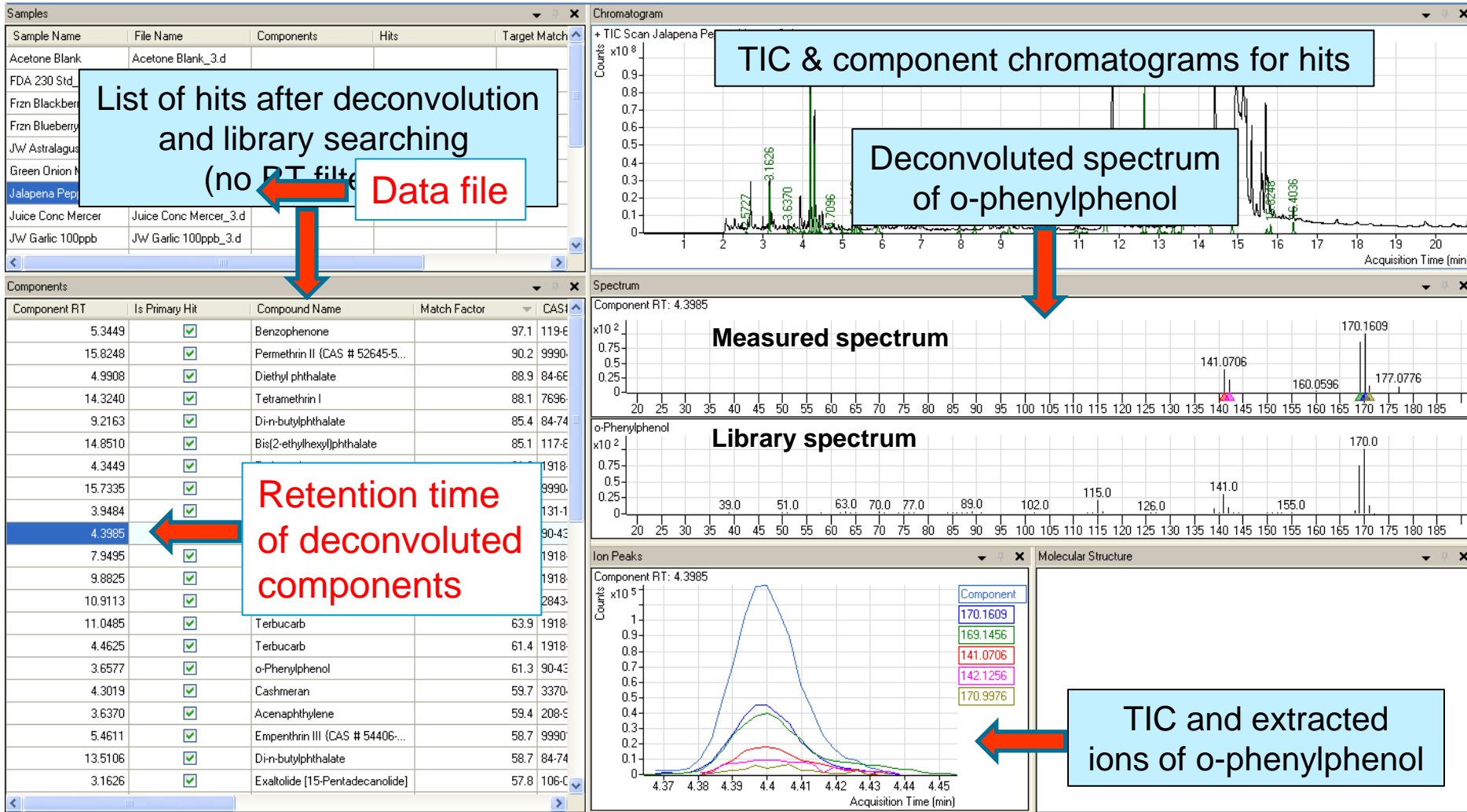
- Fast scan rates >50 spectra/second facilitate deconvolution of peaks with small retention time differences
- Using high mass accuracy spectra with fractional masses rather than integer masses provides:
 - more masses that are distinct between the two co-eluting compounds
 - better deconvolution particularly at low levels
 - better sensitivity

Deconvolution – identify unknown residues

Jalapeno Pepper Extract – GC/TOF Analysis



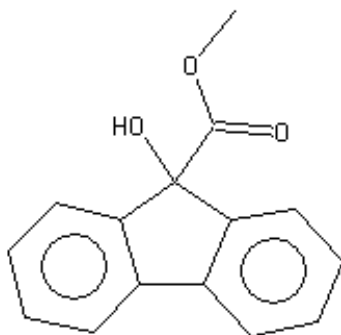
Unknowns analysis – one page for everything



How much resolution is enough?

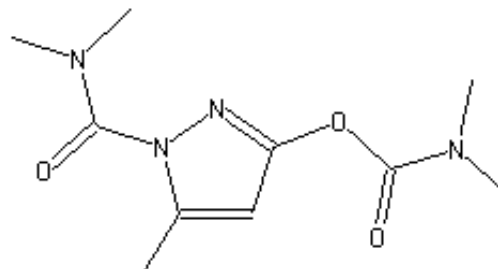
Relative abundance of co-eluting compounds affects the result

Flurenol methyl ester
 $m/z = 240.0781$
Mass error = 1.7 ppm



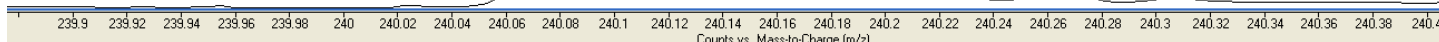
240.0785
240.1218

Dimetilan
 $m/z = 240.1217$
Mass error = 0.4 ppm



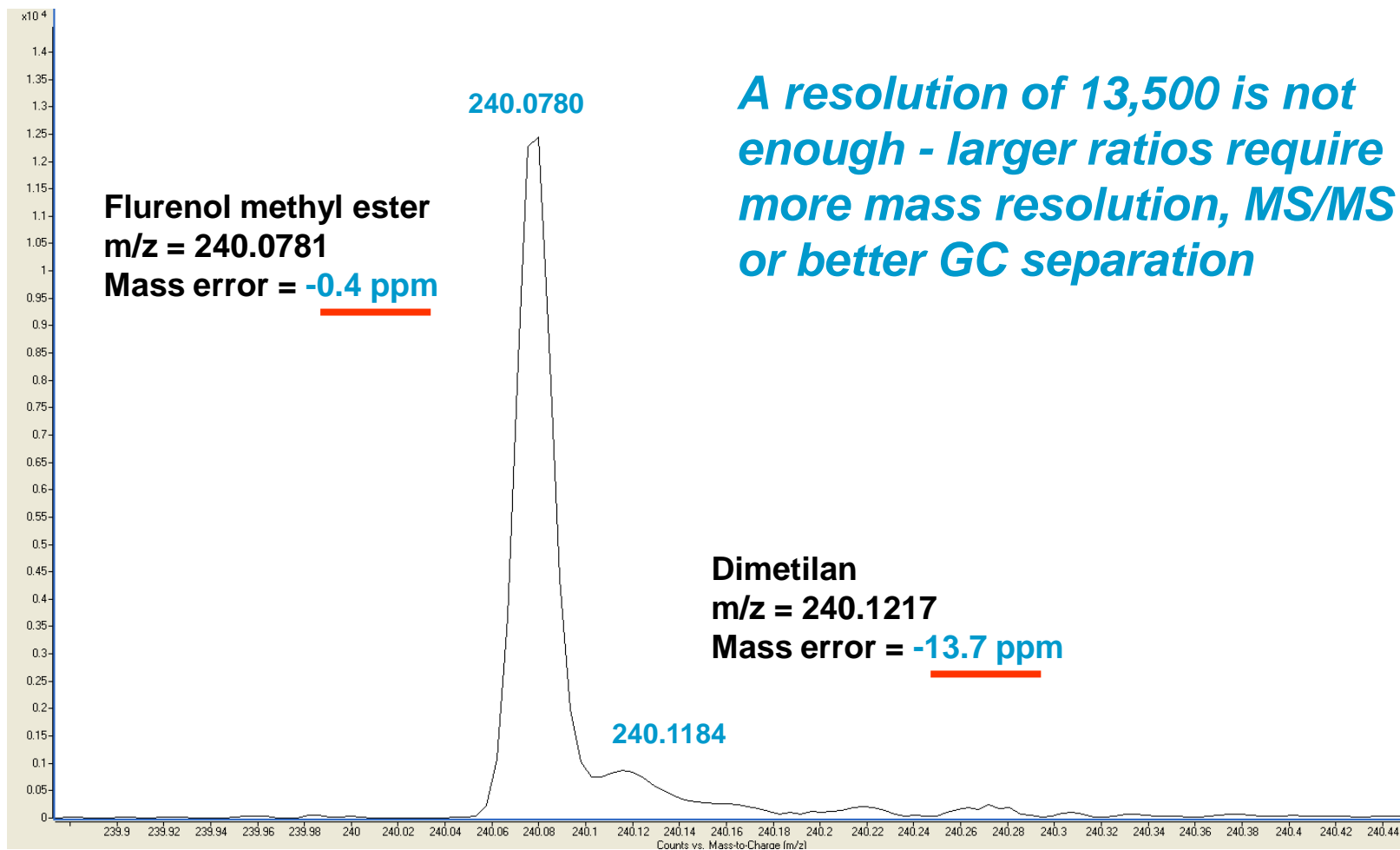
$\Delta m = 0.043$ Da.

~13,500 resolution FWHM



How much resolution is enough?

Abundance Ratio - flurenol methyl ester:Dimetilan = 10:1



A resolution of 13,500 is not enough - larger ratios require more mass resolution, MS/MS or better GC separation

What occurs when the measured mass shifts

- Large shifts in measured mass reduces value for confirmation – necessitates larger extraction windows
- Large shifts in measured mass changes EIC as peak elutes because the ratio of analyte to matrix changes and therefore the mass shift changes



What to do when you run out of resolution?

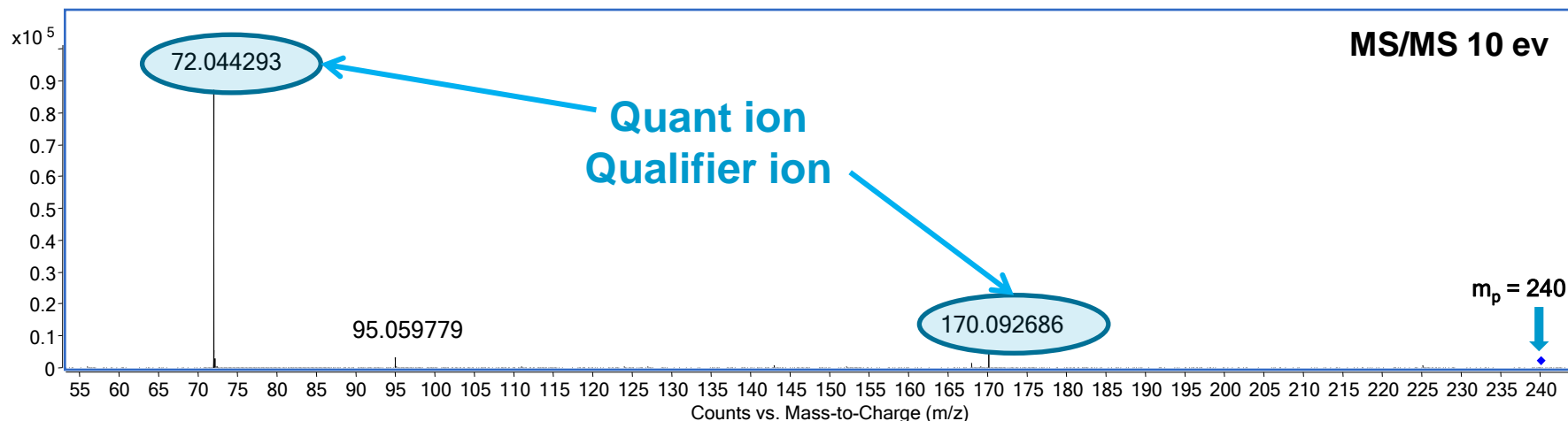
Use MS/MS

- MS/MS product ions are generally well separated in mass
- Therefore - product ion masses do not shift due to lack of resolution
- Why use Q-TOF rather than triple quadrupole?
 - Sometimes qualifier ions have low abundance
 - Accurate mass measurement of quantitation ion offers more confirmatory information than using a **low abundance qualifier ion**

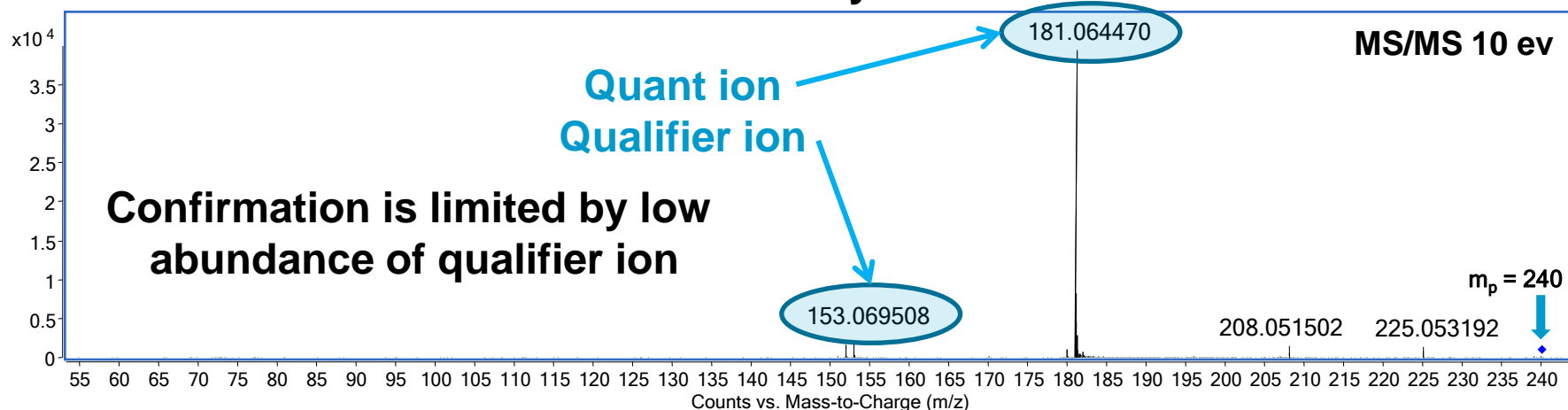
What to do when you run out of resolution?

Use MS/MS

Dimetilan



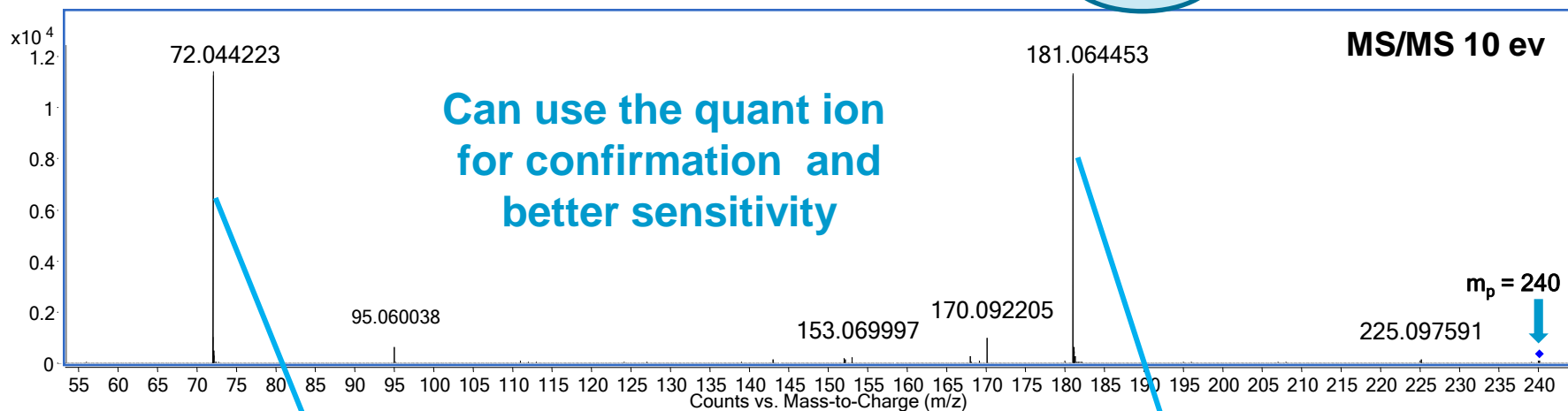
Flurenol methyl ester



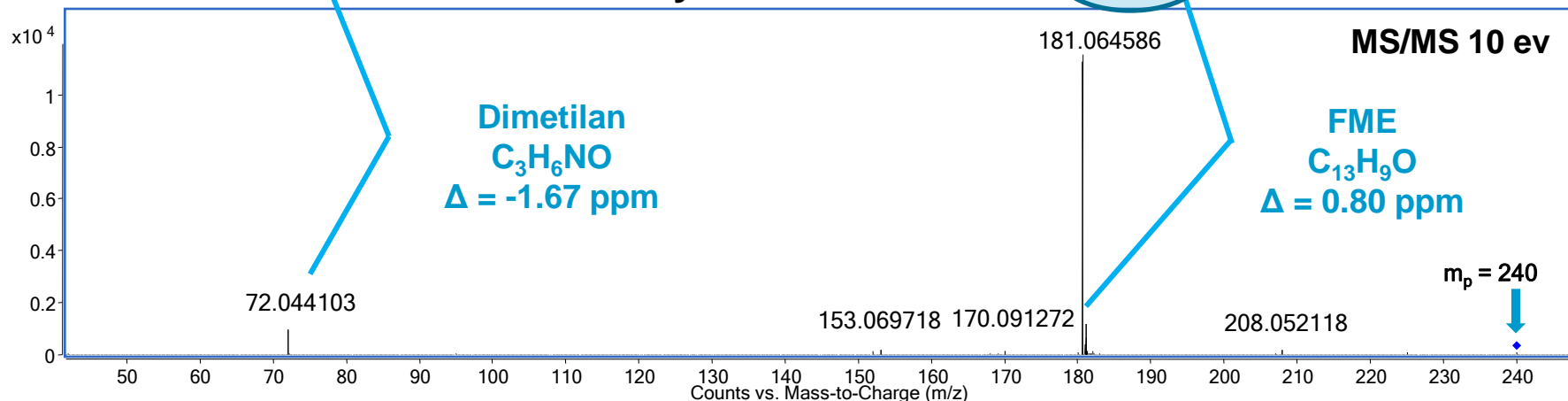
What to do when you run out of resolution?

Use MS/MS

Flurenol methyl ester + Dimetilan 1:1

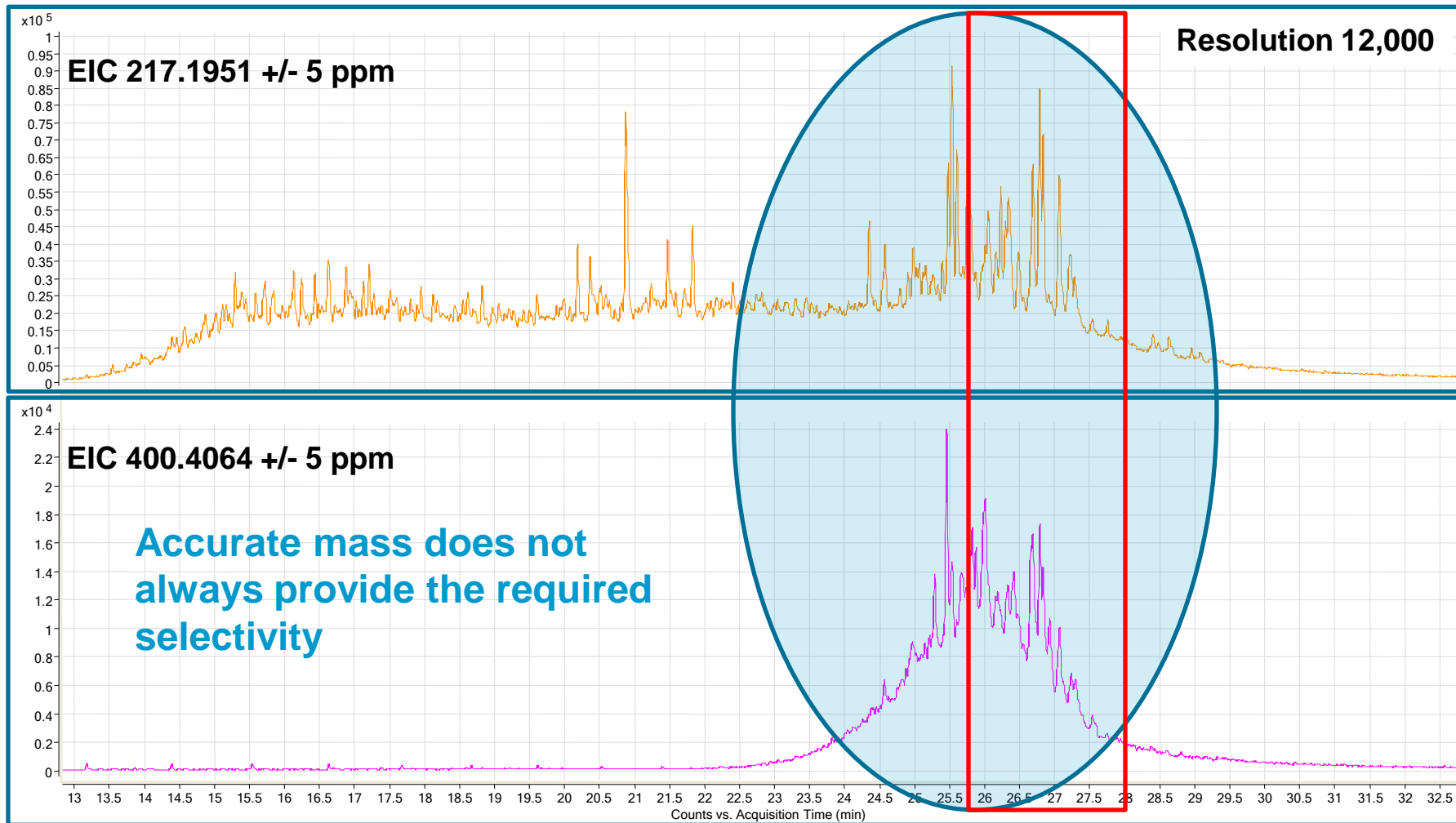


Flurenol methyl ester + Dimetilan 10:1



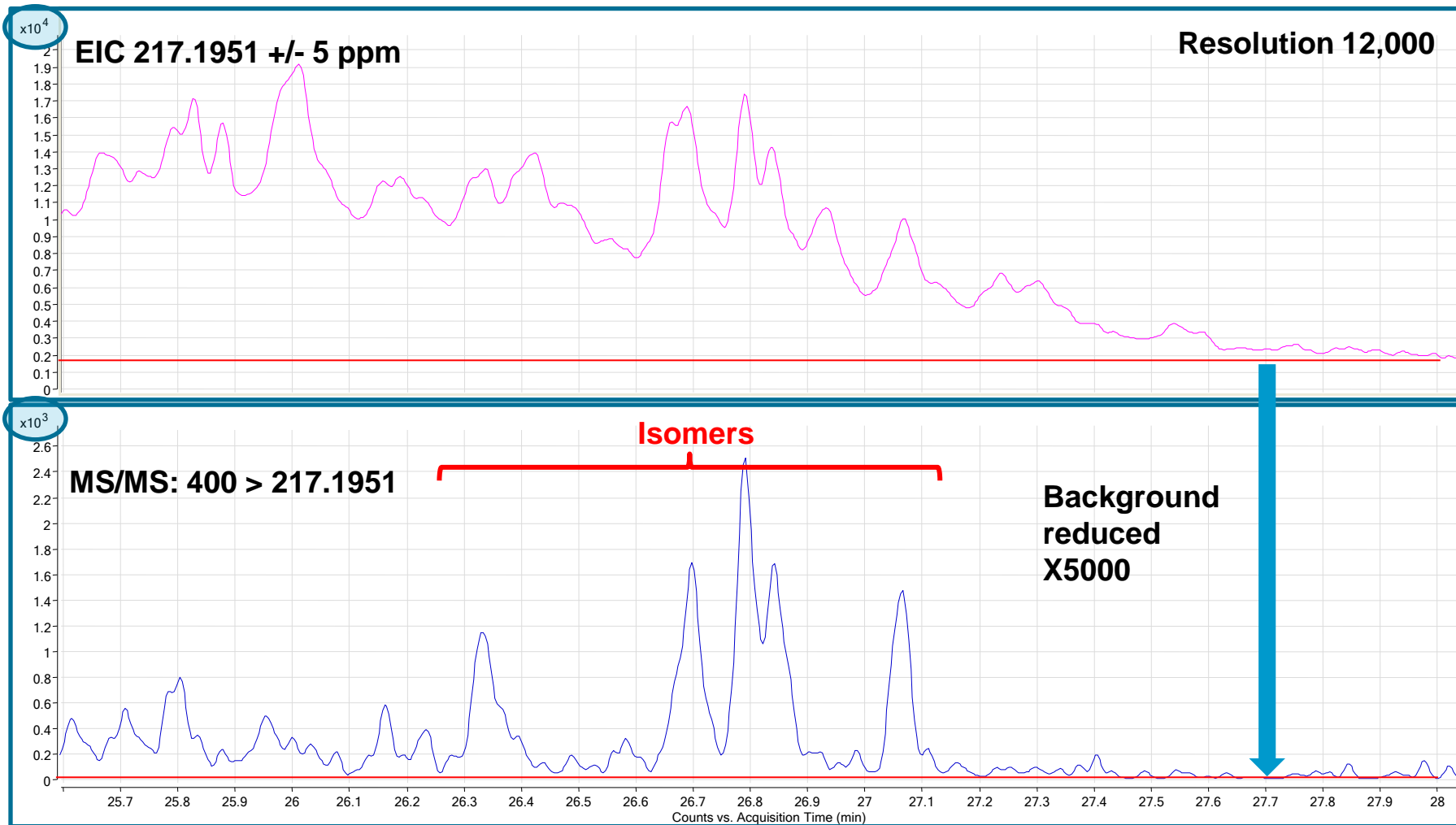
Crude oil: steranes – MS

Problem: identify and quantify isomers



Crude oil: steranes – MS/MS

Problem: identify and quantify isomers



MS/MS accurate mass EIC

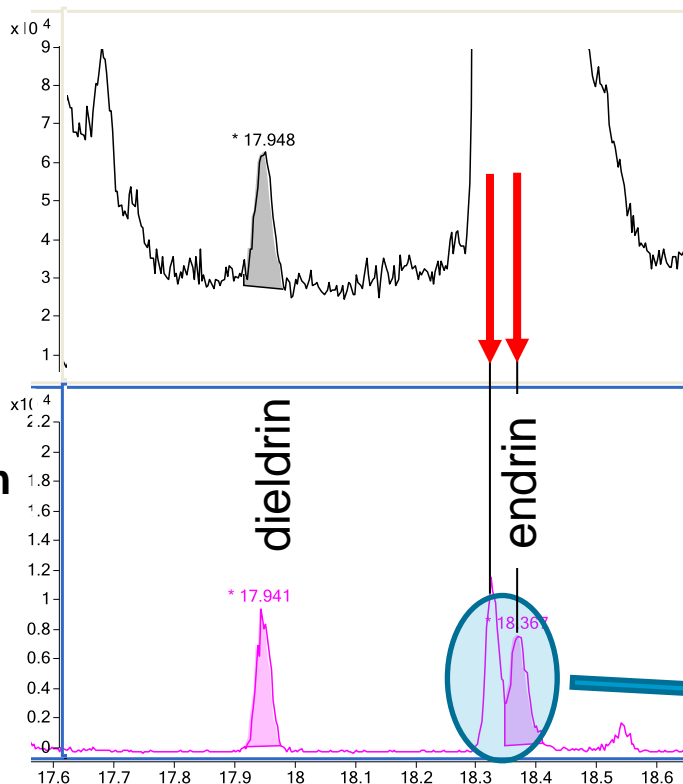
10 pg sample – resolution 13,000

Full scan MS
High resolution/accurate mass

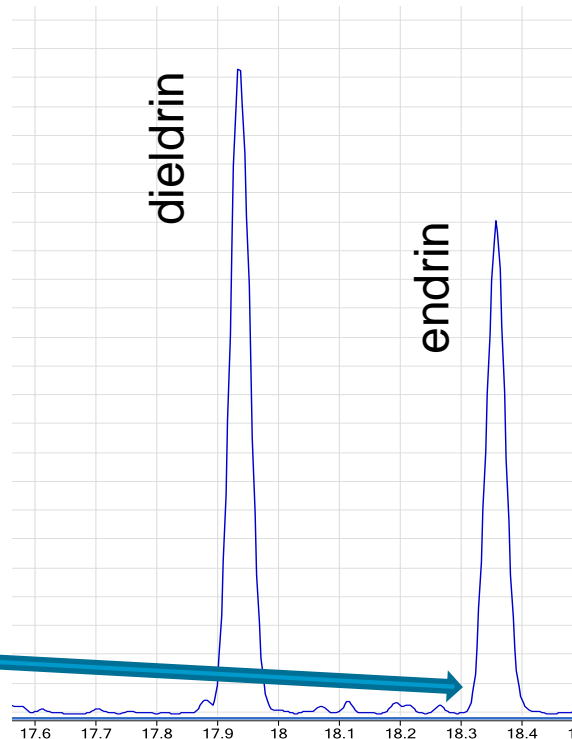
MS/MS 263>192.9150
High resolution/accurate mass

EIC 192.9150 +/- 20ppm

EIC
262.8564 ±0.5 Da.



EIC
262.8564 ±20 ppm



Resolution and accurate mass are insufficient – MS/MS solves the problem

Summary

- The resolution and mass accuracy available from a TOF instrument can solve many analytical problems
- For some analyte and matrix combinations there is not enough mass resolution to confirm and quantify over the required analyte/matrix concentration range
- Q-TOF MS/MS provides constant measured product ion mass values over a larger analyte concentration range than TOF alone when there is not enough resolution

Thank you for your attention