

Accurate Mass and Spectral Accuracy of Agilent LC/MS Single Quadrupole Systems with Cerno MassWorks

Application Note

Author

Don Kuehl, PhD
Cerno Bioscience

Introduction

Single quadrupole LC/MS instruments are the workhorses of analytical labs but fall short for applications that require unknown compound formula identification. On a single quadrupole MS system operating at unit mass resolution, the typical manufacturer specification is for ~0.1–0.5 Da mass accuracy, which is insufficient for confident elemental composition determination (formula ID). However, with the advanced calibration techniques available in Cerno MassWorks software, LC/MS single quad instruments can provide up to 100x improvement in mass accuracy, and perhaps more importantly, provide accurate modeling of the compound isotope profile to 99.0 % spectral accuracy [1]. The enabling combination of the Agilent LC/MS and MassWorks can provide a cost-effective solution for formula ID, for pure compounds. This application note evaluates the performance of an Agilent LC/MS in terms of achievable mass accuracy and spectral accuracy to enable unknown formula ID.



Agilent Technologies

Experimental

The Agilent standard tune solution (G2421-60001) was used as the calibration sample for the performance test. This solution contains a number of calibration ions at mass values from 117 Da to 2,722 Da. The instrument parameters of the Agilent 6120B Single Quadrupole LC/MS were set to acquire data in raw (profile) mode with zero ion threshold. The instrument scan range was set to 100–1,000 Da under the Scan mode with Full Scan Data Storage [2]. The test sample for evaluating the performance was reserpine, a common benchmark standard used in LC/MS, with a molecular weight of 608.68, and a molecular formula of $C_{33}H_{40}N_2O_9$. Ten microliters of both the tune solution and the reserpine were loop-injected into the LC/MS. The concentration of the reserpine was 100 ppm for a total injected amount of 1 μ g, the equivalent amount on the column in LC analysis. The reserpine was injected 10 times to obtain meaningful statistics.

The acquired Agilent MS data of the tune solution were opened directly in the MassWorks software under Calibration mode. The calibration ions were automatically located by simply loading a predefined ion list for the Agilent tune mixture. Figure 1 shows some software screenshots of the calibration process. Upon reviewing the results, the calibration file was saved and applied to all the sample runs producing fully calibrated mass spectra.

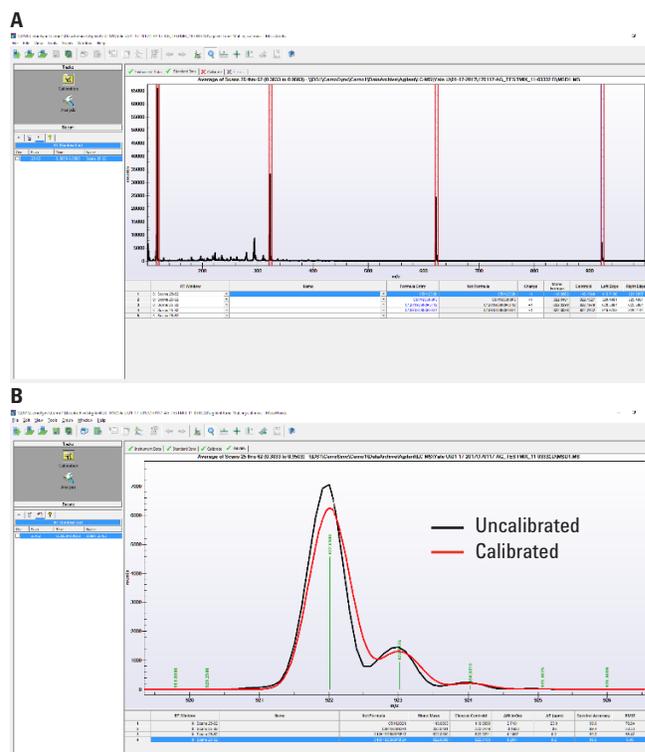


Figure 1. Screen captures show automatic selection of calibration ions (A), and a review of calibrated and uncalibrated mass spectra (B).

Results and Discussion

Figure 2 shows the formula search results under the Analysis mode from the calibrated data for one of the repeat measurements. The search was performed to find all formulae containing C, H, N, and O within a 10 mDa mass tolerance window (that is, the measured mass after MassWorks calibration is accurate to 0.00x Da), producing 38 possible elemental compositions. Since the whole MS profile including all the major and minor isotopes has also been calibrated to a known lineshape, it is a simple matter to calculate the true mass spectrum based on first principles for each possible elemental composition, and compute a corresponding spectral

accuracy by comparing it to the calibrated mass spectrum. The spectral accuracy is a simple mathematical metric that quantitates how well the true and calibrated mass spectrum match. The formula search candidates can then be sorted by spectral accuracy to provide extremely high confidence in the formula ID results. For the example shown in Figure 2, the spectral accuracy is over 99.1 % for the correct reserpine formula, with the next best match falling below 98.7 %. This advanced formula ID algorithm, Calibrated Lineshape Isotope Profile Search (CLIPS), combines both mass accuracy and spectral accuracy to enable formula ID on these otherwise conventional workhorse LC/MS instruments.

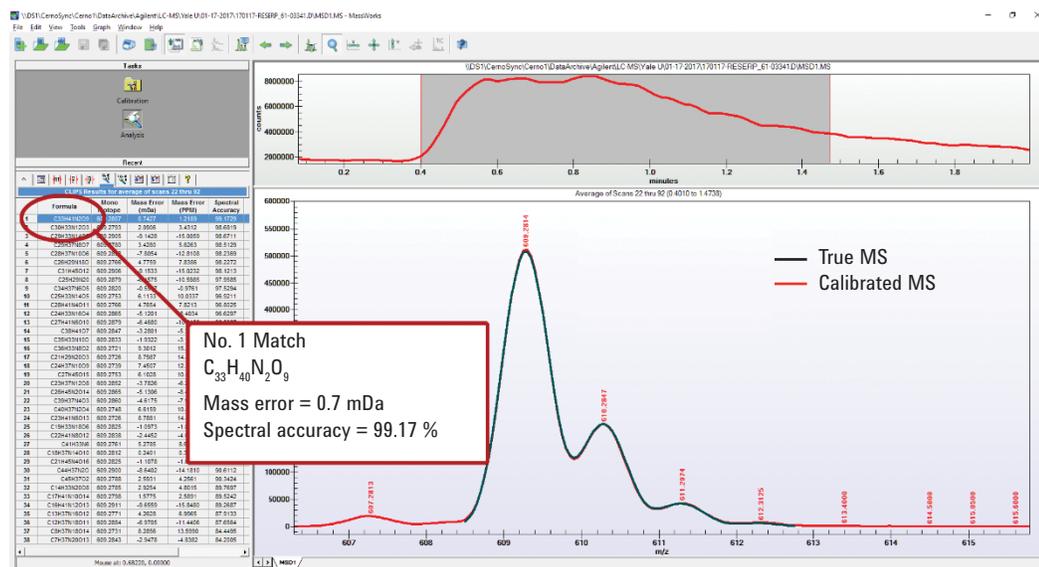


Figure 2. Formula search results for reserpine, showing the no. 1 match, with less than 1 mDa mass error and greater than 99 % spectral accuracy. The plot shows the spectral overlay of the true mass spectrum (black) and the calibrated (red) mass spectrum, and how accurate the spectral fit is.

Table 1 shows the statistical summary of the 10 repeat measurements. The repeatability of the measurements shows that the mass error attainable was within ± 0.005 Da (5 mDa), which is approximately 100x better than the uncalibrated instrument specification. In addition, the calibrated mass spectral lineshape provides an additional, and more powerful, metric for formula ID, with a spectral accuracy of 99.0 %. The well known stability of the Agilent single quadrupole instrument family has been shown to be excellent, and can hold calibration for many days [3], making frequent recalibrations unnecessary and improving the ease of use of MassWorks software for formula ID.

Table 1. Statistical Summary of the Mass Error and Spectral Accuracy from 10 Consecutive Injections of Reserpine

Injection	Accurate mass	Mass error (mDa)	Mass error (ppm)	Spectral accuracy (%)
1	609.2781	-2.6	-4.2	99.06
2	609.2822	1.5	2.5	99.05
3	609.2824	1.7	2.9	99.01
4	609.2840	3.3	5.5	98.71
5	609.2781	0.4	0.7	98.98
6	609.2814	0.7	1.2	99.17
7	609.2802	-0.5	-0.8	98.82
8	609.2800	-0.7	-1.1	99.11
9	609.2781	-1.5	-2.4	98.82
10	609.2822	-0.2	-0.3	99.00
Average	609.2807	0.3	0.4	98.97
Std dev	0.0021	1.7	2.8	0.15

Conclusion

The combination of MassWorks with the Agilent LC/MS single quadrupole is shown to provide mass accuracy of better than 5 mDa and a spectral accuracy of 99.0 % in repeat measurements of reserpine. The novel MS calibration and analysis software can transform the workhorse single quadrupole LC/MS into a powerful and cost-effective tool for formula ID. This enhanced capability can save both time and money by delivering capabilities usually restricted to more expensive and specialized high-resolution instruments to a routine analytical lab.

Acknowledgements

Special thanks to Mr. Terence Wu, Director of the Yale University West Campus Core MS Facility for his help with the experimentation and the use of the facility and the Agilent LC/MS system.

References

1. Y. Wang, M. Gu. "The Concept of spectral accuracy for MS" *Anal. Chem.* **82**, 7055-7062 (2010).
2. For a detailed guide on setting up the instrument parameters for MassWorks, see the Cerno support notes [here](#).
3. J. Mullis, F. Qiu, Y. Wang. "The Robustness of Formula Determination on a Single Quadrupole GC/MS" *ASMS* (2008).

For More Information

These data represent typical results. For more information on our products and services, visit our Web site at www.agilent.com/chem.

www.agilent.com/chem

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc., 2017
 Printed in the USA
 May 18, 2017
 5991-8065EN



Agilent Technologies