TECHNICAL NOTE

Smart MID Mode: increased instrument robustness for complex samples using the DFS Magnetic Sector GC-HRMS

Authors: Heinz Mehlmann, Dirk Krumwiede; Thermo Fisher Scientific, Bremen, Germany

Keywords: Dioxins and Furans, Magnetic Sector, mass accuracy, robustness, Smart MID

Goal

Demonstrate how the new Smart MID feature of the DFS SmartTune Operating Software for Windows 10 enables to reduce the potential negative effects of complex matrices in the analysis of Dioxins and POPs.

Introduction

The analysis of Dioxins and Furans can be challenging due to the matrix remaining in the extract, also after sample clean-up. This can affect the robustness of the analytical process.

To solve this issue, the Thermo Scientific[™] DFS SmartTune[™] Operating Software offers now an optional MID analysis mode for the Thermo Scientific[™] DFS[™] Magnetic Sector GC-HRMS: the Smart MID Mode. This mode was specifically developed for difficult samples.



The large majority of samples can be successfully analyzed with the Default MID Mode, which offers an easy-to-use approach and continues to be the default mode for MID analysis on the DFS Magnetic Sector GC-HRMS.

The Smart MID Mode should be employed for especially challenging samples, which cause analytical problems when using the default MID mode.



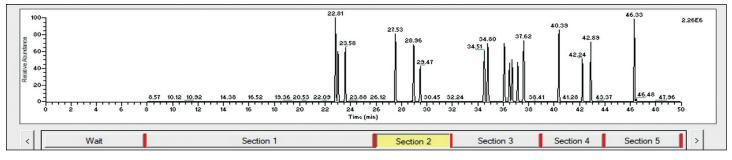
Generally the analysis of Dioxins and Furans and related persistent organic pollutants (POPs) is a targeted quantitative analysis of known contaminants with typically very low levels. For this challenging analysis the mass spectrometric approach of so-called multiple ion detection (MID), sometimes also called selected ion monitoring (SIM), is employed. In this analysis mode the data acquisition covers specifically the mass traces of the analytes only rather than full scan data. The benefit using this approach is that the instrument can concentrate analysis time to detect a higher number of target ions by using higher dwell times. The dwell time is the assigned measuring time for every single target ion according to the MID setup. Due to longer dwell times (measuring times) for the critical target analyte masses better ion statistics are achieved which is beneficial in terms of sensitivity.

For Dioxin and Furan analysis based on the isotope dilution technique, typically the two most abundant isotope masses for each native analyte and ¹³C labeled internal standard are simultaneously monitored. Isomers of the same chlorination degree and therefore with same masses, however, are chromatographically separated in time on the GC.

In this technical note we will see the advancements of the Smart MID Mode, highlighting its potential for mitigating the negative effects of challenging dioxin samples on a series of important factors for routine laboratories: robustness of the system, but also cost and time because the need for repeated analysis of difficult samples can be avoided.

As mentioned above low level target compound analysis such as Dioxin/Furan analysis is performed with the Multi lon Detection (MID) Mode. The analysis is generally divided into multiple time sections each containing different specific target masses.

In the section dialog all specific target masses for each MID time section are defined plus two reference gas masses, the Lock- and Cali-mass used for the constant automatic mass recalibration of the mass spectrometer. This approach ensures maximum mass accuracy and stability for the high resolution analysis of low level contaminants.





egment Settings Times Start [min]: 29.00 + End [min]: 37.50 + Measure [min]: 8.50 + Cycle [sec]: 1.20 +							
Section 2	L/C	Mass	Gr.	Int.	Time[ms]	Compound	Comment A
1	Lock	313.98336	1	20	11	FC43	
2		339.85915	1	1	223	pcdf	QM native pcdf
3		341.85620	1	1	223	pcdf	RM native pcdf
4		351.89941	1	5	44	PCDF	QM[13]C-PCDF
5		353.85702	1	1	223	pcdd	QM native pcdd
6		353.89646	1	5	44	PCDF	RM[13]C-PCDF
7		355.85407	1	1	223	pcdd	RM native pcdd
8	Cali	363.98017	1	20	11	FC43	
9		367.89433	1	5	44	PCDD	QM[13]C-PCDD
10		369.89138	1	5	44	PCDD	RM[13]C-PCDD

Figure 2. Example of the penta-chlorinated dibenzo dioxin/furans mass table (MID section 2).

Within each MID analysis time section the traces of target masses and reference masses are acquired. During the acquisition a permanent automatic mass re-calibration is carried out. This automatic mass calibration process includes the following steps: first step is the first so-called locking of the reference mass signal specified as lock mass, secondly, an electric calibration using a second reference mass peak, the so-called cali mass, is carried out, and in the third and final step the successive monitoring of the target masses takes place. While step 2 and 3 describe a so-called MID mass scanning cycle which is constantly repeated, the locking process is executed only once at the start of each of the MID analysis time sections.

During the lock process at the start of each MID time section the system identifies the reference mass labeled as lock mass using a small mass scan over a defined narrow mass range, called lock window. It corrects its position to the exact theoretical value which is the basis for the following calibration.

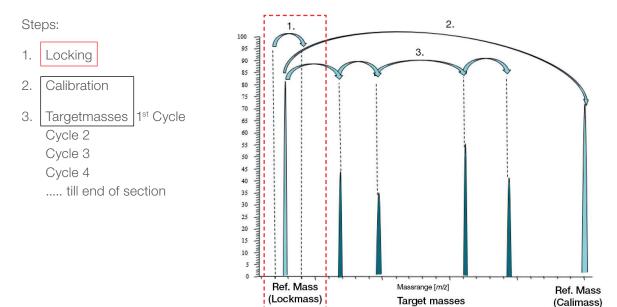


Figure 3. Steps carried out in each MID analysis section.

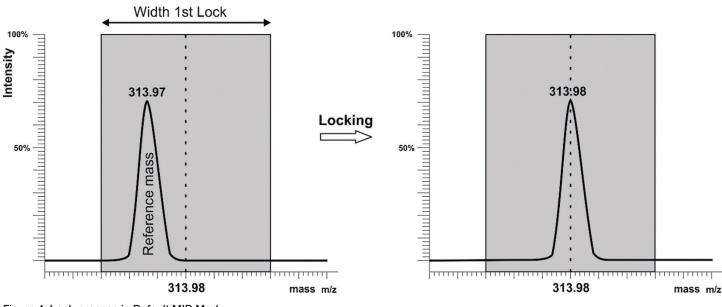


Figure 4. Lock process in Default MID Mode.

In the Default MID Mode the locking is done during the analysis of the sample at the start of each MID time section. At that point, the sample with its matrix is already present and the temperature in the GC is high so that matrix contaminants are eluting from the GC column into the ion source of the mass spectrometer together with the target analytes of interest.

With the presence of high matrix during this locking process, there is a risk in the Default MID Mode that a

mass signal from a matrix compound is close to the lock mass peak, which can disturb the locking process. An intensive matrix peak within the mass scanning range of the lock window can be misinterpreted as lock mass, which would result in a wrong locking (Figure 6). As a consequence, the complete mass calibration for the specific MID time section would be false and the MID section would fail, which means chromatograhic peaks of the target compounds are not or not correctly detected.

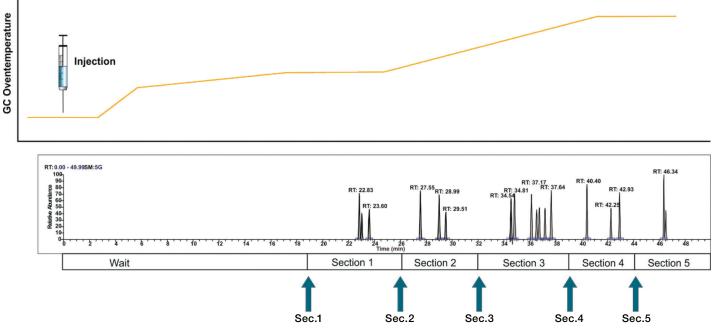


Figure 5. Default MID Mode: The locking for each section is done during the analysis when sample matrix is present, and the GC is hot.

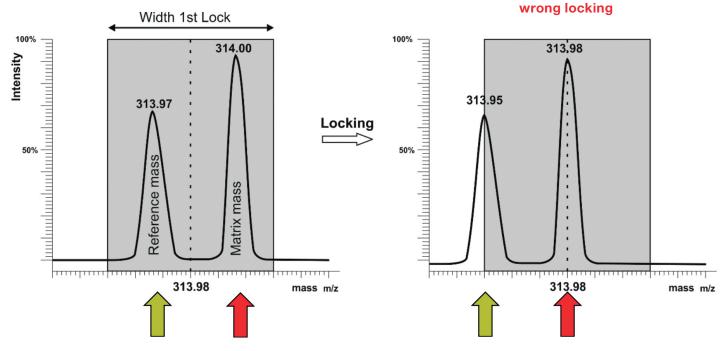


Figure 6. Matrix interference can disturb the lock process which leads to a wrong locking.

In the **Smart MID Mode**, an MID calibration table is created prior to the injection of the sample extract when no matrix is present, and the GC temperature is still low. During the analysis, this calibration table is applied.

That allows minimizing the scanning width of the lock window, as the mass position of the lock mass is already very close to its theoretical value by using the pre-calibration parameters (Figure 7). Matrix peaks cannot affect the lock process any more as they will mostly be outside of the narrow scanning mass range of the locking window. In the unlikely case where interference matrix peaks still fall within the narrow Smart MID lock window they practically overlap with the lock mass peak and thus show the same mass assignment.

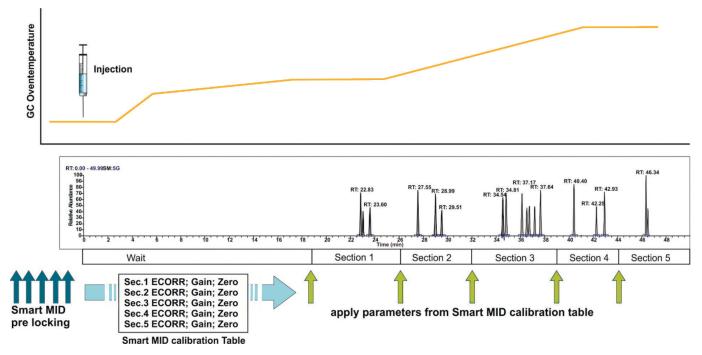


Figure 7. Principle of the Smart MID locking and calibration.

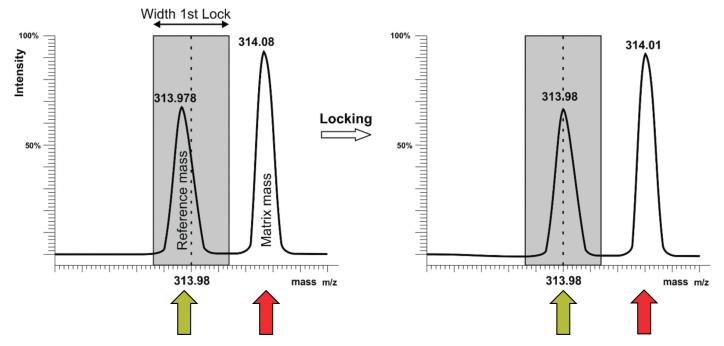


Figure 8. In the Smart MID Mode, the Lock window is much narrower so that no matrix peak can interfere.

Demonstration of the principle of the Smart MID Mode

A Matrix Simulation Experiment based on column bleed can demonstrate the advantages of the Smart MID Mode by setting up an MID analysis experiment with a reference (lock) mass close to one of the GC column bleed peaks, which is used to simulate the matrix peak. At a sufficiently high GC temperature, the column bleed peak will be more intensive than the reference peak and will accordingly interfere with a correct locking when using a standard locking window as in the Default MID Analysis Mode. Using the Smart MID mode in contrast and thus applying the pre-locking (see Figure 7) or pre-calibration, which in this matrix simulation experiment is executed before the GC is heated up. The reference (lock) mass peak is already almost in the center of the lock window. Therefore in the following MID analysis run when the GC is heated up the mass range scanning width of the lock window can be set almost as narrow as the peak base of the reference (lock) mass. The column bleed peak is still present, but far outside of this window and can not interfere with the locking process.

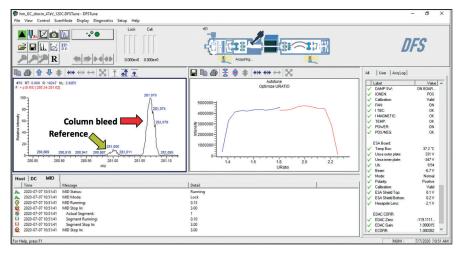


Figure 9. The column bleed peak (right) in the lock window is more intensive than the reference peak (left).

The system misinterprets the column bleed peak as lock mass. As a consequence the whole following MID analysis section fails due to a failed mass calibration.

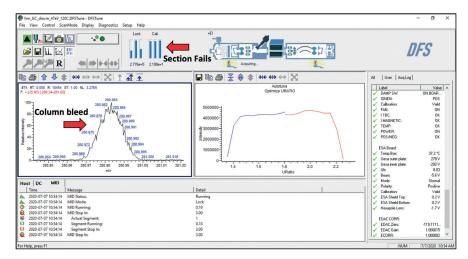


Figure 10. The system locks on the more intensive column bleed peak and not on the reference peak.

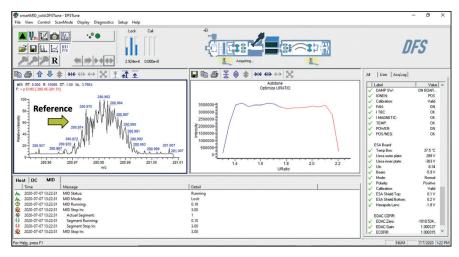


Figure 11. The system locks correctly on the reference mass, due to the pre-determined parameters.

Now, using the Smart MID mode, the system locks correctly and the MID section is acquiring analysis data correctly.

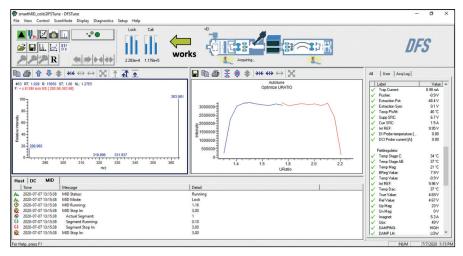


Figure 12. By using the Smart MID Mode, the MID section works correctly, and it is not affected by the column bleed peak.

Conclusion

This technical note illustrates that with the optional Smart MID Mode of the Thermo Scientific DFS SmartTune Operating Software the potential negative effects of complex matrices can be overcome for their analysis with the DFS Magnetic Sector GC-HRMS.

By determining all MID mass calibration parameters before the injection of the sample extract (when no sample matrix is present in the instrument), a very narrow lock mass scan width can be applied in the following sample analysis run, excluding potential matrix interference signals. The Smart MID Mode is offered as an option for difficult samples alongside the Default MID Analysis Mode, which continues to be the default mode, due to it's ease-of-use operation.

No matter the analysis, the DFS Magnetic Sector GC-HRMS performs compliant Dioxins and POPs Analysis, supported by constantly improved software features that are aimed at productivity and ease-of-use of your lab.

Find out more at thermofisher.com/MagneticSector

© 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representatives for details. **TN30716-EN 1020C**

