

Poster Reprint

**ASMS 2025**  
**Poster number TP 315**

# Extractable Leachable Analysis Workflow using MassHunter Explorer 2.0 with NIST and Sirius MS/MS Searching for Enhanced Compound Identification

David A. Weil\*, Sierra D. Durham, James Pyke and Matt  
Curtis

Agilent Technologies, Santa Clara, CA 95051

## Introduction

### New Software for Non-targeted HRMS Extractable and Leachable Data Analysis

One of the biggest challenges of doing extractable leachable (E&L) analysis is converting non-targeted high-mass resolution data into the confident identification of suspect, target and unknown compounds. Chromatographic deconvolution is one method utilized extensively for high-resolution GCMS data where fragment ions are aligned and searched against EI spectral libraries. Similar analysis for high-resolution LCMS data is more complicated since compounds can generate multiple adducts, charge states, and in-source fragment ions prior to generating any MSMS fragment ions.

Herein, we demonstrate the use of a new project-based software program, MassHunter Explorer 2.0, that extracts MS and data dependent acquisition (DDA) high-resolution MSMS data. MassHunter Explorer 2.0 uses a simple user-interface to import and data files and do chromatographic retention time (RT) correction to adjust for run-to-run variation. Unsupervised non-targeted data mining provided a means to generated chromatographically significant features with grouping of adducts ( $H^+$ ,  $NH_4^+$ ,  $Na^+$ ,  $K^+$ ), dimers, trimers, simple in-source fragments (e.g.  $H_2O$ ,  $H_2$ ,  $NH_3$ ,  $C_4H_8$ ,  $CO_2$ , etc) and isotopes. Non-targeted MSMS spectra are extracted along with the MS data by adduct and average collision energy. Statistical analysis such as pair-wise fold change, volcano plots, unique features and hierarchical clustering provides a means to quickly determine differences between groups.

Compound identification uses accurate mass and isotope peak abundances to search E&L application-focused commercial and user-generated databases and MS/MS libraries. For compounds not observed in these databases, MassHunter Explorer supports direct export of MSMS spectra for searching either NIST MS Search 3.0 or SIRIUS MSMS software.



Figure 1: Infinity II 1290 LC with Revident LC/Q-TOF.

## Experimental

### Extractables & Leachables from Catheters

Catheters from three suppliers were used in this study:

- Catheter A: Made from PVC, no DEHP, DINP, BPA or natural rubber latex
- Catheter B: Made with natural rubber latex base and silicone elastomer coating, no claim to be phthalate or BPA free.
- Catheter C: Made from silicone-coated natural rubber latex, no DEHP or BPA; contained in secondary plastic packaging bag

The catheters were extracted with 20 mL of pH 2.5 water, pH 10 water, or 50:50 ethanol: water at 37 °C for 24, 48 and 162 hours. The additional bag from catheter C was extracted using 50:50 ethanol: water to check for secondary migration. The extracts were all analyzed using a 1290 Infinity II HPLC system coupled to the Revident LC/Q-TOF (Figure 1, Table 1).

Non-targeted DDA MSMS data was collected using multiple collision energies (10, 20, 40 V); and both medium ion and narrow ion isolations mode. Triplicate Injections using iterative active exclusion was utilized to enhance the collection of MSMS over wide concentration range.

Table 1. Revident LC/Q-TOF Method Parameters

Parameter	Value
Column	Agilent Poroshell AQ-C18 2.1 x 100 mm
Column temperature	40 °C
Injection volume	1 µL
Flow rate	0.35 mL/min
Mobile Phase	A: Water, 2.5 mM $NH_4$ Formate, 0.05% FA B: Methanol, 2.5 mM $NH_4$ Formate, 0.05% FA
LC Gradient	2% B for 1.0 min; to 100% B at 16.0; hold 100% B to 28.0 min; to 2% at 28.1 min post time 5 min
MS Source	Dual AJS ESI
MS Mode	Auto MS/MS, positive polarity
Collision Energy	20; 40
MS Range	40-1700 $m/z$
Acquisition Rate	MS: 4 spectra/s, MSMS: 6 spectra/s
Drying Gas	250 °C, 11 L/min
Nebulizer Pressure	35 psi
Sheath Gas	300 °C, 11 L/min
Capillary Voltage	3500 V

# Results and Discussion

## MassHunter Explorer 2.0 Streamlines E&L Analysis Workflows for LC/Q-TOF

An overview of the LC/Q-TOF workflow for E&L analysis is shown in Figure 2. The LC/Q-TOF workflow for E&L analysis consists of untargeted analysis, identification using MassHunter Explorer 2.0 software, and optional exporting of remaining unknowns to either NIST MS Search 3.0 or SIRIUS for further identification (Figure 2).

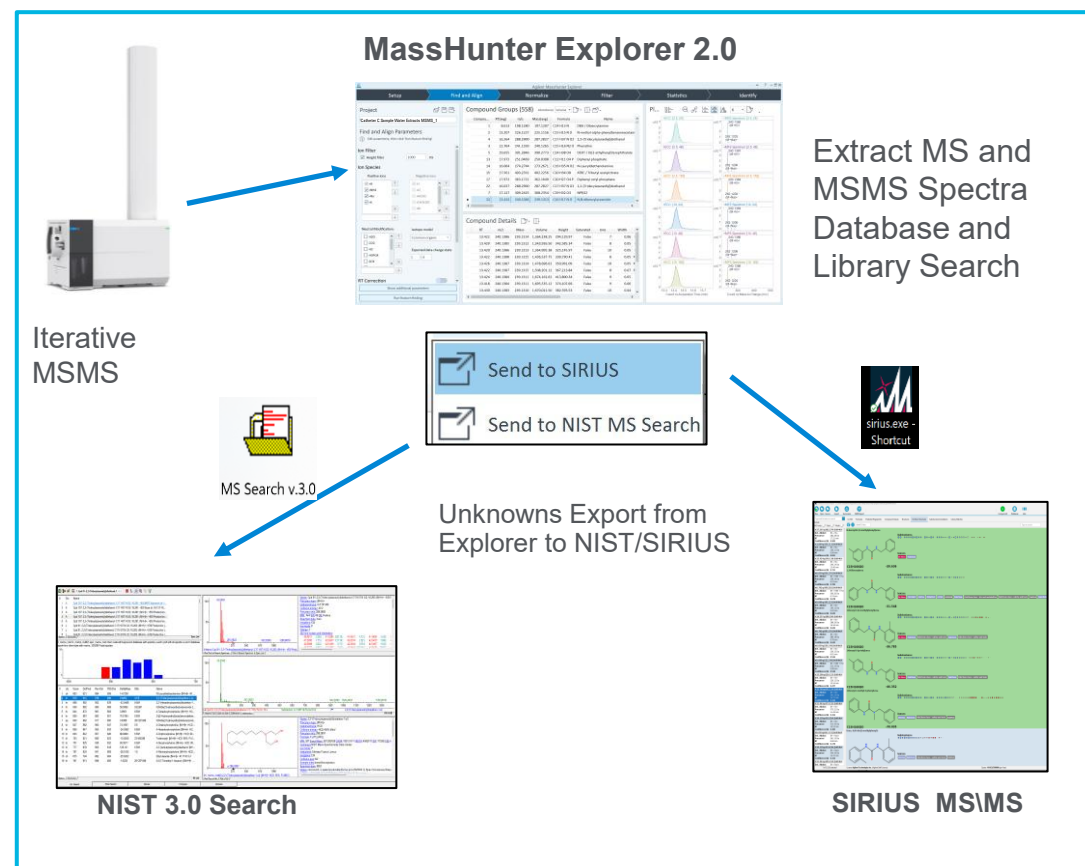


Figure 2 Data Analysis Workflow

## LC/Q-TOF MS and MS/MS Data Processing in MassHunter Explorer 2.0

MassHunter Explorer 2.0 can now extract both MS and MS/MS high-resolution mass spectra by RT, accurate mass, and abundance. The MSMS spectra are extracted by precursor mass (adducts) as average of collision energies with the option for deisotoping to simplify database searching. Results from the "Find and Align" step is shown below in Figure 3.

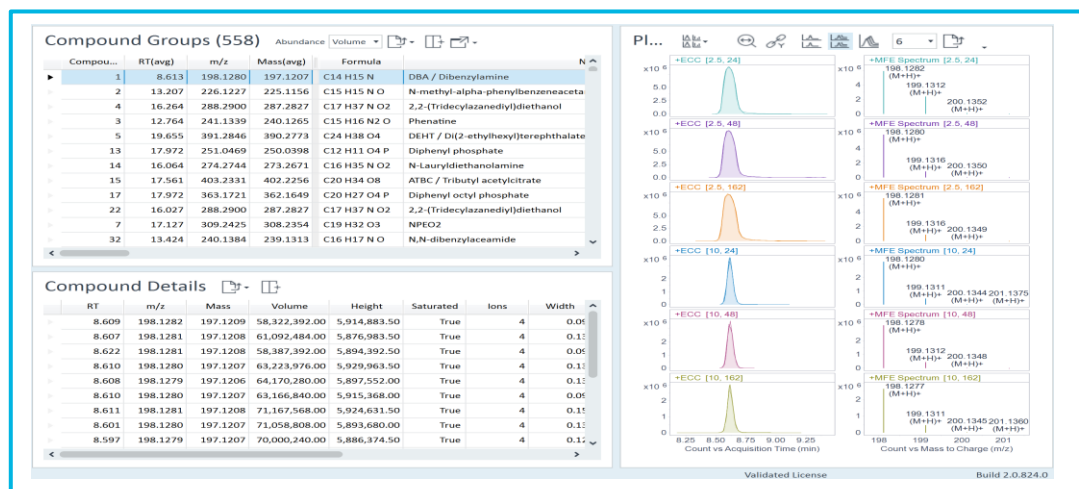


Figure 3. Results from "Find and Align" Function in MassHunter Explorer 2.0.

## Compound Identification in MassHunter Explorer 2.0 Utilizes MS and MSMS Data

Batch compound identification in MassHunter Explorer 2.0 leverages both the MS and MSMS spectra, when available, to match MS-level results based on accurate mass, isotope fidelity, and optional retention time. For compounds with MSMS data, the extracted MSMS data is then searched against the MSMS library spectra based on MSMS tolerance, and forward and reverse scoring, with example library search results shown in Figure 4.

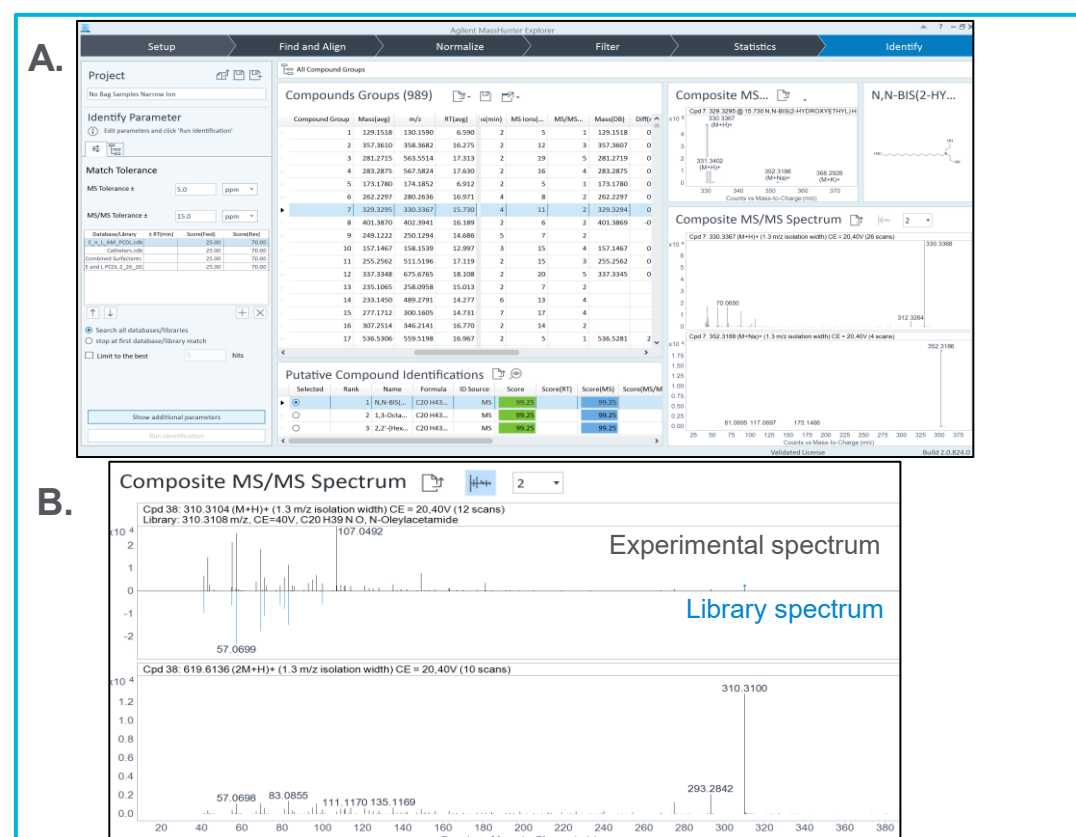


Figure 4A. MassHunter Explorer 2.0 Identification lists library search results, composite MS spectrum, proposed structure, putative compound ID, and the composite MSMS spectra. B) For the MSMS spectra with a library match, mirror plots comparing experimental and library spectra are shown.

## Development of E&L LC/Q-TOF Focused Database

Compound identification in E&L is hindered by the lack of commercial E&L focused databases. Agilent released the first commercial E&L based database that contained 1000+ compounds including over 360 with corresponding MS/MS spectra in 2017. Even large LC-MSMS search libraries like NIST, only contain a very small number of E&L focused compounds compared to metabolomics and environmental application space (Figure 5).

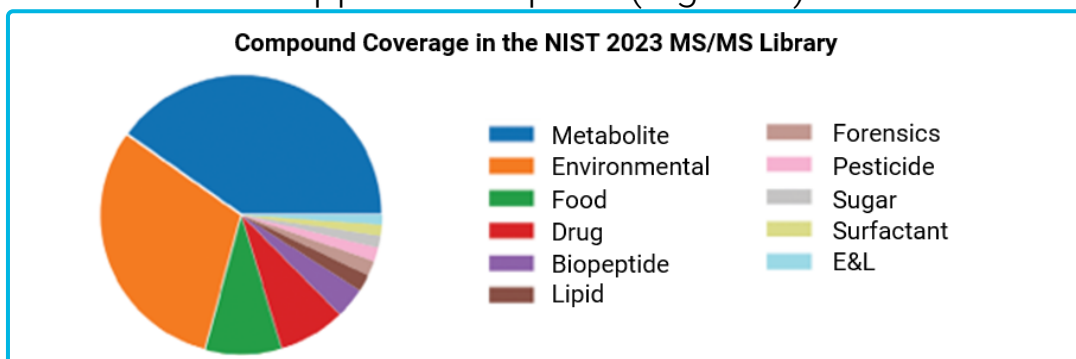


Figure 5. Application Coverage of NIST 2023 Search

# Results and Discussion

## Agilent ChemVista Combines Existing 3<sup>rd</sup> Party Data for Easy Library Building

An alternative approach is to build E&L focused database using information available from resources such as the EPA Comptox dashboard, ELSIE, SciFinder™ and the MassBank of North America (MONA). The Agilent ChemVista database and library management software provides a means to easily import, organize and export databases with chemical identifiers (CAS, SMILES, InChI), MSMS spectra, and RT (Figure 5). Compound lists with MSMS spectra can be exported as \*.SDF files and then converted and used as custom database in NIST MS Search 3.0 software. One can also export a list of compounds with SMILES structures in \*.SDF format that can be used by the SIRIUS MSMS search software.

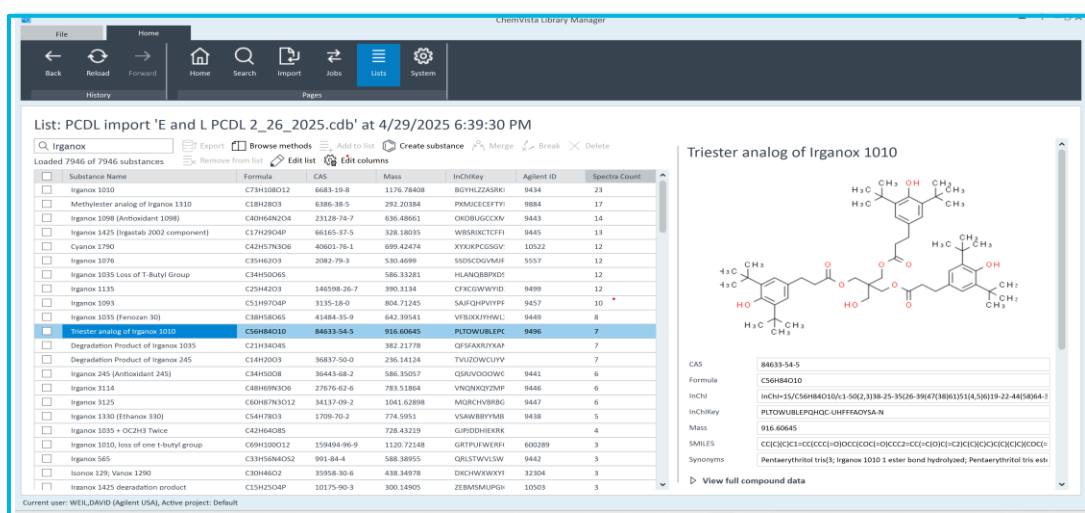


Figure 6. ChemVista database and library software.

## Streamlined Unknowns Identification with Direct Export from Explorer 2.0 to NIST MS Search 3.0

Exporting of compounds directly from Explorer 2.0 into NIST 2023 program provides a quick means to identify compounds using Hybrid MSMS Impurity Tolerant or Partial Spectrum searching. In Figure 7A, we show the MSMS search of  $m/z$  280.2636 correctly identifies adduct as  $[M+H]^+$  corresponding Linoleamide  $C_{18}H_{33}NO$ . Second example, in Figure 7B, has  $m/z$  310.2378 corresponding to  $[M+NH_4]^+$  of  $C_{18}H_{28}O_3$  Methyl 3-(3,5-di-tert-butyl-4-hydroxyphenyl) propanoate, a possible breakdown product.



Figure 7. Results from NIST MSMS Search

<https://www.agilent.com/en/promotions/asms>

This information is subject to change without notice.

DE-007086

© Agilent Technologies, Inc. 2025  
Published in USA, May 15, 2025

## Unknowns Identification using Sirius MSMS Search

A major problem in E&L compound identification based on formulas alone is that results are possibly biased towards compounds present in a database but not confirmed with RT or MSMS library matching. An issue compounded by the limited number of MSMS spectra, including very few from structural isomers, available in commercial databases.

A major advantage of SIRIUS over NIST searching is that SIRIUS will recalculate molecular formulas using MS and MSMS data, correlate MSMS data with predicted molecular fingerprints, generate fragmentation trees, and propose structures from online and custom databases or generate De Novo structures. Using SIRIUS, the presence of several sulfur containing compounds were identified in the EtOH:Water 7-day extract (Catheter C) that were not detected using the NIST 2023 search. Elemental compositions and tentative structures (some de novo generated) are shown below in Figure 8.

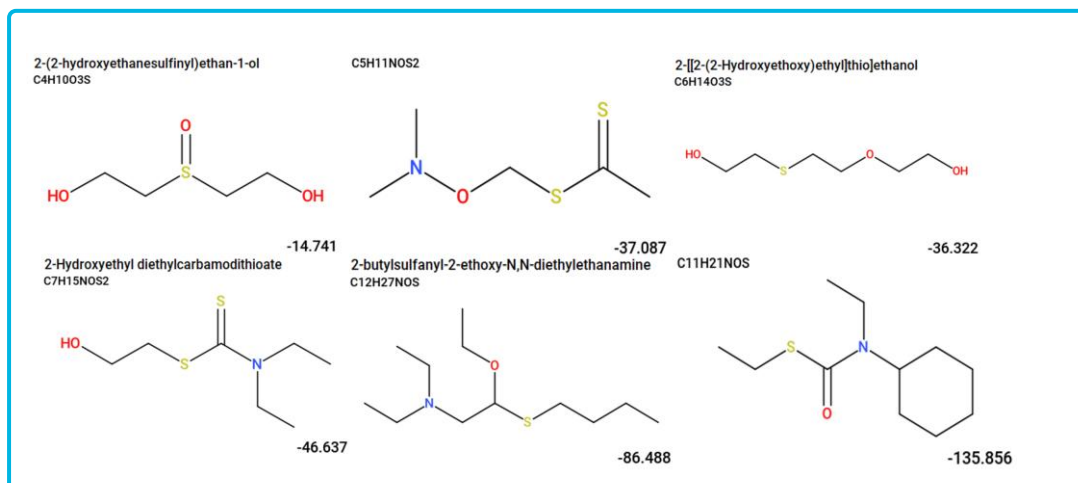


Figure 8. Proposed structures for sulfur compounds in Catheter C EtOH:Water extract

## Conclusions

MassHunter Explorer 2.0 provides a streamlined workflow that improves productivity by:

- Processing of both MS and MSMS mass spectral data
- MSMS searching using custom databases/libraries
- Streamlining export of MSMS spectra to both NIST2023 MSMS Search and SIRIUS MSMS
- Importing results using ChemVista Software

## References

1. Nature BioTechnology, 40, 411-421 (2022)
2. Bright Giant, Sirius Version 6.1.1 (Jan. 2025)

COI: David A. Weil, Sierra D. Durham, James Pyke, and Matt Curtis are employees of Agilent Technologies.