

Build Curated and Annotated HRAM MSⁿ Spectral Libraries To Aid In Unknown Structure Elucidation

Caroline Ding¹, Kate Comstock¹, Seema Sharma¹, Mark Sanders¹, Michal Raab²,
¹Thermo Fisher Scientific, San Jose, CA, USA 95134, ²HighChem, Bratislava, Slovakia



ABSTRACT

Purpose: Validate new library searching algorithms and structure ranking algorithm mzLogic™ for local proprietary MSⁿ spectral libraries to aid in structure elucidation of unknown drug metabolites, impurities or degradants.

Methods: Ten Sildenafil drug standards were infused onto the Thermo Scientific™ Orbitrap ID-X™ Tribrid™ mass spectrometer to obtain MSⁿ spectral trees for building a spectral library. LCMS data were collected with a mixture of four sildenafil standards to test library searching, mzLogic and unknown structure elucidation. Mass Frontier™ 8.0 software was used for the complete workflow including library building, data processing, identification and structure elucidation.

Results: Curated HRAM MSⁿ spectral library which contains substructure annotations helps to elucidate the structures of true unknowns. mzLogic utilizes the spectra and structure information in the local library for unknown structure ranking.

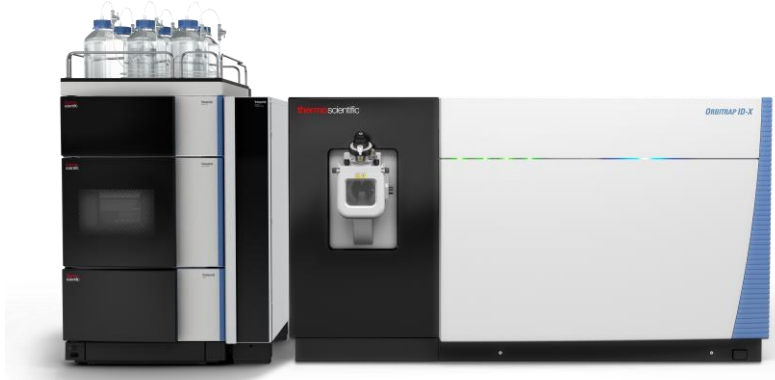
INTRODUCTION

Small molecule structure elucidation is a very challenging and time consuming task. A mass spectral library with extensive MSⁿ spectral tree and substructure information is a valuable tool for rapid identification of small molecule unknowns and unknown unknown structure characterization. The objective of this work is to demonstrate a complete workflow from building local version of “mzCloud” - HRAM MSⁿ spectral library using the Orbitrap ID-X Tribrid mass spectrometer for in-house proprietary compounds with automated curation and structure annotations in Thermo Scientific™ Mass Frontier™ structure identification software 8.0. The MSⁿ spectral library with structure annotations not only enables library searching to quickly identify unknown compounds in the library, but also enables substructure matching. The new mzLogic structure ranking algorithm is able to use the spectral knowledge in the library, rank possible structure candidates based on a combination of spectral similarity and common substructure overlapping, which results in quick and confident structure identification of unknown drug metabolites, impurities or degradants.

MATERIALS AND METHODS

Data Acquisition

An Orbitrap ID-X Tribrid mass spectrometer coupled with a Thermo Scientific™ Vanquish™ UHPLC system were used for data acquisition. As proof of concept, a library was created for ten sildenafil drug analog standards. Compounds were infused by electrospray ionization and MSⁿ data collected with varying collision energies for multiple fragmentation activation types. Four sildenafil standards with same mass (two of them were in the library, and two of them were not in the library) were mixed and analyzed by LC/MSⁿ.



The ten sildenafil drug standards for the library:

