

Poster Reprint

ASMS 2023
Poster number TP 242

Accurate Mass Library for PFAS Analysis in Environmental Samples Using High Resolution GC/Q-TOF

Sofia Nieto¹; Matthew Giardina¹; Matthew Curtis¹; Luann Wong²; Gabrielle Black²; Thomas Young², Charles Yang¹

¹Agilent Technologies, Santa Clara, CA

²Department of Civil and Environmental Engineering,
University of California, Davis, CA

Introduction

Per- and polyfluoroalkyl substances (PFAS) are emerging contaminants of increasing concern due to their environmental persistence, toxicity, and capability of bioaccumulation. There are currently thought to be over 6,000 PFAS that have been commercially produced and recent studies have shown that many emerging PFAS that have been detected in the environment can be volatile or semi-volatile in nature. Therefore, a variety of analytical techniques are necessary for their detection. GC/MS is typically used for detecting volatile and non-polar PFAS compounds. In this study we used GC/Q-TOF system to take advantage of high resolution for detecting compounds with mass defect that is different from that of complex environmental matrices. For specific and sensitive PFAS detection in soil and drinking water, we have created an accurate mass GC/MS library for PFAS. We have also identified other contaminants in drinking water such as disinfection byproducts, industrial chemicals originated from personal care products, drugs, and pesticide residues.



Figure 1. Agilent 7250 GC/Q-TOF

Experimental

GC-amenable PFAS standards have been used to obtain accurate mass spectra. Soil was sampled from two fields in California that have historically received biosolids and extracted with methylene chloride. The drinking water samples were collected at two different locations in California and represented two different water source categories: a small surface water (Weaverville) and a mixed surface and ground water (Irvine). Water samples (2.4 L) were extracted on a multi-mode SPE (HLB, WAX, WCS, Isoelut ENV) and eluted with 5% MTBE in MeOH, DCM, 0.5% NH_4OH in 1:1 EtAc:MeOH, and 1.7% formic acid in 1:1 EtAc:MeOH. The combined extracts were concentrated, solvent exchanged to EtAc and diluted 10x. GC/MS analysis was performed using an Agilent 8890 GC coupled to an Agilent 7250 high resolution Q-TOF (Figure 1) using the following the data acquisition parameters (Table 1).

Table 1. GC/Q-TOF Acquisition Parameters

GC and MS Conditions	DB-5MS	DB-624
MS	7250 Q-TOF	
GC	7890	
Inlet	MMI, 4-mm UI liner single taper with wool	
Inlet temperature	70 °C for 0.01 min; 300 °C/min to 250 °C	
Injection volume	1 μL	
Columns	DB-5MS UI, 30 m x 0.25 mm x 0.25 μm	DB-624 UI, 30 m x 0.25 mm x 1.4 μm
Oven temperature program	35 °C for 2 min; 7 °C/min to 210 °C, 20 °C/min to 300 °C, 4 min hold	30 °C for 2 min; 3 °C/min to 75 °C, 2 °C/min to 110 °C, 10 °C/min to 210 °C, 20 °C/min to 240 °C, 2 min hold
Column flow	1.2 mL/min constant flow	1 mL/min constant flow
Carrier gas	Helium	
Transfer line temperature	250 °C	
Quadrupole temperature	150 °C	
Source temperature	200 °C	
Electron energy	70 eV	
Emission current	Variable by time segment, 0.01 to 5 μA	
Spectral acquisition rate	5 Hz	
Mass range (Tune)	50 to 1200 m/z	

The chromatographic deconvolution and library search were performed in the MassHunter Unknowns Analysis 11.1. Accurate mass EI fragments were converted to the theoretical m/z using MassHunter Qualitative Analysis software version 10.0, and the spectra were exported into the accurate mass Personal Compound Database and Library (PCDL) Manager version 8.0. The PCDL for PFAS, PCDL for Pesticides and Environmental contaminants, as well as NIST 20 were used to perform initial compound identification. Retention Indices and accurate mass information were utilized to confirm the compound ID. Statistical analysis was performed in Mass Profiler Professional (MPP) 15.1.

Accurate Mass Library for PFAS

In order to create an accurate mass GC/MS PCDL, the spectra have been collected for over a hundred PFAS compounds. Accurate mass fragment ions have been automatically annotated with formulas based on accurate mass information and isotope ratios using MassHunter Qualitative Analysis software (Figure 2). The fragment formula annotations were manually verified, corrected when necessary and automatically converted to the theoretical m/z.

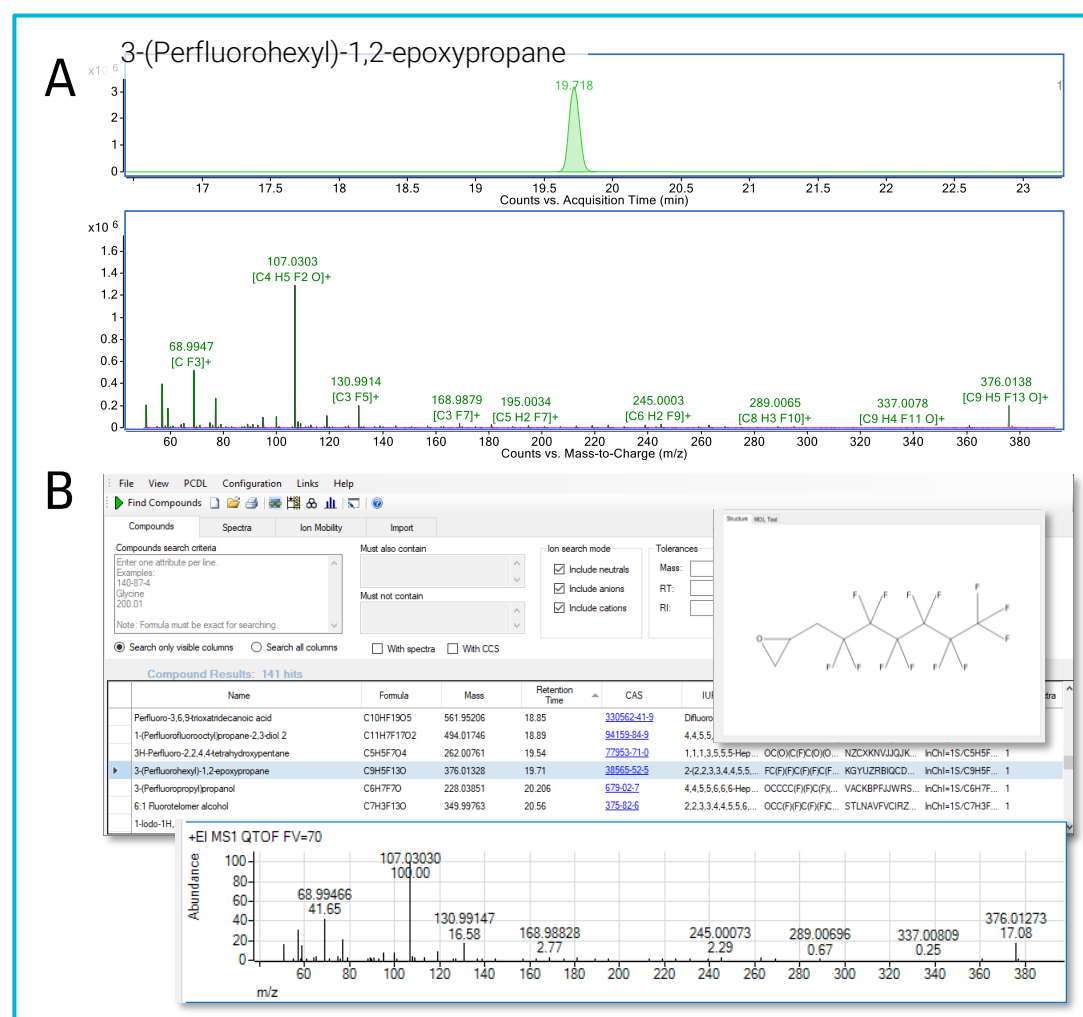


Figure 2. (A) EIC of the molecular ion and fragment formula annotation of spectrum for one of the PFAS compounds in MassHunter Qualitative Analysis software. (B) The PFAS PCDL contains EI spectra as well as the metadata including molecular structure and database identifiers.

The PFAS compound classes include perfluoroalkyl iodides (PFAIs), fluorotelomer iodides (FTIs), fluorotelomer alcohols (FTOHs), fluorotelomer olefins (FTOs), fluorotelomer acrylates (FTACs), fluorotelomer methacrylates (FTMACs) and perfluoroalkyl carboxylic acids (PFCAs) among others (Figure 3).

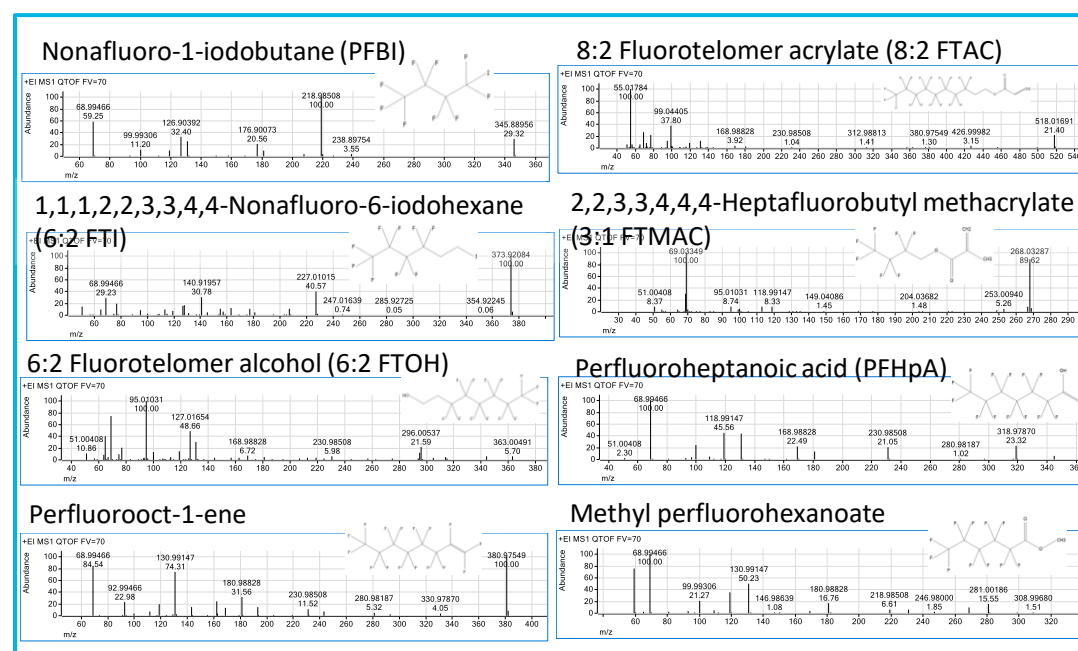


Figure 3. Examples of spectra in PFAS PCDL from different PFAS compound classes.

PFAS in Environmental Samples

For PFAS detection, the extracts of drinking water and soil were separated on DB-624 column analyzed by the GC/Q-TOF MS. The chromatographic deconvolution was performed in the Unknowns Analysis software using a SureMass algorithm that is optimized for complex high resolution EI data. The PFAS PCDL was used to search the deconvoluted spectra with RT matching. Figure 4 shows PFAS compounds identified in soil and drinking water (one in each matrix). PFAS (a derivative of PFCAs) were detected in most drinking water samples and the soil extract from Field 1.

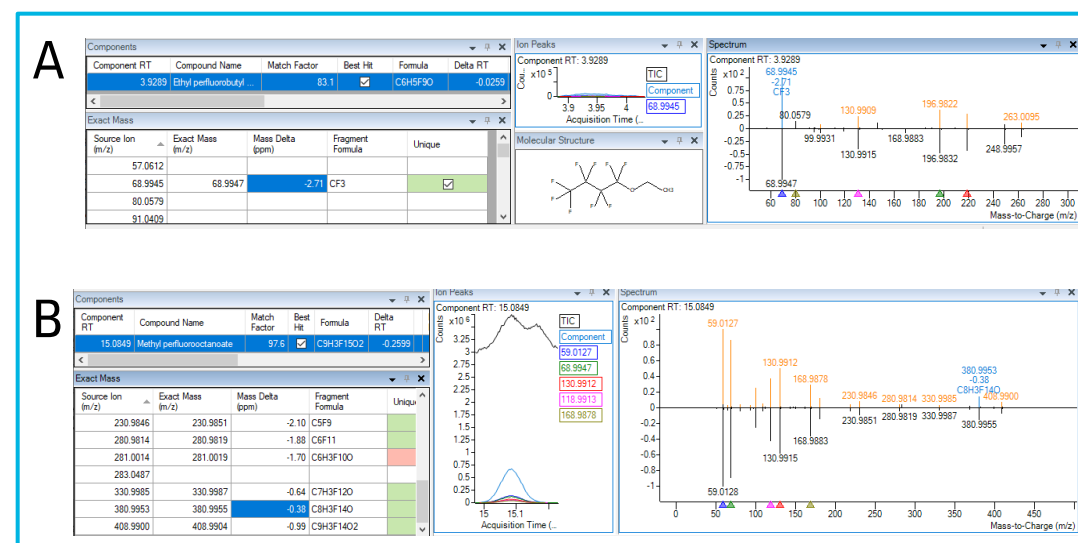


Figure 4. PFAS identified in soil Field 1 (A, ethyl perfluorobutyl ether) and drinking water Irvine (B, Methyl perfluorooctanoate) samples.

Identification of the Additional Contaminants in Drinking Water Samples

To identify other contaminants in drinking water samples the GC/Q-TOF Pesticide PCDL as well as NIST 20 library were used. Over hundred contaminants have been identified and confirmed using accurate mass information (Figure 5 and Tables 2 and 3) from sample without reinjection.

