

The effect of particle size reduction techniques on extraction and recovery of 16 PFAS in food-contact paper packaging matrices

Authors: Paulo Silva, Keith Vorst, Greg Curtzwiler, Iowa State University, Ames, IA, USA; Charles Yang, Thermo Fisher Scientific, San Jose, CA, USA

Keywords: Polyfluoroalkyl and Perfluoroalkyl Substances, PFAS, PFAA, PFOS, PFOA, GenX, PFCs, Extraction, Sample Preparation, Spike Recovery, Food Contact Materials, Orbitrap Exploris 120 mass spectrometer, TraceFinder software, Vanquish Flex UHPLC system

Goal

To investigate the effect of two particle size reduction techniques, ball milling and blade cutting-grinding, on the Perfluoroalkyl and Polyfluoroalkyl substances (PFAS) recovery of three types of food contact materials (microwave popcorn bags, molded fiber bowl, and wrappers) using Focused Ultrasound Solid-Liquid Extraction (FUSLE) technique for extraction of 16 targeted fluorinated compounds.

Introduction

Paper and board (P&B) food contact materials (FCMs) have gained special attention lately due to the broad use of commercial additive blends and raw materials of known and unknown nature that have raised health and regulatory concerns.



Natural fibers of bleached or unbleached cellulose are used to make P&B, moreover, it can also be recycled from recovered materials. P&B used as FCMs can be noted as chemically complex matrices, partly due to the naturally occurring substances in P&B, but also due to chemical treatments used to make these materials suitable for food contact (Bengtström et al., 2014). Since P&B are used in a variety of applications, one of the challenges in paper production is to achieve specific technical functionalities. Therefore, the use of chemical additives is widely employed in the manufacturing process to achieve various performance requirements. Among these additives, processing aids and functional additives are some of the main categories. Processing aids are used to improve the efficiency of the paper making processes and are not intended to be transferred into the final product, although traces can be found. Some typical examples are defoamers, biocides, felt cleaners, and deposit control agents.

Commercial fluorochemicals are very useful in P&B production as they impart water, oil, and stain repellency onto these materials and act as dispersion and levelling agents. Polyfluorinated surfactants (PFS) are the main group of chemicals in commercial blends used to improve paper technical performance. Proper quantification of these substances in extracts of P&B are not always possible, as appropriate analytical standards are not commercially available. Although they can be relatively stabilized in paper matrices, they are still of major concern as they can be precursors of poly- and per-fluorinated alkyl substances (PFAS). PFAS can function as monomers or be attached to a polymer backbone in these matrices. PFAS have been used in paper and P&B packaging since the 1950s, mostly as coatings to prevent the paper material from soaking up fats and water, but also in printing inks and as moisture barriers (Trier et al., 2017). Some examples include fast food paper wrappers, microwave popcorn bags, cake forms, sandwich and butter paper, chocolate paper, paper for dry foods and pet foods. PFAS have been linked to a variety of human health issues, including cancers, decreases in fertility, and reduced immune system function, raising concerns for health and toxicology studies as well as regulations (E.P. et al., 2009; Fair et al., 2011; Hines et al., 2009; Macon et al., 2011; Pelch et al., 2019; Rosenmai et al., 2016; Tucker et al., 2015). Long-chain PFAS (8 carbons or higher) have been phased out in the United States and European Union due to these health concerns (United States Environmental Protection Agency, 2006), but they are often replaced in manufacturing processes by short-chain PFAS (Wang et al., 2015). More studies revealed that short-chain PFAS may carry similar health concerns as long-chain PFAS, despite the short-chain compounds reduced bioaccumulation (Scheringer et al., 2014; The Danish Environmental Protection Agency, 2015).

PFAS in general have been found in surface water, groundwater, finished drinking water, rainwater, and air emissions in some areas. Currently there are no maximum contaminant levels established for PFAS in food packaging, US-EPA has established drinking water health advisories for PFOA and PFOS at 70 parts per trillion (States & Protection, 2009). PFAS as non-intentionally added substances (NIAS) can also find their way into P&B matrices. Since PFAS are ubiquitous in the environment, it can be present in processing water in paper mills.

Given the health and migration concerns associated with PFAS in FCM, it is imperative that adequate and good performance analytical methods be developed to quantify different PFAS accurately and efficiently in a variety of matrices.

Method sensitivity for detection of PFAS have been improved dramatically in the last few decades by the use of advanced analytical technologies such as triple quadrupole tandem mass spectrometry coupled to liquid chromatography (LC-MS/MS) and liquid chromatography quadrupole high resolution accurate mass spectrometry (LC-HRAM). However, efforts to improve extraction and clean-up of solid-matrix samples for subsequent analysis on those technologies are still needed in order to minimize uncertainties and assure instrumental reproducibility and accuracy in workflows. Several extraction methods have been reported for the extraction of PFAS in different matrices (Nakayama et al., 2019). Examples of methods used on solid matrices include: solid-liquid extraction (SLE), pressurized liquid extraction (PLE), ultrasound assisted extraction (UAE), microwave assisted extraction (MAE), and focused ultrasound solid-liquid extraction (FUSLE) (Martínez-Moral & Tena, 2013; Monge Brenes et al., 2019; Zabaleta et al., 2014). Among all these methods, FUSLE has been validated and shown to be a low-cost, fast, simple and safe extraction technique with PFAS recoveries on food matrices and popcorn bags of nearly 100% (Moreta & Tena, 2013, 2014). Although it is known that a reduction in particle size could always lead to increased extraction efficiency, none of these studies address this variable, which seem to pose a challenge whenever sampling solid matrices.

This application note describes a direct analysis workflow for the determination of 16 targeted PFAS in three types of food contact matrices (microwave popcorn bags, molded fiber bowl, and food wrappers), evaluating their recovery through extractions employing two particle size reduction techniques.

Experimental

Sample preparation

Sample cutting was performed on each packaging material using a trimmer (Swingline™ ClassicCut™ Lite Guillotine) with a stainless-steel blade to obtain a sample material consisting of rectangles of 25–50 mm². Sections of popcorn (no susceptor was used as part of the sample) and sandwich bags that contained adhesives were removed prior to cutting. Compressed house air was used to remove excess

sample from cutting surfaces, and all cutting, and sample preparation surfaces were cleaned with methanol between matrix types to avoid cross-contamination.

Cutting grinding—each pre-cut packaging material was pulverized into uniform particle powder using an IKA™ A11 Analytical Mill with a stainless-steel fiber cutting blade attached (30 seconds for popcorn bags; 1 minute for sandwich bags; 1 minute for the paper bowls). The mill and its components were cleaned completely with methanol between samples of different materials to avoid cross-contamination.

Ball milling—each pre-cut packaging material was pulverized into uniform particle powder using a Retsch™ MM 400 ball mill with a 25 mm diameter stainless-steel ball in each 50 cm³ stainless steel jars and stainless-steel balls, with each type requiring two 1-minute intervals to be ground into a fine powder at 30 Hz. Alconox® detergent solution was used to clean the ball/jar followed by methanol rinsing between samples of different materials to avoid cross-contamination.

Spiked samples were prepared by adding a methanolic PFAS native standard solution to each powdered packaging material (20 ppb) dispersed in ethyl acetate (Monge Brenes et al., 2019). The suspended and spiked samples were mixed thoroughly, then evaporated to dryness using a water bath set at 45 °C, and ground again to ensure homogeneity. Spiked and non-spiked samples were stored in polyethylene bags wrapped in aluminum foil and refrigerated (Frigidaire™, FFTR1814TWO) at 4 °C for subsequent analyses.

Extraction

Extractions of the different PFAS-spiked samples were carried out by following a validated method (Moreta & Tena, 2014) with few modifications. A focused ultrasonic liquid extraction (FUSLE) procedure using a Misonix™ S-4000 Ultrasonic Sonicator with a power of 600 W and an operating frequency of 20 kHz, equipped with a 3 mm titanium tip, was utilized to extract the PFAS from the samples. Each different sample material spiked with the native PFAS standard cocktail had undergone three extractions. A known amount of ground paper (~1.000 ± 0.001 g of homogenized sample) was placed into a 50 mL (34 × 100 mm) glass centrifuge tube and 24 mL of HPLC grade ethanol was added to each sample. Before each extraction, 100 µL of 300 ng mL⁻¹ of mass labeled PFAS standard solution was added to each tube. The weight of sample used in each extraction was accurately recorded

and used to normalize the concentration of PFAS obtained per gram of paper. The sonicator probe was inserted in the mixture to a depth of 2 cm from the bottom of the test tube. Each individual tube was then secured in an ice bath and subsequently sonicated. Samples were exposed to 30% amplitude at 50% pulsed cycle for 10 s. Extracts were filtered through a 60 mL Pyrex® Buchner funnel with fritted disc and porosity 10–15 µm using a vacuum pump at 550 in Hg vacuum. The probe, glassware, and extracted samples were washed twice with 2.5 mL of ethanol each rinse. The total amount of filtered extract with rinses was transferred to a 40 mL Pyrex scintillation vial without cap and immediately evaporated to dryness under a nitrogen stream using a nitrogen evaporator (N-EVAP™ 111) equipped with water bath set at 45 °C. The dry residue was reconstituted with 1 mL of LC-MS grade methanol and filtered into a 300 µL polypropylene LC vial using a disposable polypropylene medical sterile syringe equipped with a 13 mm diameter, 0.22-µm nylon filter.

System configuration

A Thermo Scientific™ Vanquish™ Flex Binary UHPLC system coupled to a Thermo Scientific™ Orbitrap Exploris™ 120 mass spectrometer was employed for the separation and quantification of the PFAS. The setup for the UHPLC and gradient profile used to elute and separate the PFAS is shown in Table 1.

Table 1. LC parameters for chromatographic separation of the PFAS

| | | |
|------------------------------------|--|-------------|
| Analytical column | Thermo Scientific™ Hypersil GOLD™ aQ column, 100 × 2.1 mm, 1.9 µm (P/N 25302-102130) | |
| PFAS Upgrade Kit for Vanquish Flex | PN 80100-62144 | |
| Trap column | Hypersil GOLD column, 50 × 3 mm, 1.9 µm (P/N 25002-053030) | |
| Column temperature | 30 °C | |
| Autosampler temperature | 7 °C | |
| Flow rate | 300 µL/min | |
| Solvent A | LC-MS grade water acidified with formic acid to 0.1% (v/v) with ammonium formate to obtain 10 mM | |
| Solvent B | LC-MS grade methanol | |
| Injection volume | 1.0 µL | |
| Gradient | Time (min) | % Solvent B |
| | 0 | 60 |
| | 4 | 100 |
| | 6 | 100 |
| | 7 | 20 |
| | 8 | 100 |
| | 9 | 20 |
| | 10 | 100 |
| | 11 | 20 |
| | 12 | 60 |
| | 13.5 | 60 |

Table 2. List of targeted PFAS included in this method.

| Native Analyte | Native Acronym (CAS) | Formula (native) | Surrogate Analyte |
|---|-----------------------------|---|-------------------|
| Perfluorobutanoic acid | PFBA (375-22-4) | C ₄ F ₇ O ₂ | M3PFBA |
| Perfluoropentanoic acid | PFPeA (2706-90-3) | C ₅ HF ₉ O ₂ | M3PFPeA |
| Perfluorobutanesulfonic acid | PFBS (375-73-5) | C ₄ HF ₉ O ₃ S | M3PFBS |
| Perfluorohexanoic acid | PFHxA (2706-90-3) | C ₆ HF ₁₁ O ₂ | MPFHxA |
| Perfluoropentanesulfonic acid | PFPeS (2706-91-4) | C ₅ HF ₁₁ O ₃ S | MPFHxS |
| 2,3,3,3-Tetrafluoro-2-(1,1,2,2,3,3,3- heptafluoropropoxy) propanoic acid (GenX) | HPFO-DA (62037-80- 3) | C ₆ HF ₁₁ O ₃ | M3HFPO-DA |
| Perfluoroheptanoic acid | PFHpA (375-85-9) | C ₇ HF ₁₃ O ₂ | MPFHxA |
| Perfluorohexanesulfonic acid | PFHxS (355-46-4) | C ₆ HF ₁₃ O ₃ S | MPFHxS |
| Sodium dodecafluoro-3H-4, 8- dioxananoate | NaDONA (958445- 44-8) | C ₇ H ₅ F ₁₂ NO ₄ | M8PFOA |
| Perfluoroheptanesulfonic acid | PFHpS (375-92-8) | C ₇ HF ₁₅ O ₃ S | MPFHxS |
| Perfluorooctanoic Acid | PFOA (335-67-1) | C ₈ HF ₁₅ O ₂ | M8PFOA |
| Perfluorooctanesulfonic acid | PFOS (1763-23-1) | C ₈ HF ₁₇ O ₃ S | M8PFOS |
| Perfluorononanoic acid | PFNA (375-95-1) | C ₉ HF ₁₇ O ₂ | M8PFOA |
| Potassium 9-chlorohexadecafluoro-3- oxanonane-1-sulfonate | 9Cl-PF3ONS (73606-19- 6) | C ₉ ClF ₁₆ KO ₄ S | MPFHxS |
| Perfluorodecanoic acid | PFDA (335-76-2) | C ₁₀ HF ₁₉ O ₂ | M8PFOA |
| 11-chloroeicosafuoro-3-oxaundecane-1- sulfonic acid | 11Cl-PF3OUdS (763051- 92-9) | C ₁₀ HClF ₂₀ O ₄ S | MPFHxS |

Table 3. Mass spectrometry conditions.

| Source Parameter | Value |
|--|---------------------|
| Ionization | H-ESI Negative mode |
| Spray voltage | 2.5 kV |
| Ion transfer tube temperature | 256 °C |
| Sheath gas | 48 |
| Aux gas | 11 |
| Vaporizer temperature | 413 °C |
| Sweep gas | 1 |
| RF lens | 70 |
| Acquisition Type | Full MS |
| Resolution | 60,000 |
| AGC target | Standard (1e6) |
| Maximum injection time | Auto |
| Mass range | 65–650 m/z |
| Acquisition Type | DIA |
| Resolution | 30,000 |
| AGC target | Standard (1e6) |
| Maximum injection time mode | Auto |
| Loop Count | 1 |
| Isolation window | 200 m/z |
| Fragmentation mode | HCD |
| HCD collision energy (%) / stepped NCE | 10, 30 |
| Number of Scan Events | 5 |

Data analysis

Identification and quantitation of PFAS targets

TraceFinder software was used to process the targeted screening quantitative data. Identification of the PFAS compounds was done by matching the retention time of the native standards, and confirmation identity was accomplished using accurate-mass measurements with spectral library matching. Confirmation of native PFAS compounds and their respective surrogate analytes are shown in Table 4. The concentration of each PFAS was determined using the response ratio of the PFAS quantitation (abundance of the precursor using MS¹ filtering mode) from the inclusion list (Table 4) to that of the relevant labeled surrogate standard.

Statistical analysis

Statistical analyses were performed using JMP Pro™ software (v. 14). Student t-test was used to compare whether the means of the two sets of data were statistically significantly different from each other.

Table 4. Inclusion list used for targeted PFAS analysis in data processing.

| Analyte | Retention Time (min) | Precursor Ion (m/z) | Product Ion (m/z) |
|--------------|----------------------|---------------------|---|
| PFBA | 1.18 | 212.9792 | 168.98937 |
| M3PFBA | 1.18 | 215.98926 | 171.9997 |
| PFPeA | 1.52 | 262.97601 | 218.9862 |
| M3PFPeA | 1.52 | 265.98636 | 222.9900 |
| PFBS | 1.54 | 298.94299 | 79.9574 98.9558 298.9430 |
| M3PFBS | 1.54 | 301.95306 | 79.9574 98.9558 |
| PFHxA | 2.12 | 312.97281 | 268.9830 |
| MPFHxA | 2.12 | 314.97952 | 268.9835 269.9867 |
| PFPeS | 2.14 | 348.9398 | 79.9574 98.9558 118.9925 |
| HPFO-DA | 2.34 | 328.96773 | 168.9895 284.9783 |
| M3HFPO-DA | 2.34 | 331.97779 | 286.9849 |
| PFHpA | 3.07 | 362.96962 | 318.9798 |
| PFHxS | 3.09 | 398.9366 | 79.9574 98.9558 118.9925 |
| MPFHxS | 3.09 | 402.945 | 84.9907 169.9891 250.9761 376.9688 398.9358 |
| NaDONA | 3.13 | 376.96887 | 84.9907 250.9761 |
| PFHpS | 3.77 | 448.93341 | 79.9574 98.9558 168.9894 |
| PFOA | 3.77 | 412.96643 | 168.9894 368.9766 |
| M8PFOA | 3.77 | 420.99326 | 118.9926 171.9995 375.9997 |
| PFOS | 4.23 | 498.93022 | 79.9574 98.9558 |
| M8PFOS | 4.23 | 506.95706 | 79.9573 98.9557 418.9731 498.9295 |
| PFNA | 4.25 | 462.96323 | 168.9894 268.983 418.9734 |
| 9CI-PF3ONS | 4.41 | 530.89558 | 82.9609 98.9557 198.9492 350.9442 |
| PFDA | 4.62 | 512.9606 | 168.9894 268.9835 318.97979 468.9701 |
| 11CI-PF3OUdS | 4.99 | 630.8902 | 82.9609 198.9493 450.9386 |
| d5-N-EtFOSAA | 5.59 | 531.0093 | 168.9896 218.9864 268.9835 330.0905 |

Results and discussion

Recovery of the 16 PFAS compounds spiked into the three different food contact matrices using two different particle-size reduction techniques is summarized in Table 5. The recovery values were higher or lower for each analyte within each food contact material between the two particle-size reduction techniques used in this study.

The great majority of compounds analyzed in these methods were within the recovery range of 70–150%, except for three compounds (HPFO-DA, PFHpA, PFHxA) that resulted in lower recoveries depending on the type of matrix that were analyzed. The lower recovery observed might be related to co-eluting matrix components that may have caused signal suppression since minimal sample clean-up was used in these methods.

It can be observed that very low coefficient of variation values were achieved through independent extractions. The values for % recoveries of the compounds PFBA and L-PFBS analyzed in the microwave popcorn bag are not available since copious amounts of these analytes were already present in this matrix. Those abundances were out of the concentration range of the analyte's responses used in the standard calibration curves.

An evaluation of the total mass of all 16 PFAS recovered (spiked + already present) per mass of the food contact matrix using both particle-size reduction techniques can be found in Table 6. The use of the ball analytical mill and blade analytical mill for particle-reduction of samples from the molded fiber bowl and brown sandwich bags show no statistical differences between the two techniques for total recovery of PFAS and also from microwave popcorn bag samples. The amounts of PFBA and L-PFBS analyzed in the microwave popcorn bag were not taken into account for this evaluation for the reason explained earlier in this section.

Table 5. PFAS recoveries (%) and (%CV) of three different food contact matrices using two particle-size reduction techniques for extraction.

| Analyte | Molded Fiber Bowl | | | |
|--------------|----------------------------------|---------|-----------------------------------|---------|
| | Ball Analytical Mill (%Recovery) | %CV>10% | Blade Analytical Mill (%Recovery) | %CV>10% |
| PFBA | 84.60 | 1.04 | 75.78 | 13.59 |
| PFPeA | 88.04 | 0.25 | 84.72 | 4.56 |
| L-PFBS | 85.82 | 3.32 | 85.76 | 1.32 |
| PFHxA | 117.81 | 3.34 | 129.02 | 6.22 |
| L-PFPeS | 82.56 | 0.83 | 81.93 | 2.02 |
| HPFO-DA | 73.08 | 0.65 | 49.32 | 2.72 |
| PFHpA | 94.09 | 3.30 | 99.53 | 2.85 |
| L-PFHxS | 90.55 | 2.00 | 89.33 | 2.10 |
| NaDONA | 89.71 | 4.22 | 87.78 | 5.14 |
| L-PFHpS | 96.75 | 1.73 | 97.38 | 0.42 |
| PFOA | 88.80 | 4.95 | 93.25 | 4.23 |
| PFOS | 91.09 | 2.31 | 90.38 | 2.87 |
| PFNA | 92.80 | 5.30 | 91.72 | 6.16 |
| 9CI-PF3ONS | 98.81 | 0.06 | 96.55 | 3.56 |
| PFDA | 88.85 | 3.69 | 93.41 | 6.07 |
| 11CI-PF3OUdS | 111.30 | 0.35 | 111.33 | 1.66 |

Table 5. Continued.

| Brown Sandwich Bag | | | | |
|-----------------------|----------------------------------|---------|-----------------------------------|---------|
| Analyte | Ball Analytical Mill (%Recovery) | | Blade Analytical Mill (%Recovery) | |
| | | %CV>10% | | %CV>10% |
| PFBA | 81.75 | 3.84 | 69.63 | 5.79 |
| PFPeA | 85.82 | 3.13 | 75.94 | 1.78 |
| L-PFBS | 81.91 | 4.36 | 79.42 | 0.14 |
| PFHxA | 110.62 | 27.14 | 47.60 | 4.54 |
| L-PFPeS | 76.10 | 3.05 | 69.94 | 2.56 |
| HPFO-DA | 88.10 | 1.57 | 70.50 | 3.06 |
| PFHpA | 106.67 | 0.49 | 101.74 | 4.27 |
| L-PFHxS | 89.97 | 2.22 | 86.32 | 2.80 |
| NaDONA | 82.61 | 0.11 | 80.44 | 4.35 |
| L-PFHpS | 94.52 | 0.51 | 87.72 | 4.03 |
| PFOA | 97.18 | 3.13 | 90.01 | 5.67 |
| PFOS | 93.22 | 0.58 | 87.93 | 4.52 |
| PFNA | 119.64 | 0.30 | 117.18 | 3.83 |
| 9CI-PF3ONS | 95.18 | 0.20 | 89.36 | 5.84 |
| PFDA | 130.49 | 1.24 | 120.92 | 3.85 |
| 11CI-PF3OUdS | 99.14 | 3.22 | 93.57 | 6.46 |
| Microwave Popcorn Bag | | | | |
| Analyte | Ball Analytical Mill (%Recovery) | | Blade Analytical Mill (%Recovery) | |
| | | %CV>10% | | %CV>10% |
| PFBA | NA | NA | NA | NA |
| PFPeA | 156.52 | 33.99 | 78.10 | 5.30 |
| L-PFBS | 88.11 | 2.41 | 87.07 | 0.74 |
| PFHxA | NA | NA | NA | NA |
| L-PFPeS | 90.36 | 3.65 | 93.89 | 1.20 |
| HPFO-DA | 87.03 | 10.98 | 80.61 | 0.60 |
| PFHpA | 123.48 | 11.56 | 130.21 | 4.28 |
| L-PFHxS | 94.14 | 2.39 | 92.63 | 1.77 |
| NaDONA | 75.70 | 1.52 | 70.99 | 7.78 |
| L-PFHpS | 109.42 | 5.75 | 112.72 | 1.44 |
| PFOA | 89.74 | 8.88 | 81.02 | 4.56 |
| PFOS | 91.60 | 0.27 | 92.17 | 1.14 |
| PFNA | 116.91 | 4.72 | 108.22 | 6.26 |
| 9CI-PF3ONS | 114.97 | 5.36 | 117.61 | 2.07 |
| PFDA | 159.79 | 1.50 | 151.41 | 10.76 |
| 11CI-PF3OUdS | 115.13 | 2.22 | 116.21 | 0.88 |

Table 6. Total PFAS recovered (ng/g) and (%CV) in three different food contact matrices using two particle-size reduction techniques for extraction.

| Molded Fiber Bowl | | | | Brown Sandwich Bag | | | | Microwave Popcorn Bag | | | |
|-----------------------------|------|------------------------------|------|-----------------------------|------|------------------------------|------|-----------------------------|------|------------------------------|------|
| Ball Analytical Mill (ng/g) | %CV | Blade Analytical Mill (ng/g) | %CV | Ball Analytical Mill (ng/g) | %CV | Blade Analytical Mill (ng/g) | %CV | Ball Analytical Mill (ng/g) | %CV | Blade Analytical Mill (ng/g) | %CV |
| 325.7 | 1.56 | 316.6 | 0.11 | 322.9 | 3.52 | 305.3 | 1.92 | 314.7 | 1.32 | 313.1 | 2.06 |

Conclusion

Particle size reduction techniques for extraction and spike recovery of 16 Perfluoroalkyl and Polyfluoroalkyl substances (PFAS) including seven perfluorocarboxylic acids (PFCAs), five perfluoroalkylsulfonates (PFASs), two chloroperfluoroether sulfonates (Cl-PFESAs), one polyfluoroether carboxylate (PFECAs), and one hexafluoropropylene oxide dimer acid (GenX) on sampling of different cellulosic-based food contact materials were developed in the present study.

No statistical differences were found between the two milling techniques for recovery of the total amount of PFAS spiked in the microwave popcorn bag, molded fiber bowl, and sandwich wrapper matrices. The ultra-high performance liquid chromatography and electrospray ionization (UHPLC/ESI) quadrupole Orbitrap Exploris 120 high-resolution mass spectrometry was reliably able to detect and quantify the analytes in relatively heavy matrices that used minimal clean up during extraction with the surrogate dilution method to correct for matrix effects. The applicability of this methodology in PFAS analysis can be improved in the future by developing strategies for more comprehensive sample clean up depending on the type of matrix used. Among those, solid phase chromatography can be suited for extraction of analytes of different polarities and matrix-interactions. As laboratories are often focused on method detection of analytes, sample preparation procedure studies are of great value to obtain reliable data and to streamline the workflow and reduce turnover time.

Acknowledgments

Research funded by the Packaging and Food Protection Consortium (PFPC) at the Iowa State University, a food industry-sponsored consortium of over 20 food processing and packaging companies and their suppliers. This research is a part of a collaboration between PFPC and Thermo Fisher Scientific.

Find out more at

thermofisher.com/FoodPackaging

References

1. Bengtström, L., Trier, X., Granby, K., Rosenmai, A. K., & Petersen, J. H. (2014). Fractionation of extracts from paper and board food contact materials for in vitro screening of toxicity. *Food Additives and Contaminants - Part A Chemistry, Analysis, Control, Exposure and Risk Assessment*. <https://doi.org/10.1080/19440049.2014.912357>
2. Martínez-Moral, M. P., & Tena, M. T. (2013). Focused ultrasound solid-liquid extraction of perfluorinated compounds from sewage sludge. *Talanta*. <https://doi.org/10.1016/j.talanta.2013.02.020>
3. Monge Brenes, A. L., Curtzwiler, G., Dixon, P., Harrata, K., Talbert, J., & Vorst, K. (2019). PFOA and PFOS levels in microwave paper packaging between 2005 and 2018. *Food Additives and Contaminants: Part B Surveillance*. <https://doi.org/10.1080/19393210.2019.1592238>
4. Moreta, C., & Tena, M. T. (2013). Fast determination of perfluorocompounds in packaging by focused ultrasound solid-liquid extraction and liquid chromatography coupled to quadrupole-time of flight mass spectrometry. *Journal of Chromatography A*. <https://doi.org/10.1016/j.chroma.2013.06.024>
5. Moreta, C., & Tena, M. T. (2014). Determination of perfluorinated alkyl acids in corn, popcorn and popcorn bags before and after cooking by focused ultrasound solid-liquid extraction, liquid chromatography and quadrupole-time of flight mass spectrometry. *Journal of Chromatography A*. <https://doi.org/10.1016/j.chroma.2014.06.018>
6. Nakayama, S. F., Yoshikane, M., Onoda, Y., Nishihama, Y., Iwai-Shimada, M., Takagi, M., Kobayashi, Y., & Isobe, T. (2019). Worldwide trends in tracing poly- and perfluoroalkyl substances (PFAS) in the environment. In *TrAC - Trends in Analytical Chemistry*. <https://doi.org/10.1016/j.trac.2019.02.011>
7. Pelch, K. E., Reade, A., Wolffe, T. A. M., & Kwiatkowski, C. F. (2019). PFAS health effects database: Protocol for a systematic evidence map. In *Environment International*. <https://doi.org/10.1016/j.envint.2019.05.045>
8. Rosenmai, A. K., Taxvig, C., Svengen, T., Trier, X., van Vugt-Lussenburg, B. M. A., Pedersen, M., Lesné, L., Jégou, B., & Vinggaard, A. M. (2016). Fluorinated alkyl substances and technical mixtures used in food paper-packaging exhibit endocrine-related activity in vitro. *Andrology*. <https://doi.org/10.1111/andr.12190>
9. Scheringer, M., Trier, X., Cousins, I. T., de Voogt, P., Fletcher, T., Wang, Z., & Webster, T. F. (2014). Helsingør Statement on poly- and perfluorinated alkyl substances (PFASs). *Chemosphere*. <https://doi.org/10.1016/j.chemosphere.2014.05.044>
10. States, U., & Protection, E. (2009). Provisional Health Advisories for Perfluorooctanoic Acid (PFOA) and Perfluorooctane Sulfonate (PFOS). Environmental Protection.
11. The Danish Environmental Protection Agency. (2015). Short-chain Polyfluoroalkyl Substances (PFAS). In Environmental project No. 1707.
12. Trier, X., Taxvig, C., Rosenmai, A. K., & Pedersen, G. A. (2017). PFAS in paper and board for food contact - options for risk management of poly- and perfluorinated substances. In Nordic Council of Ministers.
13. Tucker, D. K., Macon, M. B., Strynar, M. J., Dagnino, S., Andersen, E., & Fenton, S. E. (2015). The mammary gland is a sensitive pubertal target in CD-1 and C57Bl/6 mice following perinatal perfluorooctanoic acid (PFOA) exposure. *Reproductive Toxicology*. <https://doi.org/10.1016/j.reprotox.2014.12.002>
14. United States Environmental Protection Agency. (2006). 2010/15 PFOA Stewardship Program. United States Environmental Protection Agency.
15. Wang, Z., Cousins, I. T., Scheringer, M., & Hungerbuehler, K. (2015). Hazard assessment of fluorinated alternatives to long-chain perfluoroalkyl acids (PFAAs) and their precursors: Status quo, ongoing challenges and possible solutions. *Environment International*. <https://doi.org/10.1016/j.envint.2014.11.013>
16. Zabaleta, I., Bizkarguenaga, E., Iparragirre, A., Navarro, P., Prieto, A., Fernández, L. Á., & Zuloaga, O. (2014). Focused ultrasound solid-liquid extraction for the determination of perfluorinated compounds in fish, vegetables and amended soil. *Journal of Chromatography A*. <https://doi.org/10.1016/j.chroma.2014.01.025>
17. Zafeiraki, E., Costopoulou, D., Vassiliadou, I., Bakeas, E., & Leondiadis, L. (2014). Determination of perfluorinated compounds (PFCs) in various foodstuff packaging materials used in the Greek market. *Chemosphere*. <https://doi.org/10.1016/j.chemosphere.2013.09.092>

For Research Use Only. Not for use in diagnostic procedures. © 2021 Thermo Fisher Scientific Inc. All rights reserved. Swingline and ClassicCut are trademarks of Swingline, a division of ACCO Brands, IKA is a trademark of IKA-Werke GmbH & Co, Alconox is a trademark of Alconox Inc., Frigidaire is a trademark of Electrolux AB, Misonix is a trademark of Misonix Inc., Pyrex is a trademark of Corning Inc., N-EVAP is a trademark of Organomation Associates, Inc., and JMP Pro is a trademark of SAS Institute Inc. All other 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 Inc. 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 representative for details. **AN65960-EN 0421M**

ThermoFisher
SCIENTIFIC