Food Testing & Agriculture



# A Study of Method Limit of Quantitation for 30 PFAS in Food

Using Captiva EMR PFAS food passthrough cleanup by LC/MS/MS detection

#### **Author**

Limian Zhao Agilent Technologies, Inc.

#### **Abstract**

This application note presents a study for the method limit of quantitation (LOQ) determination for 30 per- and polyfluoroalkyl substances (PFAS) targets in foods using a newly developed method of QuEChERS extraction followed by Agilent Captiva EMR PFAS Food cartridge passthrough cleanup, then LC/MS/MS detection. The LOQ $_{\rm spiking}$  was determined as the lowest validated spiking level in food matrices, with acceptable method accuracy and repeatability results. The LOQ $_{\rm cal}$  was calculated based on the standard deviation of LOQ $_{\rm spiking-A}$  following the procedure published in 40 CFR Appendix B to Part 136.¹ Both LOQ $_{\rm spiking-A}$  and LOQ $_{\rm cal-A}$  from this method were compared against the required LOQ $_{\rm AOACr}$  from the AOAC SMPR 2023.003 guideline.² The LOQ $_{\rm cal-A}$  results from this method were also compared to LOQ $_{\rm cal-F}$  from the U.S. Food and Drug Administration (FDA) method³, which uses the same calculation method to determine method LOQ $_{\rm cal-F}$  The results showed that the newly developed method provided acceptable LOQs meeting the AOAC SMPR guideline requirements. The new method LOQ $_{\rm cal-F}$  results were lower than the LOQ $_{\rm cal-F}$  from the FDA method for foods in the same category.

#### Introduction

Determination of PFAS residues in food is an emerging topic that has gained significant attention over the last several years. In April 2023, the European Commission enforced regulations for four PFAS compounds, namely perfluorooctane sulfonic acid (PFOS), perfluorooctanoic acid (PFOA), perfluorononanoic acid (PFNA), and perfluorohexane sulfonic acid (PFHxS) in eggs, fish, seafood, meat, and edible offal. <sup>4,5</sup> In November 2023, AOAC released the Standard Method Performance Requirements (SMPR) 2023-003 for the analysis of 30 PFAS in produce, beverages, dairy products, eggs, seafood, meat products, and feed. <sup>2</sup> The FDA has released their updated method (version 3.0) for the determination of 30 PFAS in food and feed. <sup>3</sup>

A new method, using QuEChERS extraction followed by Enhanced Matrix Removal (EMR) mixed-mode passthrough cleanup on Captiva EMR PFAS Food cartridges for sample preparation and LC/MS/MS detection for sample analysis, was developed and validated in multiple food matrices. Method validation was strictly implemented based on the AOAC SMPR guideline in 15 food matrices, including infant formula, milk, eggs<sup>6</sup>, baby food<sup>7</sup>, beef, tuna, shrimp<sup>8</sup>, carrots, grapes, mushrooms, lettuce, tomatoes, orange juice<sup>9</sup>, bovine kidney<sup>10</sup>, and dry soybeans<sup>11</sup>.

The most challenging LOQ requirements are on the four critical PFAS targets – PFHxS, PFOA, PFNA, and PFOS – where the required LOQs are 10x lower than those for other PFAS targets. Considering the various requirements on method LOQs for different PFAS targets in different food matrices by the AOAC SMPR guideline and EU regulation, multiple levels of quality control (QC) samples were prepared by pre-spiking the food matrix blank with PFAS standards at different concentrations. The method  $\text{LOQ}_{\text{spiking}}$  was decided based on the acceptance criteria through the actual spiking test. However, the method  $\text{LOQ}_{\text{spiking}}$  may not reflect the true method LOQ due to multiple contributions arising from the complexity of PFAS analysis in food.

In this work, a case study of method LOQs for PFAS analysis in food was implemented based on raw data generated in previous work. The calculation of method  $LOQ_{cal}$  and method detection limit (MDL) were based on the procedure published in 40 CFR Appendix B to Part 136. The calculated results were also compared to the AOAC SMPR 2023.003 guideline requirements and results reported by FDA method C-010.03 for the corresponding food category.

## **Experimental**

Method MDL and LOQ $_{cal-A}$  for 30 PFAS in food using this method were calculated using Equations 1 and 2 based upon data obtained from LOQ $_{\rm spiking}$  measurements in previous studies.  $^{6-11}$ 

#### Equation 1.

$$MDL_A = SD_{LOQspiking-A} \times 3.14$$

#### Equation 2.

$$LOQ_{Cal-A} = SD_{LOQspiking-A} \times 10$$

where

- MDL<sub>A</sub> is the method detection limit using the Aglient method
- LOQ<sub>cal-A</sub> is the method calculated LOQ using the Agilent method
- ${\rm SD_{LOQspiking-A}}$  is the standard deviation at the validated  ${\rm LOQ_{spiking-A}}$  level using the Agilent method

The determination of LOQ<sub>spiking-A</sub> was based on the lowest spiking QC level samples that met the AOAC SMPR guideline criteria on target recovery, repeatability, and selectivity.<sup>2</sup>

A total of 15 food matrices were included in this study, where three groups were classified based on the required LOQ level for the PFAS targets. These food matrices are listed in Table 1 with the AOAC SMPR guideline and EU regulation requirements on method LOQs.

In addition, the method  $LOQ_{cal-A}$  also needs to be within the established calibration range (Table 2). The method calibration range in food matrix (µg/kg) is determined by the neat standard calibration range (ng/L) corrected by the concentration or dilution factor introduced during sample preparation. For the calculated  $LOQ_{cal-A}$  below the lowest calibration range, the lowest calibration level was reported as  $LOQ_{cal-A}$ .

**Table 1.** Evaluation food matrices and their requirements on method LOQs based on AOAC SMPR guideline and EU regulation.

			Requirements on Meth	ug/kg)		
Group	Evaluated Food Matrix	Food Category	PFOS, PFOA, PFNA, and PFHxS	PFBA and PFPeA	Other PFAS	
	Baby food	Food for				
	Infant formula	infants and young children	≤ 0.01 for all (AOAC)			
	Carrot, grape,		≤ 0.015 for PFHxS (EU)	_	0.1	
1	mushroom,	Produce	≤ 0.005 for PFNA (EU)	≤ 1	≤ 0.1	
	lettuce, and tomato		≤ 0.01 for PFOA and PFOS (EU)			
	Orange juice	Beverage				
			≤ 0.01 for all (AOAC)			
			≤ 0.02 for PFOS (EU)			
I-A	Milk	Milk	≤ 0.01 for PFOA (EU)	≤ 1	≤ 0.1	
I-A			≤ 0.005 for PFNA (EU)			
			≤ 0.015 for PFHxS (EU)			
	Beef	Fish and meat				
II	Tuna	of terrestrial animals	≤ 0.1 for all (AOAC, EU)	≤1	≤1	
	Eggs	Eggs	≤ 0.3 for all (AOAC, EU)	< 3	≤ 3	
	Shrimp	Seafood	≥ 0.3 IOI all (AUAU, EU)	≥ 3	≥ 3	
III	Bovine kidney	Edible offal	≤ 0.4 for all (AOAC)	< <b>4</b>	< 4	
	boville klaney	Edible Offal	≤ 0.5 for all (EU)	≤ 4	≤ 4	
	Dry soybean	Feed	≤ 0.5 for all (AOAC)	≤ 5	≤ 5	

Table 2. Method calibration range for 30 PFAS in food.

	Concentration	Method Calibration Range (μg/kg)							
Evaluated Matrix	or Dilution Factor	PFBA	PFPeA	Other 28 Targets					
Carrot, Grape, Mushroom, Lettuce, Tomato, Orange Juice, and Baby Food	10x Concentration	0.01 to 5	0.005 to 2.5	0.001 to 1					
Infant Formula	5x Concentration	0.02 to 10	0.004 to 2	0.002 to 1					
Milk and Eggs	10x Concentration	0.02 to 10	0.004 to 2	0.002 to 1					
Beef, Tuna, Shrimp, and Dry Soybean	5x Concentration	0.04 to 20	0.008 to 4	0.004 to 2					
Bovine Kidney	10x Dilution	2 to 1, 000	0.4 to 200	0.2 to 100					

## **Results and discussion**

## ${\sf Method}\; {\sf LOQ}_{\sf spiking}$

The validated method  $LOQ_{spiking}$  in the representative food matrices shown in Table 3 was determined based on the lowest spiking level with acceptable criteria, including method accuracy (recovery), repeatability (RSD%), and selectivity. For food matrices with the positive detection of targets, the  $LOQ_{spiking}$  has to be higher than the calculated LOQ based on the target quantitation in replicated matrix blank samples using Equation 3 based on AOAC SMPR.<sup>4</sup>

#### Equation 3.

 $LOQ = 10 \times S_s$ 

where  $\mathbf{S}_{\mathrm{s}}$  is the sample standard deviation of the replicate matrix blank samples.

Figure 1 shows the method accuracy (recovery %) at  $LOQ_{spiking-A}$  in all tested food matrices, and Figure 2 shows the method repeatability (RSD%) at  $LOQ_{spiking-A}$  in all tested food matrices.

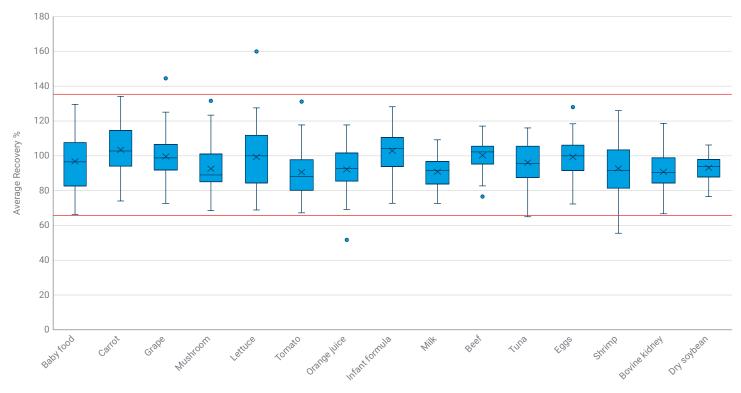
Results show that all the targets in all matrices meet the required LOQs with acceptable method accuracy, repeatability, and selectivity, except 4:2 FTS, 6:2 FTS, PFOS in carrot, PFOA in infant formula, 6:2 FTS in milk, and PFOS in beef. The significant positive detection of these targets in the specific food matrices resulted in a raised method LOQ<sub>spiking-A</sub>.

**Table 3.** The validated method  $LOQ_{spiking-A}$  for 30 PFAS targets in food.

	Method LOQ <sub>spiking-A</sub> in Food (μg/kg)														
Target	Baby Food	Carrot	Grape	Mushroom	Lettuce	Tomato	Orange Juice	Infant Formula	Milk	Beef	Tuna	Eggs	Shrimp	Bovine Kidney	Dry Soybean
PFBA	1.0	0.2	0.1	1.0	0.2	1.0	0.2	0.1	0.1	0.4	0.4	0.2	0.4	2.0	5.0
PFPeA	0.01	0.01	0.01	0.005	0.01	0.005	0.01	0.02	0.02	0.04	0.04	0.02	0.04	0.4	0.1
PFBS	0.1	0.01	0.01	0.004	0.002	0.01	0.004	0.01	0.01	0.02	0.02	0.01	0.02	0.2	0.05
4:2 FTS	0.002	0.2	0.1	0.001	0.1	0.1	0.004	0.01	0.01	0.02	0.02	0.01	0.02	0.4	0.05
PFHxA	0.004	0.004	0.01	0.004	0.02	0.02	0.02	0.02	0.02	0.02	0.02	0.02	0.02	0.4	0.05
PFPeS	0.002	0.02	0.01	0.004	0.01	0.002	0.002	0.01	0.01	0.02	0.02	0.01	0.02	0.4	0.05
HFPO-DA	0.001	0.004	0.004	0.004	0.1	0.004	0.002	0.01	0.01	0.02	0.02	0.01	0.02	0.2	0.05
PFHpA	0.002	0.01	0.004	0.004	0.002	0.002	0.004	0.01	0.01	0.02	0.02	0.01	0.02	0.2	0.05
PFHxS	0.002	0.01	0.01	0.004	0.01	0.004	0.004	0.01	0.01	0.02	0.02	0.01	0.02	0.2	0.05
DONA	0.001	0.1	0.002	0.002	0.002	0.002	0.001	0.01	0.1	0.02	0.02	0.01	0.02	0.2	0.5
6:2 FTS	0.002	0.2	0.004	0.01	0.1	0.004	0.02	0.01	0.2	0.02	0.02	0.01	0.02	0.4	0.1
PFOA	0.002	0.01	0.002	0.01	0.004	0.004	0.01	0.02	0.01	0.02	0.02	0.01	0.04	0.4	0.05
PFHpS	0.004	0.002	0.002	0.01	0.002	0.002	0.001	0.01	0.01	0.02	0.02	0.01	0.02	0.2	0.05
PFNA	0.002	0.004	0.002	0.01	0.002	0.004	0.004	0.01	0.01	0.4	0.02	0.01	0.1	0.4	0.05
PFOS	0.002	0.02	0.002	0.001	0.002	0.001	0.004	0.01	0.01	0.02	0.04	0.1	0.1	0.4	0.05
9CI-PF3ONS	0.002	0.001	0.002	0.001	0.002	0.001	0.001	0.01	0.01	0.02	0.02	0.01	0.02	0.2	0.05
8:2 FTS	0.001	0.002	0.001	0.002	0.001	0.002	0.001	0.01	0.01	0.02	0.02	0.01	0.04	0.4	0.05
PFDA	0.001	0.002	0.004	0.002	0.001	0.002	0.002	0.01	0.01	0.02	0.02	0.01	0.1	0.4	0.05
PFNS	0.001	0.002	0.002	0.01	0.001	0.004	0.001	0.01	0.01	0.02	0.02	0.01	0.02	0.2	0.05
PFDS	0.01	0.004	0.004	0.004	0.01	0.004	0.002	0.01	0.01	0.02	0.02	0.01	0.02	0.2	0.05
PFUnDA	0.002	0.004	0.002	0.01	0.002	0.002	0.001	0.01	0.01	0.02	0.04	0.01	0.4	0.4	0.05
PFOSA	0.004	0.02	0.004	0.002	0.1	0.004	0.002	0.01	0.01	0.02	0.02	0.01	0.02	0.2	0.05
11Cl-PF30UdS	0.004	0.002	0.001	0.001	0.001	0.001	0.02	0.01	0.01	0.02	0.02	0.01	0.02	0.2	0.05
PFUnDS	0.01	0.004	0.01	0.004	0.002	0.004	0.01	0.01	0.01	0.02	0.02	0.01	0.02	0.2	0.05
PFDoDA	0.004	0.004	0.002	0.01	0.002	0.004	0.001	0.01	0.01	0.02	0.04	0.01	0.04	0.4	0.05
10:2 FTS	0.002	0.002	0.004	0.002	0.004	0.001	0.001	0.01	0.01	0.02	0.02	0.01	0.02	1.0	0.05
PFDoS	0.01	0.02	0.002	0.004	0.004	0.002	0.01	0.01	0.01	0.02	0.1	0.01	0.02	0.2	0.05
PFTrDA	0.002	0.001	0.001	0.004	0.004	0.002	0.001	0.01	0.01	0.02	0.02	0.01	0.4	0.2	0.05
PFTrDS	0.02	0.01	0.004	0.01	0.02	0.02	0.01	0.01	0.01	0.02	0.4	0.01	0.02	0.2	0.05
PFTeDA	0.004	0.002	0.002	0.004	0.001	0.001	0.001	0.01	0.01	0.02	0.02	0.01	0.1	0.4	0.05

**Bold**: Highlights for four critical PFAS targets, i.e. PFHxS, PFOA, PFNA, and PFOS, and their results.

Red: Outlier due to significant positive detection of target in sample matrix blank.



 $\textbf{Figure 1.} \ \, \textbf{Method recovery at LOQ}_{\text{spiking-A}} \ \, \textbf{level for 30 PFAS targets in food (refer to Table 3 for LOQ}_{\text{spiking-A}} \ \, \textbf{level details)}.$ 

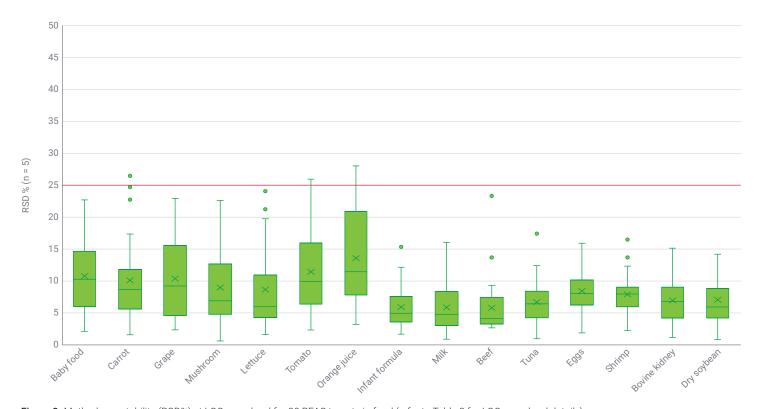


Figure 2. Method repeatability (RSD%) at LOQ<sub>spiking-A</sub> level for 30 PFAS targets in food (refer to Table 3 for LOQ<sub>spiking-A</sub> level details).

## Method MDL and LOQ<sub>cal</sub>

Agilent methods  $\mathrm{MDL_A}$  and  $\mathrm{LOQ}_{\mathrm{cal-A}}$  were determined based on validated method  $\mathrm{LOQ}_{\mathrm{spiking-A'}}$  using the aforementioned equations. Table 4 shows Agilent method  $\mathrm{MDL_A}$ , and Table 5 shows Agilent method  $\mathrm{LOQ}_{\mathrm{cal-A}}$ . The results were compared to those from FDA method C-010.03³ for targets in the representative food matrices from the same category, including lettuce, milk, eggs, fish, and feed. Since data for PFBA and PFPeA are not available for FDA method C-010.03, the comparison is limited to the remaining 28 PFAS targets. The results are shown in Figure 3.

The calculated LOQ $_{cal-A}$  were comparable to the LOQ $_{spiking-A}$  levels, with many of LOQ $_{cal-A}$  below the LOQ $_{spiking-A}$ . When the LOQ and was calculated below the lowest calibration level, the lowest calibration level was considered as method LOQ<sub>cal</sub> instead and differentiated in italics. Based on calculation, LOQ and for all targets in all matrices were shown to be below the required LOQ that met both the AOAC SMPR guideline and EU regulation for required LOQs. The only two exceptions, 4:2 FTS in carrot and PFNA in beef (highlighted in red), were caused by the significant positive detection of targets in the sample matrices. For four critical targets - PFHxS, PFOA, PFNA, and PFOS - the results are highlighted in bold. Comparing the targeted LOQ suggested in the EU regulation<sup>4,5</sup> for the four critical PFAS targets, the Agilent method LOQ sal-4 satisfied the EU-targeted LOQs in milk, which are  $\leq 0.01 \,\mu\text{g/kg}$  for PFOS and PFOA,  $\leq 0.02 \,\mu\text{g/kg}$  for PFNA, and ≤ 0.04 µg/kg for PFHxS. For these four targets, the ultralow EU-targeted LOQs in fruits, vegetables, and food for infants and young children are significantly challenging, i.e.,  $\leq 0.001 \,\mu g/kg$  for PFOA and PFNA,  $\leq 0.002 \,\mu g/kg$  for PFOS. and ≤ 0.004 µg/kg for PFHxS. The matrix background was the biggest barrier for the method to achieve these ultra-low LOQ levels in real food matrices. However, the Agilent method still met these ultra-low LOQs for the critical four targets in some food matrices with cleaner matrix background, which included 0.001  $\mu$ g/kg LOQ<sub>cal.</sub> for PFOA in grape and tomato, 0.001  $\mu g/kg$  LOQ<sub>cal-A</sub> for PFNA in grape, 0.002  $\mu g/kg$ (and below)  $\mathsf{LOQ}_{\mathsf{cal-A}}$  for PFOS in grape, mushroom, lettuce, and tomato, and 0.004 µg/kg (and below) LOQ and for PFHxS in baby food, mushroom, and lettuce. These results demonstrate that the Agilent method can reach ultra-low LOQ levels, given an acceptably clean food matrix.

The LOQ<sub>cal</sub> comparison between the Agilent method (LOQ<sub>cal-A</sub>) and FDA method C-010.03 (LOQ<sub>cal-F</sub>) is based on careful selection of food matrices, covering produce, milk, eggs, fish, and feed categories. For the produce category, lettuce was used in both method validations. For the eggs category, egg was used in both methods. Therefore, the comparison in these two matrices was guite straightforward. For the milk category, whole milk was used in the Agilent method validation, while chocolate milk was used in the FDA method. For the fish category, salmon was used in the FDA method, while canned tuna was used in the Agilent method. For the feed category, corn was used in the FDA method, while soybean was used in the Agilent method. Even with the slight differences in these food matrices, they are considered comparable sample matrices, and thus were also chosen for method comparison. Including these food matrices also expands the method comparison for more complicated and challenging food categories. The calculation of LOQ<sub>cal-A</sub> was based on validated  $\mathrm{LOQ}_{\mathrm{spiking-A'}}$  while  $\mathrm{LOQ}_{\mathrm{cal-F}}$  was based on results from a 0.05 µg/kg spiking level using the FDA method.3

The results summarized in Figure 3 show that Agilent method LOQ<sub>cal-A</sub> levels are overall below LOQ<sub>cal-E</sub> using the FDA C-010.03 method, with few exceptions due to the matrix positive detection impacts. The significantly lower  $LOQ_{spiking-A}$ is attributed to the much more simplified sample preparation method procedure, where improved matrix removal efficiency allows the collection of validated data at low LOQ sniking-A levels. The higher matrix removal efficiency also allows the use of larger sample sizes for extraction, results in a higher concentration factor, and reduces matrix effect for target determination. This also contributes to the lower LOQ and A using the Agilent method. Even with the equivalent spiking level for method LOQ $_{\text{spiking}}$  (0.05  $\mu$ g/kg) in the feed matrix comparison, the LOQ by the Agilent method still ranges lower than that of the FDA method. This confirms that the reduced matrix effect and larger sample size can improve method LOQs overall.

**Table 4.** Method  $MDL_A$ s for 30 PFAS targets in food.

	Method MDL <sub><sub>A</sub></sub> (μg/kg)														
Target	Baby Food	Carrot	Grape	Mushroom	Lettuce	Tomato	Orange Juice	Infant Formula	Milk	Beef	Tuna	Eggs	Shrimp	Bovine Kidney	Dry Soybean
PFBA	0.282	0.078	0.018	0.044	0.027	0.163	0.149	0.019	0.008	0.042	0.099	0.026	0.041	0.159	0.974
PFPeA	0.002	0.001	0.002	0.002	0.002	0.002	0.006	0.004	0.002	0.006	0.011	0.004	0.015	0.081	0.010
PFBS	0.011	0.002	0.004	0.001	0.001	0.003	0.002	0.004	0.001	0.003	0.005	0.001	0.005	0.047	0.010
4:2 FTS	0.001	0.065	0.014	0.001	0.008	0.013	0.002	0.002	0.001	0.006	0.006	0.001	0.005	0.070	0.007
PFHxA	0.001	0.001	0.003	0.002	0.002	0.012	0.002	0.005	0.002	0.005	0.009	0.001	0.004	0.124	0.015
PFPeS	0.001	0.001	0.003	0.002	0.001	0.002	0.002	0.001	0.001	0.004	0.008	0.001	0.005	0.086	0.007
HFPO-DA	0.001	0.005	0.002	0.002	0.047	0.002	0.002	0.001	0.001	0.007	0.005	0.003	0.003	0.088	0.015
PFHpA	0.001	0.005	0.002	0.001	0.001	0.001	0.001	0.002	0.002	0.005	0.005	0.002	0.010	0.031	0.021
PFHxS	0.001	0.003	0.003	0.001	0.001	0.002	0.003	0.002	0.002	0.005	0.004	0.001	0.006	0.024	0.007
DONA	0.001	0.016	0.001	0.001	0.001	0.001	0.001	0.001	0.022	0.004	0.005	0.001	0.005	0.018	0.065
6:2 FTS	0.002	0.011	0.001	0.001	0.014	0.001	0.002	0.007	0.006	0.005	0.004	0.002	0.006	0.111	0.029
PFOA	0.002	0.003	0.001	0.001	0.001	0.001	0.001	0.005	0.003	0.001	0.005	0.003	0.008	0.072	0.016
PFHpS	0.001	0.001	0.001	0.003	0.001	0.001	0.001	0.003	0.002	0.005	0.002	0.002	0.005	0.076	0.008
PFNA	0.001	0.002	0.001	0.002	0.003	0.002	0.001	0.002	0.002	0.059	0.005	0.003	0.031	0.113	0.015
PFOS	0.002	0.004	0.001	0.001	0.001	0.001	0.002	0.001	0.001	0.005	0.008	0.011	0.031	0.122	0.009
9CI-PF3ONS	0.001	0.001	0.001	0.001	0.001	0.001	0.001	0.001	0.001	0.006	0.006	0.001	0.005	0.036	0.007
8:2 FTS	0.001	0.001	0.001	0.001	0.001	0.001	0.001	0.002	0.001	0.003	0.005	0.003	0.009	0.103	0.004
PFDA	0.001	0.001	0.002	0.001	0.001	0.001	0.001	0.002	0.003	0.008	0.013	0.003	0.052	0.097	0.008
PFNS	0.001	0.001	0.002	0.001	0.002	0.001	0.001	0.003	0.001	0.002	0.005	0.001	0.005	0.058	0.004
PFDS	0.001	0.003	0.002	0.002	0.003	0.003	0.001	0.001	0.003	0.003	0.005	0.002	0.004	0.024	0.001
PFUnDA	0.001	0.001	0.001	0.001	0.001	0.001	0.001	0.001	0.001	0.008	0.008	0.001	0.007	0.178	0.020
PFOSA	0.003	0.012	0.002	0.001	0.019	0.001	0.001	0.004	0.004	0.003	0.005	0.005	0.008	0.024	0.008
11Cl-PF30UdS	0.001	0.001	0.001	0.001	0.001	0.001	0.014	0.001	0.002	0.004	0.005	0.001	0.004	0.053	0.003
PFUnDS	0.002	0.004	0.003	0.003	0.002	0.003	0.006	0.005	0.001	0.004	0.008	0.002	0.008	0.061	0.013
PFDoDA	0.001	0.002	0.001	0.001	0.001	0.001	0.001	0.003	0.003	0.004	0.011	0.003	0.012	0.115	0.014
10:2 FTS	0.001	0.001	0.003	0.001	0.001	0.001	0.001	0.001	0.001	0.003	0.005	0.001	0.006	0.066	0.008
PFDoS	0.006	0.005	0.001	0.003	0.002	0.001	0.005	0.002	0.003	0.003	0.005	0.001	0.004	0.071	0.021
PFTrDA	0.001	0.001	0.001	0.001	0.001	0.001	0.001	0.002	0.003	0.012	0.173	0.008	0.082	0.065	0.022
PFTrDS	0.006	0.007	0.003	0.01	0.015	0.005	0.007	0.004	0.003	0.007	0.005	0.001	0.007	0.030	0.010
PFTeDA	0.004	0.001	0.001	0.001	0.001	0.001	0.001	0.001	0.003	0.008	0.009	0.002	0.029	0.084	0.025

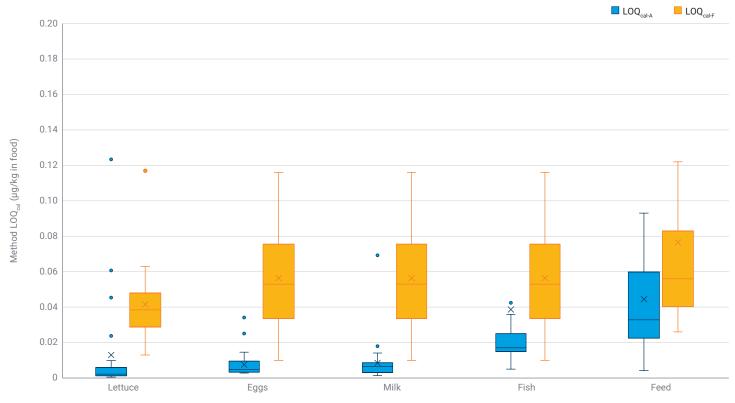
**Bold**: Highlights four critical PFAS targets, i.e. PFHxS, PFOA, PFNA, and PFOS, and their results.

**Table 5.** Method LOQ  $_{\mbox{\tiny cal-A}}$  for 30 PFAS targets in food.

		Method LOQ <sub>cal-A</sub> (µg/kg)														
Target	Baby Food	Carrot	Grape	Mushroom	Lettuce	Tomato	Orange Juice	Infant Formula	Milk	Beef	Tuna	Eggs	Shrimp	Bovine Kidney	Dry Soybean	
PFBA	0.899	0.247	0.056	0.139	0.085	0.519	0.474	0.061	0.025	0.134	0.316	0.082	0.130	0.506	3.103	
PFPeA	0.008	0.005	0.006	0.005	0.006	0.007	0.018	0.011	0.005	0.019	0.036	0.014	0.047	0.258	0.033	
PFBS	0.035	0.007	0.013	0.001	0.002	0.008	0.006	0.014	0.002	0.008	0.016	0.004	0.017	0.200	0.033	
4:2 FTS	0.003	0.208	0.046	0.001	0.024	0.041	0.005	0.006	0.003	0.019	0.019	0.003	0.017	0.223	0.021	
PFHxA	0.004	0.004	0.010	0.005	0.006	0.039	0.007	0.016	0.007	0.016	0.030	0.003	0.014	0.394	0.047	
PFPeS	0.002	0.017	0.010	0.006	0.004	0.005	0.007	0.004	0.002	0.014	0.024	0.005	0.016	0.272	0.022	
HFPO-DA	0.002	0.002	0.007	0.006	0.123	0.005	0.005	0.003	0.003	0.023	0.015	0.009	0.010	0.281	0.047	
PFHpA	0.002	0.015	0.005	0.004	0.001	0.003	0.003	0.006	0.005	0.014	0.017	0.006	0.033	0.200	0.067	
PFHxS	0.002	0.010	0.009	0.004	0.002	0.006	0.010	0.005	0.005	0.016	0.013	0.004	0.018	0.200	0.022	
DONA	0.001	0.049	0.001	0.001	0.001	0.002	0.001	0.004	0.069	0.012	0.016	0.004	0.014	0.200	0.208	
6:2 FTS	0.005	0.034	0.002	0.005	0.045	0.004	0.007	0.021	0.018	0.016	0.012	0.005	0.018	0.353	0.093	
PFOA	0.005	0.009	0.001	0.002	0.002	0.001	0.004	0.014	0.009	0.004	0.015	0.010	0.025	0.230	0.049	
PFHpS	0.004	0.004	0.003	0.009	0.002	0.001	0.001	0.008	0.007	0.016	0.005	0.007	0.016	0.243	0.025	
PFNA	0.002	0.006	0.001	0.006	0.005	0.007	0.003	0.007	0.007	0.189	0.017	0.008	0.100	0.359	0.047	
PFOS	0.006	0.014	0.002	0.002	0.002	0.001	0.006	0.003	0.003	0.014	0.025	0.034	0.100	0.388	0.029	
9CI-PF30NS	0.002	0.001	0.001	0.001	0.001	0.002	0.001	0.003	0.003	0.018	0.019	0.003	0.017	0.200	0.023	
8:2 FTS	0.002	0.001	0.002	0.003	0.002	0.004	0.002	0.006	0.004	0.009	0.017	0.010	0.030	0.327	0.011	
PFDA	0.002	0.005	0.006	0.004	0.001	0.003	0.002	0.006	0.004	0.024	0.042	0.010	0.167	0.309	0.026	
PFNS	0.002	0.003	0.005	0.003	0.005	0.003	0.002	0.009	0.003	0.006	0.017	0.004	0.015	0.200	0.014	
PFDS	0.002	0.008	0.006	0.005	0.010	0.010	0.003	0.004	0.008	0.009	0.016	0.005	0.013	0.200	0.004	
PFUnDA	0.002	0.001	0.001	0.004	0.003	0.002	0.001	0.002	0.004	0.025	0.025	0.003	0.023	0.566	0.063	
PFOSA	0.008	0.038	0.006	0.003	0.061	0.004	0.001	0.014	0.014	0.008	0.017	0.015	0.027	0.200	0.026	
11Cl-PF30UdS	0.004	0.001	0.002	0.001	0.001	0.001	0.045	0.004	0.006	0.014	0.015	0.004	0.013	0.200	0.011	
PFUnDS	0.007	0.012	0.012	0.008	0.005	0.010	0.019	0.015	0.003	0.012	0.026	0.005	0.025	0.200	0.043	
PFDoDA	0.005	0.005	0.005	0.004	0.002	0.002	0.002	0.010	0.009	0.013	0.036	0.010	0.039	0.365	0.045	
10:2 FTS	0.003	0.003	0.003	0.002	0.001	0.001	0.002	0.002	0.003	0.009	0.015	0.003	0.018	0.210	0.025	
PFDoS	0.018	0.015	0.015	0.009	0.005	0.001	0.016	0.008	0.010	0.011	0.015	0.003	0.013	0.225	0.066	
PFTrDA	0.002	0.001	0.001	0.003	0.001	0.002	0.001	0.006	0.009	0.039	0.551	0.025	0.262	0.205	0.069	
PFTrDS	0.019	0.021	0.021	0.031	0.047	0.015	0.021	0.013	0.008	0.024	0.017	0.003	0.022	0.200	0.033	
PFTeDA	0.012	0.002	0.002	0.003	0.001	0.002	0.001	0.003	0.008	0.026	0.030	0.007	0.092	0.268	0.080	

**Bold**: Highlights four critical PFAS targets, i.e. PFHxS, PFOA, PFNA, and PFOS, and their results.

*Italic blue*: Highlights for LOQ<sub>cal-A</sub> where the calibration range lowest level was used, when calculated LOQ<sub>cal</sub> was below the method calibration range. Red: Outlier due to significant positive detection of target in sample matrix blank.



**Figure 3.** Method  $LOQ_{cal}$  comparison of 28 PFAS determinations in representative foods;  $LOQ_{cal-A}$  using the Agilent method versus  $LOQ_{cal-F}$  using the FDA C-010.03 method.

## Conclusion

A case study on the method limits of quantitation was conducted for the validated Agilent method for 30 PFAS in food. Method  $\mathrm{LOQ}_{\mathrm{spiking-A}}$ ,  $\mathrm{MDL}_{\mathrm{A}}$ , and  $\mathrm{LOQ}_{\mathrm{cal-A}}$  were summarized for 30 PFAS targets in all 15 food matrices. Both method  $\mathrm{LOQ}_{\mathrm{spiking-A}}$  and  $\mathrm{LOQ}_{\mathrm{cal-A}}$  were compared to the LOQ levels required by the AOAC SMPR 2023.003 guideline and EU regulation; all met the accepted criteria, with few exceptions due to the significant positive detection of target in food matrices. Method  $\mathrm{LOQ}_{\mathrm{cal-A}}$  results were then compared between the Agilent and FDA C-010.03 methods, demonstrating a much lower method  $\mathrm{LOQ}_{\mathrm{cal}}$  achieved by the Agilent method, which is attributed to the simplified method, higher matrix removal efficiency, and reduced matrix effect.

#### References

- Definition and Procedure for the Determination of the Method Detection Limit, Revision 2. 40 CFR Appendix B to Part 136. Washington (DC). Definition and Procedure for the Determination of the Method Detection Limit, Revision 2; December 2016.
- AOAC (2023) Standard Method Performance Requirements (SMPRs) for Per- and Polyfluoroalkyl Substances (PFAS) in Produce, Beverages, Dairy Products, Eggs, Seafood, Meat Products, and Feed (AOAC SMPR 2023.003).
- 3. Determination of 30 Per- and Polyfluoroalkyl Substances (PFAS) in Food and Feed using Liquid Chromatography-Tandem Mass Spectrometry (LC-MS/MS), FDA Foods Program Compendium of Analytical Laboratory Methods: Chemical Analytical Manual (CAM) **2024**, C-010.03.

- EUR-Lex (2023) Consolidated text: Commission Regulation (EU) 2023/915 of 25 April 2023 on Maximum Levels for Certain Contaminants in Food and Repealing Regulation (EC) No 1881/2006.
- 5. Guidance Document on Analytical Parameters for the Determination of Per- and Polyfluoroalkyl Substances (PFAS) in Food and Feed, Version 1.2, May **2022**.
- 6. Zhao, L.; Giardina, M.; Parry, E. Agilent Technologies, Inc. Determination of 30 Per- and Polyfluoroalkyl Substances in Infant Formula, Milk, and Eggs, *Agilent Technologies application note*, publication number **5994-7366EN**, **2024**.
- Zhao, L.; Giardina, M.; Parry, E. Determination of 30 Perand Polyfluoroalkyl Substances in Baby Food, Agilent Technologies application note, publication number 5994-7367EN, 2024.
- 8. Zhao, L.; Giardina, M.; Parry, E. Determination of 30 Perand Polyfluoroalkyl Substances in Beef, Tuna, and Shrimp, *Agilent Technologies application note*, publication number **5994-7368EN**, **2024**.
- 9. Zhao, L.; Giardina, M. Determination of 30 Per- and Polyfluoroalkyl Substances in Fruits, Vegetables, and Juices, *Agilent Technologies application note*, publication number **5994-7369EN**, **2024**.
- Zhao, L.; Giardina, M.; Parry, E. Determination of 30 Per- and Polyfluoroalkyl Substances in Bovine Kidney, Agilent Technologies application note, publication number 5994-7370EN, 2024.
- 11. Zhao, L.; Giardina, M.; Parry, E. Determination of 30 Per- and Polyfluoroalkyl Substances in Dry Soybeans, *Agilent Technologies application note*, publication number 5994-7371EN. 2024.

#### www.agilent.com

DE-003306

This information is subject to change without notice.

