

## PFAS – Analysis Arium® Mini Plus Extend

Sample	Detection threshold	Detected Concentration	Unit	Method
PFBA	50	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFPeA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFHxA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFHpA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFOA branched	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFOA linear	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFOA total	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFNA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFDA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFUnDA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFDoDA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFTTrDA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFTeDA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFHxDA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFBS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFPeS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFHxS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFHpS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFOS linear	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFOS branched	0.6	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFOS total	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFNS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFDS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFUnDS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFDoDS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
PFTTrDS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
4:2 FTS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
6:2 FTS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
8:2 FTS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
10:2 FTS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>
N-MeFOSAA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:A</sup>

**Execution and Analysis Procedure**

The water analysis was executed by TÜV Rheinland Energy GmbH, an internationally recognized testing laboratory for special analytics, based on following measurement method: MS-0047387 Rev. 0, in accordance with DIN 38407-42, 2011-03. Relative expanded measurement uncertainty (k=2): 50 %.

<sup>A</sup>=The method has been accredited.

<sup>PV</sup>=The method has been partially validated.

The tests were performed with the Arium® Mini Plus Extend, without final filter, fed with tap water.

Sample	Detection threshold	Detected Concentration	Unit	Method
N-EtFOSAA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:PV</sup>
8:2diPAP	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:PV</sup>
PFECHS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 <sup>1:PV</sup>

#### Execution and Analysis Procedure

The water analysis was executed by TÜV Rheinland Energy GmbH, an internationally recognized testing laboratory for special analytics, based on following measurement method: MS-0047387 Rev. 0, in accordance with DIN 38407-42, 2011-03. Relative expanded measurement uncertainty (k=2): 50 %.

<sup>^</sup>=The method has been accredited.

<sup>PV</sup>=The method has been partially validated.

The tests were performed with the Arium® Mini Plus Extend, without final filter, fed with tap water.

#### Germany

Sartorius Lab Instruments  
GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Goettingen  
Phone +49 551 308 0

#### USA

Sartorius Corporation  
565 Johnson Avenue  
Bohemia, NY 11716  
Phone +1 631 254 4249  
Toll-free +1 800 635 2906

 For further information, visit  
[sartorius.com](https://www.sartorius.com)