

PFAS – Analysis Arium® Sterile Plus Filter

Sample	Detection threshold	Detected Concentration	Unit	Method
PFBA	50	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFPeA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFHxA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFHpA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFNA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFDA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFUnDA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFDoDA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFTTrDA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFTeDA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFHxDA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFBS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFPeS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFHxS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFHpS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFNS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFDS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFUnDS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFDoDS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFTTrDS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
4:2 FTS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
6:2 FTS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
8:2 FTS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
10:2 FTS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
N-MeFOSAA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
N-EtFOSAA	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:PV}
8:2diPAP	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:PV}
PFECHS	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:PV}

Execution and Analysis Procedure

The water analysis was executed by TÜV Rheinland Energy & Environment 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 %.

^A = The method has been accredited.

^{PV} = The method has been partially validated.

The tests were performed with the Arium® Sterile Plus Filter attached to an Arium® Mini Plus, fed with tap water. The Sterile Plus Filter was flushed with 5 liters of ultrapure water before the sample was collected.

Sample	Detection threshold	Detected Concentration	Unit	Method
PFOA (linear)	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1,A}
PFOA (branched)	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1,A}
PFOA (total)	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1,A}
PFOS (linear)	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1,A}
PFOS (branched)	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1,A}
PFOS (total)	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1,A}

Execution and Analysis Procedure

The water analysis was executed by TÜV Rheinland Energy & Environment 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 %.

^A = The method has been accredited.

^{PV} = The method has been partially validated.

The tests were performed with the Arium® Sterile Plus Filter attached to an Arium® Mini Plus, fed with tap water. The Sterile Plus Filter was flushed with 5 liters of ultrapure water before the sample was collected.

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