

PFAS – Analysis Arium® CellPlus Ultrafilter

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® CellPlus Ultrafilter attached to an Arium® Mini Essential, fed with tap water. The CellPlus Ultrafilter was flushed with 20 liters of ultrapure water before the sample was collected.

Sample	Detection threshold	Detected Concentration	Unit	Method
PFOA (lin)	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFOA (br)	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFOA (sum)	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFOS (lin)	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFOS (br)	0.5	Under detection threshold	ng/L (ppt)	MS-0047387 ^{1:A}
PFOS (sum)	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® CellPlus Ultrafilter attached to an Arium® Mini Essential, fed with tap water. The CellPlus Ultrafilter was flushed with 20 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