

PFAS – Analysis Arium® Comfort I

Sample	Detection threshold	Detected Concentration	Unit	Method
PFBA	5	Under detection threshold	ng/l (ppt)	MS-0047387
PFPeA	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFHxA	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFHpA	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFOA total	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFNA	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFDA	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFUnDA	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFDODA	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFTTrDA	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFTeDA	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFHxDA	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFBS	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFPeS	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFHxS	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFHpS	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFOS total	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFNS	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFDS	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFUnDS	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFDODS	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFTTrDS	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
4:2 FTS	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
6:2 FTS	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
8:2 FTS	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
10:2 FTS	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
N-MeFOSAA	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
N-EtFOSAA	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
8:2diPAP	0.5	Under detection threshold	ng/l (ppt)	MS-0047387
PFECHS	0.5	Under detection threshold	ng/l (ppt)	MS-0047387

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 %.

^A=The method has been accredited.

^{PV}=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 Göttingen
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)