

PFAS – Analysis Arium® Pro

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}
PFOA total	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}
PFOS total	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:PV}
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 GmbH, an internationally recognized testing laboratory for special analytics, based on following measurement method: MS-0047387. The method have been partially validated. The tests were performed with the Arium® Pro VF, without final filter, fed with DI water.

¹ MS-0047387 Rev. 0, following DIN 38407-42, 2011-03. Relative expanded uncertainty of measurement (k=2): 50%.

^A Accredited method

^{PV} Partially validated method

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)