

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 linear	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}
PFTrDA	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 linear	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}
PFTrDS	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

For further information, visit www.sartorius.com

USA

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