

PFAS – Analysis Arium® Comfort I

Sample	Detection threshold	Detected Concentration	Unit	Method
PFBA	5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFPeA	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFHxA	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFHpA	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFOA linear	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFOA total	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFNA	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFDA	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFUnDA	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFDoDA	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFTTrDA	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFTeDA	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFHxDA	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFBS	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFPeS	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFHxS	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFHpS	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFOS linear	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFOS total	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFNS	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFDS	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFUnDS	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFDoDS	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFTTrDS	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
4:2 FTS	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
6:2 FTS	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
8:2 FTS	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
10:2 FTS	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
N-MeFOSAA	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
N-EtFOSAA	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
8:2diPAP	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A
PFECHS	0.5	Under detection threshold	ng/l (ppt)	QMA-504-197 ^A

Execution and Analysis Procedure


The water analysis was executed by ProChem GmbH, an internationally recognized testing laboratory for special analytics, based on following measurement method: QMA-504-197^A. The method have been partially validated. The tests were performed with the Arium® Comfort I, 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
www.sartorius.com