

PFAS – Analysis Arium<sup>®</sup> Comfort II

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

**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<sup>A</sup>. The method have been partially validated. The tests were performed with the Arium<sup>®</sup> Comfort II, 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](http://www.sartorius.com)