

VOC – Analysis Arium® Mini Plus Extend

| Sample | Detection Threshold | Detected Concentration | Unit | Method |
|-------------------------------------|---------------------|---------------------------|------------|----------|
| 1, 1 Dichloropropene | 0.3 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 1 Dichlorethane | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 1 Dichlorethene | 0.2 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 1, 1, 2 Tetrachlorethane | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 1, 1, 2, 2 Tetrachlorethane | 0.2 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 1, 1, Trichlorethane | 0.25 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 1, 2 Trichlorethane | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 2 Dibromo-3-chloropropane (DBCP) | 0.5 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 2 Dibromoethane | 0.3 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 2 Dichlorethane | 0.4 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 2 Dichlorobenzene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 2 Dichloropropane | 0.5 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 2, 3 Trichlorobenzene | 0.05 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 2, 3 Trichloropropane | 0.2 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 2, 4 Trichlorobenzene | 0.05 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 2, 4 Trimethylbenzene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 3 Dichlorobenzene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 3 Dichloropropane | 0.2 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 3, 5 Trimethylbenzene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 1, 4 Dichlorobenzene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 2 Chlorotoluene | 0.2 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 2, 2 Dichloropropane | 0.4 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 4 Chlorotoluene | 0.2 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| 4 Isopropyltoluene | 0.05 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Benzene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Bromobenzene | 0.2 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Bromochloromethane | 0.4 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Bromodichloromethane | 0.2 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Bromoform | 0.5 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Bromomethane | 0.25 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Chlorobenzene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Chloroethane | 0.25 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Chloroform | 0.5 | Under detection threshold | µg/L (ppb) | HS-GC/MS |

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 methods: HS-GC/MS, GC/MS. The methods have been partially validated. The tests were performed with the Arium® Mini Plus Extend, without final filter, fed with tap water.

| Sample | Detection Threshold | Detected Concentration | Unit | Method |
|---------------------------|---------------------|---------------------------|------------|----------|
| cis-1,2-Dichloroethene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| cis-1,3-Dichloropropene | 0.3 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Dibromochloromethane | 0.5 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Dibromomethane | 0.25 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Dichloromethane | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Ethylbenzene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Hexachlorobutadiene | 0.005 | 0.010 | µg/L (ppb) | GC/MS |
| Isopropylbenzene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| m-Xylene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Naphthaline | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| n-Butylbenzene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| n-Propylbenzene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| o-Xylene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| p-Xylene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| sec-Butylbenzene | 0.05 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Styrene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| tert-Butylbenzene | 0.05 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Tetrachloroethene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Tetrachloromethane | 0.4 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Toluene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| trans-1,2-Dichloroethene | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| trans-1,3-Dichloropropene | 0.5 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Trichloroethene | 0.2 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Trichlorofluoromethane | 0.2 | Under detection threshold | µg/L (ppb) | HS-GC/MS |
| Vinyl chloride | 0.1 | Under detection threshold | µg/L (ppb) | HS-GC/MS |

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 methods: HS-GC/MS, GC/MS. The methods have 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 Goettingen
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)