

Trace Analysis Arium® Advance EDI Series

	Detection threshold	Unit	Calculated concentration Arium® Advance EDI	Procedure
Elements				
Antimony Sb	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Arsenic As	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Bismuth Bi	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Bromide Br ⁻	0.005	µg/L (ppb)	Under detection threshold	IC
Cadmium Cd	0.0005	µg/L (ppb)	Under detection threshold	ICP-MS
Cesium Cs	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Cerium Ce	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Chromium Cr	0.005	µg/L (ppb)	Under detection threshold	ICP-MS
Dysprosium Dy	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Erbium Er	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Europium Eu	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Gadolinium Gd	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Gallium Ga	0.005	µg/L (ppb)	Under detection threshold	ICP-MS
Germanium Ge	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Gold Au	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Hafnium Hf	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Holmium Ho	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Indium In	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Iridium Ir	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Lanthanum La	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Lead Pb	0.005	µg/L (ppb)	Under detection threshold	ICP-MS
Lutetium Lu	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Mercury Hg	0.005	µg/L (ppb)	Under detection threshold	ICP-MS
Molybdenum Mo	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Neodymium Nd	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Niobium Nb	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Osmium Os	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Palladium Pd	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Platinum Pt	0.002	µg/L (ppb)	Under detection threshold	ICP-MS
Praseodymium Pr	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Rhenium Re	0.001	µg/L (ppb)	Under detection threshold	ICP-MS

Execution and Analysis Procedure

The purified water analysis was executed by ATU GmbH, Analytics for Technology and Environment of the Arium® Advance EDI. ATU GmbH is an internationally recognized testing laboratory for the analysis of purified water. The following procedures were used for the analysis: "IC" (Dionex), "GF-AAS" (Dionex, Perkin Elmer), "ICP-MS 7500 cs" (Agilent) and "GC-MS." The purified water was taken directly from the product water outlet of the Arium® Advance EDI.

	Detection threshold	Unit	Calculated concentration Arium® Advance EDI	Procedure
Rhodium Rh	0.002	µg/L (ppb)	Under detection threshold	ICP-MS
Rubidium Rb	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Ruthenium Ru	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Samarium Sm	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Scandium Sc	0.002	µg/L (ppb)	Under detection threshold	ICP-MS
Selenium Se	0.01	µg/L (ppb)	Under detection threshold	ICP-MS
Silver Ag	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Tantalum Ta	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Tellurium Te	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Terbium Tb	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Thallium Tl	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Thorium Th	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Thulium Tm	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Tin Sn	0.002	µg/L (ppb)	Under detection threshold	ICP-MS
Tungsten W	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Uranium U	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Vanadium V	0.0005	µg/L (ppb)	Under detection threshold	ICP-MS
Ytterbium Yb	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Yttrium Y	0.001	µg/L (ppb)	Under detection threshold	ICP-MS
Zirconium Zr	0.002	µg/L (ppb)	Under detection threshold	ICP-MS
Anorganic				
Trimethylamine TMA	0.005	µg/L (ppb)	Under detection threshold	ICP-MS
Phosphate PO ₄ ³⁻	0.01	µg/L (ppb)	Under detection threshold	ICP-MS

Execution and Analysis Procedure


The purified water analysis was executed by ATU GmbH, Analytics for Technology and Environment of the Arium® Advance EDI. ATU GmbH is an internationally recognized testing laboratory for the analysis of purified water. The following procedures were used for the analysis: "IC" (Dionex), "GF-AAS" (Dionex, Perkin Elmer), "ICP-MS 7500 cs" (Agilent) and "GC-MS." The purified water was taken directly from the product water outlet of the Arium® Advance EDI.

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