

## Lab Water Applications

### Arium® Lab Water System

Lab Water Quality by System	Advanced RO	Advanced EDI	Mini	Mini UV	Mini Essential	Mini Essential UV	Mini Plus	Mini Plus UV	Pro	Pro DI	Pro UV	Pro UF	Pro VF	Comfort I	Comfort I UV	Comfort II	Comfort II UV
Type 1 Water			■	■	■	■	■	■	■	■	■	■	■	■	■	■	■
Type 2 Water		■														■	■
Type 3 Water	■							■	■					■	■		

Daily Water Consumption by System <sup>1</sup>	Advanced RO	Advanced EDI	Mini	Mini UV	Mini Essential	Mini Essential UV	Mini Plus	Mini Plus UV	Pro	Pro DI	Pro UV	Pro UF	Pro VF	Comfort I	Comfort I UV	Comfort II	Comfort II UV
Type 1 ultrapure water < 10 liter/day			■	■	■	■	■	■									
Type 1 ultrapure water 10 - 40 liter/day														■	■	■	■
Type 1 ultrapure water 40- 100 liter/day									■	■	■	■	■				
Type 2 pure water < 120 liter/day (5 l/h)		■														■	■
Type 2 pure water < 150 liter/day (10 l/h)		■														■	■
Type 3 pure water < 140 liter/day (8 l/h)	■													■	■		
Type 3 pure water < 200 liter/day (16 l/h)	■													■	■		
Type 3 pure water < 270 liter/day (24 l/h)	■													■	■		

### Lab Water Application Overview | System Requirements by Application

Feed Application	Advanced RO	Advanced EDI	Mini	Mini UV	Mini Essential	Mini Essential UV	Mini Plus	Mini Plus UV	Pro	Pro DI	Pro UV	Pro UF	Pro VF	Comfort I	Comfort I UV	Comfort II	Comfort II UV
Feed ultrapure water systems	■	■															
Feed distilled systems		■															
Water for Laboratory devices (Autoclaves   Washing Machine etc.)	■	■												■	■	■	■

General Laboratory & Analytical Application	Advanced RO	Advanced EDI	Mini	Mini UV	Mini Essential	Mini Essential UV	Mini Plus	Mini Plus UV	Pro	Pro DI	Pro UV	Pro UF	Pro VF	Comfort I	Comfort I UV	Comfort II	Comfort II UV
Buffer, media and pH solutions	■	■	■	■	■	■	■	■	■	■	■	■	■	■	■	■	■
Histology		■	■	■	■	■	■	■	■	■	■	■	■	■	■	■	■
ELISA (Enzyme-Linked Immunosorbent Assay)		■	■	■	■	■	■	■	■	■	■	■	■	■	■	■	■
AAS (Atomic Absorption Spectroscopy)		■	■	■	■	■	■	■	■	■	■	■	■	■	■	■	■
Solutions for chemical analysis and synthesis		■	■	■	■	■	■	■	■	■	■	■	■	■	■	■	■
GF-AAS (Graphite Furnace Atomic Absorption Spectrometry)			■	■	■	■	■	■	■	■	■	■	■	■	■	■	■
Preparation of reagents, blank samples, etc.			■	■	■	■	■	■	■	■	■	■	■	■	■	■	■
Photometry			■	■	■	■	■	■	■	■	■	■	■	■	■	■	■

Molecular Biology   Lifescience Application	Advanced RO	Advanced EDI	Mini <sup>2</sup>	Mini UV	Mini Essential <sup>2</sup>	Mini Essential UV	Mini Plus <sup>2</sup>	Mini Plus UV	Pro	Pro DI	Pro UV	Pro UF	Pro VF	Comfort I <sup>2</sup>	Comfort I UV	Comfort II <sup>2</sup>	Comfort II UV
Electrophoresis			■		■		■					■	■	■		■	
Northern Blot			■		■		■					■	■	■		■	
Southern Blot			■		■		■					■	■	■		■	
Western Blot			■		■		■					■	■	■		■	
Endotoxin analysis			■		■		■					■	■	■		■	
Immunocytochemistry			■		■		■					■	■	■		■	
Production of monoclonal antibodies			■		■		■					■	■	■		■	
PCR (Polymerase Chain Reaction)			■		■		■					■	■	■		■	
DNA Sequencing			■		■		■					■	■	■		■	
Nutrient media for cell culture (Mammalia & plant)			■		■		■					■	■	■		■	
Chromatography)			■		■		■					■	■	■		■	

Analytical Application	Advanced RO	Advanced EDI	Mini	Mini UV	Mini Essential	Mini Essential UV	Mini Plus	Mini Plus UV	Pro	Pro DI	Pro UV	Pro UF	Pro VF	Comfort I	Comfort I UV	Comfort II	Comfort II UV
SPE (Solid phase extraction)				■		■		■			■		■	■	■		■
Trace metal analysis				■		■		■			■		■	■	■		■
IC (Ion chromatography)				■		■		■			■		■	■	■		■
ICP-MS (Inductively Coupled Plasma Mass Spectrometry)				■		■		■			■		■	■	■		■
GC-MS (Gas Chromatography-Mass Spectrometry)				■		■		■			■		■	■	■		■
HPLC (High-Performance Liquid Chromatography)				■		■		■			■		■	■	■		■
TOC analysis				■		■		■			■		■	■	■		■

All displayed applicable systems starting with the minimal requested water quality criteria

1. Average indication depending on incoming feed water
2. Only in combination with an Arium® Cell Plus ultrafilter.