

Find the Right Arium® Mini for Your Laboratory

Feed Water	Arium® Mini	Arium® Mini Plus	Arium® Mini Essential	Arium® Mini Plus Extend	Arium® Mini Essential Extend
Tap Water		✓		✓	
Pretreated water (RO DI line)			✓		✓
Pretreated water (from supply container)	✓				
Flexible Removable Dispenser				✓	✓
Bagtank Technology	✓	✓		✓	
Purified Water					
Type 1	✓	✓	✓	✓	✓
Type 3		✓		✓	