

Arium® Bag

The Most Innovative Bagtank System

- Fast and easy replacement of the Arium® Bag
- High user safety as the Bagtank eliminates the need for cleaning chemicals

Description

Pure water is stored inside the laboratory water system, which reliably protects pretreated pure water from secondary contamination.

Sartorius Bagtank technology enables consistent water quality over a prolonged period, ensuring continuously reproducible results.

Unlike conventional water tanks, the Arium® Bag ensures high user safety and saves time as it eliminates the need for a complicated cleaning procedure with chemicals.



Technical Specifications | Ordering Information

Materials	
Bag	S71 proprietary film
Tubing	TuFlux®

Bag dimensions* [H × W]	
5-Liter bag	40 × 33 cm (15.7" × 12.9")
20-Liter bag	86.5 × 43.0 cm (34.1" × 16.9")
50-Liter bag	90.0 × 58.1 cm (35.4" × 22.9")

*When empty

Order Number	Description
H2O-CBS-5-S	Arium® 5-Liter bag for Arium® Mini and Mini Plus; qty. per pkg.: 1
H2O-CBS-20	Arium® 5-Liter bag for Arium® Bagtank 20; qty. per pkg.: 2
H2O-CBS-50	Arium® 50-Liter bag for Arium® Bagtank 50 and 100; qty. per pkg.: 2

Intended Use

Device type:

- Arium® Mini and Mini Plus
- Arium® Bagtank