

Arium® Mini Plus Pretreatment Cartridge



Reliable Protection for the Pretreatment of Feed Water

- Fast and effective adsorption of impurities by high-grade activated carbon
- Highly efficient catalyzer for removing oxidating agents such as chlorine
- Highly efficient reverse osmosis membranes; optimized water usage
- Low-energy membranes for ecological and economic operation

Description

Efficient purification is performed by a combination of activated carbon, a catalyzer and a downstream reverse osmosis membrane.

The spherical, catalytic activated carbon and an additional catalyst reliably remove oxidants, such as chlorine and ozone, heavy metal ions and particulate contaminants, from feed water.

In addition, due to the downstream reverse osmosis membrane, up to 98% of all salts, as well as bacteria and particles, are retained.

Technical Specifications | Ordering Information

Materials	
Housing	High-grade polypropylene
Filter media	Spherical, catalytic activated, carbon
Dimensions [W × H × D]	18 × 26 × 11 cm (7 × 10.2 × 4.3")
Operating weight	3.5 kg (7.7 lbs.)
Feed water requirements	See Technical Specifications on page 3

Order Number	Description
H2O-CPR	Arium® Mini Plus pretreatment cartridge; qty. per pkg.: 1

Intended Use

- Device type:
- Arium® Mini Plus