

# Arium® Softener Cartridge

## For maximum service life of the EDI module

- Consistently high water quality
- Long service life
- Effective CaCO<sub>3</sub> elimination

## Description

It is sensible to soften the feed water to improve protection of the EDI module.

The cartridge reliably removes traces of alkaline earth ions from the water, thereby guaranteeing consistently high water quality and a long service life of the EDI module.



## Technical Specifications | Ordering Information

| Materials        |                              |
|------------------|------------------------------|
| Housing          | High-purity polypropylene    |
| Filling material | Ultrapure ion exchange resin |

| Order number | Description                     |
|--------------|---------------------------------|
| H2O-CSO-1    | Arium® Softener Cartridge, 1 pc |

### Intended Use

Device type:

- Arium® Advance EDI
- Arium® Comfort II