

Arium® Cleaning Solution

Effective cleaning for a long lifetime

- Highly effective cleaning
 - Free of TOC (organic components)
- Surfactant-free
- Environmentally safe and gentle on the materials

Description

With this cleaning agent, the regular removal of impurities that develop during water purification is easy and effective, this results in higher flow rates and longer service lives.

This highly effective agent is already filled in 50 ml syringes and ready for direct connection.



Technical Specifications | Ordering Information

Ingredients

Sodium hypochlorite

Order number

611CDS1

Description

Arium® Cleaning Solution,
50 ml cleaning solution
filled into syringes, 1 pc

Intended Use

Device type:

- Arium® Pro DI, Pro UF, Pro UV and Pro VF