

Arium® UV Lamp (185 | 254 nm)

Ultrapure water, free of TOC

- Horizontal installation, optimized temperature gradient
- Effectively destroys organic compounds
- Easy replacement

Description

The horizontally arranged UV lamp provides particularly reliable results. In contrast to vertical devices, the temperature gradient is less pronounced and thus prevents the influence on the activity of UV waves.

The two different wavelengths reliably remove organic substances up to a TOC (total organic carbon) content of ≤ 5 ppb*.



Technical Specifications | Ordering Information

Materials

TOC value for product water* ≤ 5 ppb

* Depends on the type of system and on the feed water

Order number

H2O-CEL1

Description

Arium® UV Lamp (185 | 254 nm), 1 pc

Intended Use

Device type:

- Arium® Mini
- Arium® Mini Essential
- Arium® Mini Plus