

# 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  $\leq 2$  ppb\*.



## Technical Specifications | Ordering Information

---

### Materials

---

TOC value for product water\*  $\leq 2$  ppb

---

\* Feed water < 50 ppb TOC

---

### Order number

---

611CEL1

---

### Description

---

Arium® UV Lamp (185 | 254 nm), 1 pc

---

### Intended Use

Device type:

- Arium® Comfort (UV & TOC version)
- Arium® Pro UV & VF (also TOC version)