## S/RICH SILS

## Arium® Mini

## Compact Ultrapure Water Systems for 10L per Day

- Constant low conductivity for reliable results
- Integrated Bagtank technology for best economic water storage
- Easy-to-use touch display for intuitively menu navigation
- Space-saving unit with a width of only 28 cm (11 inches)
- Highest quality standards made in Germany

## Simplifying Progress

