

## arium® Pressure Tanks

Product Water Storage Tanks for Reverse Osmosis System 61316 and Type 2 Water System 61215



### Description

The arium® pressure tanks are designed to store water produced by the arium® 61316 reverse osmosis system and the 61215 Type 2 water system.

Each of these vessels is a completely enclosed pressure tank incorporating an internal diaphragm, which expands and contracts relative to the preset tank pressure and the volume of water in storage. An FDA-approved double butyl diaphragm isolates the product water within the pressure tank, ensuring water purity is maintained during storage.

Several benefits are gained by the use of a closed pressure tank:

- Contamination by airborne bacteria and gases is reduced.
- Tanks can be mounted in any position for delivery of water to optimize space and storage location.
- The pressurized reservoir is capable of delivering RO water directly to ultrapure water systems and other laboratory equipment without the need of an additional pump.

### Specifications

#### Materials of Construction

|            |  |
|------------|--|
| Exterior:  | Two part polyurethane   epoxy primed paint |
| Reservoir: | Patented double butyl diaphragm            |
| Plumbing:  | PE and PP                                  |

#### Operating Pressure

|                     |   |
|---------------------|---|
| Tank:               | 0–10 bar   0–146 psi max.                 |
| Standard precharge: | 0–2.6 bar   0–38 psi<br>± 0.3 bar   5 psi |

#### Operating Temperature

5°C–90°C (41°F–200°F)  
Not to exceed 60°C (140°F) ambient

#### Connectors

|         |  |
|---------|--|
| Inlet:  | 6 mm (1/4") John Guest   |
| Outlet: | 9 mm (3/8") quick connect,<br>6 mm (1/4") NPT for optional pressure switch or outlet |

#### Pressure Gauge

0–700 kPa | 0–100 psi

#### Shutoff Valves

Manual ball valve and quick-disconnect shutoff valve included

#### Optional Accessories

Pressure switch, order no. 611APS1, to connect pressure tank to arium® 611 systems

#### Certification

The arium® tanks meet CE and NSF [58] standards

## Laboratory Water Systems

| <b>Order Number</b> | <b>Capacity<br/>@ 2.5 bar</b> | <b>Height</b> | <b>Diameter</b> | <b>Weight<br/>Empty</b> | <b>Weight<br/>Full</b> |
|---------------------|-------------------------------|---------------|-----------------|-------------------------|------------------------|
| 613APV31            | 30 litres                     | 57 cm         | 41 cm           | 11 kg                   | 41 kg                  |
|                     | 8 gallons                     | 23"           | 16"             | 24 lbs                  | 90 lbs                 |
| 613APV70            | 70 litres                     | 111 cm        | 41 cm           | 21 kg                   | 91 kg                  |
|                     | 19 gallons                    | 44"           | 16"             | 46 lbs                  | 201 lbs                |
| 613APV100           | 100 litres                    | 94 cm         | 53 cm           | 26 kg                   | 126 kg                 |
|                     | 26 gallons                    | 37"           | 21"             | 57 lbs                  | 278 lbs                |

Sartorius AG  
Weender Landstrasse 94–108  
37075 Goettingen, Germany  
Phone +49.551.308.0  
Fax +49.551.308.3289  
[www.sartorius.com/arium](http://www.sartorius.com/arium)

USA +1.631.2544249  
UK +44.1372.737100  
France +33.1.69192100  
Italy +39.055.505671  
Spain +34.91.3588566  
Japan +81.3.33295533

Specifications subject to change without notice.  
Printed in Germany on paper that has been  
bleached without any use of chlorine.  
W/sart-254a - G  
Publication No.: SLG2032-e06094  
Order No.: 85030-517-67