



## arium<sup>®</sup> Dispenser Gun

The solution for greater flexibility in the laboratory.



The arium dispenser gun is an ergonomic, easy-to-handle dispensing unit which, thanks to its different designs, can be used for dispensing both pure and ultrapure water. The variants are available as bench top units with a stand or as a wall-mounted system depending on space requirements.

Thanks to the extended tube guide, the work area for dispensing pure or ultrapure water is extended to up to 2.5 meters. Furthermore, the height-adjustable stand enables optimal adjustment to different sampling chamber sizes. (The stand is also optionally available for the display | dispensing unit.)

A sterile Sartopore<sup>®</sup> 2 150 membrane final filter capsule with a 0.2 µm membrane (validated as a sterile filter according to HIMA and ASTM F-838-83 guidelines) guarantees the sterile dispensing of pure and ultrapure water (EU > 1 CFU/1000 ml).



### Areas of Application:

- Dispensing ultrapure water
- Dispensing pure water\*

\* In conjunction with an arium 613 APV pressure tank

### Features

- Suitable for pure and ultrapure water dispensing
- Sterile water dispensing through a 0.2 µm membrane final filter capsule
- Available with a stand or wall mounting bracket
- Ergonomically-designed dispenser gun
- Height-adjustable stand
- Range up to 2.5 meters
- Easy-to-use
- Light weight

### Specifications

#### Material

Stand	Aluminum (gray anodized)
Dispenser gun	Plastic (white finish)
Tubing	PE   PVDF

#### Dimensions (W × H × D)

Dispenser gun with stand (without tubing)	185 × 595 × 510 mm
Dispenser gun with wall mounting bracket (without tubing)	90 × 100 × 285 mm

#### Weight

Dispenser gun with stand (without tubing)	5.60 kg
Dispenser gun with wall mounting bracket (without tubing)	0.46 kg

#### Feed Water Requirements

Input Pressure	0 – 3 bar
Temperature	5 – 35 °C

#### Product Water Quality

Particles*	< 1/ml
Bacteria*	< 1 CFU/1000 ml

\* When using a Sartopore<sup>®</sup> 2 150 membrane final filter capsule

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen, Germany  
Phone +49.551.308.0  
Fax +49.551.308.3289  
www.sartorius-stedim.com  
  
USA Toll-Free +1.800.368.7178  
UK +44.1372.737159  
France +33.442.845600  
Italy +39.055.63.40.41  
Spain +34.90.2110935  
Japan +81.3.3740.5407

Specifications subject to change without notice. Printed and copyrighted by Sartorius Stedim Biotech GmbH. | · G  
Publication No.: SLG2047-e11051  
Order No.: 85032-540-13  
Ver. 05 | 2011

Order Number	Description
H2Opro-AMDG1	Dispenser gun including height-adjustable stand
H2Opro-AMDG2	Dispenser gun including wall mounting kit
613-AMDG1	Dispenser gun including height-adjustable stand for 613APV pressure tank
613-AMDG2	Dispenser gun including wall-mounting kit for 613APV pressure tank
5441307H4--CE--B	Sartopore <sup>®</sup> 2 150 sterile final filter, 0.2 µm pore size (5 units)