

# Certificate of Conformity

## **Cartridge Kit H2O-CRO-H-2, H2O-CRO-H-1, 613CPM4 and 613CPM4-----V for arium® advance, comfort, 613 & 612 laboratory water systems**


This document certifies that the designated product was manufactured according to Sartorius Weighing Technology acceptance criteria.

The manufacturer hereby declares that the products below are manufactured to meet his published specifications for these products. The manufacturer also declares that the product H2O-CRO-H-2, H2O-CRO-H-1, 613CPM4 and 613CPM4-----V have been manufactured from approved materials of construction and comply with design and manufacturing documentation.

When used unmodified and according to good manufacturing practices, all H2O-CRO-H-2, H2O-CRO-H-1, 613CPM4 and 613CPM4-----V membranes will comply with the U.S. Food, Drug and Cosmetic Act as amended under Food Additive Regulation 21 CFR 177-2550. The product may be used in processing liquid food and to purify water for food manufacturing compliance with the FDA 21 CFR 177-2550 regulatory citation. In use, the products should be maintained in a sanitary manner in accordance with good manufacturing practice. The product mentioned should be rinsed prior to first use according to the corresponding user manual of the related water purification device.

01.01.2013

Date



---

Dr. Susan Gerlghausen  
Director of Quality

Sartorius Lab Instruments GmbH & Co. KG  
Weender Landstrasse 94-108  
37075 Goettingen, Germany

[www.sartorius.com](http://www.sartorius.com)