

# Certificate of Conformity

## Material Origin Certificate **arium<sup>®</sup> mini H2O-CPR Pre-Treatment Cartridge**

This product has passed Sartorius Stedim Biotech in-house tests and thus meets Sartorius Stedim Biotech stringent quality control standards.

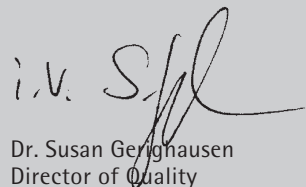
We hereby declare that this product are manufactured to meet the published specifications for these products. The manufacturer also declares that the product have been manufactured from approved materials of construction and comply with design and manufacturing documentation.

Cartridge housings are constructed of 100% virgin polypropylene.

Furthermore, we declare, to the best of his knowledge, that the product does not contain any hazardous material as supplied.

June 2015

Date



Dr. Susan Gerigshausen  
Director of Quality