

# Certificate of Approval

This is to certify that the Management System of:

## Promega GmbH Euro Hub

Pfingstweidstr. 23, 68199 Mannheim, Germany

has been approved by LRQA to the following standards:

ISO 13485:2016



P.G. Cornelissen - Area Operations Manager - Western and Central Europe

Issued By: Lloyd's Register Deutschland GmbH

for and on behalf of: Lloyd's Register Quality Assurance Limited

Current Issue Date: 5 December 2017

Expiry Date: 13 January 2020

Certificate Issue Number: 10054676

Original Approvals:

ISO 13485 14 January 2011

Approval Certificate Number: ISO 13485 – 0020402

The scope of this approval is applicable to:

Warehousing and distribution of reagents for basic and applied research for life science.



001