

# 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 Lloyd's Register to the following standards:

ISO 13485:2016



P.G. Cornelissen - Area Manager North Europe

Issued by: Lloyd's Register Nederland B.V.

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

Current issue date: 14 January 2020  
Expiry date: 13 January 2023  
Certificate identity number: 10233689

Original approval(s):  
ISO 13485 – 14 January 2011

Approval number(s): 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