

# Certificate of Approval

This is to certify that the Management System of:

## Promega GmbH

Schildkrötstr. 15, 68199 Mannheim, Germany

has been approved by LRQA to the following standards:

ISO 9001:2015



P.G. Cornelissen - Area Manager North Europe

Issued by: Lloyd's Register Deutschland GmbH

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

Current issue date: 21 January 2019  
Expiry date: 20 January 2022  
Certificate identity number: 10153531

Original approval(s):  
ISO 9001 – 21 January 2004

Approval number(s): ISO 9001 – 0018332

The scope of this approval is applicable to:

Distribution and service of reagents and (instrument-) system solutions for basic and applied research with life sciences.



001