

Certificate of Approval

This is to certify that the Management System of:

Promega GmbH

Gutenbergring 1, 69190 Walldorf, Germany

has been approved by Lloyd's Register to the following standards:

ISO 9001:2015

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.



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



001