

Certificate of Approval

This is to certify that the Management System of:

Promega GmbH

Gutenbergring 10, 69190 Walldorf, Germany

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

ISO 13485:2016

Approval number(s): ISO 13485 – 0025822

The scope of this approval is applicable to:

Distribution and service of reagents and (instrument-) system solutions for basic and applied research within life sciences and In Vitro Diagnostics.



Paul Graaf

Area Operations Manager North Europe

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

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



001