18 December 2023 17 December 2026 10563341

Original approval(s): ISO 13485 - 18 December 2020

LRQA

## **Certificate of Approval**

This is to certify that the Management System of:

## Promega GmbH

Gutenbergring 10, 69190 Walldorf, Germany

has been approved by LRQA 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, Europe

Issued by: LRQA Limited



LRQA Group Limited, its affiliates and subsidiaries and their respective officers, employees or agents are, individually and collectively, referred to in this clause as 'LRQA'. LRQA assumes no responsibility and shall not be liable to any person for any loss, damage or expense caused by reliance on the information or advice in this document or howsoever provided, unless that person has signed a contract with the relevant LRQA entity for the provision of this information or advice and in that case any responsibility or liability is exclusively on the terms and conditions set out in that contract. Issued by: LRQA Limited, 1 Trinity Park, Bickenhill Lane, Birmingham B37 7ES, United Kingdom