

Certificate of Approval

This is to certify that the Management System of:

Promega GmbH Euro Hub

Gutenbergring 10, 69190 Walldorf, Germany

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

ISO 13485:2016

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.



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