

Certificate of Approval

This is to certify that the Management System of:

Promega GmbH Euro Hub

Gutenbergring 10, 69190 Walldorf, Germany

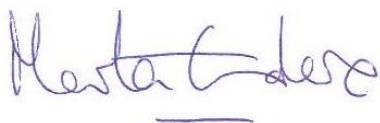
has been approved by LRQA to the following standards:

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.



Marta Escudero

Regional Director, Europe

Issued by: LRQA Limited



0001