

Certificate



This is to confirm that the quality management system and its application, as employed by

Certificate number: EBN.ISO.QMS.4592
Issue date: 21-03-2024
Valid from: 21-03-2024
Valid until: 21-03-2027

Promega Benelux B.V.

located at Leiden

meets with the requirements of the quality management system

ISO 9001:2015

This certificate is valid for the following scope of activities:

Distribution of molecular and cellular biology reagents and related instrumentation for life science research

Ing. C.A.N. van Zwienen,
General director

