

Certificate of Approval

This is to certify that the Management System of:

Promega Benelux B.V.

Schipholweg 1 (6e verdieping), 2316 XB Leiden, Netherlands

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

ISO 9001:2015

Approval number(s): ISO 9001 – 00023197

The scope of this approval is applicable to:

Supplier of Biological Reagents and Instruments for life science Research.



P.G. Cornelissen

Area Manager North Europe

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

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



001