

---

# Certificate of Approval

This is to certify that the Management System of:

## Promega Benelux B.V.

Schipholweg 103 (3e verdieping), 2316 XC Leiden, The Netherlands

has been approved by LRQA 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 Scientific and Medical Research.



**Paul Graaf**

---

Area Operations Manager, Europe

Issued by: LRQA Limited

