



Lloyd's Register  
LRQA

## CERTIFICATE OF APPROVAL

This is to certify that the Quality Management System of:

**Promega Benelux B.V.**  
**Schipholweg 1 (6e verdieping)**  
**2316 XB Leiden**  
**The Netherlands**

has been approved by Lloyd's Register Quality Assurance  
to the following Quality Management System Standard:

**ISO 9001 : 2015**

The Quality Management System is applicable to:

**The supply of biological reagents and  
instruments for scientific research, human identification,  
molecular diagnostics and quality control.**

Approval Certificate No: RQA653404	Original Approval	:	15 January 1999
	Current Certificate	:	1 January 2016
	Certificate Expiry	:	31 December 2019

Issued by: Lloyd's Register Nederland B.V. for and on behalf of  
Lloyd's Register Quality Assurance Limited



001

K.P. van der Mandelelaan 41a, 3062 MB Rotterdam, Nederland

For and on behalf of 1 Trinity Park, Bickenhill Lane, Birmingham, B37 7ES, United Kingdom

This approval is carried out in accordance with the LRQA assessment and certification procedures and monitored by LRQA

The use of the UKAS Accreditation Mark indicates Accreditation in respect of those activities covered by the Accreditation Certificate Number 001.