



## CERTIFICATE OF APPROVAL

This is to certify that the Quality Management System of:

**Promega Benelux B.V.  
Leiden, The Netherlands**

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

**ISO 9001 : 2008**

The Quality Management System is applicable to:

**Supplier of Biological Reagents and Instruments  
for Scientific and Medical Research.**

Approval  
Certificate No: 653404

Original Approval: 15 January 1999  
Current Certificate: 1 January 2008  
Certificate Expiry: 31 December 2010

Issued by: LRQA (Rotterdam)



This document is subject to the provision on the reverse  
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  
LRQA 10-12-1