



CERTIFICATE OF APPROVAL

This is to certify that the Quality Management System of:

Promega GmbH
Schildkrötstr. 15, 68199 Mannheim
Germany

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:

**Distribution and service of reagents and (instrument-)
system solutions for basic and applied research
within life sciences.**

Approval
Certificate No: KLN 4000095

Original Approval: 21 January 2004

Current Certificate: 21 January 2010

Certificate Expiry: 20 January 2013

Issued by: Lloyd's Register Quality Assurance GmbH



001

This document is subject to the provision on the reverse.

Innere Kanalstr. 15, 50823 Germany, Deutschland, Handelsregister Nr. B 34587

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

Macro Revision 13