

# Certificate of Registration



This is to certify that the Quality Management System of:

**Promega UK Limited**

Delta House, Enterprise Road, Southampton Science Park, Chilworth, Southampton, SO16 7NS, United Kingdom

applicable to:

**The stockholding and distribution of molecular and cellular biology reagents and related instrumentation for life science research. Provision of technical support for all supplied products and instrument service and repair**

has been assessed and registered by NQA against the provisions of:

**BS EN ISO 9001:2015**

This registration is subject to the company maintaining a quality management system, to the above standard, which will be monitored by NQA

Managing Director

Certificate No.	9668
ISO Approval Date	22 June 1998
Reissued	13 April 2017
Valid Until:	12 April 2020
EAC Code:	29



015