

# Certificate of Registration



This is to certify that the Quality Management System of:

**Promega UK Limited**

2 Benham Road, Southampton Science Park, Chilworth, Hampshire, SO16 7QJ, 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:

**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



0015

Certificate No.	9668
ISO Approval Date:	22 June 1998
Reissued:	6 March 2026
Valid Until:	11 April 2029
EAC Code:	29, 31, 29, 35