

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



A handwritten signature in black ink, appearing to read 'N. Wray', is positioned above the title 'Managing Director'.

Managing Director



015

Certificate No.	9668
ISO Approval Date:	22 June 1998
Reissued:	3 April 2020
Valid Until:	12 April 2023
EAC Code:	29