

Certificate of Registration



This is to certify that the Occupational Health & Safety Management System of:

Promega UK Limited

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

applicable to:

All activities associated with our stockholding and distribution of molecular and cellular biology reagents and related instrumentation for the life sciences sector. The use of a life sciences laboratory for teaching and demonstration purposes

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

ISO 45001:2018

This registration is subject to the company maintaining an occupational health & safety management system, to the above standard, which will be monitored by NQA.



Managing Director

| | |
|--------------------|-------------|
| Certificate No. | 187387 |
| ISO Approval Date: | 15 May 2023 |
| Valid Until: | 14 May 2026 |
| EAC Code: | 29, 31 |



0015