

Certificate of Registration

This is to certify that the Management System of:

Promega UK Ltd
2 Benham Road, Southampton Science Park, Southampton,
SO167QJ

has been approved by Alcumus ISOQAR and is compliant
with the requirements of:

ISO 45001: 2018

SCOPE OF REGISTRATION

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.

CERTIFICATE NUMBER: 23544-OHS-001

Initial Registration Date: 15/05/2023
Re-issue Date: 29/05/2025
Current Expiry Date: 14/05/2026

SIGNED



Jim Anderson, Chief Executive Officer
(on behalf of Alcumus ISOQAR)

This certificate will remain current subject to the company maintaining its system to the required standard. This will be monitored regularly by Alcumus ISOQAR. Further clarification regarding the scope of this certificate and the applicability of the relevant standards' requirement may be obtained by consulting Alcumus ISOQAR.

Alcumus ISOQAR Limited, Cobra Court, 1 Blackmore Road, Stretford, Manchester M32 0QY

T: 0161 865 3699 **E:** isoqarenquires@alcumus.com **W:** isoqar.com

This certificate is the property of Alcumus ISOQAR and be returned of request.

