



ISOQAR



Certificate of Registration

This is to certify that the Management System of:

Promega UK Ltd

**2 Benham Road, Southampton Science Park, Southampton, SO16
7QJ**

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

ISO 14001: 2015

SCOPE OF REGISTRATION

All activities associated with our stockholding and managed 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-EMS-001

SIGNED

Initial Registration Date: 05/02/2024
Previous Expiry Date: 05/02/2027
Recertification Audit Date: 02/04/2026
Re-issue Date: 17/04/2026
Current Expiry Date: 30/04/2029

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

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

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 ISOQAR and be returned of request.

