



# 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 45001: 2018

### SCOPE OF REGISTRATION

All activities associated with the provision of stockholding and managed distribution of molecular and cellular reagents and related instrumentation, including servicing & maintenance, for the life sciences sector. The use of a life sciences laboratory for teaching and demonstration purposes.

**CERTIFICATE NUMBER: 23544-OHS-001**

**SIGNED**

**Initial Registration Date:** 15/05/2023  
**Previous Expiry Date:** 14/05/2026  
**Recertification Audit Date:** 05/02/2026  
**Re-issue Date:** 26/02/2026  
**Current Expiry Date:** 14/05/2029

**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.

