

# 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 14001: 2015



**Certificate Number:** 23544-EMS-001  
**Initial Registration Date:** 05/02/2024  
**Expiry Date:** 05/02/2027

## 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 life science laboratory for teaching and demonstrating purpose.

Signed:  
Alyn Franklin, 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:** [isoqarenquiries@alcumus.com](mailto:isoqarenquiries@alcumus.com) **W:** [alcumus.com/isoqar](http://alcumus.com/isoqar)

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