



# ENVIRONMENTAL MANAGEMENT SYSTEM CERTIFICATE

Certificate No.: ZHJT25E11513R0S

Promega (Beijing) Biotech Co., Ltd

Unified social credit code: 91110101717856885T

The management system is up to standard

GB/T24001-2016/ISO14001:2015

The certificate is valid to the following product(s)/service

Sales and related management activities of specialty chemical products (biological reagents) (excluding hazardous chemicals), instruments, and electronic products.

Registration Address: Global Trade Center B907-909, No. 36 Eastern Road, 3rd North Ring Road, Dong Cheng District, Beijing

Audit Address: Global Trade Center B907-909, No. 36 Eastern Road, 3rd North Ring Road, Dong Cheng District, Beijing

Date of Issue: December 09th, 2025

Issued By:

Valid until: December 08th, 2028

**Beijing Zhongheng Zhengtong Certification Co., Ltd.**

Company address: Room 602, unit 2, floor 6, building 11, Chaoyang North Road, Chaoyang District, Beijing

The effectiveness of the Certificate is subject to QR Code in the lower left corner.

Meanwhile, you can search the website of certification body: [www.bjzhzt.com](http://www.bjzhzt.com)

or search the CNCA website: [www.cnca.gov.cn](http://www.cnca.gov.cn)

