

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 9001:2015

This is to certify that: Shanghai Promega Biological Products Ltd.
Building 12B
No. 88 Xinjun Ring Road
Caohejing HiTech Park of Pujiang
Minhang
Shanghai
201114
China

上海普洛麦格生物产品有限公司
91310000607216836N
中国
上海
闵行区
漕河泾开发区浦江高科技园
新骏环路88号
12B
邮编: 201114

Holds Certificate No: **FM 603098**

and operates a Quality Management System which complies with the requirements of ISO 9001:2015 for the following scope:

The design and development, production and relevant activities for S1 Nuclease, M-MLV reverse Transcriptase H⁻, RNasin Ribonuclease Inhibitor, Taq DNA Polymerase, Pfu DNA Polymerase.

S1核酸酶、M-MLV反转录酶H⁻、核糖核酸酶抑制剂、Taq DNA聚合酶、Pfu DNA聚合酶的设计开发、生产及相关活动。

For and on behalf of BSI:

Chris Cheung, Head of Compliance & Risk - Asia Pacific

Original Registration Date: 2010-11-01

Latest Revision Date: 2017-11-10

Effective Date: 2016-10-31

Expiry Date: 2019-10-30

Page: 1 of 1



...making excellence a habit.™

This certificate was issued electronically and remains the property of BSI and is bound by the conditions of contract.

An electronic certificate can be authenticated [online](#).

Printed copies can be validated at www.bsi-global.com/ClientDirectory or telephone +86 10 8507 3000, as well as at www.cnca.gov.cn, the official website of CNCA – Certification and Accreditation Administration of the People's Republic of China.

Further clarifications regarding the scope of this certificate and the applicability of ISO 9001:2015 requirements may be obtained by consulting the organization. The certified organization shall be subject to surveillance audit periodically with acceptable results for maintaining the validity of this certificate. This certificate is valid only if provided original copies are in complete set.