



Certificate of Registration

OUALITY MANAGEMENT SYSTEM - ISO 13485:2016

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: MD 603096

and operates a Quality Management System which complies with the requirements of ISO 13485:2016 for the following scope:

The design & manufacture of CT/ NG/ UU DNA qPCR Detection Kit, CT DNA qPCR Detection Kit, HR-HPV (14 types) DNA qPCR detection Kit, HSV 1 and 2 DNA qPCR Detection Kit. 沙眼衣原体/淋球菌/解脲脲原体核酸检测试剂盒、沙眼衣原体核酸检测试剂盒、高危型人乳头瘤病毒(14个型)核酸检测试剂盒、 单纯疱疹病毒I和II型核酸分型检测试剂盒的研发和生产。

For and on behalf of BSI:

Graeme Tunbridge, Senior Vice President Medical Devices

Original Registration Date: 2013-11-25

Latest Revision Date: 2022-11-09

bsi.



Effective Date: 2022-11-25 Expiry Date: 2025-11-24

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.