

# Certificate of Registration

## QUALITY MANAGEMENT SYSTEM - ISO 13485: 2003

*This is to certify that:*

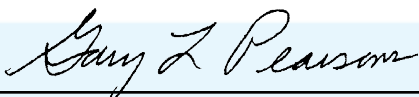
**Promega Corporation**  
**2800 Woods Hollow Road**  
**Madison**  
**Wisconsin**  
**53711-5399**  
**USA**

*Holds Certificate No:* **FM 505517**

*and operates a Quality Management System which complies with the requirements of ISO 13485: 2003 for the following scope:*

The design and manufacture of products, service and information applicable to the delivery of reagents, reagent support systems, instrumentation and medical devices for the biotechnology, DNA identification and in vitro diagnostic markets.

*For and on behalf of BSI:*



*President, BSI Management Systems (Americas)*

Originally registered: **08/24/2006**

Latest Issue: **08/24/2006**

Expiry Date: **08/23/2009**



Page: 1 of 2

This certificate remains the property of BSI, Inc. and shall be returned immediately upon the request.  
An electronic certificate can be authenticated [online](#). Printed copies can be validated at [www.bsi-global.com/ClientDirectory](http://www.bsi-global.com/ClientDirectory)  
To be read in conjunction with the scope above or the attached appendix.  
Americas Headquarters: 12110 Sunset Hills Road, Suite 200, Reston, VA 20190, USA.  
Group Headquarters: 389 Chiswick High Road, London, W4 4AL, UK.



Certificate No: **FM 505517**

**Location**

**Registered Activities**

Promega Corporation  
2800 Woods Hollow Road  
Madison  
Wisconsin  
53711-5399  
USA

The design and manufacture of products, service and information applicable to the delivery of reagents, reagent support systems, instrumentation and medical devices for the biotechnology, DNA identification and in vitro diagnostic markets.

Originally registered: **08/24/2006**

Latest Issue: **08/24/2006**

Expiry Date: **08/23/2009**

Page: 2 of 2

This certificate remains the property of BSI, Inc. and shall be returned immediately upon the request.  
An electronic certificate can be authenticated [online](http://www.bsi-global.com/ClientDirectory). Printed copies can be validated at [www.bsi-global.com/ClientDirectory](http://www.bsi-global.com/ClientDirectory)  
To be read in conjunction with the scope above or the attached appendix.  
Americas Headquarters: 12110 Sunset Hills Road, Suite 200, Reston, VA 20190, USA.

Group Headquarters: 389 Chiswick High Road, London, W4 4AL, UK.