



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 America, Inc.

Originally Registered: **08/24/2006**

Latest Issue: **07/28/2009**

Expiry Date: **08/23/2012**



Page: 1 of 1

This certificate remains the property of BSI and shall be returned immediately upon request. An electronic certificate can be authenticated [online](http://www.bsigroup.com/ClientDirectory). Printed copies can be validated at www.bsigroup.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.

