


Ion PGM™ Controls Kit v2

Catalog Number 4482010

Pub. No. MAN0013492 Rev. A.0

Contents	Cap color	Quantity	Volume	Shipping and Storage
E. coli DH10B Genomic DNA Control (100 ng/μL)	Red	1 tube	30 μL	-30°C to -10°C
Ion Sphere™ Test Fragments (150,000/μL)	Green	1 tube	18 μL	
E. coli DH10B Control 200 Library	Blue	1 tube	12 μL	
E. coli DH10B Control 400 Library	Orange	1 tube	16 μL	

 **WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from www.lifetechnologies.com/support.

Description

The Ion PGM™ Controls Kit v2 (Cat. no. 4482010) includes the following components:

- E. coli DH10B Genomic DNA Control (100 ng/μL) sufficient for preparing 3 libraries (1 μg control DNA per library)
- Ion Sphere™ Test Fragments sufficient for 3 sequencing runs (750,000 ISPs per chip)
- E. coli DH10B Control 200 Library sufficient for 10+ template preparation reactions (1 μL per reaction)
- E. coli DH10B Control 400 Library sufficient for 10+ template preparation reactions (1 μL per reaction)

Store the kit components as directed in the table.

The Ion PGM™ Controls Kit v2 is used in conjunction with the following kits:

- Ion Xpress™ Plus Fragment Library Kit (Cat. no. 4471269)
- Ion Plus Fragment Library Kit (Cat. no. 4471252)
- Ion PGM™ Template OT2 200 Kit (Cat. no. 4480974)
- Ion PGM™ Template OT2 400 Kit (Cat. no. 4479878)
- Ion PGM™ IC 200 Kit (Cat. no. 4484080)
- Ion PGM™ Hi-Q™ Chef Kit (Cat. no. A25948)
- Ion PGM™ Sequencing 200 Kit (Cat. no. 4474004)
- Ion PGM™ Sequencing 400 Kit (Cat. no. 4482002)
- Ion PGM™ Hi-Q™ Sequencing Kit (Cat. no. A25592)

See the appropriate user guide for your kit, available on the Ion Community at <http://ioncommunity.lifetechnologies.com/welcome>, for more information.

Precautions

Handle biological samples and all other materials according to local laboratory guidelines and regulations. Avoid cross-contamination of samples and reagents to ensure quality results.

If equipment or reagents are used in a manner not specified by the manufacturer, the protection provided by the equipment may be impaired.

Note: Refer to the expiration date on the package labels. Do not use kits after the expiration date.

Customer and technical support

Visit www.lifetechnologies.com/support for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
 - Product FAQs
 - Software, patches, and updates
- Order and web support
- Product documentation, including:
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

The information in this guide is subject to change without notice.

DISCLAIMER

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

Important Licensing Information: This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2015 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For support visit lifetechnologies.com/support or email techsupport@lifetech.com

lifetechnologies.com

16 February 2015

