

Ion S5™ Controls Kit Plus

Catalog Number A30729

Pub. No. MAN0016206 Rev. A.0

 **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 thermofisher.com/support.

Product description

The Ion S5™ Controls Kit Plus includes the following components:

- Human CEPH Genomic DNA Control (100 ng/μL) sufficient for preparing 3 libraries (1 μg control DNA per library)
- Human CEPH 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 at 1 μL per reaction, or 3 template preparation reactions at 5 μL per reaction

The Ion S5™ Controls Kit Plus can be used with the following kits:

- Ion 520™ & Ion 530™ Kit – Chef
- Ion 540™ Kit – Chef
- Ion 520™ & Ion 530™ Kit – OT2
- Ion 540™ Kit – OT2
- Ion 520™ & Ion 530™ ExT Kit – Chef

Kit contents and storage

Contents	Amount	Storage
Human CEPH Genomic DNA Control (red cap)	30 μL	-30°C to -10°C
Human CEPH Control 200 Library (yellow cap)	12 μL	
<i>E. coli</i> DH10B Control 400 Library (orange cap)	16 μL	

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. Store kit components as directed.

IMPORTANT! Do not substitute components in this kit with components from other Ion Torrent™ controls kits. This protocol has been verified using these specific materials. Substitution may adversely affect system performance.

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

Customer and technical support

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

- Worldwide contact telephone numbers
- Product support, including:
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- 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.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

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

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

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.

Corporate entity: Life Technologies Corporation | Carlsbad, CA 92008 USA | Toll Free in USA 1 800 955 6288

©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For Research Use Only. Not for use in diagnostic procedures.