


Ion 510™ Chip Kit

Catalog Numbers A34292

Pub. No. MAN0016856 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](https://www.thermofisher.com/support).

Description

The Ion 510™ Chip Kit contains barcoded chips for sample tracking and sequencing runs using the Ion S5™ and Ion S5™ XL Sequencer. During sequencing, the Ion 510™ Chip detects polymerase-driven base incorporation and translates this information into digital form.

The Ion 510™ Chip offers:

- eFuse technology providing sample tracking functionality.
- Compatibility with all current library preparation methods.

Note: This product is compatible with the Ion Chef™ Instrument and Ion S5™ or Ion S5™ XL Sequencer workflows.

Contents and storage

Contents	Amount	Storage
Ion 510™ Chip	2 x 4 chips	15°C–30°C

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.

IMPORTANT! Do not substitute components in this kit with components from other Ion Torrent™ chip 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](https://www.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.



Manufacturer: Life Technologies Corporation | 5781 Van Allen Way | Carlsbad, CA 92008

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.

Revision history: Pub. No. MAN0016856

Revision	Date	Description
A.0	12 June 2017	New document

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

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.