

# Ion 540™ Kit – Chef

Catalog Number A27759

Pub. No. MAN0010848 Rev. B.0

**Note:** For safety and biohazard guidelines, see the “Safety” appendix in the following product documentation: *Ion 540™ Kit - Chef User Guide* (Pub. no. MAN0010851). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

## Description

The Ion 540™ Kit – Chef contains 8 boxes (Cat. No. [A34461](#)) or 10 boxes (Cat. No. [A34019](#)) total (see “Kit summary” on page 1 for details).

The Ion 540™ Kit – Chef contains template preparation reagents sufficient for 8 reactions for automated template preparation and chip loading using the Ion Chef™ System, and reagents for sequencing using the Ion S5™ or Ion S5™ XL System. Store the kit components as directed. For more information, see the *Ion 540™ Kit – Chef User Guide* (Pub. No. MAN0010851).

## Kit summary

| Component                    | Part No. | Quantity per kit |
|------------------------------|----------|------------------|
| Ion S5™ Chef Supplies        | A27755   | 4 boxes          |
| Ion S5™ Chef Solutions       | A27754   | 1 box            |
| Ion 540™ Chef Reagents       | A27758   | 1 box            |
| Ion S5™ Sequencing Solutions | A27767   | 1 box            |
| Ion S5™ Sequencing Reagents  | A27768   | 1 box            |

## Ion Chef™ reagents and materials

| Contents  | Amount / box | Storage      |
|---|--------------|--------------|
| <b>Ion S5™ Chef Supplies (Part No. A27755)</b>  |              |              |
| Chip Adapter                                    | 2            | 15°C to 30°C |
| Enrichment Cartridge v2                         | 1            |              |
| Tip Cartridge v2                                | 1            |              |
| PCR Plate                                       | 1            |              |
| PCR Plate Frame                                 | 1            |              |
| Frame Seal v2                                   | 1            |              |
| Recovery Station Disposable Lid v2              | 2            |              |
| Recovery Tube v2                                | 12           |              |
| <b>Ion S5™ Chef Solutions (Part No. A27754)</b> |              |              |
| Ion S5™ Chef Solutions                          | 4 cartridges | 15°C to 30°C |

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

| Contents  | Amount / box | Storage        |
|---|--------------|----------------|
| <b>Ion 540™ Chef Reagents (Part No. A27758)</b> |              |                |
| Ion 540™ Chef Reagents                          | 4 cartridges | -30°C to -10°C |

## Ion S5™ sequencing reagents and materials

| Contents  | Amount / box | Storage                       |
|---|--------------|-------------------------------|
| <b>Ion S5™ Sequencing Solutions (Part No. A27767)</b> |              |                               |
| W2 Solution   | 4 × 1.5 L    | 15°C to 30°C                  |
| Ion S5™ Cleaning Solution                             | 250 mL       |                               |
| <b>Ion S5™ Sequencing Reagents (Part No. A27768)</b>  |              |                               |
| Ion S5™ Sequencing Reagents                           | 4 cartridges | -30°C to -10°C <sup>[1]</sup> |

<sup>[1]</sup> Cartridges ship at 2°C to 8°C. Store as indicated, do not store on dry ice.

**IMPORTANT!** Do not store the Ion S5™ Sequencing Reagents (Part No. A27768) on dry ice or in a closed environment containing dry ice.

## 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™ template or sequencing 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.

## 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 at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://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](http://www.thermofisher.com/support).

## Customer and technical support

Visit [thermofisher.com/support](http://thermofisher.com/support) for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
  - Product FAQs
  - Software, patches, and updates
  - Training for many applications and instruments
- Order and web support
- Product documentation
  - 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.



**Revision history:** Pub. No. MAN0010848 B.0

| Revision | Date           | Description   |
|----------|----------------|---|
| B.0      | 11 August 2023 | Updated Ion S5™ Chef Supplies (Part No. A27755) kit to include the PCR Plate Frame. |
| A.0      | 12 August 2015 | New product information sheet for Ion 540™ Kit – Chef.                              |

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

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. 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:** 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.

©2015-2023 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.