

Ion 510™ & Ion 520™ & Ion 530™ Kit – Chef

Catalog Numbers A34461, A34019

Pub. No. MAN0016955 Rev. C.0

Note: For safety and biohazard guidelines, see the “Safety” appendix in the following product documentation: *Ion 510™ & Ion 520™ & Ion 530™ Kit – Chef User Guide* (Pub. no. MAN0016854). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Description

The Ion 510™ & Ion 520™ & Ion 530™ 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 510™ & Ion 520™ & Ion 530™ 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 510™ & Ion 520™ & Ion 530™ Kit – Chef User Guide* (Pub. No. MAN0016854).

Kit summary

Component	Part No.	Quantity per kit	
		A34461	A34019
Ion S5™ Chef Supplies	A27755	4 boxes	4 boxes
Ion S5™ Chef Solutions	A27754	1 box	1 box
Ion 510™ & Ion 520™ & Ion 530™ Chef Reagents	A34018	1 box	1 box
Ion S5™ Sequencing Solutions	A27767	1 box	2 boxes
Ion S5™ Sequencing Reagents	A27768	1 box	2 boxes

Ion Chef™ reagents and materials

Contents	Amount / box	Storage
Ion S5™ Chef Supplies (Part No. A27755)		
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	
Ion S5™ Chef Solutions (Part No. A27754)		
Ion S5™ Chef Solutions	4 cartridges	15°C to 30°C

Contents	Amount / box	Storage
Ion 510™ & Ion 520™ & Ion 530™ Chef Reagents (Part No. A34018)		
Ion 510™ & Ion 520™ & Ion 530™ Chef Reagents	4 cartridges	–30°C to –10°C

Ion S5™ sequencing reagents and materials

Contents	Amount / box	Storage
Ion S5™ Sequencing Solutions (Part No. A27767)		
Ion S5™ Wash Solution	4 × 1.5 L	15°C to 30°C
Ion S5™ Cleaning Solution	250 mL	
Ion S5™ Sequencing Reagents (Part No. A27768)		
Ion S5™ Sequencing Reagents	4 cartridges	–30°C to –10°C ^[1]

^[1] 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.

Customer and technical support

Visit [thermofisher.com/support](https://www.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.

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. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Revision history: Pub. No. MAN0016955 C.0

Revision	Date	Description
C.0	11 August 2023	Updated Ion S5™ Chef Supplies (Part No. A27755) kit to include the PCR Plate Frame.
B.0	22 August 2017	Corrected the number of boxes listed in the kit summary.
A.0	12 June 2017	New product information sheet for Ion 510™ & Ion 520™ & Ion 530™ 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.

©2017-2023 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.