

Ion 520™ & 530™ Kit – Chef

Catalog Number A27757

Pub. No. MAN0010847 Rev. A.0

Ion S5™ Chef Supplies (Part no. A27755; 4 boxes supplied per kit)		
Component	Quantity per box	Shipping and storage
Chip Adapter	2 adapters	15°C to 30°C
Enrichment Cartridge v2	1 cartridge	
Tip Cartridge v2	1 cartridge	
PCR Plate and Frame Seal v2	1 each	
Recovery Station Disposable Lid v2	2 lids	
Recovery Tube v2	12 tubes	

Ion S5™ Chef Solutions (Part no. A27754)		
Component	Quantity per kit	Shipping and storage
Ion S5™ Chef Solutions	4 cartridges	15°C to 30°C

Ion 520™ & 530™ Chef Reagents (Part no. A27756)		
Component	Quantity per kit	Shipping and storage
Ion 520™ & 530™ Chef Reagents	4 cartridges	-30°C to -10°C

Ion S5™ Sequencing Solutions (Part no. A27767; 2 boxes supplied per kit)			
Component	Quantity per box	Volume	Shipping and storage
Ion S5™ Wash Solution	4 bottles	1.5 L	15°C to 30°C
Ion S5™ Cleaning Solution	1 bottle	250 mL	

Ion S5™ Sequencing Reagents (Part no. A27768; 2 boxes supplied per kit)		
Component	Quantity per box	Shipping and storage
Ion S5™ Sequencing Reagents	4 cartridges	-30°C to -10°C

Note: For safety and biohazard guidelines, refer to the “Safety” appendix in the *Ion 520™ & 530™ Kit – Chef User Guide* (Pub. no. MAN0010846). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Description

The Ion 520™ & 530™ Kit – Chef (Cat. no. A27757) contains 10 boxes total:

- Ion S5™ Chef Supplies (Part no. A27755; 4 boxes per kit)
- Ion S5™ Chef Solutions (Part no. A27754)
- Ion 520™ & 530™ Chef Reagents (Part no. A27756)
- Ion S5™ Sequencing Solutions (Part no. A27767; 2 boxes per kit)
- Ion S5™ Sequencing Reagents (Part no. A27768; 2 boxes per kit)

The Ion 520™ & 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, refer to the *Ion 520™ & 530™ Kit – Chef User Guide* (Pub. no. MAN0010846).

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 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 found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

Customer and technical support

Visit thermofisher.com/techresources 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.

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: 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 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 thermofisher.com/techresources or email techsupport@lifetech.com

thermofisher.com/lifescience