

Ion S5™ and Ion S5™ XL Instrument

Catalog Numbers A27212, A27214

Pub. No. MAN0010842 Rev. C.0

Ion S5™ System (Catalog No. A27212)			
Components			Part No.
Ion S5™ Sequencer			A27211
Ion S5™ Installation Kit			A27215
Ion S5™ XL System (Catalog No. A27214)			
Contents			Part No.
Ion S5™ XL Sequencer			A27213
Ion Torrent™ Server			A28563
Ion S5™ Installation Kit			A27215
Ion S5™ Installation Kit (Part No. A27215; Not available for separate purchase.)			
Contents	Part No.	Quantity	Storage
Ion 540™ Chip Kit	A27765	4 pack	15°C to 30°C
Ion S5™ Sequencing Solutions	A27767	1	
Ion S5™ Sequencing Reagents	A27768	1	-30°C to -10°C
Ion S5™ Cartridge Tool	A28308	1	15°C to 30°C

Note: For safety and biohazard guidelines, see the “Safety” appendix in the *Ion S5™ and Ion S5™ XL Instrument User Guide* (Pub. No. MAN0010811). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Description

Each Ion S5™ and Ion S5™ XL System ships with an Ion S5™ Installation Kit (Part No. A27215) containing all the reagents and materials necessary to install an Ion S5™ or Ion S5™ XL Sequencer.

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

Customer and technical support

Visit 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.



Life Technologies Holdings Pte Ltd | Block 33 | Marsiling Industrial Estate Road 3 | #07-06, Singapore 739256

Revision history: Pub. No. MAN0010842

Revision	Date	Description
C.0	1 February 2023	Removed the following skus from the Ion S5™ Installation Kit contents table. <ul style="list-style-type: none">• A28195• A27897• A29022
B.0	21 February 2022	Update the quantity of the Ion S5™ Cartridge Tool in the Ion S5™ Installation Kit.
A.0	1 September 2015	New product information sheet for the Ion S5™ Sequencer and Ion S5™ XL Sequencer instruments.

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.

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