

Ion AmpliSeq™ Exome RDY Kits

Catalog Numbers A38262, A38264

Pub. No. MAN0010195 Rev. D

Note: For safety and biohazard guidelines, see the “Safety” appendix in the following product documentation: *Ion AmpliSeq™ Exome RDY Library Preparation User Guide* (Pub. No. [MAN0010084](#)). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Product description

The Ion AmpliSeq™ Exome RDY Kits (Cat. Nos. [A38262](#) and [A38264](#)) provide reagents and materials for preparing eight Ion AmpliSeq™ exome libraries.

The kits include the following:

- Ion AmpliSeq™ Library Kit Plus
- Ion AmpliSeq™ Exome RDY Panel
 - 1×8 format (one plate with eight rows filled; Cat. No. [A38262](#)) or
 - 4×2 format (four plates with two rows filled per plate; Cat. No. [A38264](#))

Store the kit components as directed.

Ion AmpliSeq™ Exome RDY Kit configurations

Ion AmpliSeq™ Exome RDY Kit	Cat. No.	Ion AmpliSeq™ Exome RDY Panel plate format
Ion AmpliSeq™ Exome RDY Kit 1×8	A38262	1 plate × 8 rows
Ion AmpliSeq™ Exome RDY Kit 4×2	A38264	4 plates × 2 rows/plate

Kit contents and storage

Ion AmpliSeq™ Library Kit Plus (Cat. No. [4488990](#); sufficient for preparing eight exome libraries)

Component	Amount	Storage
5X Ion AmpliSeq™ HiFi Mix (red cap)	120 µL	–30°C to –10°C
FuPa Reagent (brown cap)	48 µL	
Switch Solution (yellow cap)	96 µL	
DNA Ligase (blue cap)	48 µL	
25X Library Amp Primers (pink cap)	48 µL	
1X Library Amp Mix (black cap)	1.2 mL	
Low TE	6 mL	15°C to 30°C ^[1]

^[1] Can be stored at –30°C to –10°C.

Ion AmpliSeq™ Exome RDY Panel (one format provided per kit)

Component, format	Amount	Storage
Ion AmpliSeq™ Exome RDY Panel, 1×8 (Part No. 4489838)	1 plate for 8 libraries	2°C to 8°C
Ion AmpliSeq™ Exome RDY Panel, 4×2 (Part No. 4489840)	4 plates for 2 libraries each	

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 can be impaired.

Refer to the expiration date on the package labels. 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 its affiliates 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 questions, contact Life Technologies at www.thermofisher.com/support.



Life Technologies Corporation | 5781 Van Allen Way | Carlsbad, California 92008 USA

For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](https://www.thermofisher.com/symbols-definition).

Revision history: Pub. No. MAN0010195 D

Revision	Date	Description
D	4 December 2024	Updated storage temperature for the Ion AmpliSeq™ Exome RDY Panel.
C.0	17 January 2018	New catalog numbers added for the Ion AmpliSeq™ Exome RDY Kits, and sequencing chips removed from the kits.
B.0	19 September 2017	<ul style="list-style-type: none">• Ion AmpliSeq™ Exome RDY S5 Kits included• Ion AmpliSeq™ Exome RDY Kit catalog numbers updated• Reagent volumes in the Ion AmpliSeq™ Library Kit Plus updated• Rebranding and other general updates
A.0	28 March 2014	New Product Information Sheet for the Ion AmpliSeq™ Exome RDY Kits.

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.

©2014-2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.