

Ion AmpliSeq™ Transcriptome Human Gene Expression Kit

Catalog Number A26325, A26326, A26327

Pub. No. MAN0010744 Rev. B.0

Note: For safety and biohazard guidelines, see the “Safety” appendix in the *Ion AmpliSeq™ Transcriptome Human Gene Expression Kit User Guide* (Pub. No. MAN0010742). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Product description

The Ion AmpliSeq™ Human Gene Expression Core Panel provides gene-level expression information from a single multiplexed panel targeting over 20,000 genes (> 95% of the RefSeq gene database). The small amplicon designs enable the use of the panel with RNA isolated from fixed tissues, such as formalin-fixed paraffin-embedded (FFPE) samples. Library preparation with the kit requires 10 ng of total RNA input, without the need for poly (A) selection or ribosomal RNA depletion.

Note: For preparing automated libraries with the Ion Chef System, use the Ion AmpliSeq™ Transcriptome Human Gene Expression Panel: Chef-ready Kit (Cat. No. A31446, ordered separately). For more information, see the *Ion AmpliSeq™ Library Preparation on the Ion Chef™ System User Guide* (Pub. No. MAN0013432).

Contents and storage

The Ion AmpliSeq™ Transcriptome Human Gene Expression Kit (Cat. Nos. A26325, A26326, and A26327) provides reagents for 24, 96, and 384 libraries, respectively.

Component	Number of kits		
	Cat. No. A26325 (24 reactions)	Cat. No. A26326 (96 reactions)	Cat. No. A26327 (384 reactions)
Ion AmpliSeq™ Library Kit Plus	1	4	16
SuperScript™ VIL0™ cDNA Synthesis Kit	1	2	8
Ion AmpliSeq™ Transcriptome Human Gene Expression Core Panel	1	4	16

The kit components for each kit are shown in the following table. The amount of each component is indicated for a single kit.

Contents	Amount	Storage
Ion AmpliSeq™ Library Kit Plus [1]		
5X Ion AmpliSeq™ HiFi Mix (red cap)	130 µL	-30°C to -10°C
FuPa Reagent (brown cap)	60 µL	
Switch Solution (yellow cap)	130 µL	
DNA Ligase (blue cap)	60 µL	
25X Library Amp Primers (white cap)	60 µL	
1X Library Amp Mix (black cap)	1.5 mL	
Low TE (clear cap)	1.7 mL	Room temp (15–30°C)
SuperScript™ VIL0™ cDNA Synthesis Kit [2, 3]		
5X VIL0™ Reaction Mix	200 µL	-30°C to -10°C
10X SuperScript™ III Enzyme Mix	100 µL	

Contents	Amount	Storage
Ion AmpliSeq™ Transcriptome Human Gene Expression Core Panel^[3]		
Ion AmpliSeq™ Transcriptome Human Gene Expression Core Panel	192 µL	-30°C to -10°C

^[1] Kit is shipped on frozen gel packs. Store as indicated.

^[2] May also be ordered separately (Cat. No. 11754-050).

^[3] Shipped on dry ice. Store as indicated.

Customer and technical support

Visit thermofisher.com/support for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- 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.

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.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

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.

Corporate entity: Life Technologies Corporation | Carlsbad, CA 92008 USA | Toll Free in USA 1 800 955 6288

©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.