

Genexus™ Barcodes 1–32 HD

Catalog Number A40261

Pub. No. MAN0018964 Rev. A.0

Note: For safety and biohazard guidelines, see the “Safety” appendix in the *Genexus™ Integrated Sequencer User Guide* (Pub. No. MAN0017910). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Product description

The Ion Torrent™ Genexus™ Barcodes 1–32 HD provides one plate that is filled with 32 dried-down, single-use barcodes, and is covered with a pierceable foil seal that eliminates cross-contamination between wells. The barcodes can be used in increments of one to 32, until all barcodes are exhausted. When used with Genexus™ Library Strips 1 and 2-HD (Cat. No. A40255), the kit supports flexible, fully automated preparation and normalization of up to 32 Ion AmpliSeq™ HD libraries on the Ion Torrent™ Genexus™ Integrated Sequencer. For Ion AmpliSeq™-based assays, select from Ion Torrent™ Genexus™ Barcodes 1–32 AS (Cat. No. A40258), Genexus™ Barcodes 33–64 AS (Cat. No. A40259), Genexus™ Barcodes 65–96 AS (Cat. No. A40260), and Genexus™ Barcodes 1–96 AS (Cat. No. A40257), which are ordered separately.

Contents and storage

Component	Label color	Quantity	Storage
Genexus™ Barcodes 1–32 HD	Purple	1 plate	15°C to 30°C

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.

IMPORTANT! Do not substitute components in this kit with components from other Ion Torrent™ templating or sequencing kits. This protocol has been verified using these specific materials. Substitution can adversely affect system performance.

Note: Refer to the expiration date on the package label. Do not use products 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.



Life Technologies Corporation | 7335 Executive Way | Frederick, MD 21704 | USA
For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

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.

Revision history: Pub. No. MAN0018964

Revision	Date	Description
A.0	2 December 2019	New product information sheet for Genexus™ Barcodes 1–32 HD

Important Licensing Information: This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.