

Genexus™ Pipette Tips

Catalog Number A40266

Pub. No. MAN0018966 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™ Pipette Tips contain 12 individually wrapped racks of 96 tips each, that are specially designed to fit into designated spaces on the deck of the Ion Torrent™ Genexus™ Integrated Sequencer. The lid and wrapping keeps tips free of dust and contaminants before use, and the yellow base protects tips from contacting surfaces during handling. Each run utilizes two to six racks of pipette tips, depending on configuration.

Contents and storage

Component	Quantity	Storage
Genexus™ Pipette Tips	12 racks	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 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. MAN0018966

Revision	Date	Description
A.0	2 December 2019	New product information sheet for Genexus™ Pipette Tips

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.