Capillary arrays for the 3500/3500xL Genetic Analyzer

Catalog Numbers 4404683, 4404685, 4404687, 4404689

Pub. No. 100070886 Rev. A



WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from **thermofisher.com/support**.

Product description

The capillary array for the 3500/3500xL Genetic Analyzer is available in 8-capillaries or 24-capillaries.



WARNING! SHARP The load-end of the capillary array has small, blunt ends that can lead to piercing injury.

Contents and storage

Cat. No.	Size	Storage
4404683	8-capillary, 36-cm	Room temperature
4404685	8-capillary, 50-cm	
4404687	24-capillary, 36-cm	
4404689	24-capillary, 50-cm	

Expiration date/on-instrument use limits

The on-instrument life is determined by the limit that is reached first—number of days after first installation, samples run, injections performed, or expiry date. Usage is tracked by the system software.

For the limits set by the software, refer to the user guide provided with the software or instrument.

Instrument	On-instrument use limits Lower of:	Guidelines
8-capillary	160 injections or expiration date listed on packaging and	Store capillary arrays with the loading-end of the
24-capillary	RFID label when used with Thermo Fisher Scientific reagents.	capillary array in distilled water to prevent the polymer from drying in the capillaries.

Install the capillary array

See the instrument user guide for instructions on how to install the capillary array.

Important notice regarding use of consumables that exceed supported limits

BEFORE DISMISSING THE WARNING THAT THE CONSUMABLES HAVE REACHED SUPPORTED LIMITS AND CONTINUING WITH OPERATION OF THE INSTRUMENT, PLEASE READ AND UNDERSTAND THE FOLLOWING IMPORTANT NOTICE AND INFORMATION:

Life Technologies does not recommend the use of consumables that exceed supported limits. The recommended limits are designed to promote the production of high quality data and minimize instrument downtime. Reagent and consumable lifetime minimum performance are based on testing and studies that use reagents and consumables that have not exceeded supported limits.

The use of consumables beyond the supported limits may impact data quality or cause damage to the instrument or capillary array. The cost of repairing such damage is *NOT* covered by any Life Technologies product warranty or service plan. Customer use of expired consumables is at customer's own risk and without recourse to Life Technologies. For example, product warranties do not apply to defects



resulting from or repairs required due to misuse, neglect, or accident including, without limitation, operation outside of the environmental or use specifications or not in conformance with Life Technologies instructions for the instrument system, software, or accessories.

Please see your specific service contract or limited product warranty for exact language regarding coverage and ask your Life Technologies representative if you have further questions.

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.



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

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.

Revision history: Pub. No. 100070886

Revision	Date	Description
A.0	11 April 2018	New document.

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.

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.



11 April 2018