POP-7[™] Polymer for 3730/3730xl DNA Analyzer

Catalog Numbers 4363929, 4335615, 4363935, and 4335611

Pub. No. 100032432 Rev. B

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 polymer is the separation matrix for capillary electrophoresis on 3730/3730*xl* DNA Analyzers. It can be used for sequencing analysis and fragment analysis applications. The polymer is pre-formulated.

Contents and storage

Cat. No.	Amount	Storage	
4363929	1 x 28-mL bottle (8,000–12,000 samples/bottle) ^[1]	Store at 2–8°C.	
4335615	5 x 28-mL bottles (8,000–12,000 samples/bottle) ^[1]		
4363935 10 x 28-mL bottles (8,000–12,000 samples/bottle) ^[1]			
4335611	30 x 28-mL bottles (8,000–12,000 samples/bottle) ^[1]		

^[1] The polymer has been validated for HID applications.

Expiration date/on-instrument supported limits

The on-instrument life is determined by the limit that is reached first—number of days after first installation, samples run, or expiration date.

Bottle size	Array size	On-instrument supported limits ^[1] Lower of:	Guidelines
28-mL	36-capillary	7 days, 8,000–12,000 samples, or expiration date	The polymer has been verified for use for up to 7 days on the instrument. See "Important notice regarding use of consumables that exceed supported limits" on page 2.
	50-capillary	7 days, 8,000–12,000 samples, or expiration date	

[1] The bottle has adequate polymer to support the stated number of samples, plus additional volume to accommodate installation and changing the polymer type.

Replenish or change the polymer

1. Check the expiration date on the label to ensure that the polymer is not expired and will not expire during intended use.

IMPORTANT! Do not use if the product is expired, if the bottle or label is damaged, or if the cap is missing or damaged.

2. Allow the refrigerated polymer to equilibrate to ambient temperature before use.



3. In the software, click Wizards to replenish or change the polymer.

Option	Select	
To replenish	Wizards > Replenish Polymer Wizard, then select Same Lot or Different Lot Wizards > Change Polymer Wizard	
To change		

4. Follow the prompts in the Wizard window.

See the instrument user guide for instructions on starting the runs.

Important notice regarding use of consumables that exceed supported limits

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 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 | 35 Wiggins Avenue | Bedford, MA 01730

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. 100032432

Revision	Date	Description
В	31 October 2018	Updates to template and Wizard options.
A	26 March 2015	New document.

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.

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

