

Conditioning Reagent Kit 3500/Flex Series

SeqStudio™ Flex and 3500 series instruments

Catalog Number 4393718

Pub. No. 4408237 Rev. E

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](https://www.thermofisher.com/support).

Product description

The conditioning reagent is used to prime the polymer delivery pump, wash the polymer delivery pump between polymer type changes, perform instrument shutdown procedures, and perform routine maintenance. It is supplied as a ready-to-use pouch with a radio frequency identification (RFID) tag incorporated into the label. The instrument uses the RFID tag to track the conditioning reagent usage and expiration.

Contents and storage

Cat. No.	Description	Storage
4393718	Conditioning Reagent Kit 3500/Flex Series, 1 pouch	2–8°C After removing from storage, use the pouch within 24 hours.

Expiration date/on-instrument supported limits

On-instrument supported limits	Guidelines
After it is installed on the instrument, the pouch is good for a one-time use.	See the expiration date on the label. See "Important notice regarding use of consumables that exceed supported limits" on page 2.

Install the conditioning reagent

The instrument wizards dictate use of the conditioning reagent. Install the pouch only when requested to do so by the wizard.

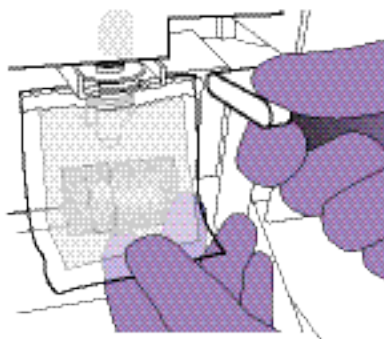
1. Check the expiration date on the label to ensure that the conditioning reagent is not expired and will not expire during intended use.

IMPORTANT! Do not use if the product is expired, if the pouch or label is damaged, or if the top seal is missing or damaged.

2. Allow the refrigerated conditioning reagent to equilibrate to ambient temperature (15–30°C) before use.
3. Peel off the seal at the top of the pouch fitment.

Note: You may notice a tiny droplet of conditioning reagent inside the fitment (residual from the pouch filling process). This is **not** expected to cause any performance issues.

4. With the RFID label facing the instrument, slide the pouch fitment onto the slot of the lever assembly. Push the lever up to snap the pouch into the connector end of the instrument pump.



Note: The RFID label must face the instrument (away from you) to ensure that the RFID information is read accurately by the instrument.

5. Follow the wizard instructions or refer to the instrument user guide to complete the necessary procedure.

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:

Thermo Fisher Scientific 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 Thermo Fisher Scientific product warranty or service plan. Customer use of expired consumables is at customer's own risk and without recourse to Thermo Fisher Scientific. 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 Thermo Fisher Scientific 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 Thermo Fisher Scientific 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 | 1455 Concord Street | Framingham, Massachusetts 01701 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. 4408237

Revision	Date	Description
E	31 January 2022	Added the SeqStudio™ Flex Series Genetic Analyzer. In "Product description", added routine maintenance to the list of uses for the conditioning reagent. Updated branding and legal information.
D	31 March 2015	Baseline for this revision.

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.

©2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.