

# HT ExoSAP-IT™ High-Throughput PCR Product Cleanup

Brief Protocol

Catalog Number 78395

Pub. No. MAN0016839 Rev. A.0 (02/2017)

**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](http://thermofisher.com/support).

## Product description

HT ExoSAP-IT™ reagent treats PCR products ranging in size from less than 100 bp to over 20 kb with absolutely no sample loss by degrading unused primers and nucleotides. This HT formulation and packaging is designed for use with liquid handling robots or a multichannel pipette. The 8-tube strip format holds 60 reactions in each tube for a total of 480 reactions per strip. The plate format holds 12 of the 8-tube strips for a total of 5,760 reactions per plate. The reaction volume is based on a 2 µL quantity of HT ExoSAP-IT™ reagent for 5 µL of post-PCR reaction product. Simply remove the strip cap and begin pipetting.

For 96-well plates, all 12 of the 8-tube strip caps can be removed for robotic pipetting. Add HT ExoSAP-IT™ reagent directly to the reaction products following PCR. HT ExoSAP-IT™ reagent is active in commonly used PCR buffers, so no buffer exchange is required. After treatment, it is inactivated by heating to 80°C for 15 minutes. The treated PCR products are now ready for subsequent analysis in applications that require DNA to be free of excess primers and nucleotides.

## PCR cleanup protocol

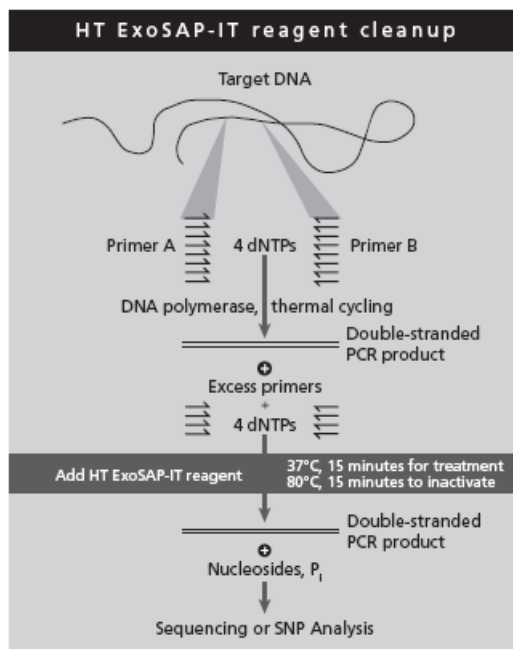
**Note:** Store HT ExoSAP-IT™ reagent in a non-frost-free freezer.

1. Remove HT ExoSAP-IT™ reagent strips or plate from -20°C freezer and thaw on ice. Keep on ice throughout this procedure.
2. Mix 5 µL of a post-PCR reaction product with 2 µL of HT ExoSAP-IT™ reagent for a combined 7 µL reaction volume.

When treating PCR product volumes greater than 5 µL, simply increase the amount of HT ExoSAP-IT™ reagent proportionally.

3. Incubate at 37°C for 15 minutes to degrade excess primers and nucleotides.
4. Incubate at 80°C for 15 minutes to inactivate HT ExoSAP-IT™ reagent.
5. The PCR product is now ready for use in DNA sequencing, SNP analyses, or other primer-extension applications. Treated PCR products may be stored at -20°C until required.

## HT ExoSAP-IT™ reagent cleanup overview



## Customer and technical support

Visit [thermofisher.com/support](http://thermofisher.com/support) for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
  - Product FAQs
  - Software, patches, and updates
  - Training for many applications and instruments
- Order and web support
- Product documentation, including:
  - 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 found on Life Technologies' website at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://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](http://www.thermofisher.com/support).

---

**Manufacturer's address:** Affymetrix Inc. | 3450 Central Expressway | Santa Clara, CA 95051 | USA

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.

**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.

**Corporate entity:** Life Technologies Corporation | Carlsbad, CA 92008 USA | Toll Free in USA 1 800 955 6288

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.