

USB® HT ExoSAP-IT® High-Throughput PCR Product Cleanup

Product number 78395

Brief protocol

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.

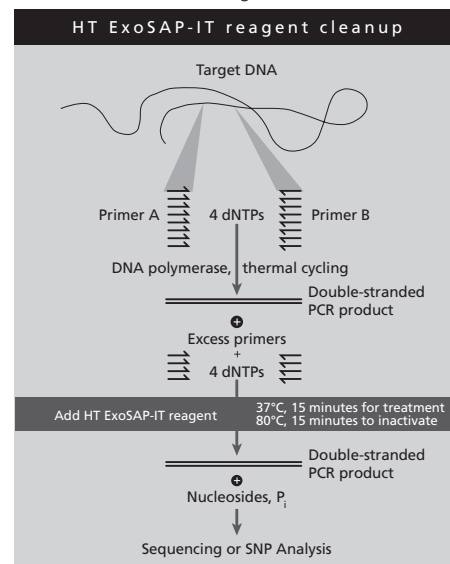
Add HT ExoSAP-IT reagent directly to the reaction products following PCR. HT ExoSAP-IT reagent is active in commonly used PCR buffers, so no buffer exchange is required. After treatment, it is inactivated by heating to 80°C for 15 minutes. The treated PCR products are now ready for subsequent analysis in applications that require DNA to be free of excess primers and nucleotides.

PCR cleanup protocol:

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



For research use only. Not for use in diagnostic procedures.

USB, Affymetrix and ExoSAP-IT are registered trademarks of Affymetrix, Inc.

ExoSAP-IT PCR Product Cleanup is covered by patents owned by Affymetrix, including US

Patent Nos. 6,379,940 and 6,387,634 and corresponding patents in Canada and Japan.

ExoSAP-IT PCR Product Cleanup—This product or portions thereof are sold under license from GE Healthcare under US Patent Nos. 5,741,676 and 5,756,285.

© 2015 Affymetrix, Inc. All rights reserved.

Affymetrix, Inc.

USA: USBtechsupport@affymetrix.com

Europe: USBtechsupporteurope@affymetrix.com

usb.affymetrix.com

P 78395B

rev 06/15