

GeneChip™ T7-Oligo(dT) Promoter Primer Kit

Catalog Number 900375

Pub. No. 701182 Rev. 3

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 use

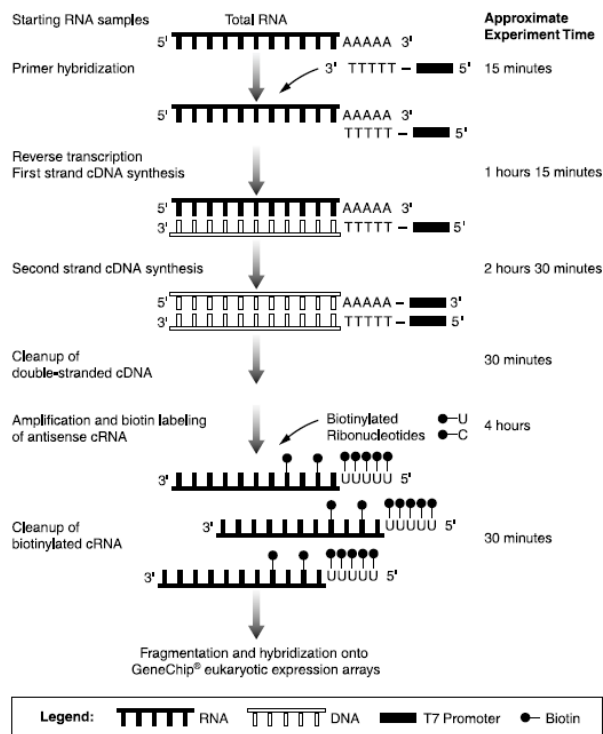
The Applied Biosystems™ GeneChip™ T7-Oligo(dT) Promoter Primer Kit provides reagents essential for the primer hybridization step in the first strand cDNA synthesis reaction of eukaryotic target preparation.

Introduction

In situ synthesized oligonucleotide probes on GeneChip™ eukaryotic expression arrays are designed in the sense orientation of corresponding transcripts. In GeneChip™ eukaryotic expression analysis, sample amplification and labeling is used to produce biotinylated antisense targets for hybridization to probes on the arrays. Instructions for the primer hybridization step are described below. The remaining procedure for target preparation is in the *GeneChip™ Expression Analysis Technical Manual* (Technical Manual), available on our website and will not be duplicated here.

The complete procedure involves hybridizing a chimeric T7-oligo(dT) primer to the sample mRNA in the reverse transcription reaction for first strand cDNA synthesis, followed by second strand cDNA synthesis. The purified double-stranded cDNA containing a T7 promoter at its 5' end can now serve as a template in the *in vitro* transcription reaction that produces many copies of complementary RNA (cRNA) that has incorporated biotinylated nucleotides. The biotinylated cRNA targets are then fragmented and hybridized to GeneChip™ expression arrays.

Eukaryotic Target Labeling GeneChip® Probe Arrays



The GeneChip™ T7-Oligo(dT) Promoter Primer Kit provides the essential reagents that are used in the primer hybridization step of the first strand cDNA synthesis reaction. This kit has modified the stock concentration of the T7-oligo(dT) primer to 50 μM from 100 μM as previously recommended. Use the following procedure for the primer hybridization step, then follow the protocols that are described in the Technical Manual for the subsequent cDNA synthesis and cRNA amplification.

Kit components and storage

Unless otherwise indicated, all materials are available through thermofisher.com.

MLS: Fisher Scientific (fisherscientific.com) or other major laboratory supplier.

Cat. No.	Size	Components
900375	150 reactions	50 μM T7-oligo(dT) primer, HPLC purified 300 μL × one vial (sequence: 5' - GGCCAGTGAATTGTAATACG ACTCACTATAGGGAGGCGG -(dT) ₂₄ -3') RNase-free water 825 μL × 2 vials

IMPORTANT! The primer should be stored at -20°C. Performance has been shown to be unaffected for up to 25 freeze-thaw cycles.

Materials not provided with the kit

See the Technical Manual for a complete list of reagents that are needed to complete cDNA synthesis, cleanup, and cRNA amplification and labeling.

Before starting

The quality of the RNA samples is essential to the overall success of the analysis. We recommend routine assessment of RNA quality and integrity for all samples.

Either eukaryotic total RNA or poly-A RNA can be used as starting material for GeneChip™ target labeling. A minimum amount of 5 µg of total RNA or 0.2 µg of poly-A RNA is recommended.

Follow routine laboratory procedures for handling RNA samples and components of this kit to prevent RNase contamination.

Primer hybridization procedure

1. Remove the T7-oligo(dT) primer from the freezer. Ensure that the primer is thawed before proceeding.
2. Lightly vortex the vial of primer for complete mixing. Briefly spin down the reagent in a microcentrifuge.
3. In a microfuge tube, combine sample RNA with 2 µL of 50 µM T7-oligo(dT) primer.
4. Calculate the amount of RNase-free water that is needed to make up a final volume of 20 µL for the entire first strand cDNA synthesis reaction. See the Technical Manual for the volume calculation. Add the appropriate amount of RNase-free water to the reaction.
5. Heat the reaction mix at 70°C for 10 minutes to denature the sample RNA/promoter primer mix.

6. Quickly spin the microfuge tube and immediately cool on ice.
7. The RNA/primer complex is now ready for cDNA synthesis. See the Technical Manual for complete GeneChip™ target amplification instructions.

Customer and technical support

Visit 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. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Thermo Fisher Scientific Baltics UAB | V.A. Graiciuno 8, LT-02241 | Vilnius, Lithuania

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. 701182

Revision	Date	Description
3	06 September 2017	Update document to current template.
2	07 November 2002	Baseline for revision history.

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.

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