

GeneChip™ WT PLUS Reagent Kit and GeneChip™ HT WT PLUS Reagent Kit

Catalog Numbers 902280, 902281, and 902414

Doc. Part No. 703172 Pub. No. MAN0017939 Rev. A.0

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™ WT PLUS Reagent Kit (manual version) and GeneChip™ HT WT PLUS Reagent Kit (high-throughput version) generate amplified and biotinylated sense-strand DNA targets from total RNA without the need for an up-front selection or enrichment step for mRNA. The kits are optimized for use with our Whole Transcript (WT) Expression Arrays.

The GeneChip™ WT PLUS Reagent Kit is offered in three configurations.

- 10-reaction kit for generating ten target preps using a manual assay protocol.
- 30-reaction kit for generating 30 target preps using a manual assay protocol, OR 24 target preps using an automated target prep assay on the Beckman Coulter™ Biomek™ FX^P.
- 96-reaction kit for generating 96 target preps using an automated target prep assay on the Beckman Coulter™ Biomek™ FX^P.

Introduction

Both the manual and high-throughput (HT) versions of the Applied Biosystems™ GeneChip™ WT PLUS Reagent Kit use a reverse transcription priming method that primes the entire length of RNA, including both poly(A) and nonpoly(A) mRNA to provide complete and unbiased coverage of the transcriptome. Each kit consists of reagents and a protocol for preparing hybridization-ready targets from 50–500 ng of total RNA. WT PLUS Reagent is optimized to work with a wide range of samples including tissues, cells, cell lines, and whole blood. The total RNA samples can be used directly without removal of ribosomal or globin RNA before target preparation with WT PLUS Reagent.

Kit components

Module	Component	Storage
WT Amplification Kit Module 1	First-Strand Enzyme	-20°C
	First-Strand Buffer	-20°C
	Second-Strand Enzyme	-20°C
	Second-Strand Buffer	-20°C
	IVT Enzyme	-20°C
	IVT Buffer	-20°C
	Control RNA (1 mg/mL HeLa total RNA)	-20°C
	2nd-Cycle Primers	-20°C
	2nd-Cycle ss-cDNA Enzyme	-20°C
	2nd-Cycle ss-cDNA Buffer	-20°C
	RNase H	-20°C
Nuclease-free Water	any temp ^[1]	
WT Amplification Kit Module 2	Purification Beads	4°C ^[2]
GeneChip™ Eukaryotic Poly-A RNA Control Kit	Poly-A Control Stock	-20°C
	Poly-A Control Dil Buffer	-20°C
GeneChip™ WT Terminal Labeling Kit	10X cDNA Fragmentation Buffer	-20°C
	UDG, 10 U/μL	-20°C
	APE 1, 1,000 U/μL	-20°C
	5X TdT Buffer	-20°C
	TdT, 30 U/μL	-20°C
	DNA Labeling Reagent, 5 mM	-20°C
	RNase-free Water	any temp ^[1]
GeneChip™ Hybridization Control Kit	20X Hybridization Controls	-20°C
	3 nM Control Oligo B2	-20°C

^[1] Store the Nuclease-free Water at -20°C, 4°C, or room temperature. Not included in Cat. No. 902414.

^[2] Do not freeze.

Shipping information

The five components of the GeneChip™ WT PLUS Reagent Kit and the GeneChip™ HT WT PLUS Reagent Kit are shipped at different temperatures.

- WT Amplification Kit Module 1 (–20°C component)
- WT Amplification Kit Module 2 (4°C component)
- GeneChip™ Eukaryotic Poly-A RNA Control Kit (–20°C component)
- GeneChip™ WT Terminal Labeling Kit (–20°C component)
- GeneChip™ Hybridization Control Kit (–20°C component)

Ordering information

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

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

Cat. No.	Product name	Description
902280	GeneChip™ WT PLUS Reagent Kit	10 reactions
902281	GeneChip™ WT PLUS Reagent Kit	30 reactions
902414	GeneChip™ HT WT PLUS Reagent Kit	96 reactions
Supporting products		
900720	GeneChip™ Hybridization, Wash, and Stain Kit	Sufficient for 30 reactions
901667	GeneAtlas™ Hybridization, Wash, and Stain Kit for WT Array Strips	Sufficient for 60 reactions
901622	GeneTitan™ Hybridization, Wash, and Stain Kit for WT Array Plates	Sufficient for 96 reactions

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.



Manufacturer: 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, 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. MAN0017939

Revision	Date	Description
A.0	10 July 2018	Initial release in Thermo Fisher Scientific document control system. Supersedes legacy Affymetrix publication number 703172. Updated to correct component in shipping information section.

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.

©2018 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.

thermofisher.com/support | thermofisher.com/askaquestion

thermofisher.com

10 July 2018

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.

ThermoFisher
SCIENTIFIC