

GeneChip™ Pico Kit and GeneChip™ HT Pico Kit

Catalog Numbers 902789, 902790, and 902791

Pub. No. 703361 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.

Intended use

This product information sheet provides general instruction for the Applied Biosystems™ GeneChip™ Pico Kit and the Applied Biosystems™ GeneChip™ HT Pico Kit. Each kit generates amplified and biotinylated double-stranded cDNA (ds-cDNA) from total RNA sample and is optimized for use with gene-level expression arrays such as 3' IVT, Gene ST, among many others.

The GeneChip™ Pico Kit is offered in three configurations:

1. 12-reaction kit for generating 12 target preparations using a manual assay protocol.
2. 30-reaction kit for generating 30 target preparations using a manual assay protocol; or 24 target preparations using an automated target preparation assay on the Beckman Coulter Biomek FX^P.
3. 96-reaction kit for generating 96 target preparations using an automated target preparation assay on the Beckman Coulter Biomek FX^P.

Introduction

Both the manual and high throughput versions of the Pico Reagent Kit use a reverse transcription priming method that primes at the poly-A tail as well as throughout the entire length of RNA to ensure unbiased coverage of the transcriptome. Each kit is comprised of reagents and a protocol for preparing hybridization ready targets from 100 pg to 10 ng of purified total RNA from cells or tissues and 500 pg to 50 ng of purified total RNA from FFPE tissues. The kit is ideal for amplification of intact, partially degraded, and compromised RNA samples, including cell lines, fresh tissues, whole blood, and formalin-fixed paraffin-embedded (FFPE) tissues. The total RNA samples can be used directly without removal of ribosomal or globin RNA prior to target preparation with Pico Reagent.

Instructions for use

Please refer to the *GeneChip™ Pico Kit User Guide* (Pub. No. 703308 for manual assay and Pub. No. 703398 for automation assay) for procedures on DNA target preparation, target hybridization, fluidics setup, array scanning, and data analysis instructions.

Kit components

Module	Component	Storage
Pico Amplification Kit, Module 1	Pico First-Strand Enzyme	-20°C
	Pico First-Strand Buffer	-20°C
	Pico Cleanup Reagent	-20°C
	Pico 3' Adaptor Enzyme	-20°C
	Pico 3' Adaptor Buffer	-20°C
	Pico PCR Enzyme	-20°C
	Pico PCR Buffer	-20°C
	Pico IVT Enzyme	-20°C
	Pico IVT Buffer	-20°C
	Poly-A Control Stock	-20°C
Pico Amplification Kit, Module 2	Nuclease-Free Water	Any temp ^[1]
	Pico 2nd-Cycle Primers	-20°C
	Pico 2nd-Cycle ds-cDNA Enzyme	-20°C
	Pico 2nd-Cycle ds-cDNA Buffer	-20°C
	Pico Frag. & Label Enzyme	-20°C
Pico Amplification Kit, Module 3	Pico Frag. & Label Buffer	-20°C
	Purification Beads	4°C ^[2]
	Pico Hydrolysis Buffer	4°C
Control HeLa RNA	Pico Neutralization Buffer	4°C
	Control RNA (100 ng/μL HeLa total RNA)	-20°C
GeneChip™ Hybridization Control Kit	20X Hybridization Controls	-20°C
	3 nM Control Oligo B2	-20°C
Tubes Organizer	Plastic vinyl template for organization and storage of components in 9 x 9 array, 81-places square wells, 5 1/4 in. x 5 1/4 in (e.g., Nalgene™ CryoBox Cat. No. 5026-0909, or equivalent).	

^[1] Store the Nuclease-Free Water at -20°C, 4°C, or room temperature.

^[2] Do not freeze.

Shipping information

The five components of the GeneChip™ Pico Kit are shipped at different temperatures:

- Pico Amplification Kit Module 1 (–20°C component)
- Pico Amplification Kit Module 2 (–20°C component)
- Pico Amplification Kit Module 3 (4°C component)
- Control HeLa RNA (–20°C component)
- GeneChip™ Hybridization Control Kit (–20°C component)

Functional testing








Every lot is rigorously tested for performance.

Ordering information

Cat. No.	Product name	Description
902789	GeneChip™ Pico Kit	Sufficient for 12 reactions
902790	GeneChip™ Pico Kit	Sufficient for 30 reactions
902791	GeneChip™ HT Pico Kit	Sufficient for 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
901531	GeneAtlas™ Hybridization, Wash, and Stain Kit for 3' IVT Arrays	Sufficient for 60 reactions
901622	GeneTitan™ Hybridization, Wash, and Stain Kit for WT Array Plates	Sufficient for 96 reactions
901530	GeneTitan™ Hybridization, Wash, and Stain Kit for 3' IVT Arrays	Sufficient for 96 reactions

Symbol table

Table below shows a legend of the graphic symbols used for product label, product information sheets and user guides.

Symbol/Letters	Statement
	CATALOG NUMBER
	BATCH CODE
	USE BY YYYY-MM-DD
	TEMPERATURE LIMITATION
	CONTAINS SUFFICIENT FOR <n> TESTS
	CONSULT INSTRUCTIONS FOR USE
	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.



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, 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. 703361

Revision	Date	Description
3	20 July 2017	Updated document to current template.
2	10 March 2016	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.