

# GeneChip™ Eukaryotic Poly-A RNA Control Kit

Catalog Numbers 900433

Pub. No. 701379 Rev. 7



**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).

## Introduction

The Applied Biosystems™ GeneChip™ Eukaryotic Poly-A RNA Control Kit is designed specifically to provide exogenous positive controls to monitor the entire GeneChip™ eukaryotic target labeling process.

Each eukaryotic GeneChip™ expression array contains probe sets for several *B. subtilis* genes that are absent in the samples analyzed (*lys*, *phe*, *thr*, and *dap*). This Poly-A RNA Control Kit contains *in vitro* synthesized, polyadenylated transcripts for these *B. subtilis* genes that are pre-mixed at staggered concentrations to allow GeneChip™ expression array users to evaluate the overall success of the assay. The concentrated Poly-A Control Stock can be diluted with the Poly-A Control Dil Buffer (supplied as part of the kit) and spiked directly into the RNA samples to achieve the final concentrations (known as a ratio of copy number) summarized below:

Poly-A RNA Controls	Final concentration
<i>lys</i>	1:100,000
<i>phe</i>	1:50,000
<i>thr</i>	1:25,000
<i>dap</i>	1:6,667

The controls are then amplified and labeled together with the samples. Examining the hybridization intensities of these controls on GeneChip™ arrays helps you monitor the labeling process independently from the quality of the starting RNA samples.

## Instructions for use

Detailed procedures for using this kit are provided in the *GeneChip™ Expression Analysis Technical Manual*, *GeneChip™ WT PLUS Reagent Kit User Guide*, and the *GeneChip™ 3' IVT PLUS Reagent Kit User Guide*, available at [thermofisher.com](http://thermofisher.com).

## Kit components

Component	Volume
Poly-A Control Stock	16 µL
Poly-A Control Dil Buffer	3.8 mL

## Storage

The GeneChip™ Eukaryotic Poly-A RNA Control Kit must be kept at -20°C in a non-frost-free freezer. Performance of the Poly-A Control Stock has been shown to be unaffected for up to eight freeze-thaw cycles.

## Functional testing

**GeneChip™ array functional testing:** Each lot is functionally tested in the GeneChip™ expression assay following the recommended protocols.

**Nuclease testing:** Each component is tested in rigorous nuclease assays.

## 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).

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. 701379

Revision	Date	Description
7	01 August 2017	Updated document to current template.
6	23 May 2014	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.