


FlashTag™ Biotin HSR RNA Labeling Kit

Catalog Numbers 901910 and 901911

Doc. Part No. 703098 Pub. No. MAN0017699 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

This document provides information for the Applied Biosystems™ FlashTag™ Biotin HSR RNA Labeling Kit. This kit is available in two configurations: 10 or 30 reactions. The kit includes all the necessary reagents for labeling any RNA sample for analysis on GeneChip™ miRNA Arrays.

See the *FlashTag™ Biotin HSR RNA Labeling Kit User Guide* (Pub. No. 703095) for target preparation, target hybridization, and array processing procedures.

Kit components

Vial number	Component
1	10X Reaction Buffer
2	25mM MnCl ₂
3	ATP Mix
4	PAP Enzyme
5	5X FlashTag™ Biotin HSR Ligation Mix
6	T4 DNA Ligase
7	HSR Stop Solution
8	RNA Spike Control Oligos
9	ELOSA Spotting Oligos
10	ELOSA Positive Control
11	Nuclease-Free Water
12	27.5% Formamide

Storage

The Applied Biosystems™ FlashTag™ Biotin HSR RNA Labeling Kit should be stored at -20°C.

Ordering information

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

Product name	Description	Source
FlashTag™ Biotin HSR RNA Labeling Kit	10 reactions	901910
	30 reactions	901911
Supporting products		
GeneChip™ Hybridization, Wash, and Stain Kit ^[1]	Sufficient for 30 reactions	900720
GeneChip™ Hybridization Control Kit	30 reactions	900454
	150 reactions	900457
GeneChip™ Control Oligo B2, 3 nM	30 reactions	900301
GeneAtlas™ Hybridization, Wash, and Stain Kit for miRNA Array Strips	Sufficient for 60 reactions	902134
GeneTitan™ Hybridization, Wash, and Stain Kit for miRNA Array Plates	Sufficient for 96 reactions	902276

^[1] Each kit contains 1 Hybridization Module, 1 Stain Module, 3 bottles of Wash Buffer A, and 1 bottle of Wash Buffer B sufficient for 30-reactions. Individual kit components may be ordered separately.

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. MAN0017699

Revision	Date	Description
A.0	13 May 2018	Initial release in Thermo Fisher Scientific document control system. Supersedes legacy Affymetrix publication number 703098. Updated to the current document template, with associated updates to trademarks, logos, licensing, and warranty.

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.