

GeneTitan™ Hybridization, Wash, and Stain Kit for WT Array Plates

Catalog Number 901622

Pub. No. 702899 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

This sheet provides information for the Applied Biosystems™ GeneTitan™ Hybridization, Wash, and Stain Kit for WT Array Plates (Cat. No. 901622). The kit is composed of two boxes:

- Part No. 901621: GeneTitan™ Hybridization Module for WT Array Plates—sufficient for 12 × 8 reactions, 6 × 16 reactions, 4 × 24 reactions, or 1 × 96 reactions

Note: The GeneTitan™ Hybridization, Wash, and Stain Kit for WT Array Plates (Cat. No. 901622) consists of Part No. 901621 and Cat. No. 901583. Part No. 901621 is not available for purchase separately.

- Cat. No. 901583: GeneTitan™ Wash Buffers A and B Module—sufficient for one array plate.

The GeneTitan™ Hybridization, Wash, and Stain Kit for WT Array Plates includes all the necessary reagents for processing Whole Transcript (WT) Array Plates.

Module components

Component	Volume	Storage
GeneTitan™ Hybridization Module for WT Array Plates (Part No. 901621)		
5X WT Hyb Add 1	4 × 925 µL	2°C to 8°C
15X WT Hyb Add 4	4 × 355 µL	2°C to 8°C
2.5X WT Hyb Add 6	4 × 1,525 µL	2°C to 8°C
Stain Cocktail 1 & 3	22 mL	2°C to 8°C
Stain Cocktail 2	13 mL	2°C to 8°C
Array Holding Buffer	110 mL	2°C to 8°C
GeneTitan™ Wash Buffers A & B Module (Cat. No. 901583)		
Wash Buffer A	1,045 mL	2°C to 8°C
Wash Buffer B	450 mL	2°C to 8°C

Note: The WT Hyb Add reagent names match the order in which reagents are added. For example, WT Hyb Add 4 is the fourth component added during preparation of the Hybridization Mix. WT Hyb Add 2, 3, and 5 are not used and are not part of the Hybridization Module.

Shipping information

The two components of the GeneTitan™ Hybridization, Wash, and Stain Kit for WT Array Plates are shipped at different temperatures. The GeneTitan™ Hybridization Module for WT Array Plates (Part No. 901621) is sent at 5°C. The GeneTitan™ Wash Buffers A & B Module (Cat. No. 901583) is sent at ambient temperature.

Functional testing

Every lot is rigorously tested for performance.

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/size
901622	GeneTitan™ Hybridization, Wash, and Stain Kit for WT Array Plates Contains the GeneTitan™ Hybridization Module for WT Array Plates (Part No. 901621) and the GeneTitan™ Wash Buffers A and B Module (Cat. No. 901583).	Sufficient for 96 reactions
Related products		
902414	GeneChip™ HT WT PLUS Reagent Kit	Sufficient for 96 reactions
902281	GeneChip™ WT PLUS Reagent Kit	Sufficient for 30 reactions (manual) or 24 reactions (automated)
902508	HT WT PLUS Consumables for Biomek™ FX ^P Target Prep Express	Sufficient 4 × 24 or 4 × 96 reactions

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.



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. 702899

Revision	Date	Description
3	06 September 2017	Update document to current template.
2	24 April 2015	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.