

# HT WT PLUS Consumables for Biomek™ FX<sup>P</sup> Target Prep Express

Catalog Numbers 902508

Pub. No. 703236 Rev. 2

**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](https://www.thermofisher.com/support).

## Product use

The Applied Biosystems™ HT WT PLUS Consumables for Biomek™ FX<sup>P</sup> Target Prep Express contains plastic consumables that are needed for target preparation on the Biomek™ FX<sup>P</sup> using the GeneChip™ HT WT PLUS Reagent Kit. The accessory kit contains consumables sufficient for 4x24 or 4x96 reaction runs. See the *GeneChip™ HT WT PLUS Reagent Kit User Guide* (Pub. No. 703208) for instructions on using the consumables.

## Kit components

Component	Number per run	Number per kit	Figure
96-well deep well microplate, polypropylene, 2 mL, non-sterile	2	8	Figure 1
Greiner U-bottom (round) plate, polypropylene	1	4	Figure 2
Bio-Rad™ Hard-Shell™ PCR plate, blue shell/clear well	18	72	Figure 3
Greiner UV-Star™ plate, nonsterile, clear, flat bottom	2	8	Figure 4
Costar™ 96-well U-bottom plate, polystyrene	4	16	Figure 5
24-well plate, 10 mL, square well, pyramid bottom, polypropylene	1	4	Figure 6
FrameStar™ Break-A-Way plate	1	4	Figure 7
'P' Pad for lid, autosealing microplates, arched with Microseal™	1	2 <sup>[1]</sup>	Figure 8

<sup>[1]</sup> Reusable. See the *GeneChip™ HT WT PLUS Reagent Kit User Guide* (Pub. No. 703208) for instructions.

## Storage

The HT WT PLUS Consumables for Biomek™ FX<sup>P</sup> Target Prep Express must be stored at room temperature.

## Ordering information

Cat. No.	Product name	Description
902508	HT WT PLUS Consumables for Biomek™ FX <sup>P</sup> Target Prep Express	Labware sufficient for 4x24 or 4x96 reaction runs

## Component images



Figure 1 96-well deep well microplate



Figure 2 Greiner U-bottom (round) plate

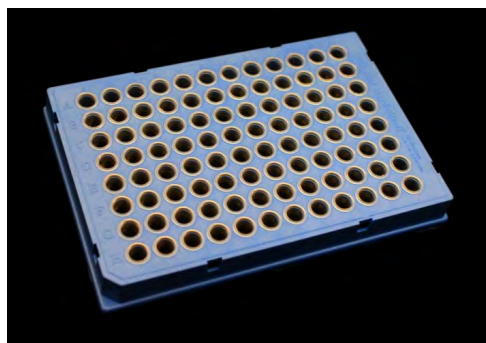


Figure 3 Bio-Rad™ Hard-Shell™ PCR plate



Figure 4 Greiner UV-Star™ plate



Figure 5 Costar™ 96-well U-bottom plate



Figure 6 24-well plate

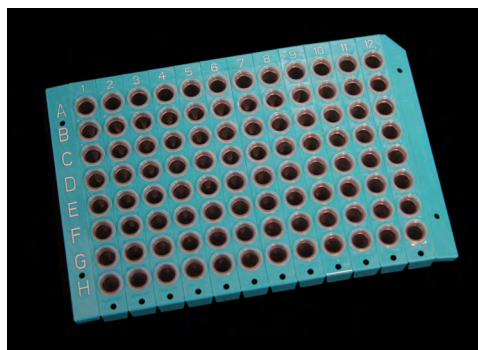


Figure 7 FrameStar™ Break-A-Way plate

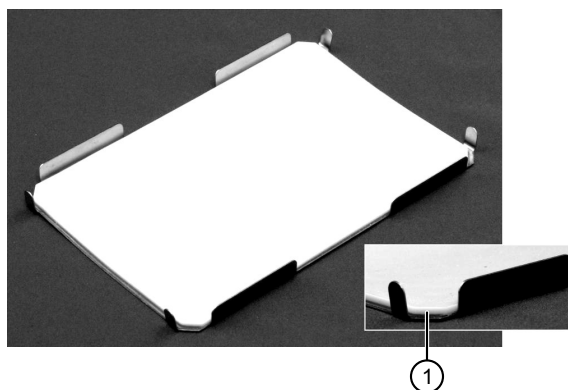


Figure 8 'P' Pad for lid, arched with Microseal™

① 'P' Pad

### 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. 703236

Revision	Date	Description
2	15 August 2017	Update document to current template.
1	18 March 2014	Initial release of document.

**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. Biomek is a registered trademark of Beckman Coulter, Inc. Bio-Rad, Hard-Shell, and Microseal are registered trademark of Bio-Rad Laboratories, Inc. UV-Star is a registered trademark of Greiner Bio-One GmbH. Costar is a registered trademark of Corning, Inc. FrameStar is a registered trademark of 4titude Ltd.