A large, vertical, metallic DNA double helix structure that runs down the left side of the page, starting from the top and ending at the bottom.

Site Preparation Guide

GeneChip® HT Methods for
Expression on the Beckman
Coulter® Biomek® FX^P Target
Prep Instrument

**For research use only.
Not for use in diagnostic procedures.**

Trademarks

GeneChip®, NetAffx®, Command Console®, Powered by Affymetrix™, GeneChip-compatible™, Genotyping Console™, DMET™, GeneTitan®, Axiom™, and GeneAtlas™ are trademarks or registered trademarks of Affymetrix, Inc. All other trademarks are the property of their respective owners.

Limited License

Subject to the Affymetrix terms and conditions that govern your use of Affymetrix products, Affymetrix grants you a non-exclusive, non-transferable, non-sublicensable license to use this Affymetrix product only in accordance with the manual and written instructions provided by Affymetrix. You understand and agree that, except as expressly set forth in the Affymetrix terms and conditions, no right or license to any patent or other intellectual property owned or licensable by Affymetrix is conveyed or implied by this Affymetrix product. In particular, no right or license is conveyed or implied to use this Affymetrix product in combination with a product not provided, licensed, or specifically recommended by Affymetrix for such use.

Patents

Cartridge Arrays: Products may be covered by one or more of the following patents: U.S. Patent Nos. 5,445,934; 5,744,305; 5,945,334; 6,140,044; 6,261,776; 6,291,183; 6,346,413; 6,399,365; 6,420,169; 6,551,817; 6,610,482; 6,733,977; 6,955,915 and D430,024 and other U.S. or foreign patents. Products are manufactured and sold under license from OGT under 5,700,637 and 6,054,270

Array Plates: Products may be covered by one or more of the following patents: U.S. Patent Nos. 5,445,934; 5,744,305; 6,261,776; 6,291,183; 6,346,413; 6,399,365; 6,610,482; 6,733,977; and other U.S. or foreign patents. Products are manufactured and sold under license from OGT under 5,700,637 and 6,054,270.

Software: Products may be covered by one or more of the following patents: U.S. Patent Nos. 5,733,729; 5,795,716; 5,974,164; 6,066,454; 6,090,555; 6,185,561; 6,188,783; 6,223,127; 6,228,593; 6,229,911; 6,242,180; 6,308,170; 6,361,937; 6,420,108; 6,484,183; 6,505,125; 6,510,391; 6,532,462; 6,546,340; 6,687,692; 6,607,887; 7,062,092 ; 7,451,047; 6,090,555; 6,611,767; 6,687,692; 6,829,376; 7,130,458; 7,451,047 and other U.S. or foreign patents.

Reagents: Products may be protected by one or more of the following patents: U.S. Patent Nos. 6,864,059; 7,468,243; 7,491,818.

Copyright

© 2010 Affymetrix Inc. All rights reserved.

Contents

Chapter 1	Beckman Coulter Biomek FX^P Target Prep Express Configuration	1
	Equipment	1
	Related Documentation	2
	Other Equipment Required	3
	Oven	3
	Centrifuge	3
	Thermal Cyclers	3
	Microplate Reader	3
	Miscellaneous Equipment Required	3
Chapter 2	Site Preparation Guide — GeneChip[®] HT 3' IVT Express Kit Assay	5
	Equipment and Software Required	5
	Reagents and Arrays Required	5
	Labware Required for the GeneChip HT 3' IVT Express Assay	6
	Photographs of the Labware Required	7
	Labware Templates	11
	Tube Block / Cold Block Template	11
	Modular Reservoir Frame Template Stickers	11
	Related Affymetrix Documentation	12
Chapter 3	Site Preparation Guide — GeneChip[®] HT WT Assay	13
	Equipment and Software Required	13
	Consumables Required	13
	Reagents Required	14
	Labware Required for the GeneChip HT WT Assay	15
	Labware Templates	21
	Reagent Cold Block Template	21
	Modular Reservoir Frame Template Stickers	21
	Related Affymetrix Documentation	22

Beckman Coulter Biomek FX^P Target Prep Express Configuration

This chapter details the configuration required for the Beckman Coulter Biomek[®] FX^P TPE System for expression assays.

Equipment

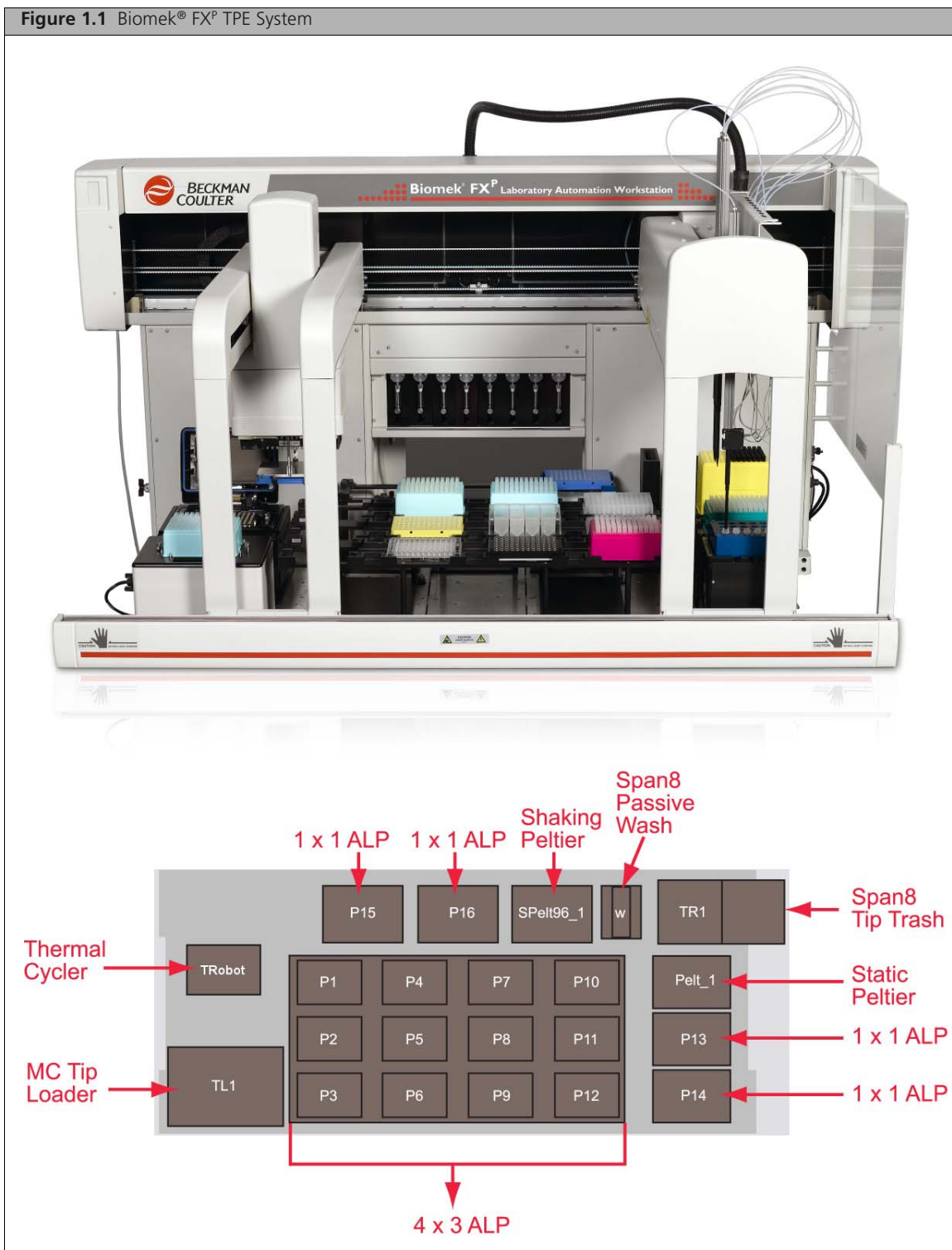
The Biomek FX^P Target Prep Express (P/N A83103) is required to run GeneChip HT expression assays. This workstation includes the accessories, software and deck configuration listed in [Table 1.1](#). For information on upgrading a Biomek liquid handler already at your site, contact Beckman Coulter. In addition to this workstation, also required are the components listed in [Chapter 2](#), if running the GeneChip HT 3' IVT Express Assay, or [Chapter 3](#) if running the GeneChip HT WT Assay.

Table 1.1 Biomek FX^P Target Prep Express required for running the GeneChip[®] HT Methods for Expression

✓	Biomek FX ^P Target Prep Express	Quantity	Beckman Coulter Part Number
□	Biomek FX^P Target Prep Express The part number for this workstation includes the components listed below.	1	A83103
	■ Biomek FX ^P Dual Multichannel Span-8	1	A31844
	■ Automation Controller XP, Biomek system software and monitor	1	A16170
	■ Packaged Tip Loader	1	719356
	■ Span-8 Disposal ALP	1	719590
	■ Disposable Tip Hardware Kit - 8 pk	1	A39380
	■ BFX 96-channel Disp, Tip Head, 200U	1	719368
	■ Pkg Static ALP Platform	4	719357
	■ 4 X 3 ALP Kit	1	719948
	■ Adaptor, Deep Well Plate	1	A83050
	■ Tube Block	1	A83054
	■ Integrated Solutions Components Ref Quote 14202A	1	
	■ Integrated Solution – Class 5 Peltier ALP Biomek FX ^P v3.3 Integration - 1 unit	1	969125
	■ Integrated Solution – Class 7 Peltier Shaking ALP FX ^P v3.3 Integration - 1 unit	1	969126
	■ Integrated Solution – Class 2 Special Lid Covers - 2 units	1	
	■ Required: Integrated Solution – Class 5 Biometra TRobot FX ^P 3.3 integration - 1 unit	1	969125

Related Documentation

- *Biomek® Liquid Handler User's Manual*, Beckman Coulter P/N 987834
- *Biomek® FX Laboratory Automation Workstation Preinstallation Manual*, Beckman Coulter P/N 719450, rev AH
- *Biomek® Software User's Manual*, Beckman Coulter P/N 987835
- *TRobot Thermocycler Instruction Manual*, biometra



Other Equipment Required

Oven

GeneChip[®] Hybridization Oven 645: Affymetrix, P/N 00-0331

Centrifuge

Eppendorf centrifuge 5810R to spin 96 well PCR plates

Eppendorf centrifuge 5415C to spin 1.5 mL to 2 mL microcentrifuge tubes

Thermal Cyclers

A Biomek FX^P Target Prep Express robot with an on-deck Biometra TRobot 96 Thermal cycler (P/N 050-991) is required. The Biometra Thermal cycler should be purchased directly from the local Biometra distributor (<http://www.biometra.com/distributors.0.html>). The Biometra TRobot 96 integration kit must be purchased from Beckman Coulter Inc.

Microplate Reader

We recommend using the microplate reader listed in [Table 1.2](#).

Table 1.2 Microplate Reader

✓	Item	Supplier	Part Number
☐	DTX 880 Multimode Detector with Genomic Filter Slide	Beckman Coulter	987921 – detector A30184 – filter slide

Miscellaneous Equipment Required

The remaining equipment required is listed in [Table 1.3](#).

Table 1.3 Other Equipment Required

✓	Item	Supplier	Part Number
☐	P1000 Rainin pipetor with correct pipette tips		
☐	Benchtop Cooler, –20°C	Stratagene	400012
☐	Freezer, –20°C	Your choice	—
☐	Ice bucket, 4 to 9 liters	Any vendor	—
☐	Microfuge (for tubes and strip tubes)	Any vendor	—
☐	Refrigerator, 2 to 8°C	Any vendor	—
☐	Vortex-Genie 2 (or equivalent; for plates and microtubes)	Scientific Industries	SI-0236 (120 V/60 Hz) SI-0246 (230 V/50 Hz)

Site Preparation Guide — GeneChip® HT 3' IVT Express Kit Assay

This site preparation guide includes the supplier and part number information for the equipment, software, reagents, arrays, labware and other consumables that have been validated for use with the GeneChip® HT 3' IVT Express Kit used on the Beckman Coulter Biomek® FX^P TPE System.

Equipment and Software Required

Table 2.1 Affymetrix equipment, software, reagents and arrays required

✓	Item	Part Number
Equipment		
	Beckman Coulter Biomek FX ^P Target Prep Instrument (see Chapter 1)	Beckman Coulter, A83103
□	GeneTitan® Instrument*	Contact Affymetrix, Inc.
Software		
□	GeneChip® Command Console (AGCC)	version 3.0 or later

*Required if running the GeneTitan setup for array plates.

Reagents and Arrays Required

Table 2.2 Affymetrix equipment, software, reagents and arrays required

✓	Item	Part Number
Reagents (array plates)		
□	GeneChip® HT 3' IVT Express Kit Kit packaging options: <ul style="list-style-type: none"> ■ 1 x 96 reaction Kit ■ 4 x 24 reactions Kit 	901253 901225
□	GeneChip® HT Hybridization, Wash and Stain Kit for GeneTitan™ and 3' IVT Express	901530
Reagents (cartridge arrays)		
□	GeneChip® 3' IVT Express Kit Kit packaging options: <ul style="list-style-type: none"> ■ Sufficient for 10 Reactions ■ Sufficient for 30 Reactions 	901228 901229
□	GeneChip® Hybridization, Wash and Stain Kit	900720
Arrays		
□	Any GeneChip 3' IVT cartridge array or Affymetrix 3' IVT array plate may be used with this assay.	see website

Labware Required for the GeneChip HT 3' IVT Express Assay

The labware required to run the GeneChip HT 3' IVT Express Assay on the Biomek FX^P Target Prep Express is listed [Table 2.3](#). Photographs of the labware are provided in [Table 2.5 on page 7](#).

Table 2.3 Labware for the Biomek FX^P Target Prep Express

✓	Item	Quantity Required 24-Array Plate Run	Quantity Required 96-Array Plate Run	Supplier	Part Number
Pipette Tips, barrier, 96 tips/rack					
☐	Biomek Span P50, pre-sterile, barrier	365 tips	576 tips		A21586
☐	Biomek AP96, P250, pre-sterile, barrier	480 tips 576 tips*	480 tips 576 tips*	Beckman Coulter	717253
☐	Biomek Span P250, pre-sterile, barrier	85 tips	59 tips		379503
☐	Biomek Span P1000, pre-sterile, barrier, conductive	47 tips 71 tips*	69 tips 93 tips*		987925
Labware					
Affymetrix Labware Kit for IVT Express Method on Beckman Biomek FX ^P Target Prep Instrument [†] , consists of:				Affymetrix	901561
☐	Plate, 1.2 mL Square Well Storage, Low Profile (4)	1	1		
☐	Plate, Bio-Rad Hardshell PCR (16)	2	2		
☐	Plate, Costar Brand Serocluster U-bottom (8)	2	2		
☐	Plate, UV Star Greiner (8)	2	2		
☐	Plate, 96 Well, Conical, PP, 340 µL/well (4)	1	1		
☐	Tube, Conical Bottom, 2.0mL, Screwcap, Sterile (24)				
☐	Lid, Autosealing Microplates, Arched (1)	1	1		
☐	Pad, Microseal, P (1)	1	1		
Reservoirs, modular for reagents					
☐	Frame for Modular Reservoirs	2	2		372795
☐	Quarter Reservoir	8 11*	8 11*	Beckman Coulter	372790
☐	Quarter Reservoir, Divided by Width	3	3		372792

*Quantity needed if running GeneTitan setup.

[†]The labware kit contains consumables sufficient for 4x24 or 4x96 rxn runs. Kit quantities shown in parenthesis.

Table 2.4 Miscellaneous Reagents and Materials for Target Preparation

✓	Item	Supplier	Part Number
☐	DNAZap	Ambion	9890
☐	RNaseZap	Ambion	9786
☐	100% ETOH	various	
☐	Cold Reagent Block template: <ul style="list-style-type: none"> ■ 24 samples ■ 96 samples 	Affymetrix	15-0431 15-0432
☐	Affymetrix® GeneChip® HT 3' IVT Reservoir Template	Affymetrix	90-0905
☐	SPRIPlate 96R Super Magnet	Agencourt	A32782

Photographs of the Labware Required

The photographs in [Table 2.5](#) are provided to help visualize the labware required on the Biomek FX^P Target Prep Express to run the target preparation portion of the GeneChip HT 3' IVT Express Assay.

Table 2.5 Labware used on the Biomek Workstation deck

Labware	Supplier and Part Number	Labware Image
Biomek AP96 – P250 Pipette Tips (aqua box; pre-sterile, barrier)	Beckman Coulter P/N 717253	
Biomek Span P1000 Pipette Tips (yellow box; pre-sterile, barrier, conductive)	Beckman Coulter P/N 987925	

Table 2.5 Labware used on the Biomek Workstation deck

Labware	Supplier and Part Number	Labware Image
<p>Biomek Span P250 Pipette Tips (green box; pre-sterile, barrier)</p>	<p>Beckman Coulter P/N 379503</p>	
<p>Biomek Span P50 Pipette Tips (magenta box; pre-sterile, barrier)</p>	<p>Beckman Coulter P/N A21586</p>	
<p>Plate, Bio-Rad Hardshell PCR (available in multiple colors)</p>	<p>Affymetrix P/N 901561 (Part of the Labware Kit for IVT Express Method on Beckman Biomek FX^P Target Prep Instrument)</p>	
<p>Plate, 1.2 mL Square Well Storage, Low Profile</p>	<p>Affymetrix P/N 901561 (Part of the Labware Kit for IVT Express Method on Beckman Biomek FX^P Target Prep Instrument)</p>	

Table 2.5 Labware used on the Biomek Workstation deck

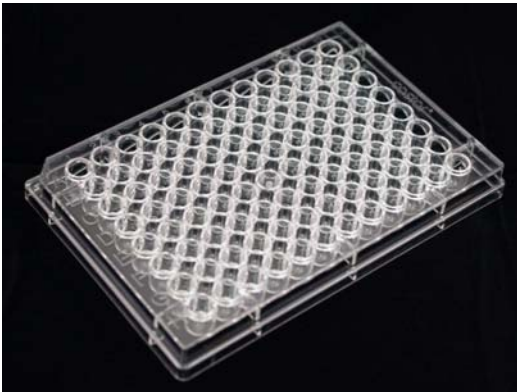



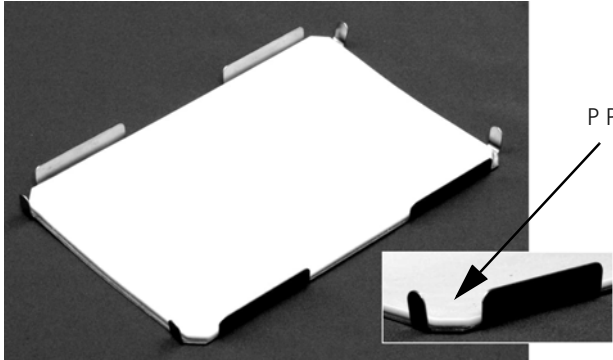

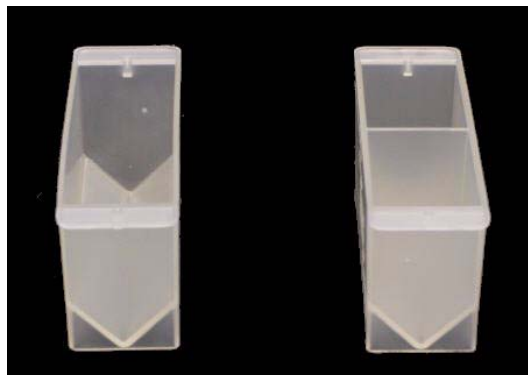
Labware	Supplier and Part Number	Labware Image
Plate, Costar Brand Serocluster U-bottom	Affymetrix P/N 901561 (Part of the Labware Kit for IVT Express Method on Beckman Biomek FX [®] Target Prep Instrument)	
Plate, UV Star Greiner	Affymetrix P/N 901561 (Part of the Labware Kit for IVT Express Method on Beckman Biomek FX [®] Target Prep Instrument)	
Plate, 96 Well, Conical Bottom, Polypropylene	Affymetrix P/N 901561 (Part of the Labware Kit for IVT Express Method on Beckman Biomek FX [®] Target Prep Instrument)	
Tube, Conical Bottom, 2.0mL, Screwcap, Sterile	Affymetrix P/N 901561 (Part of the Labware Kit for IVT Express Method on Beckman Biomek FX [®] Target Prep Instrument)	

Table 2.5 Labware used on the Biomek Workstation deck

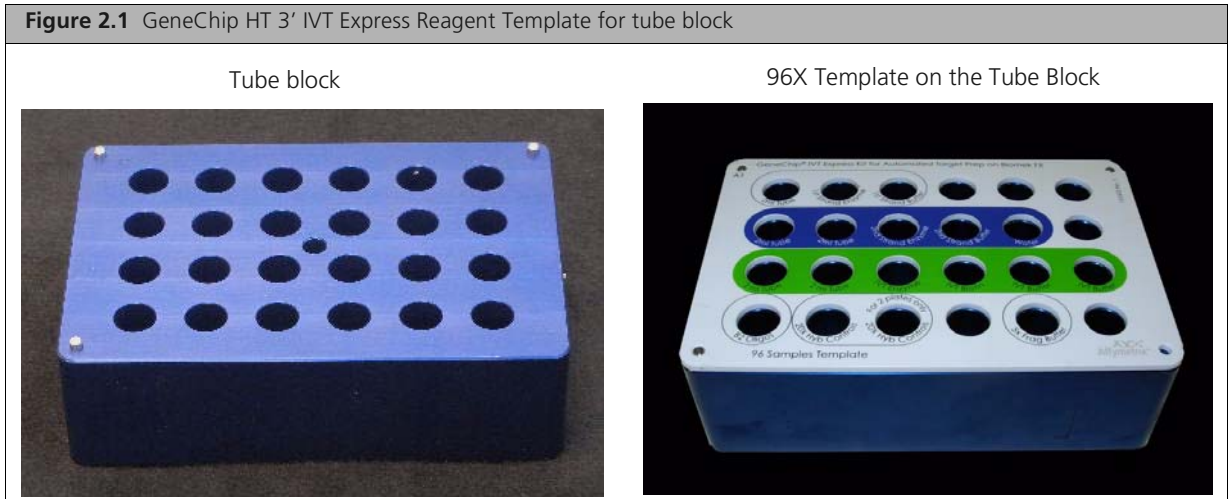
Labware	Supplier and Part Number	Labware Image	
<p>Lid, Autosealing Microplates, Arched with Microseal, P Pad</p>	<p>Affymetrix P/N 901561 (Part of the Labware Kit for IVT Express Method on Beckman Biomek FX^P Target Prep Instrument)</p>		
<p>Quarter Reservoir Frame</p>	<p>Beckman Coulter P/N 372795</p>		
<p>Quarter Reservoirs</p> <ul style="list-style-type: none"> ■ Quarter module, 40 mL capacity ■ Quarter module divided by width, 19 mL capacity each receptacle 	<p>Beckman Coulter</p> <p>P/N 372790 (40 mL) P/N 372792 (19 mL)</p>	<p>Undivided 40 mL capacity</p>	<p>Divided by width 19 mL capacity</p>



Labware Templates

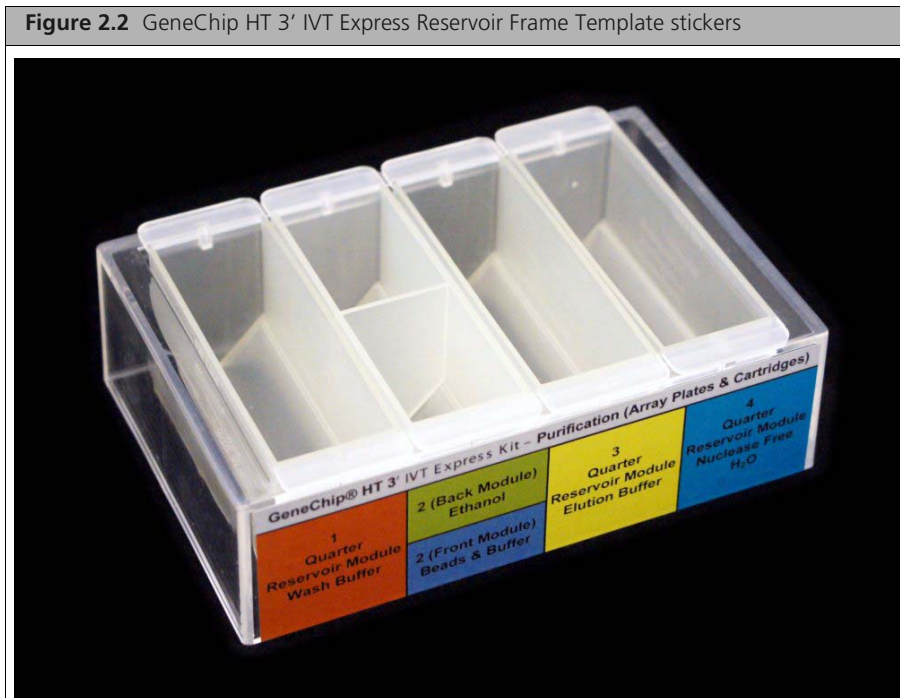
Tube Block / Cold Block Template

The GeneChip HT 3' IVT Express Kit cold block template was designed to fit precisely onto the top of the Beckman Coulter Tube Block (P/N A83054). The template is held in place by the metal posts on the block (Figure 2.1). Use template P/N 15-0431 for 24X runs and template P/N 15-0432 for 96X runs to ensure the proper placement of reagent tubes onto the block for each method.



Modular Reservoir Frame Template Stickers

Template stickers are used with the modular reservoir frame to ensure proper placement of reagents in the reservoirs. Template stickers are available for the (1) Purification, (2) Normalization & Hybridization, and (3) GeneTitan Setup methods.



Related Affymetrix Documentation

- *GeneChip® HT 3' IVT Express Kit User Manual*, P/N 702662
- *GeneChip® Expression Wash, Stain and Scan User Manual for Cartridge Arrays*, P/N 702731
- *GeneTitan® Instrument User Guide for Expression Array Plates*, P/N 702933
- *Affymetrix GeneTitan® Instrument User's Guide*, P/N 08-0296
- *Affymetrix GeneTitan® Instrument Site Preparation Guide*, P/N 08-0300
- *Affymetrix GeneChip® Command Console™ User Manual*, P/N 702569

Site Preparation Guide — GeneChip® HT WT Assay

This site preparation guide includes the supplier and part number information for the equipment, software, reagents, arrays, labware and other consumables that have been validated for use with the GeneChip® HT WT Assay used on the Beckman Biomek® FX^P TPE System.

Equipment and Software Required

Table 3.1 Affymetrix equipment, software, reagents and arrays required

✓	Item	Part Number
Equipment		
	Beckman Coulter Biomek FX ^P Target Prep Instrument (see Chapter 1)	Beckman Coulter, A83103
□	GeneTitan® Instrument*	Contact Affymetrix, Inc.
Software		
□	GeneChip® Command Console (AGCC)	version 3.0 or later

*Required if running the GeneTitan setup for array plates.

Consumables Required

Table 3.2 Affymetrix Installation Kit

✓	Component	Quantity	Source	P/N
□	Affymetrix Installation Kit consists of:		Affymetrix	90-0931
	■ HT WT Reservoir Stickers	1 set		
	■ HT WT Cold Block Template	1		
	■ Alpaqua Magnetic Plate	1		
	■ Alpillio® Plate Cushion	1		
	■ WT ATP Biomek method file (and support files)	Installed by Affymetrix FAS		
	■ Arched Auto-Sealing Lids With Wide Tabs	Pkg (Pkg of 4)		

Table 3.3 Affymetrix® and Ambion® HT WT Expression Assay Consumables for Biomek® FX^P Target Prep Express

✓	Component	Number per Kit	Source	P/N
□	Affymetrix® and Ambion® HT WT Expression Assay Consumables for Biomek® FX ^P Target Prep Express - Labware sufficient for 4x24 or 4x96 Rxn runs consists of:		Affymetrix	901646
	■ 96-Well Deep Well Microplate, polypropylene, 2 mL, non-sterile	8		
	■ Greiner U-bottom (round) Plate, polypropylene	8		
	■ Bio-Rad Hard Shell PCR Plate, blue shell/clear well	72		
	■ Greiner UV Star Plate, non-sterile, clear, flat bottom	8		
	■ Costar 96-Well U-bottom plate, polystyrene	16		
	■ 24-Well Plate, 10 mL, square well, pyramid bottom, polypropylene	4		
	■ Auto-sealing Microplate Lid with Microseal P Pad	2		

Table 3.4 Biomek® Tips and Universal labware

✓	Product	Packaging	Total 8x	Total 16x	Total 24x	Total 96x	Source	P/N
□	Biomek Span P50 Barrier Tips (Span 8 Pipettor) (magenta)	96 ea./box- 10 box/case	10 Boxes	10 Boxes	10 Boxes	10 Boxes	Beckman Coulter	A21586
□	Biomek AP96 – P250 Barrier Tips Sterile (AP 96 Multichannel Pipettor) (aqua)	96 ea./box - 10 box/case	8 Boxes	8 Boxes	8 Boxes	8 Boxes	Beckman Coulter	717253
□	Biomek Span P250 Barrier Tips (Span 8 Pipettor) (green)	96 ea./box - 10 box/case	1 Box	1 Box	1 Box	1 Box	Beckman Coulter	379503
□	Biomek Span P1000 LLS Barrier Tips (Span 8 Pipettor) (yellow)	96 ea./box - 5 box/case	2 Boxes	2 Boxes	2 Boxes	2 Boxes	Beckman Coulter	987925
□	Quarter Reservoir, Divided by Width	40 unit/case	4	4	4	4	Beckman Coulter	372792
□	Quarter Reservoir	40 unit/case	8	8	8	8	Beckman Coulter	372790
□	Frame for Reservoirs	1 ea.	reusable		reusable		Beckman Coulter	372795

Reagents Required

Table 3.5 HT WT Express Reagents, Miscellaneous Reagents, and Materials

✓	Product	Source	P/N
□	Ambion® WT Expression Kit for High Throughput Robotics Kit packaging options: <ul style="list-style-type: none"> ■ 1 x 24 reaction Kit* ■ 1 x 96 reaction Kit 	Ambion	4440537 4440536
□	Affymetrix® HT WT Terminal Labeling and Controls Kit [†] (Sufficient for 24 Rxn) Contains: <ul style="list-style-type: none"> ■ GeneChip® HT WT Terminal Labeling Kit, P/N 901555 (1 x 24 Rxn Kit) ■ GeneChip® Hybridization Control Kit, P/N 900454 (30 Rxn) ■ GeneChip® Poly-A Control Kit, P/N 900433 (100 Rxn) 	Affymetrix	901647
□	Affymetrix® HT WT Terminal Labeling and Controls Kit [†] (Sufficient for 96 Rxn) Contains: <ul style="list-style-type: none"> ■ GeneChip® HT WT Terminal Labeling Kit, P/N 901518 (1 x 96 Rxn Kit) ■ GeneChip® Hybridization Control Kit, P/N 900457 (150 Rxn) ■ GeneChip® Poly-A Control Kit, P/N 900433 (100 Rxn) 	Affymetrix	901648
□	GeneChip® Hybridization, Wash and Stain Kit – (30 Rxns, for cartridges)	Affymetrix	900720
□	GeneTitan® Hybridization, Wash and Stain Kit for WT Array Plates (96 Rxns)	Affymetrix	901622
□	DNAZap	Ambion	9890
□	RNaseZap	Ambion	9786
□	100% ETOH	various	
□	Isopropanol	various	

*One 24 rxn kit is good for (2) 8x runs, (1) 16x run or (1) 24x run.

[†]Individual components available for purchase separately.

Labware Required for the GeneChip HT WT Assay

Table 3.6 shows the types of labware used in the GeneChip HT WT Assay. For labware requirements specific to each part of the assay, refer to the GeneChip® HT WT User Guide (P/N 702923).

Table 3.6 Labware used on the Biomek Workstation Deck

Labware	Supplier & Part Number	Labware Image
Biomek AP96 – P250 Barrier Tips (aqua box; pre-sterile, barrier)	Beckman Coulter P/N 717253	
Biomek Span P1000 LLS Barrier Tips (yellow box; pre-sterile, barrier, conductive)	Beckman Coulter P/N 987925	
Biomek Span P250 Barrier Tips (green box; pre-sterile, barrier)	Beckman Coulter P/N 379503	
Biomek Span P50 Pipette Tips (magenta box; pre-sterile, barrier)	Beckman Coulter P/N A21586	

Table 3.6 Labware used on the Biomek Workstation Deck (Continued)

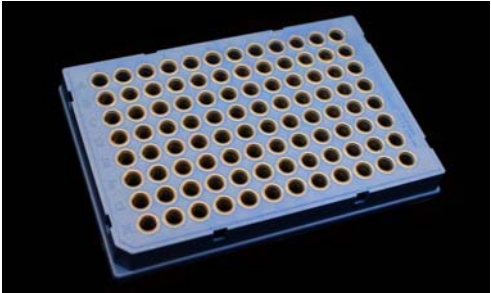
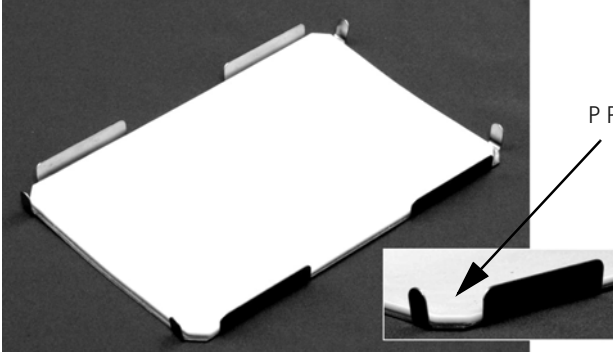


Labware	Supplier & Part Number	Labware Image
Bio-Rad Hard Shell 96-Well Plate (available in multiple colors)	Bio-Rad P/N HSP9631 (A component of the Affymetrix® and Ambion® HT WT Expression Assay Consumables for Biomek® FX ^P Target Prep Express Kit P/N 901646. Kit = 72)	
Lid, Autosealing Microplates, Arched with Microseal, P Pad	Bio-Rad P/N MSL-2032 (A component of the Affymetrix® and Ambion® HT WT Expression Assay Consumables for Biomek® FX ^P Target Prep Express Kit P/N 901646. Kit = 2)	
AbGene 96 Square Well Storage Plate (2.2 mL)	AbGene P/N AB-0932 (A component of the Affymetrix® and Ambion® HT WT Expression Assay Consumables for Biomek® FX ^P Target Prep Express Kit P/N 901646. Kit = 8)	
24 well mixing plate, 10 mL, square well, pyramid bottom	Seahorse Bioscience P/N S30024 (A component of the Affymetrix® and Ambion® HT WT Expression Assay Consumables for Biomek® FX ^P Target Prep Express Kit P/N 901646. Kit = 4)	

Table 3.6 Labware used on the Biomek Workstation Deck (Continued)





Labware	Supplier & Part Number	Labware Image
Greiner UV Star Plate	Greiner P/N 655801 (A component of the Affymetrix® and Ambion® HT WT Expression Assay Consumables for Biomek® FX ^P Target Prep Express Kit P/N 901646. Kit = 8)	
Greiner U-Bottom (Round), Polypropylene Plate	Greiner P/N 650201 (A component of the Affymetrix® and Ambion® HT WT Expression Assay Consumables for Biomek® FX ^P Target Prep Express Kit P/N 901646. Kit = 8)	
Costar Polystyrene 96-well, U-bottom Plate (non-sterile)	Costar P/N 3798 (A component of the Affymetrix® and Ambion® HT WT Expression Assay Consumables for Biomek® FX ^P Target Prep Express Kit P/N 901646. Kit = 16)	
Alpaqua Magnetic Plate	Part of the HT WT Installation Kit from Affymetrix, P/N 90-0931	

Table 3.6 Labware used on the Biomek Workstation Deck (Continued)






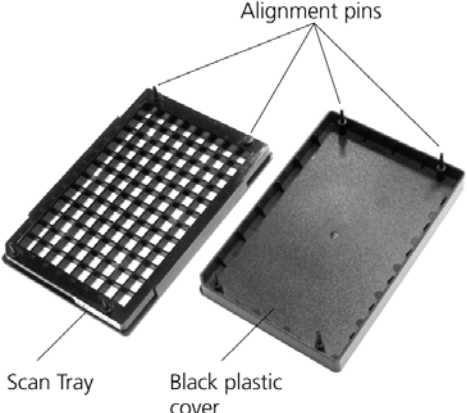
Labware	Supplier & Part Number	Labware Image																																
<p>Modular Reservoir Template Stickers</p> <p>Modular reservoir template stickers are available to place on the front of the modular reservoir frame to assist with the proper placement of modular reservoirs and reagents. See Figure 3.2 for further detail.</p>	<p>Affymetrix P/N 90-0907</p> <p>Part of the HT WT Installation kit from Affymetrix, P/N 90-0931</p>	<table border="1"> <thead> <tr> <th colspan="4">Affymetrix® and Ambion® WT Expression Modular Reservoir #1</th> </tr> </thead> <tbody> <tr> <td>1 Quarter Reservoir Module Water</td> <td colspan="3">NO Reservoir Module</td> </tr> </tbody> </table> <table border="1"> <thead> <tr> <th colspan="4">Affymetrix® and Ambion® WT Expression Modular Reservoir #2</th> </tr> </thead> <tbody> <tr> <td rowspan="2">1 Quarter Reservoir Module Water</td> <td rowspan="2">2 Quarter Reservoir Module Wash Solution</td> <td>3 (Back Module) Empty</td> <td>4 (Back Module) Elution Solution</td> </tr> <tr> <td>3 (Front Module) Binding Buffer</td> <td>4 (Front Module) IPA / EtOH</td> </tr> </tbody> </table> <table border="1"> <thead> <tr> <th colspan="3">Affymetrix® and Ambion® WT Expression Modular Reservoir #3</th> </tr> </thead> <tbody> <tr> <td>1 Quarter Reservoir Module Water</td> <td>2 Quarter Reservoir Module 2X Hybridization Mix</td> <td>NO Reservoir Module</td> </tr> </tbody> </table> <table border="1"> <thead> <tr> <th colspan="4">Affymetrix® and Ambion® WT Expression Modular Reservoir #4</th> </tr> </thead> <tbody> <tr> <td>1 Quarter Reservoir Module Holding Buffer</td> <td>2 Quarter Reservoir Module Stain 1 & 3</td> <td>3 Quarter Reservoir Module Stain 2</td> <td>NO Reservoir Module</td> </tr> </tbody> </table>	Affymetrix® and Ambion® WT Expression Modular Reservoir #1				1 Quarter Reservoir Module Water	NO Reservoir Module			Affymetrix® and Ambion® WT Expression Modular Reservoir #2				1 Quarter Reservoir Module Water	2 Quarter Reservoir Module Wash Solution	3 (Back Module) Empty	4 (Back Module) Elution Solution	3 (Front Module) Binding Buffer	4 (Front Module) IPA / EtOH	Affymetrix® and Ambion® WT Expression Modular Reservoir #3			1 Quarter Reservoir Module Water	2 Quarter Reservoir Module 2X Hybridization Mix	NO Reservoir Module	Affymetrix® and Ambion® WT Expression Modular Reservoir #4				1 Quarter Reservoir Module Holding Buffer	2 Quarter Reservoir Module Stain 1 & 3	3 Quarter Reservoir Module Stain 2	NO Reservoir Module
Affymetrix® and Ambion® WT Expression Modular Reservoir #1																																		
1 Quarter Reservoir Module Water	NO Reservoir Module																																	
Affymetrix® and Ambion® WT Expression Modular Reservoir #2																																		
1 Quarter Reservoir Module Water	2 Quarter Reservoir Module Wash Solution	3 (Back Module) Empty	4 (Back Module) Elution Solution																															
		3 (Front Module) Binding Buffer	4 (Front Module) IPA / EtOH																															
Affymetrix® and Ambion® WT Expression Modular Reservoir #3																																		
1 Quarter Reservoir Module Water	2 Quarter Reservoir Module 2X Hybridization Mix	NO Reservoir Module																																
Affymetrix® and Ambion® WT Expression Modular Reservoir #4																																		
1 Quarter Reservoir Module Holding Buffer	2 Quarter Reservoir Module Stain 1 & 3	3 Quarter Reservoir Module Stain 2	NO Reservoir Module																															
<p>24-Position Tube Rack</p>	<p>Beckman Coulter P/N 373661 (rack) P/N 373696 (insert)</p>	<p>Tube rack with 24 inserts</p> 																																
<p>Adaptor, Thin PCR (installed on the Shaking Peltier)</p> <p>This adaptor is typically installed by a Beckman Coulter field service technician during new system installation or a system upgrade. Ensure that you have one of these adaptors on the deck prior to running this assay.</p>	<p>Beckman Coulter P/N A19709</p>																																	

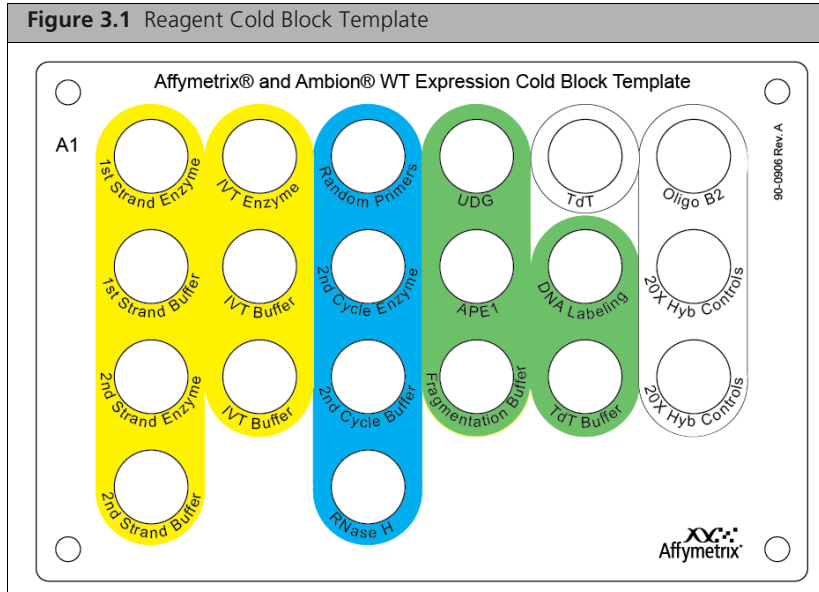
Table 3.6 Labware used on the Biomek Workstation Deck (Continued)

Labware	Supplier & Part Number	Labware Image
Alpillo Plate Cushion	Part of the HT WT Installation kit from Affymetrix, P/N 90-0931	 <p>A blue metal frame with four corner brackets. The word 'Alpillo' is printed vertically on the right side.</p>
GeneTitan Hyb Tray	Affymetrix P/N 500867 (GeneTitan trays and lids are provided with the Array Plate when ordered. The P/N provided here is for identification purposes only.)	 <p>A clear plastic tray with a grid of wells. A barcode and part number are visible on the bottom edge.</p>
GeneTitan Stain Tray	Affymetrix P/N 501025 (GeneTitan trays and lids are provided with the Array Plate when ordered. The P/N provided here is for identification purposes only.)	 <p>A clear plastic tray with a grid of wells. A barcode and part number are visible on the bottom edge.</p>
GeneTitan Scan Tray	Affymetrix P/N 501006 or 500860 (GeneTitan trays and lids are provided with the Array Plate when ordered. The P/N provided here is for identification purposes only.)	 <p>Two components are shown: a black grid-like Scan Tray and a Black plastic cover. Alignment pins are indicated by lines pointing to the cover.</p>

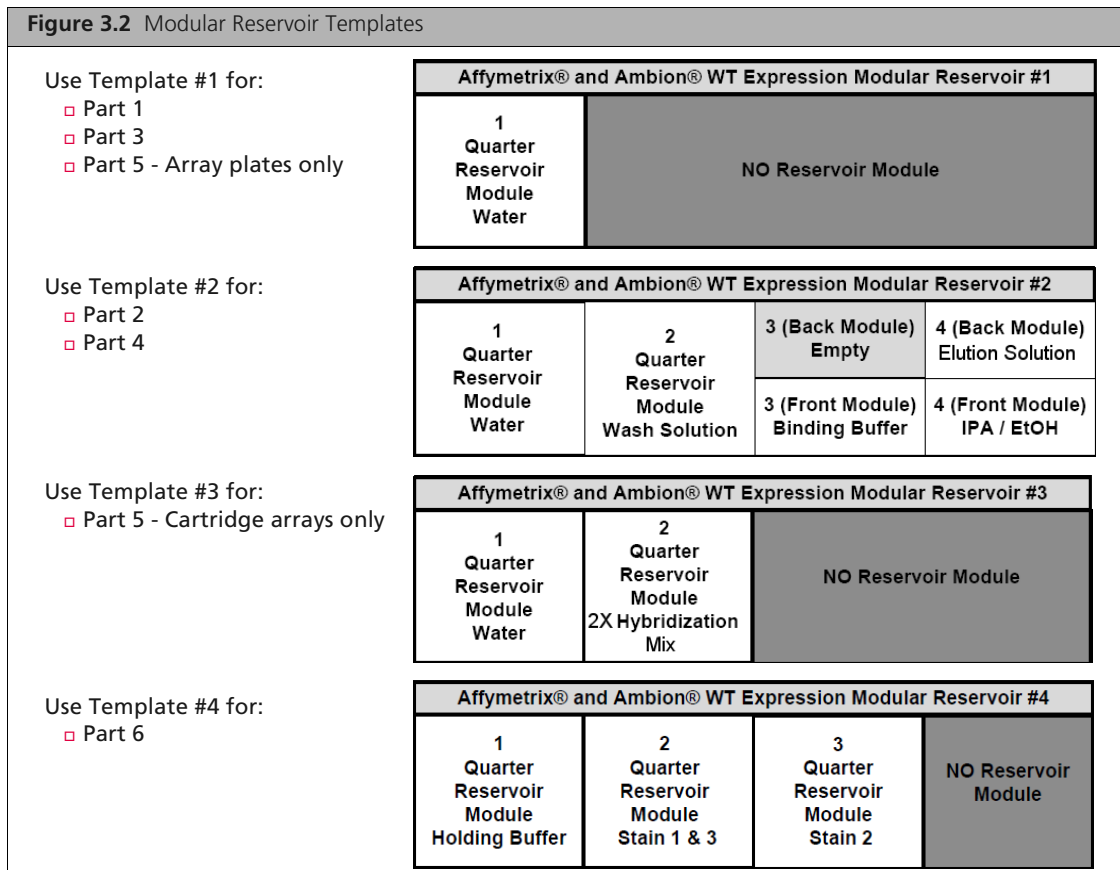
Labware Templates

The Ambion® WT Expression Kit and GeneChip® HT WT Terminal Labeling Kit provide templates to help ensure the proper placement of reagents in labware.

Reagent Cold Block Template



Modular Reservoir Frame Template Stickers



Related Affymetrix Documentation

- *GeneChip® HT WT User Guide*, P/N 702923
- *GeneChip® Expression Wash, Stain and Scan User Manual for Cartridge Arrays*, P/N 702731
- *Affymetrix GeneTitan® Instrument User's Guide*, P/N 08-0296
- *Affymetrix GeneTitan® Instrument Site Preparation Guide*, P/N 08-0300
- *Affymetrix GeneChip® Command Console™ User Manual*, P/N 702569