

# The Freedom EVO<sup>®</sup> 75 GenePaint™ Application Configuration

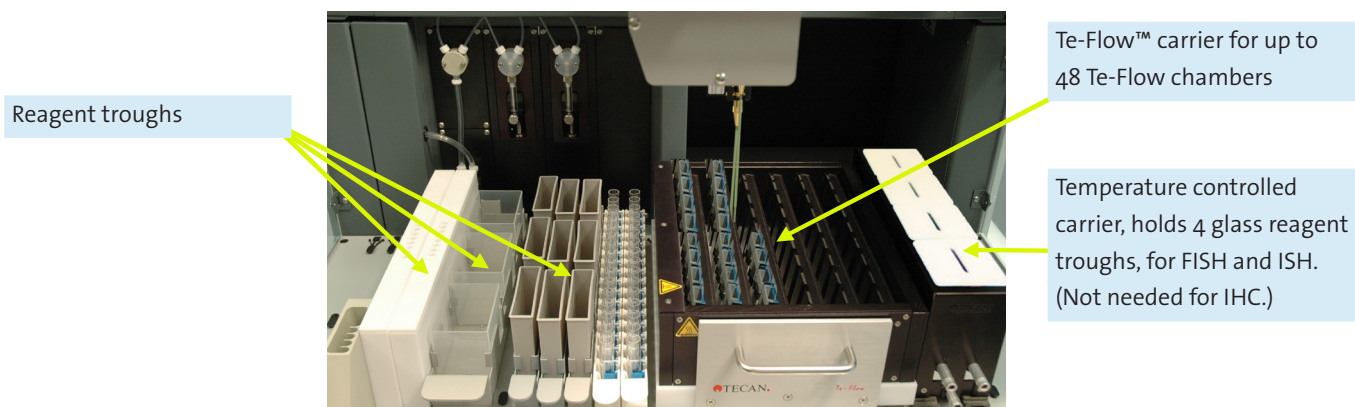
## Automated System for In Situ Hybridization (ISH), Fluorescence In Situ Hybridization (FISH), and Immunohistochemistry (IHC)

Reference 10003BF100597



Tecan's GenePaint system automates all of the steps required for performing in situ hybridization of tissues on microscope slide formats. One of the key steps in the study of gene expression is the validation of results obtained by in vitro gene expression methods (microarrays, blots, etc.). This validation typically involves the use of appropriate cell or tissue samples and the performance of in situ hybridization (ISH) to confirm the in vitro results. The unique advantage of in situ hybridization over destructive expression techniques such as microarrays is the localization of transcripts in intact cells and tissues, giving a much higher level of information.

Processing of large numbers of cell or tissue samples in these type of procedures is an extremely time consuming and laborious process. GenePaint alleviates this throughput bottleneck with the elimination of 80% of the manual steps of conventional ISH, FISH or IHC analysis. All fixing, staining, washing and hybridization processes of the slides are handled by the GenePaint system. The Freedom EVO 75 ensures reproducible and fast liquid transfer from reagent troughs to flow-through chambers.



## Suggested Configuration

Part Number	Quantity	Product Description
<b>Tecan Robotic &amp; Liquid Handling Hardware and Software</b>		
30025019	1 PC	Instrument Freedom EVO 75 base
30025021	1 PC	Right safety panel
30025026	1 PC	Liquid handling arm 2-channel
10643050	2 PC	Tips fixed washable for 96 plate
10619426	2 PC	Syringe 2.5 ml
30025868	1 PC	Wash station for 2 tips
30025024	2 PC	Fast tip wash option
30032130	1 PC	Back panel with recess (for rear routing of Te-Flow tubing)
<b>Tecan GenePaint Basics</b>		
10760810	48 PC	Chamber Flow through Te-Flow
10760800	1 PC	Rack chamber 48-position Te-Flow
10760825	2 PC	Plate mounting Te-Flow, 8 grids
10760815	1 PC	Spacers 500 pieces
10760949	1 PC	Manual Operating Te-Flow
<b>GenePaint Accessories</b>		
30014666	2 PC	Carrier trough and cover 350 ml
10613020	3 PC	Carrier additive trough 100 ml
30019985	2 PC	Carrier tubes 16-position, tube 10 mm
30016581	1 PC	Carrier trough reagent 320 ml, 3-position, incl. 3 troughs with lids
10613049	1 PC	Trough disposable 100 ml, 108 pieces
30023119	1 PC	Trough 320 ml, 5 pieces
<b>Tecan Temperature Control Items</b>		
30017432	1 PC	Carrier Thermo controlled 4 glass troughs 250 ml (Not needed for IHC)
10760831	1 PC	Thermometer Te-Flow
30017604 (EU)	1 PC	Thermostated recirculator Lauda Ecoline 205 230V
10760938	1 PC	Tubing set Te-Flow Lauda Integra T2200
10760939	1 PC	Quick coupling Te-Flow
10760035 (EU)	1 PC	Thermostated recirculator Lauda Ecoline 203 230V (not for IHC)
10626106	1 PC	Cable RS232 (for 10760035, not needed for IHC)
<b>Tecan Waste Disposal Items</b>		
30011583 (EU)	1 PC	Peristaltic pump Heidolph PD 5201
10760855	1 PC	Option head pump quick SP Heidolph
10760940	1 PC	Tubing set for Te-Flow chamber carrier waste
<b>Optional Accessories</b>		
10619677	1 PC	Bottle system liquid 10 L
10619819	1 PC	Bottle waste 20 L

Note: Waste and system liquid containers, PC and monitor may also be ordered from Tecan

Scientific instrumentation; not for use in human, clinical or diagnostic procedure.

Tecan is in major countries a registered trademark and Freedom EVO 75, Flow-Thru Technology, Freedom EVOware Standard, TouchTools Suite, Instant Pipetting, Freedom EVOware Normalization Wizard, PosID are trademarks of Tecan Group Ltd., Männedorf, Switzerland.  
© 2009, Tecan Trading AG, Switzerland, all rights reserved.

Please contact your local dealer for more information. [www.tecan.com](http://www.tecan.com), p. 2/2