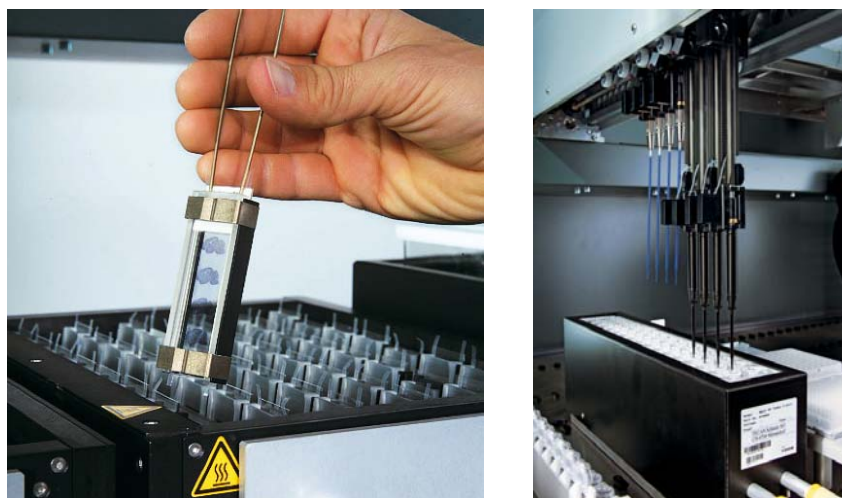


GenePaint™

An innovative solution for automated *in situ* hybridization (ISH), fluorescence *in situ* hybridization (FISH) and immunohistochemistry (IHC)

Reference 10001BO100394



GenePaint system for *in situ* hybridization (1-96 slides)

Tecan's GenePaint system, based on a new flow-through technology, automates all of the steps required for performing *in situ* hybridization (ISH) 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 performing ISH on appropriate cell or tissue samples to confirm the *in vitro* results.

The unique advantage of ISH 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 using this type of procedure, however, is an extremely time-consuming and laborious process. GenePaint alleviates this throughput bottleneck.

The advanced benefits for functional genomic studies include:

- eliminating 80 % of the manual steps in conventional ISH, FISH or IHC analysis
- extending walkaway time through unattended hybridization protocols
- allowing the use of any commercial set of reagents for hybridization studies
- assuring consistent and reproducible results of excellent quality
- saving on the amounts of reagents used

The applications described here are not available in the US outside of the research market.

Scientific instrumentation; not for use in human, clinical or diagnostic procedures. All systems or products are tailor-made one-offs that were exclusively developed on customer demand. All displayed pictures and images originate from development, test or installation phases and do not claim to represent the final installation at customer site.

Tecan is in major countries a registered trademark and GenePaint is a trademark of Tecan Group Ltd., Männedorf, Switzerland.
© 2009, Tecan Trading AG, Switzerland, all rights reserved.