

HS 4800™ Pro Hybridization Station and new PowerScanner™

Ultimate hybridization reproducibility and high resolution scanning for your microarrays

Reference 10001BO100396



HS 4800 Pro Hybridization Station



PowerScanner

The HS 4800 Pro Hybridization Station provides automated processing of microarrays on slides, from pre-hybridization and on-board denaturation to automatic slide drying, with Tecan's unique OSND™ (On Slide Nitrogen Drying) Technology. The system has been designed to avoid hands-on operation and give researchers better reproducibility, sensitivity and reliability than they can expect from traditional procedures, with even less labor involved. The system offers standardization and strict control of the entire process that are key factors in enhancing data quality and consistency.

Tecan's unique patented ABS™ (Active Bubble Suppression) system effectively minimizes the risk of bubble formation during hybridization and thus hybridization artifacts. Agitation during hybridization gives higher stringency and faster kinetics, resulting in higher sensitivity and specificity, and a wider ratiometric range of signals compared to manual methods. Precise temperature control and highly uniform washing provide consistent and reproducible results between slides.

Tecan's HS 4800 Pro hybridization station is the perfect automated solution for different array and slide formats. Easily interchangeable chambers allow one instrument to offer multiple chamber formats and sizes. The novel QuadChamber™ is the first fourfold segmented chamber for fully automated processing of microarrays. It guarantees no cross-contamination between the segments and fits perfectly on Agilent 4-Pack Microarrays and other four-segment slides.

Tecan offers a seamless integration of key functions while addressing the need for high resolution and perfect data quality. Tecan's new general purpose PowerScanner performs the following functionalities: 2 µm resolution dual color in 20 minutes at top sensitivity; < 3 % CV reproducibility intra- and inter-array; dynamic range of 10⁵; multiplex up to 24 arrays per slide, random access of up to 48 slides from a compact magazine; integrated ozone filters (dye degradation prevention); compatibility with glass slides, nitrocellulose and highly reflective coatings supporting nucleic acids, proteins and tissues; ability to support almost all main third party analysis software packages, and can be ordered together with ArrayPro™ for fully automated microarray image batch analysis.

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 HS 4800 Pro, PowerScanner, OSND Technology, ABS, QuadChamber and ArrayPro are trademarks of Tecan Group Ltd., Männedorf, Switzerland.

© 2009, Tecan Trading AG, Switzerland, all rights reserved.