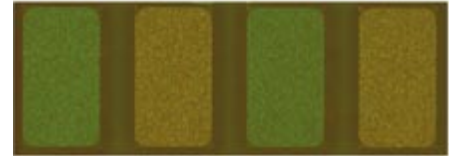


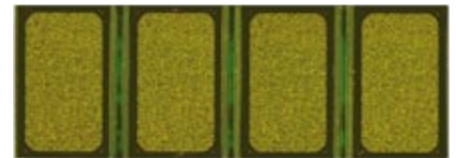
# Tecan's new QuadChamber™ is validated for use with Agilent's 4-pack Microarrays



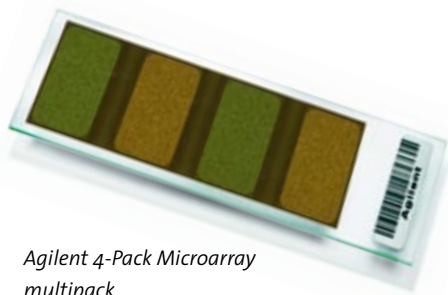
Tecan HS 4800™ with four extension units



Agilent aCGH 4-Pack Microarray



Agilent Gene Expression 4-Pack Microarray



Agilent 4-Pack Microarray multipack

Tecan has recently developed the QuadChamber for fully automated processing of four different microarrays simultaneously on one slide, using the HS Pro™ automated hybridization station. The QuadChamber was specially developed for use with Agilent's Technologies new 4 x 44k 4-Pack Gene Expression as well as CGH Microarrays, which consist of four individual, whole-genome microarrays printed on a single glass slide. This represents the first fully automated system that can independently handle four arrays on one slide with no cross-contamination between the arrays, and is the latest development in an ongoing, global collaboration between the two companies.

The QuadChamber provides a completely sealed environment around each of the

four arrays on Agilent's 4-Pack slides, with independent channels for wash buffers, independent agitation mechanisms and drying for each of the four areas. With the new chamber, the HS Pro fully automates the hybridization of the arrays from start to finish while maintaining separate compartments around each array on the slide, ensuring high quality processing with excellent reproducibility and minimal slide handling.

Agilent's 4-Pack Microarrays concentrate sample hybridization, so provide enhanced sensitivity and allow the detection of even very low abundance transcripts. Having four arrays on each slide allows researchers to generate more genomics data per experiment while considerably reducing the required volumes of samples and reagents.

Tecan and Agilent have worked together to develop a number of protocols, including automated array-based CGH, ensuring excellent quality data and results for customers straight away. "Agilent has worked closely with Tecan to develop our 4-Pack Microarrays and protocols for fully automated

hybridization," said Jeff McMillan, product manager for Genomics Automation at Agilent. "The HS Pro is the only fully automated instrument that can process our 4-Pack Microarrays."

The QuadChamber and the 4-Pack Microarrays will be presented together at a number of shows and workshops over the coming months, beginning with the Lausanne Genomics Days 2007, a symposium in Switzerland on 13-14 September, and continuing to the American Society of Human Genetics conference on 23-27 October in San Diego, USA.



Tecan's new HS Pro QuadChamber

The products described are for research use only