

PRODUCT FACT SHEET

The new Axon GenePix® 4300A and Axon GenePix® 4400A microarray scanners represent the latest in a long line of industry-leading microarray detection and analysis systems from MDS Analytical Technologies. These groundbreaking microarray instruments will enable high resolution scanning of the latest generation of high-density microscope slide-based microarrays, with the flexibility of multi-label imaging, and platform independence. The Axon GenePix 4300A and 4400A microarray scanners represent the first significant advance in microarray detection technology since the release of the GenePix Autoloader 4200AL scanner in 2004. These improvements ensure that the Axon GenePix 4300A and 4400A microarray scanners will offer the broadest applicability in the microarray industry, providing the optimal tools for all microscope slide format microarray platforms.



- **High density array scanning:** The Axon GenePix 4400A model increases the maximum optical resolution two-fold over previous microarray detection and analysis systems, with a scanning capability of 2.5 microns per pixel. This enables researchers to confidently image the upcoming generation of ultra high density microarrays from commercial providers such as Agilent Technologies and Roche-Nimblegen.
- **Reduced cost per data point:** high density microarray platforms increase the amount of data obtained from a single microarray, greatly improving the efficiency of a given experiment and significantly reducing the cost per data point by limiting the number of microarrays needed.
- **Simple upgrade path:** for researchers who do not require the high resolution, but who would benefit from the increased speed, uniformity, and sensitivity of the systems, the Axon GenePix 4300A offers 5 micron maximum resolution, which can be upgraded at any time.
- **Flexible scanner configuration:** researchers have the choice of up to four internally-installed lasers, and up to 16 emission filters for ultimate fluorophore flexibility.
- **All-new version of Axon GenePix® Pro software:** enables the acquisition and analysis of increasingly complex microarray assays.