
Sole Source Specification for Axon GenePix[®] 4200AL

From the industry-leader in performance, the Axon GenePix[®] 4200AL scanner from MDS Analytical Technologies is the only scanner available today that provides:

Features & Benefits

- Walk-Away Operation: Automatically loads and scans up to 36 slides; Batch scan and batch analysis options; Automatic PMT balancing removes need to set sensitivity parameters for each slide; E-mail notification upon completion of batch
- Flexibility in Fluorophore Selection: Houses up to 4 laser sources for excitation; Has 16-position user-accessible emission filter wheel
- Non-Confocal Technology: Provides 64 μ m depth of field; Expanded focal range is ideal for thick specimens, alternate substrates, and non-planar slide irregularities
- Sensitivity of 0.1 fluor/ μ m²: Allows detection of weak signals, while maintaining a superior signal to noise ratio
- Dynamic Range of 4 Orders of Magnitude: Provides linear results for low- and high-intensity features
- Selectable Scan Resolution: User can choose from 5 to 100 μ m resolution to optimize image quality and file size for each application
- Dynamic Laser Power Monitoring: Feedback circuit produces constant signal output throughout each scan
- Adjustable Focus: Facilitates scanning of a variety of substrate types
- Adjustable Laser Power: Allows user to optimize signal output for a variety of sample types
- Dynamic Photomultiplier Adjustment: Allows user to optimize detector voltages dynamically during image acquisition; PMTs use integrators, ensuring highest collection efficiency
- Uses Low-Noise Digitization Electronics: Provides maximum signal-to-noise ratio
- Integrated Hardware Barcode Reader: Automatically records slide ID into image file
- One License of GenePix Pro 6.0 Software with each scanner order
- Feature Viewer Provides Real-Time Results: Navigates through data while acquiring an image
- Integrated Software and Hardware: Ensures seamless communication and superior performance; Software monitors hardware (lasers and moving components) for system integrity, and provides a report on system performance
- One-Step Data Transfer to Acuity Database and Informatics Software: One platform for scanning, image analysis and data mining
- Automatic Calibration: Instantly tunes scanners to factory performance
- 16-bit TIFF Format Output (with other user-selectable formats): Images can easily be exported to, or imported from third-party programs
- Plug and Play Device: Simple SCSI connection to PC workstation
- Image Acquisition and Analysis with One Computer Operating System: Transition instantly from image acquisition to image analysis