

## Sole Source Specification for Axon GenePix<sup>®</sup> 4100A

---

From the industry-leader in performance, the Axon GenePix<sup>®</sup> 4100A scanner from MDS Analytical Technologies is the only scanner available today that provides:

### Features & Benefits

#### Features

- Greater Value: Affordable microarray scanning can be brought right into your lab.
  - Single laser version available for an even lower price.
- Smaller Size: Similar in size to a micro centrifuge or a PCR machine
- Fluorophore Flexibility: Emission filter wheel with two factory-installed filters for green and red fluorophores and up to six additional filters for greater customization.
- Powerful Software: Includes GenePix Pro, the highly popular and powerful acquisition and analysis software – widely recognized as the industry standard.
- Superior Optics and Digitization: Highly efficient non-confocal optical path ensures greater light-gathering capacity coupled with MDS AT's unparalleled low-noise digitization electronics for maximum signal-to-noise.
- Scanning Design: Sequential scan design permits flexible optical configuration at reduced cost
  - Scans standard microscope slides sequentially at two laser excitation wavelengths – full-slide scan time for two colors is approximately 13 min. at 10  $\mu$ m resolution, shorter for smaller scan regions.
- Quality You Expect from MDS Analytical Technologies: Many of the same powerful features as the popular GenePix 4000B microarray scanner:
  - Two factory-installed lasers: green (532 nm peak excitation) and red (635 nm peak excitation).
  - Automatic calibration instantly tunes scanners to factory performance.
  - Ready-to-use scripts for scan control, scanner monitoring, and batch processing.
  - Multiple pixel resolutions between 5 and 100  $\mu$ m.
  - Barcodes read automatically by software.
  - All scan events are logged automatically.
  - Scanner control is accessible through scripting for customized automation, analysis, and reports.
  - Control multiple scanners from a single computer.

#### Specifications

- Sample type: Standard microscope slides (1"  $\times$  3", 25  $\times$  75 mm, or 26  $\times$  76 mm)
- Maximum scan area: 22  $\times$  71.5 mm
- Excitation: 532 nm and 635 nm
- Scan time: 6.5 minutes per channel, full scan area, 10  $\mu$ m resolution
- Pixel resolution: Adjustable from 5 to 100  $\mu$ m
- Digital resolution: 16-bit
- Dynamic range: Four orders of magnitude
- Maximum S:N (PMT): 10,000:1
- Images: Single- or multi-image TIFF
- Connection: SCSI (adapter included)
- Power supply: 110 / 220V Universal power supply
- Dimensions: 14 x 11 x 14 inches (WxHxD)
- Weight: 40 lb.