

Sole Source Specification for GenePix® 4100A USB

From the industry-leader in performance, the GenePix® 4100A USB scanner from Molecular Devices is the only scanner available today that provides

Features & Benefits

- 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 Molecular Devices' 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 µm resolution, shorter for smaller scan regions.
- Quality You Expect from Molecular Devices: 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 µm.
 - 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" × 3", 25 × 75 mm, or 26 × 76 mm)
- Maximum scan area: 22 × 71.5 mm
- Excitation: 532 nm and 635 nm
- Scan time: 6.5 minutes per channel, full scan area, 10 µm resolution
- Pixel resolution: Adjustable from 5 to 100 µm
- Digital resolution: 16-bit
- Dynamic range: Four orders of magnitude
- Maximum S:N (PMT): 10,000:1
- Images: Single- or multi-image TIFF
- Connection: USB high-speed 2.0 (cable included)
- Power supply: 110 / 220V Universal power supply
- Dimensions: 14 x 11 x 14 inches (WxHxD)
- Weight: 40 lb.