

## Sole Source Specification for GenePix® SL50 Slide Loader

The GenePix® SL50 Slide Loader is an add-on unit for the GenePix 4300A or 4400A from Molecular Devices and the only system available today that provides:

### Features & Benefits

- Walk-Away Operation: Automatically loads up to 50 slides; Batch scan and batch analysis options;
- Light protected slide enclosure: Less than 0.25% of ambient light
- Reliable robotics: Encoder controlled close loop feedback
- High-precision gripper: Careful transfer and safe positioning in six-point slide holder
- Controlled nudging: Proper handling of a variety of different microarray slides
- Bench top size (h x w x d, mm): 620 x 807 x 870
- Slides are always accessible from the slide enclosure
- Operate in manual or in batch mode
- 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
- Image acquisition and analysis with one computer operating system: Transition instantly from image acquisition to image analysis