

MAGPIX[®]



System Specifications

Accuracy and Precision

- Sample uptake volume $\pm 5\%$
- Classification of microspheres $>80\%$
- Misclassification of microspheres $\leq 2\%$
- Distinguishes 1 to 50 unique xMAP[®] microspheres in a single sample
- Temperature control $\pm 1^\circ\text{C}$ of target
- Sample carryover $\leq 4\%$
- Soluble background fluorescence emissions at $590\text{ nm} \pm 24\text{ nm}$ automatically subtracted from fluorescence intensity values

Sensitivity

- Detect ≤ 700 fluorochromes phycoerythrin (PE) per xMAP microsphere

Optics

- Reporter channel, LED excitation: $511\text{nm} \pm 27\text{ nm}$
- Minimum reporter channel dynamic range: 3 decades of detection
 - Typical Dynamic Range: ≥ 3.5 decades of detection
- Reporter / Classification detector: CCD
- Reporter / Classification A/D resolution 16 bits
- Focus Lens: 5x magnification

Fluidics

- 96 well plate capability with additional off plate reagent area
- Piercing probe capability
- Auto adjusting probe capability
- Sample uptake volume $20\mu\text{L} - 200\mu\text{L}$.
- Internal / removable drive bottle and waste bottle
- Disposable Drive bottle volume: 650mL (8 plates)
- Waste bottle volume: 850mL

Electronics:

- USB 2.0 communications
- Input voltage range: $100 - 120$ or $200 - 240\text{ VAC}$ and $50 - 60\text{Hz}$

Setup

- Installation ≤ 2 hours
- System calibration < 10 minutes

Capacity

- Analyze multiple 96-well plates per batch
- Analyze multiple assay templates per plate
- Automatic sampling from a 96-well plate. The following microtiter plates are compatible with the MAGPIX plate holder: flat bottom, conical, round, filter bottom, [overall height no more than $0.75''$ (19mm)], any color
- Drive and waste bottles hold enough volume to run up to eight 96-well plates between refills
- Detect and distinguish surface reporter fluorescence emissions at $590\text{ nm} \pm 24\text{ nm}$ on the surface of 1-50 unique xMAP microsphere sets in a single sample

General

- Physical dimensions: 16.5cm (6.5 inches) W x 60cm (23.5 inches) D x 43cm (17 inches) H. Additional space required for the monitor/PC stand, keyboard, mouse and barcode scanner does not exceed 64.8cm (25.5 inches) W x 61cm (24 inches)
- Weight: up to 18kg (40lbs)
- Operating Environment:
 - Temperature: $15^\circ\text{C} - 35^\circ\text{C}$ ($59^\circ\text{F} - 95^\circ\text{F}$)
 - Humidity: $20\% - 80\%\text{RH}$, non-condensing

General (continued)

- Altitude: up to 2400m (7874 ft) above mean sea level
- Temperature control: Maintains samples at a constant temperature from 35 °C - 60 °C (95°F - 131 °F)
 - Warm up time: 15 minutes
- Plate throughput time: ≤ 60 minutes

Monitor Specifications

- Screen resolution and number of colors: SXGA 1280 x 1024 with 32-bit color
- Screen size: 43cm (17")

PC Specifications

- Processor: 2.8 GHz Intel® Core 2 or higher
- Main memory: 4 G RAM or higher
- Hard disk drive: 80 GB or higher (160 GB recommended)

- Communication/Ports: four USB version 2.0 compatible high speed ports
- DVD-RW drive
- Operating system: English USA Microsoft® Windows® XP Professional, SP3 or English USA Windows® 7, 32-bit (not 64-bit)
- All Luminex instrumentation is CE and Safety Agency marked (MET and/or UL and/or TUV and/or NEMKO) to electrical/safety device standards. For details on approvals and standards compliance please contact Luminex.

Order Information

- Contact your Luminex Sales Representative

For technical support, call: 877-785-2323 (U.S. and Canada) or +1 512-381-4397 (International)

Fax: 512-219-5114

E-mail: support@luminexcorp.com

Luminex Corporation
12212 Technology Blvd.
Austin, Texas 78727
www.luminexcorp.com