

# MAGPIX<sup>®</sup>

*Multiply the Power of Your Potential.*

MAGPIX is a compact fluorescent-based detection system, capable of simultaneously measuring up to 50 targets of interest in a single microtiter well. Based on Luminex's proven xMAP<sup>®</sup> Technology, the MAGPIX system uses color-coded magnetic microspheres as the substrate on which assays are performed. Ideal for low to medium-throughput life science research laboratories, MAGPIX can perform quantitative analysis of proteins and nucleic acids in a variety of sample matrices.

## Experience the benefits of MAGPIX:

**Sensitivity:** Approximately  $10^6$  copies of DNA or single-digit picogram levels of protein

**Dynamic Range (Typical):**  $\geq 3.5$  logs

**Read Time:** 96-well plate in  $\leq 60$  min

**Daily Start-Up/Shut-Down:**  $\leq 15$  min



## Features and Benefits

### Multiplex

- Capability to simultaneously measure up to 50 analytes in a single sample results in lower labor and reagent costs, lower sample requirements and improved laboratory efficiencies

### Compact Size

- MAGPIX is a compact instrument requiring less bench space

### Assays

- MAGPIX is supported with a broad menu of commercially available pre-optimized assay kits, as well as reagents for user-developed assays
- Compatible with an extensive range of sample types and requires less sample input than many other current methods

# Specifications\*

## PERFORMANCE

Multiplexing Capability	Up to 50 individual analytes
Reagent Compatibility	MagPlex® magnetic microspheres
Dynamic Range (Typical)	≥ 3.5 logs
Microplate Type	96-well plate
Reading Speed	96 wells ≤ 60 mins
Sample Temperature Control	35° – 60° C (95° – 131° F)
Sample Volume Uptake	20 – 200 µL
Probe Piercing	Yes
Auto Adjust-Probe Height	Yes
Daily Start-Up	≤ 15 minutes

## OPTICS

Detection	CCD Imager
Reporter Channel Excitation	511 nm
Classification Channel Excitation	621 nm

## PHYSICAL CHARACTERISTICS

Power	110 – 120 VAC or 200 – 240 VAC and 50 – 60 Hz
Dimensions	16.5 cm W x 60 cm D x 43 cm H (6.5" W x 23.5" D x 17" H)
Weight	17.5 kg (38.5 lbs)

## REGULATORY

Regulatory	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.
------------	---

\* For more detail see MAGPIX® system specifications (PN 89-60000-00-074) on the download tab at [www.luminexcorp.com/MAGPIX](http://www.luminexcorp.com/MAGPIX)

**UNITED STATES**  
Luminex Corporation  
Austin, Texas  
Tel: +1.512.219.8020  
Fax: +1.512.219.5195

**CANADA**  
Luminex Molecular Diagnostics  
Toronto, Ontario  
Tel: +1.416.593.4323  
Fax: +1.416.593.1066

**EUROPE**  
Luminex B.V.  
Oosterhout, The Netherlands  
Tel: +31.162.408333  
Fax: +31.162.408337

[www.luminexcorp.com](http://www.luminexcorp.com)

**CHINA**  
Luminex Shanghai Trading Co.  
Shanghai, China  
Tel: +86.21.616.50809  
Fax: +86.21.616.50811

[info@luminexcorp.com](mailto:info@luminexcorp.com)

**JAPAN**  
Luminex Japan Corporation, Ltd.  
Tokyo, Japan  
Tel: +81.3.5545.7440  
Fax: +81.3.5545.0451

**AUSTRALIA**  
BSD Robotics, A Luminex Company  
Brisbane, Queensland  
Tel: +61.7.3273.0273  
Fax: +61.7.3273.0274