

Luminex

Why Limit Yourself?

Experience more with the affordable MAGPIX[®] multiplexing system.



MAGPIX[®]

Multiply the Power of Your Potential.

www.luminexcorp.com/magpix

Less Sample Input. More Results.

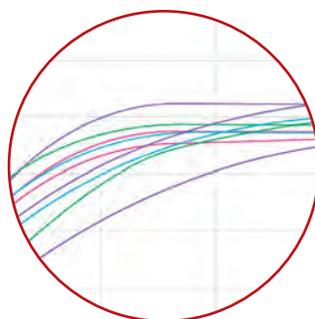
The **MAGPIX** system is a versatile multiplexing platform capable of performing qualitative and quantitative analysis of proteins and nucleic acids in a variety of sample matrices. This affordable system requires less sample input than many other current technologies and can perform up to 50 tests in a single reaction volume, greatly reducing sample input, reagents and labor while improving productivity.

The MAGPIX innovative design is based on CCD imaging technology, which allows for a compact, more robust system. Streamlined start-up and shut-down protocols and minimal maintenance requirements make the system easy to operate and maintain.

MAGPIX is supported with a broad menu of commercially available assay kits, as well as reagents to build your own assays. With MAGPIX, multiplexing is more accessible than ever.



96-well plate



Up to 50 analytes/well



Up to 4,800 results in one hour

A Simple and Intuitive Solution.

The **MAGPIX** system is designed for ease of use allowing new and experienced users alike to set up and run an assay in a minimal amount of time. ELISA-like workflow, intuitive software interface, online training and 24/7 technical support will minimize your ramp-up time to multiplexing.

MAGPIX System 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.
------------	---

Specifications subject to change.

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

A Versatile Platform to Meet Your Needs.

Functions that previously required multiple instruments can now be performed on a single platform, giving you more flexibility and a greater array of options. Whether testing for a single analyte or multiple analytes, proteins or nucleic acids, MAGPIX delivers accurate analytical performance and efficient, easy-to-follow protocols.

MAGPIX is based on xMAP® Technology, a proven platform that has supported cutting-edge research for more than a decade. Innovations in optics have resulted in a compact, robust system that is user installable in less than two hours. These advancements and low maintenance requirements make MAGPIX the ideal platform for small laboratories with limited budgets.

- Ability to perform single and multiplex analysis
- Analysis of both proteins and nucleic acids
- Broad menu of commercially available assay kits and reagents for user-developed assays

Applications

Protein Research

Cytokines and Chemokines
Cell Signaling
Inflammatory Mediators
Cancer Markers
Cardiac Markers
Protein Expression

Infectious Disease
Transcription Factors
Apoptosis
Hormone Receptors
Neurobiology
Others

Molecular Research

Gene Expression
SNP Genotyping
MicroRNA
Others

xPONENT® SOFTWARE SOLUTIONS

xPONENT offers industry-standard curve-fitting modules for quantitative analysis, allelic ratios for genotyping and qualitative algorithms for serology assays, as well as data visualization tools. The software can also be configured to automatically launch third-party data analysis programs for advanced statistical analysis.

Software Options:

- **Security Levels:** The secure xPONENT package allows system administrators to set up user access levels
- **21 CFR Part 11:** Offers multilevel user management, full audit trail, electronic records and electronic signatures
- **IQ/OQ Protocol:** Document system qualification to help laboratories meet regulatory compliance burdens



MAGPIX Features and Benefits.

Let the Luminex MAGPIX system accelerate your research by putting the power of multiplexing technology well within your reach. This groundbreaking new product gives you more data, more efficiency and more options – all at a price you won't believe. Multiply your performance, multiply your efficiency and multiply your potential with the new Luminex MAGPIX system.

Affordability

Benefit from an affordable multiplexing solution that costs up to 4x less than comparable ELISA assays.*

Versatility

Transition to multiplexing at your own pace and exponentially grow your data.

Flexibility

Achieve a wider array of goals with the ability to analyze both proteins and nucleic acids.

Compact Size

Reduce lab clutter with a smaller, more compact instrument that requires less linear bench space.

Efficiency

Measure up to 50 targets in a single reaction volume and dramatically reduce sample input.

Assay Support

Leverage a broad range of commercially available, pre-optimized assay kits, as well as reagents for user-developed assays.



* Price per result based on comparison of xMAP 42-plex kit to ELISA kit containing the same analyte.

Luminex Instrumentation.

FEATURE	INSTRUMENT		
	MAGPIX	Luminex 100/200	FLEXMAP 3D
Software	xPONENT®	xPONENT	xPONENT
Optics	LED/CCD Camera	Lasers/APDs/PMTs	Lasers/APDs/PMTs
Hardware	Fluorescent Imager	Flow Cytometry based	Flow Cytometry based
Bead Compatibility	MagPlex®	MagPlex MicroPlex® SeroMAP™ LumAvidin® xTAG®	MagPlex MicroPlex SeroMAP LumAvidin xTAG
Multiplex Capacity	50	100 (80 for MagPlex)	500
Read Time	~ 60 mins/96-well plate	~ 40 mins/96-well plate	~ 20 mins/96-well plate ~ 75 mins/384-well plate
Applications	Protein/Nucleic Acid	Protein/Nucleic Acid	Protein/Nucleic Acid
510(k) Assays	---	Over 50	---
Dynamic Range	≥ 3.5 logs (typical)*	≥ 3.5 logs	≥ 4.5 logs
Microtiter Plate	96 well	96 well	96 well and 384 well
Footprint Including PC (Linear Bench Space)	64.8 cm (24")	80.0 cm (32")	64.8 cm (24")
Weight (Analyzer)	17.5 kg (38.5 lbs)	49 kg (113 lbs)	77.1 kg (170 lbs)

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

Luminex Instrumentation: Solutions You Can Rely On

Since its incorporation in 1995, Luminex has become a global leader in multiplexing solutions. Luminex offers a full family of instruments designed to meet all your multiplexing needs. Whether it's the efficient MAGPIX system, the dynamic Luminex® 100/200™ system or the powerful FLEXMAP 3D® system, we're sure to have a solution that will help you optimize your research capabilities. Learn more about our suite of enabling multiplexing solutions. Only Luminex offers the piece of mind that comes from:

- Peer reviewed publications
- Patented xMAP technology
- Strong network of market leading partners
- Over 50 510k cleared assays
- Thousands of instruments placed worldwide

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

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

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

www.luminexcorp.com

info@luminexcorp.com