



# Luminex xMAP<sup>®</sup> Instrumentation

## Flexible Multiplexing Technology

Luminex offers three xMAP<sup>®</sup>-based instruments to run multiplexed assays for both genomic and proteomic assays, whether kit-based or lab-developed. Our platforms provide:

- Open architecture to allow development of your own unique tests
- Broad menu of commercially available assays
- Ability to multiplex up to 500 targets in one well of a microtiter plate

Feature*	Instrument		
	MAGPIX <sup>®</sup>	Luminex <sup>®</sup> 100/200 <sup>™</sup>	FLEXMAP 3D <sup>®</sup>
Software	xPONENT <sup>®</sup>	xPONENT <sup>®</sup>	xPONENT <sup>®</sup>
Optics	LED/CCD Camera	Lasers/APDs/PMTs	Lasers/ APDs/ PMTs
Hardware	Fluorescent Imager	Flow Cytometry based	Flow Cytometry based
Beads Compatibility	MagPlex <sup>®</sup> MagPlex <sup>®</sup> -TAG	MagPlex <sup>®</sup> MicroPlex <sup>®</sup> SeroMAP <sup>™</sup> LumAvidin <sup>®</sup> MagPlex <sup>®</sup> -TAG	MagPlex <sup>®</sup> MicroPlex <sup>®</sup> SeroMAP <sup>™</sup> LumAvidin <sup>®</sup> MagPlex <sup>®</sup> -TAG
	Refer to MagPlex <sup>®</sup> Ordering Guidelines for region selection.		
Multiplex Capacity	50	100 (80 for MagPlex <sup>®</sup> & MagPlex <sup>®</sup> -TAG)	500
Read Time	~60 mins/96 well plate	~40 mins/96well plate	~20 mins/96well plate
Applications	Protein/Nucleic acid	Protein/Nucleic acid	Protein/Nucleic acid
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 lb)	49 kg (113 lb)	77.1 kg (170 lb)

\*For more detail, please refer to the Market Specification documents available for each instrument on the Luminex website.



To learn more, please visit: [www.luminexcorp.com](http://www.luminexcorp.com)

Specifications and features subject to change without notice. The FLEXMAP 3D and Luminex 100/200 are class I(I) laser products.

©2016 Luminex Corporation. All rights reserved. Luminex, xMAP, FLEXMAP 3D, MAGPIX, MicroPlex, MagPlex, and LumAvidin are trademarks of Luminex Corporation, registered in the U.S. and other countries. Luminex 100/200 is a trademark of Luminex Corporation.

#### HEADQUARTERS

#### UNITED STATES

+1.512.219.8020

info@luminexcorp.com

www.luminexcorp.com

#### EUROPE

+31.73.800.1900

europe@luminexcorp.com

www.luminexcorp.com

#### CANADA

+1.416.593.4323

info@luminexcorp.com

www.luminexcorp.com

#### CHINA

+86.21.802.31150

info@luminexcorp.com

www.luminexcorp.com

#### JAPAN

+81.3.5545.7440

info@luminexcorp.com

www.luminexcorp.com

FL1503.0316