

MAGPIX[®] System Specifications

For Use With xPONENT[®] 4.3 Software

This document is intended to provide key system specifications for MAGPIX[®] instrument hardware, including the minimum specifications of the PC that will run xPONENT[®] Software.

General

Physical Dimensions	16.5 cm (6.5 in.) x 60cm (23.5 in.) x 43 cm (17 in.) (W x D x H) Additional space required for the monitor/PC stand, keyboard, mouse, and barcode scanner does not exceed 64.8 cm (25.5 in.) x 61 cm (24 in.)
Weight	18 kg (40 lbs.)
Operating Temperature	15°—35°C (59°—95°F)
Operating Humidity	20—80%, non-condensing
Altitude	Operation up to 2,400 m (7,874 ft) above mean sea level
System Warmup Time	15 min
System Initialization	<15 min. (Includes fluidics flush and weekly calibration)
System Verification	5 min
Temperature Control	Samples are maintained at a constant temperature when using the heater block (35—60°C +/- 2°C of set point; 95—131°F)
Plate Run Time	96-well plate ≤60 min

Electronics

USB 2.0-compatible communications link for fast data transfer

Input Voltage Range	100—120 V- ±10%, 1.4 Amp, and 200—240 V- ±10%, 0.8 Amp, 47—63 Hz
Installation Category	II - As defined in IEC 61010-1:2010
Pollution Degree	2 - As defined in IEC 61010-1:2010

Optics

Reporter Channel	LED excitation: 511 nm ± 27 nm
Reporter Channel Dynamic Range	3.5 decades of detection
Reporter/Classification Detector	CMOS
Reporter Channel Detection	A/D resolution 16 bits
Focus Lens	5x magnification

Fluidics

Sample Uptake Volume	20—200 µL
Well-to-well Carryover	<4%
Piercing Probe Capability	Yes
Auto Adjusting Probe Capability	Yes
Internal/Removable Drive Bottle and Waste Bottle	Yes
96-Well Plate Capability with Additional Off-Plate Reagent Area	Yes

Microspheres

- Distinguish 1 to 50 unique xMAP® MagPlex® microspheres in a single sample
- Detects a minimum of 700 fluorochromes of phycoerythrin (PE) per xMAP® microspheres
- Soluble background fluorescence emissions at 590 nm ± 24 nm automatically subtracted from fluorescence intensity values

Classification of xMAP® Microspheres	≥80%
Total System Misclassification of xMAP® Microspheres	≤2%
Temperature Control	±1°C of target

PC and Monitor Specifications

Processor	3.0 GHz Intel Core i5 (or higher)
Main Memory	8 GB RAM
Hard Drive	1 TB Hard drive space (or higher)
Ports	1x USB 3.1 Type-C 5x USB 3.1 Type-A 4x USB 2.0 Type-A 2x Display ports 1x VGA port 1x DVD R/W type drive PCI USB Adapter Card
Operating System	Microsoft® Windows® 10 Professional 64 bit, version 1709 or greater
Screen Resolution	SXGA 1,280 x 1,024, 32 bit color
Screen Size	49.3 cm (19 in.)
Display Settings	96 DPI font Default Windows® 10 theme

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.

Luminex
A DiaSorin Company

For more information, please visit: www.luminexcorp.com/MAGPIX

Products are region specific and may not be approved in some countries/regions. Please contact Luminex at support@luminexcorp.com to obtain the appropriate product information for your country of residence.

©2022 Luminex Corporation. A DiaSorin Company. All rights reserved. Luminex, MAGPIX, MagPlex, xMAP, and xPONENT are trademarks of Luminex Corporation, registered in the US. and other countries. Microsoft and Windows are registered trademarks of Microsoft Corporation.

HEADQUARTERS

UNITED STATES

+1 512 219 8020

info@luminexcorp.com

EUROPE

+31 73 800 1900

europe@luminexcorp.com

CANADA

+1 416 593 4323

info@luminexcorp.com

CHINA

+86 21 8036 9888

infocn@luminexcorp.com

JAPAN

+81 3 5545 7440

infojp@luminexcorp.com

FL416233