

Multiplex Perfected

Bio-Plex Manager MP software ensures high quality results and optimal instrument performance. Use for maintenance and data acquisition on the Bio-Plex® MAGPIX™ multiplex reader.

- Optimize performance with automated instrument management
- Integrate with powerful Bio-Plex analysis software
- Quickly adopt and master multiplex technology



Now with Automated Instrument Management

Using Bio-Plex Manager MP software is like having a technical expert monitoring your MAGPIX instrument. It suggests maintenance procedures based on current instrument state and performance. Bio-Plex Manager MP software is a dynamic, responsive member of your team. It's multiplex perfected.

Automated Instrument Management

- Automatically loads recommended start-of-day maintenance routines based on current instrument status
- Monitors performance during data acquisition, alerts the user to performance issues, and loads recommended maintenance to resolve issues
- Recommends simple step-wise troubleshooting

Software Benefits

- Detection and resolution of performance issues ensures data quality
- Confidence in proper instrument maintenance leads to confidence in results
- Highest quality results maximize the value of your experiments
- Simple integration with Bio-Plex analysis software to quickly assemble and understand your data
- Simple interface and built-in guidance enable a user of any experience level to easily run the Bio-Plex MAGPIX multiplex reader

Convert any MAGPIX Multiplex Reader into a Self-Monitoring Instrument

Even if you purchased your MAGPIX multiplex reader from another vendor you can benefit from the intelligence and simplicity of Bio-Plex Manager MP software. Contact your local Bio-Rad representative about upgrading your Bio-Plex MAGPIX multiplex reader today.

Ordering Information

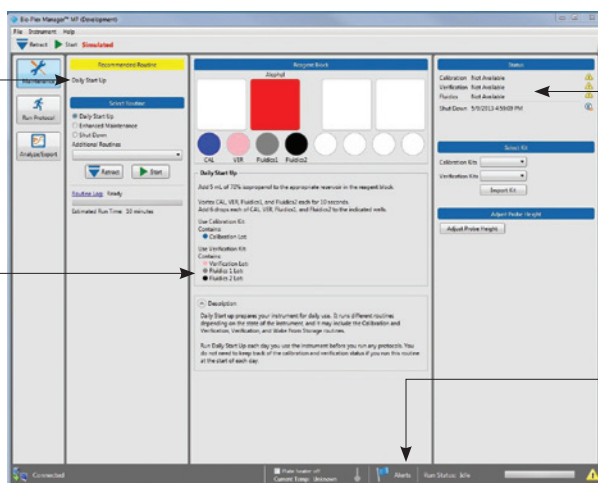
Catalog #	Description
171-051555	Bio-Plex Manager MP Software Upgrade , includes Bio-Plex Manager MP software, software activation code, Getting Started Guide, probe height adjustment plate, Upgrade Quick Guide for Bio-Plex Manager 6.1
171-015001	Bio-Plex MAGPIX Reader , includes Bio-Plex MAGPIX instrument and accessories, PC and accessories, MS Excel, Bio-Plex Manager MP software, software activation code, Getting Started Guide, and probe height adjustment plate
171-STND01	Bio-Plex Manager Desktop , 1 seat, for review and optimization of data from a single instrument run
171-001513	Bio-Plex Data Pro™ , 5 seats, for assembly of multiple instrument runs and interpretation of experimental data
171-001523	Bio-Plex Data Pro Plus , 5 seats, for assembly of multiple instrument runs and interpretation of experimental data, includes saved analysis sessions and standard curve overlays

For more information, visit www.bio-rad.com/magpix.

MAGPIX is a trademark of Luminex Corporation.

Recommended routine is presented for maintenance and performance optimization

Detailed instructions for performing routines and step-by-step troubleshooting steps



Status center tracks calibration state and whether end-of-day maintenance has been performed

Alert center tracks important messages. Complete message can be accessed at any time



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026 674 55 05
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000