



Bio-Plex® MULTIPLEX SYSTEM

Bio-Plex Pro™ Wash Station

Quick Guide

Using the Wash Station

1. Prime Channel 1.
2. Perform your Wash Programs.
3. Rinse when finished using the instrument.
 - RINSE DAY if you will use the instrument later that day
 - RINSE NIGHT at the end of the day

Which Program Do I Use?

Assay/Step	Program
Magnetic Bead–Based Assay*	
First wash	MAG x2
All other washes	MAG x3

* Magnetic bead–based assays are Bio-Plex Pro Assays or assays that use MagPlex Beads.

Manual Steps

Wash Station Protocols

Dispense 50 µl bead suspension Wash 2 cycles	MAG x2
Dispense 50 µl standards and solutions Incubate for 30–120 min* (refer to assay manual) Wash 3 cycles	MAG x3
Dispense 25 µl detection antibody Incubate for 30 min* Wash 3 cycles	MAG x3
Dispense 50 µl SA-PE Incubate for 10 min* Wash 3 cycles	MAG x3
Dispense 125 µl assay buffer Place on shaker and shake for 1 min Read on Bio-Plex System Array Reader	—

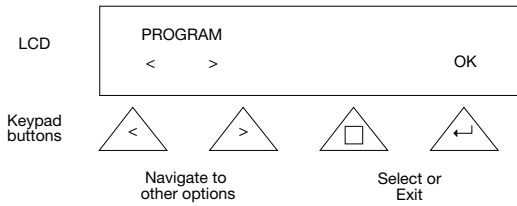
SA-PE, streptavidin-phycoerythrin.

* Incubating plates should be:

1. At room temperature.
2. Covered with aluminum foil or otherwise protected from light.
3. Shaken according to the recommended shaker settings in the product manual.

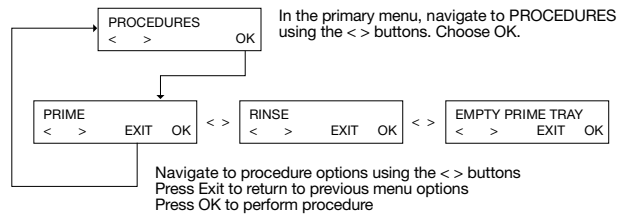


Using the Keypad



If you don't see the above screen, press Exit until you see PROGRAM.

Example Navigation



Follow the charts below to perform each of the indicated operations.

Using the Wash Station for Daily Operation

1. Prime Procedure

Caution: Do not run any programs until you have primed Channel 1.

Prime Channel 1 if:

1. Last action was a rinse with Channel 2.
2. The system has been idle for more than an hour.
3. You are changing buffer in bottle 1.

2. Wash Programs

3. Rinse Procedures

Use RINSE DAY if you will use the instrument later that day.

Use RINSE NIGHT at the end of the day.

or RINSE NIGHT

Key: <> Select option shown ↵ Use default option

Bottle Contents

- Liquid 1 — Wash buffer
- Liquid 2 — Distilled water or equivalent

MagPlex is a trademark of Luminex Corporation.

The Bio-Plex Suspension Array System includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 177 **Belgium** 32 (0)3 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 45 44 52 10 00 **Finland** 358 09 804 22 00
France 33 01 47 95 69 65 **Germany** 49 89 31 884 0 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 972 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 (0)318 540 666
New Zealand 64 9 415 2280 **Norway** 47 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 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 46 08 555 12700 **Switzerland** 41 026 674 55 05
Taiwan 886 2 2578 7189 **Thailand** 66 662 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 020 8328 2000

