

colorimetric systems

MICROPLATE WASHERS

THE 1575 IMMUNOWASH MICROPLATE WASHER OFFERS CUSTOM ASSAYS INCLUDING:

- STANDARD ELISA
- ADHERENT CELL WASHING
- BLOOD TYPING ASSAYS
- CELL GROWTH REQUIRING BOTTOM WASHING
- OVERFLOW WASH FOR WEAKLY BOUND ANTIBODIES
- CUSTOM ASSAY WASH PROCEDURES

The Model 1575 Immunowash Microplate Washer

Performance and Flexibility

The Model 1575 Immunowash microplate washer offers the total solution for laboratories requiring versatility in a microplate strip washer. It addresses the wash requirements for standard ELISA, cell-based assays, blood typing, and other custom assays.

The Model 1575 Immunowash microplate washer eliminates the problem of residual buffer in flat bottom wells with custom adjustment of needle positions for crosswise aspiration. Its residual volume of less than 2 μ l represents the best performance of any commercially available microplate washer.

Flexible Wash Programs

Sophisticated on-board software permits virtually unlimited flexibility in creation of wash programs and the capacity to save up to 75 of these programs for easy recall and use. The Model 1575 Immunowash microplate washer employs any combination of up to six single wash types or four two cycle wash types. This allows the user to customize any type of wash procedure, set soak times, and wash either individual strips or the entire microplate.

You retain control over liquid dispense speed, horizontal and vertical positions of aspirate needles, crosswise aspiration, bottom washing (Coombs washing), plate shaking, and vertical speed of the wash manifold. User-friendly on-board software greatly simplifies programming of these procedures and allows you to save up to 75 methods. The option of either 8- or 12-channel manifolds provides additional control over throughput requirements while a high power vacuum provides superior dispense and aspirate efficiency.



Simple Maintenance

The Model 1575 Immunowash microplate washer eliminates complex and time-consuming cleaning operations necessary for fixed wash manifolds with an easily accessible manifold with removable access cover. Unplugging two tubes allows the user to completely remove the manifold for additional cleaning and/or sterilization procedures. A waste bottle sensor alerts the user when the bottle becomes full and eliminates liquid backup in the system.

Laboratory Safety

A standard aerosol protection cover completely encloses the plate and manifold to both protect the user from accidental liquid splashes and eliminate interference from airborne contaminants.

BIO-RAD

Specifications

ELECTRICAL

Voltage	220/110 VAC, 50/60 Hz
Consumption	100 VA max
Fuses	2.5 AT

PHYSICAL

Width	32.5 cm
Length	40.05 cm (plate carrier in) 52.55 cm (plate carrier out)
Height	19.3 cm
Weight	9.6 Kg
Operating conditions	15–30 °C/15-85% R.H.

HARDWARE

Manifolds	8-channel or 12-channel
Vacuum	Integrated vacuum pump 9 l/min
Waste bottle volume	2,000 ml

SOFTWARE

Kits	Up to 75 Wash Kits programmable. Each kit is saved under its own 15 character name and is a combination of methods with or without soak times in strip or plate mode.
Method	6 single cycle washing methods: Wash, Aspirate, Dispense, Bottom Wash, Bottom Aspirate, Shake. 4 two cycle washing methods: Wash + Aspirate, Wash + Bottom Aspirate, Bottom Wash + Aspirate, Bottom Wash + Bottom Aspirate.
Soak time	0 to 9.9 seconds in Strip Mode. 0 to 59 minutes in Plate Mode.
Repetition	Every method can be repeated up to 9 times
Wash mode	Strip Mode or Plate Mode
Omit strip	Any strip can be omitted during kit programming
Plates	Up to 10 plate parameter sets programmable; accepts flat or round bottom microplates. Vertical and horizontal speeds, vertical and horizontal positions of aspirate needle inside well are programmable.
Other features	Priming sequence of hydraulic system during change of wash solution. Separate disinfection program Password program protect
Performance aspiration	Less than 2 µl maximum residual volume in a flat bottom well
Volume accuracy	<2.5% CV across a 96-well plate

Ordering Information

Catalog # Description

MODEL 1575 IMMUNOWASH MICROPLATE WASHER

170-7009	Model 1575 Immunowash Microplate Washer, 100/120/220/240 VAC, includes bottle and tubing set, 8-channel manifold, aerosol protection cover, power cord, and instruction manual
170-7010	Model 1575 Immunowash 12-channel manifold
170-7011	Model 1575 Immunowash Standard Maintenance Kit

BIO-RAD

**Bio-Rad
Laboratories**

Life Science
Group

Website www.bio-rad.com U.S. (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Canada (905) 712-2771
China 86-10-62051850/51 Denmark 45 39 17 99 47 Finland 358 (0)9 804 2200 France 01 43 90 46 90 Germany 089 318 84-0
Hong Kong 852-2789-3300 India (91-11) 461-0103 Israel 03 951 4127 Italy 39-02-216091 Japan 03-5811-6270 Korea 82-2-3473-4460
Latin America 305-894-5950 Mexico 514-2210 The Netherlands 0318-540666 New Zealand 64-9-4152280 Norway 22-74-18-70 Russia 7-095-4585822
Singapore 65-2729877 Spain 34-91-661-7085 Sweden 46 (0)8 627 50 00 Switzerland 01-809 55 55 United Kingdom 0800-181134