

# ImmunoWash™ 1575 Microplate Washer

## Essential Microplate Washer that Streamlines any Assay Workflow

A flexible and user-friendly instrumentation for delivering, mixing, and removing liquids from microtiter plates, the ImmunoWash 1575 microplate washer improves efficiency for many microplate applications.

### Feature

### Benefit

Up to 110 protocols can be saved and easily modified. Each protocol is a combination of 4 adjustable procedures (washing, aspiration, dispensing, and agitation) and multiple cycles	<ul style="list-style-type: none"> <li>Washing procedures are accurate and efficient at highest level of optimization</li> <li>Delivers exact reproducibility of assays and repeatability of workflows</li> </ul>
Programmable horizontal and vertical needle positions (accurate to within 0.1 mm) to support bottom washing, crosswise aspiration, and overflow washing	<ul style="list-style-type: none"> <li>Accommodates assays with special requirements (sediments, coagulants, low volumes, low liquid agitation requirements)</li> <li>Saves reagents and time</li> </ul>
Overflow washing with circular flow of liquid in each well	<ul style="list-style-type: none"> <li>Provides high wash efficiency without compromising throughput</li> <li>Replaces tedious sequences of repeated aspiration and dispensing steps</li> </ul>
Crosswise aspiration	<ul style="list-style-type: none"> <li>Minimizes residual volume to below 6 µl per well</li> </ul>
Dispenser speed control	<ul style="list-style-type: none"> <li>Delivers higher volume accuracy</li> </ul>
Adjustable plate shaking	<ul style="list-style-type: none"> <li>Minimizes bubbles and liquid adherence to well sides</li> </ul>
Easily accessible for cleaning and maintenance Removable and autoclavable plate carrier Aerosol protection cover Preprogrammed disinfection protocol	<ul style="list-style-type: none"> <li>Satisfies concerns about use of infectious agents and samples that are sensitive to the environment</li> <li>Easy decontamination and prevention of contamination</li> </ul>
Integrated and automated vacuum and dispensing pumps	<ul style="list-style-type: none"> <li>Ensures accurate and quiet washing, eliminating the need for bulky, noisy external pumps</li> <li>Improves working conditions for lab personnel</li> <li>Lowers maintenance costs</li> </ul>
Easy to install 8- and 12-channel manifolds	<ul style="list-style-type: none"> <li>Allows for different orientations of the washing procedure in plate and individual strips</li> </ul>
Multilingual (English, French, Russian) interface via LCD display	<ul style="list-style-type: none"> <li>Accommodates localized regulatory requirements and communication standards</li> </ul>

**BIO-RAD**

# ImmunoWash™ 1575 Microplate Washer

## Specifications

Residual well volume	<6 µl
Soak time	0–9.9 sec in strip mode 0–59 min in plate mode
Available manifolds	8- and 12-way
Protocol storage capacity	110 protocols
Compatible plate types	Any flat- or round-bottomed plate
Waste bottle volume	2,000 ml
Vacuum	Integrated vacuum pump, 9 L/min
Dimensions (W x D x H)	34.5 x 42.3 x 20.0 cm
Weight	10.3 kg
Operating conditions	15–40°C, 15–85% relative humidity
Voltage	200/100 VAC; 50/60 Hz
Consumption	100 VA maximum
User interface	2 x 20 character LCD screen; 5 diaphragm keys; flat keyboard

## Ordering Information

Catalog #	Description
170-7009	<b>ImmunoWash 1575 Microplate Washer</b> , 100–240 V, includes bottle and tubing set, 8-port manifold, aerosol protection cover, power cord, instructions
170-7021	<b>ImmunoWash 1575/MICROTECH™ 12-Channel Manifold</b>
170-7026	<b>ImmunoWash 1575 Standard Maintenance Kit</b>



**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://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 31 3689 6600  
**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** 30 210 777 4396 **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 **Malaysia** 60 3 2117 5260 **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 3170 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700  
**Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 6518311 **United Kingdom** 020 8328 2000