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.



Benefit
 Washing procedures are accurate and efficient at highest level of optimization Delivers exact reproducibility of assays and repeatability of workflows
 Accommodates assays with special requirements (sediments, coagulants, low volumes, low liquid agitation requirements) Saves reagents and time
 Provides high wash efficiency without compromising throughput Replaces tedious sequences of repeated aspiration and dispensing steps
■ Minimizes residual volume to below 6 µl per well
■ Delivers higher volume accuracy
Minimizes bubbles and liquid adherence to well sides
 Satisfies concerns about use of infectious agents and samples that are sensitive to the environment Easy decontamination and prevention of contamination
 Ensures accurate and quiet washing, eliminating the need for bulky, noisy external pumps Improves working conditions for lab personnel Lowers maintenance costs
 Allows for different orientations of the washing procedure in plate and individual strips
 Accommodates localized regulatory requirements and communication standards



ImmunoWash™ **1575 Microplate** Washer

Specifications

-poomounomo	
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 #

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

Description



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 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 23473 4460 Malaysia 60 3 2177 5260 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland 48 22 331 999 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

Bulletin 5786 Rev B US/EG 11-0515 0611 Sig 0211