

Specifications Sheet



Model 1575 Microplate Washer

Ordering Information

Catalog No.	Product Description
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, instruction manual
170-7021	Model 1575 Immunowash 12-Channel Manifold
170-7026	Model 1575 Immunowash Standard Maintenance Kit

Electrical voltage: 220/110 VAC, 50/60 Hz

Consumption: 100 VA maximum

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% Relative Humidity

Hardware manifolds: 8- or 12-channel

Vacuum: Integrated vacuum pump 9 L/minute

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

Methods: 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: 0-9.9 seconds in strip mode; 0-59 minutes in plate mode

Repetition: Every method can be repeated up to 9 times

Wash mode: Strip 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 protected

Performance aspiration: Less than 2 µl maximum residual volume in a flat bottom well

Distribution accuracy: 15 µl average volume variation across one plate

BIO-RAD

Bio-Rad
Laboratories



Molecular
Bioscience Group

U.S. (800) 4BIORAD • California (510) 741-1000 • Australia 02-9914-2800 • Austria (1)-877 89 01 • Belgium 09-385 55 11 • Canada (905) 712-2771 • China (01) 2046622 • Denmark 39 17 9947 • Finland 90 804 2200 • France (1) 49 60 68 34 • Germany 089 318 84-0 • India 91-11-461-0103 • Italy 02-21609 1 • Japan 03-5811-6270 • Hong Kong 7893300 • The Netherlands 0318-540666 • New Zealand 09-443 3099 • Singapore (65) 272-9877 • Spain (91) 661 70 85 • Sweden 46 (0) 8-735 83 00 • Switzerland 01-809 55 55 • United Kingdom 0800 181134