

MultiWash+ Microplate Washer

A 96- and 384-well automated washer

BENEFITS

- Walk-up usability
- Start simply with wash and rinse bottles (included)
- Flexible system washes 96- or 384-well plates
- Automated rinse reduces clogging
- Quiet and efficient

Introduction

MultiWash™+ washer is an automated microplate washer that provides maximum flexibility with a minimum footprint. The MultiWash+ washer can be configured for both 96- and 384-well plates and comes with four wash/rinse bottles and one waste bottle for out-of-the-box use.

Wash protocols & modes

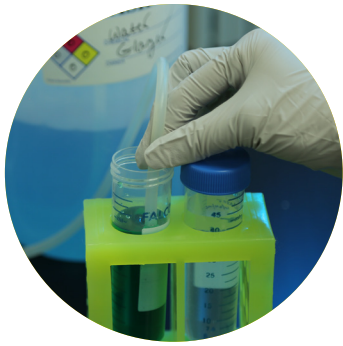
The on-board touch panel comes with the option of 20 different wash protocols, including up to 8 cycles within each protocol for walk-up useability and flexibility. Washing variations include adjustable speed and volume, adjustable aspiration speed and time, adjustable soak times and three modes of shaking are available to mix solutions.

Efficiency

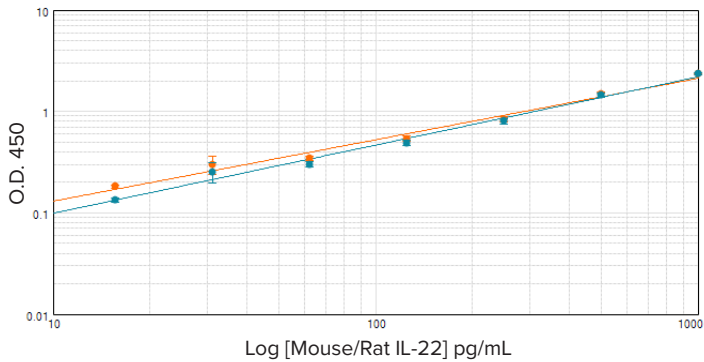
Efficient washing is achieved with cross-wise aspiration, reducing residual volume within each well. The washer has a small footprint and is vacuum and pressure free, with on-board pumps that provide a quiet wash experience.

Mouse/Rat IL-22 Quantikine ELISA using the MultiWash+ washer

A Mouse/Rat IL-22 Quantikine ELISA from R&D systems was used to evaluate the MultiWash+ washer. An IL-22 standard curve was prepared and a sandwich ELISA was performed using the MultiWash+ plate washer in strip mode to wash the wells. After reading the ELISA plate on the EMax® Plus Microplate Reader, the R2 value was 0.984. A control well was set up to verify the accuracy of the standard curve. The standard was within the range described with the kit (data not shown).



MultiWash+ washer can dispense in 50 µl increments and tubing can be inserted into a variety of containers since pressurized bottles are not required.



ELISA Assay

Technical specifications

Plate wash time (300 µl)	<ul style="list-style-type: none"> • 96-well plate, 1 wash: 1 min. 9 sec. • 96-well plate, 3 washes: 3 min. 35 sec.
Dispense volume	50 – 2000 µL in 50 µL increments
Residual volume	< 1 µL per well
Dead volume	15.62 mL
Precision	< 5% at 300 µL across the plate
User interface	2 X 16 characters backlit display keyboard with 5 function keys
Shaking	3 speeds
Microplates	<ul style="list-style-type: none"> • 96-well plates flat and round bottom (384-well plates with 16-manifold) • Fits plates with standard ANSI SBS footprint
Power requirement	90 – 250B, 60V auto-sensing, 50/60 Hz
Instrument dimensions (WxHxD)	<ul style="list-style-type: none"> • 21 x 21 x 44 cm, • 8.3" x 8.3" x 17.3"
Weight	8Kg (17.8lbs)

Ordering information

Part number

MultiWash+ washer for 96-well plates: 4 liquid lines, 8 way manifold and tubing, 4 wash/rinse bottles (2L), 1 waste bottle (2.5L)	MultiWash+
Optional 16-way manifold	5032338
Optional 12-way manifold	5032336
Washer Maintenance Kit: 2 X T 2.5A spare fuses, 2 cleaning needles, tubing sets for all bottles, cleaning tool for needles, tubing set for manifold, and spare tubing set for manifold	5032339
External Tubing Kit (4-way)	5032340
Wash bottle (2L)	5032341
Rinse bottle (2L)	5032342

Contact Us

Phone: [+1-800-635-5577](tel:+18006355577)
 Web: www.moleculardevices.com
 Email: info@moldev.com
 Check our website for a current listing of worldwide distributors.