

Bio-Rad Laboratories
Improving lives with every test.

Introducing the MedTec

AutoBlot 3000

Strip processing

Following the manual addition of test samples, the AutoBlot incubates, washes and performs subsequent reagent additions as defined by the operator during the programming phase. It permits easy setup with walk-away performance, sounding an alarm when the test is complete.

The AutoBlot is fully programmable from the front panel and will store up to ten protocols within the unit, allowing for full customization of blot assays for dispense, incubation and aspiration.

The AutoBlot dispenses and aspirates up to twenty strips in 90 seconds. This means the automated assay runs quickly, compared to performing the assay manually, with no operator intervention after sample addition. The AutoBlot 3000 dispenses up to six reagents, depending on the number of pumps.

Convenient

- Instrument, reagent and accessories package

Flexible

- Compact, light-weight, benchtop processor

Easy-to-Use

- Software and hardware programs

Reliable

- History of successful performance

*MedTec
AutoBlot 3000
Strip Processor*



BIO-RAD

AutoBlot 3000

Technical Specifications

Dimensions

22 in x 18 in x 6.5 in (556 mm x 457 mm x 165 mm)

Shipping Box Size

31 in x 23.5 in x 12 in (787 mm x 597 mm x 305 mm)

Unit Weight

30 lbs (14 kg)

Capacity

One 20-strip tray (non-heated platform)

Power

100–24 V; 50 or 60 Hz; 3.2 amp max (heated),
1.3 amp max (unheated)

Software Updates

PC download

Reagent Bottle

500 mL (wash buffer); 250 mL (DI water)

Waste Bottle

1000 µL

Dispense Volume Range

0.5 mL – 3.0 mL in 0.1 mL increments (accuracy ±10%)

Incubation Time

Up to 24 hours for all incubation periods in increments
of 1 minute

Number of Programmable Steps per Assay

15 steps

Maximum Relative Humidity

90%

Ordering Information

Catalog No.	Description	
439MT01049	Pump tubing kit.....	1
439MT01001	Tray, 20-well, AutoBlot	3-pack
439MT01008	Bottle, 1000 mL, waste with lid, AutoBlot.....	1
439MT01009	Bottle, 500 mL, wash with lid, AutoBlot.....	1
439MT01010	Bottle, 250 mL, waste with lid, AutoBlot.....	1
439MT01089	Lid, brown reagent bottle, AutoBlot.....	2 units
439MT01096	Complete tubing kit, AutoBlot 3000	1
439MT01097	Power supply, AutoBlot 3000, 24V.....	1
439MT07600	Guide to operation, manual, AutoBlot 3000	1
25224R	AutoBlot 3000, non-heated, refurbished.....	1
25224	AutoBlot 3000, non-heated, includes: 20-well tray (3-pack); 1000 mL waste bottle with lid; 500 mL wash bottle with lid; 250 mL waste bottle with lid; brown reagent bottle lid; 24 V power supply; operation manual	1 unit



**Bio-Rad
Laboratories**

For further information, please contact the Bio-Rad office
nearest you or visit our website at www.bio-rad.com.

Clinical
Diagnostics Group

Website www.bio-rad.com U.S. 1-800-2-BIO-RAD Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium 32-9-385-5511 Brazil 5521-3461-5202 Canada 1-514-334-4372 Czech Republic 420-2-41430532 China 86-21-63052255 Denmark 45-4452-1000 Finland 358-9-804-22-00 France 33-1-4795-6000 Germany 49-89-31884-0 Hong Kong 852-2789-3300 India 91-124-6398112 Israel 972-3-9514127 Italy 39-02-216091 Japan 81-3-5811-6290 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 5255-5534-2552 The Netherlands 31-318-540666 New Zealand 64-9-415-2280 Norway 47-23-38-41-30 Poland 48-22-331-99-99 Portugal 351-21-472-7700 Russia 7-095-721-14-00 Singapore 65-6415-3188 South Africa 27-11-442-85-08 Spain 34-91-590-5200 Sweden 46-8-555-127-00 Switzerland 41-61-717-95-55 Thailand 662-651-8311 United Kingdom 44-208-328-2000