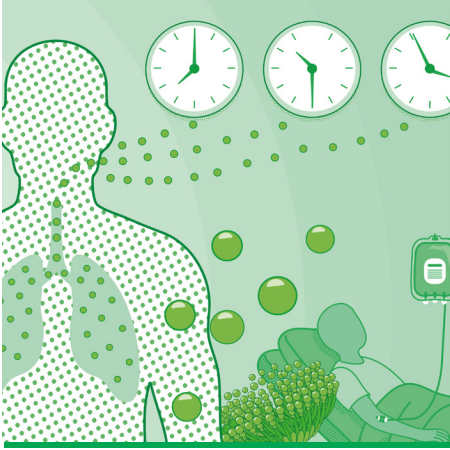




Platelia™ *Aspergillus* Ag Assay Automated Processing Quick Guide

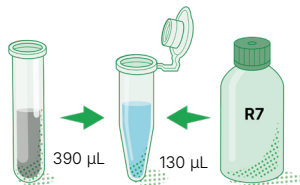
The Platelia *Aspergillus* Ag assay protocol is validated for automated processing on the Bio-Rad EVOLIS™ or EVOLIS Twin Plus™ Systems.

Important: Strict compliance with the EVOLIS and EVOLIS Twin Plus APF (Assay Protocol File) manuals is essential for successful automated processing. Bring reagents to room temperature (18–30°C) for at least 30 minutes before use.

Treatment of Samples & Controls

- 1 Mix 390 µL control sera or patient samples (sera and/or BAL fluids) with 130 µL R7 in polypropylene tubes.

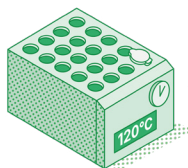
Control sera or patient sample



- 2 Tightly close tubes to prevent opening during heating.

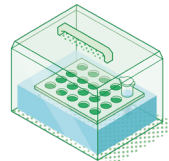
Heat Block

Vortex, then heat tubes for 6 min in a 120°C heat block.



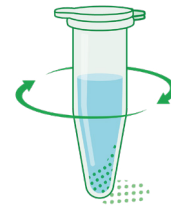
Water Bath

Vortex, then heat tubes for 3 min in a 100°C water bath.



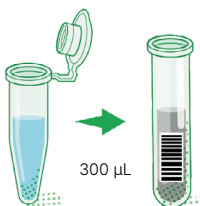
< or >

- 3 Centrifuge 10 min at 10,000 g. Supernatant will be used in step 4 and can be stored at 2–8°C for up to 48 hr prior to testing.



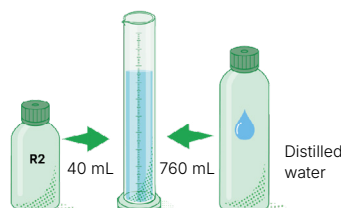
Preparation of Sample and Control Tubes

- 4 Transfer 300 µL supernatant of each treated sample or each treated control into 5 mL hemolysis tubes. Affix appropriate barcode labels onto tubes.



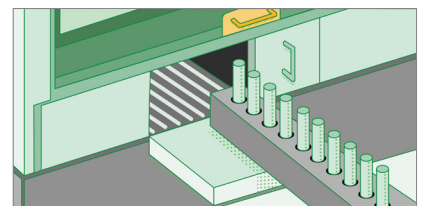
Preparation of Washing Solution

- 5 Dilute R2 to 1/20 in distilled water. The total volume will depend on your testing needs.



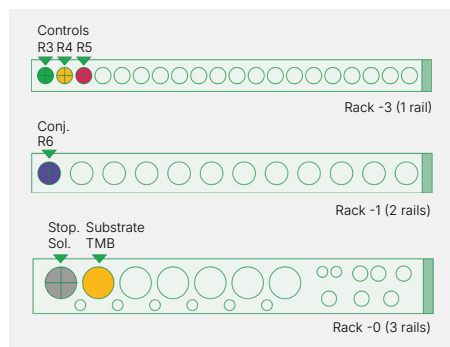
Load Samples, Controls, Reagents, and Buffers

- 6 Load barcoded treated sample tubes onto rack T of the EVOLIS System and rack \$T of the EVOLIS Twin Plus System.

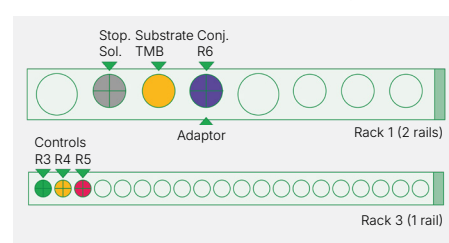


- 7** Place controls and reagents onto their appropriate racks for the EVOLIS System or EVOLIS Twin Plus System.
- Load treated control tubes (R3, R4, R5) on rack -3 of EVOLIS System; rack 3 of EVOLIS Twin Plus System
 - Load conjugate R6 on rack -1 of EVOLIS System; rack 1 of EVOLIS Twin Plus System
 - Load substrate R9 and stop solution R10 on rack -0 of EVOLIS System; rack 1 of EVOLIS Twin Plus System

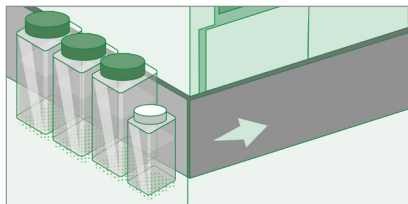
**Platelia Aspergillus Ag Assay
EVOLIS System Rack Configuration**



**Platelia Aspergillus Ag Assay
EVOLIS Twin Plus System Rack Configuration**

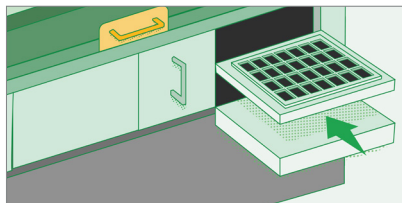


- 8** Load wash buffer, cleaning fluid, and tip racks.



Load Microplate and Start Run

- 9** Insert the plate on the plate holder and load on the EVOLIS System or EVOLIS Twin Plus System.



Automatic Reading and Interpretation

- 10** Validation criteria have been optimized for automated processing to improve the performance of the assay. Please see the full details in the user manuals: EVOLIS APF and EVOLIS Twin Plus APF.

Ordering Information

Catalog #	Description	Instruments
62794	Platelia Aspergillus Ag Assay, 96 tests	89818B EVOLIS System, 1 unit 93500B EVOLIS Twin Plus System, 1 unit
62792	Platelia Aspergillus Ag Verification Panel, 6-member panel, 3 tests	89611 EVOLIS 300 µL Tips, 18 x 960 89612 EVOLIS 1,100 µL Tips, 10 x 960

Visit [bio-rad.com/aspergillus-ag](https://www.bio-rad.com/aspergillus-ag) for more information.

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. Platelia, EVOLIS, and EVOLIS Twin Plus are trademarks of Bio-Rad Europe GmbH in certain jurisdictions. © 2026 Bio-Rad Laboratories, Inc. Supporting documents are available at downloads.bio-rad.com. EVOLIS APF—Product File Identifier 0001387 EVOLIS Twin Plus APF—Product File Identifier 0001386



**Bio-Rad
Laboratories, Inc.**

**Clinical Diagnostics
Group**

Website www.bio-rad.com/diagnostics **U.S.** 1 800 224 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 55 11 3065 7550 **Canada** 1 800 361 1808 **China** +86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23 **France** 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Greece** 30 210 7774396 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 00 800 00 24 67 23 **Italy** 00 800 00 24 67 23 **Japan** 81 3 6361 7070 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 55 5488 7670 **The Netherlands** 00 800 00 24 67 23 **New Zealand** 64 9415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23 **Republic of Korea** 82 080 007 7373 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3170 **South Africa** 27 11 442 8508 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23 **Switzerland** 00 800 00 24 67 23 **Taiwan** 886 2 2578-7189 **Thailand** 662 651 8311 **United Arab Emirates** 971 4 818 7300 **United Kingdom** 00 800 00 24 67 23