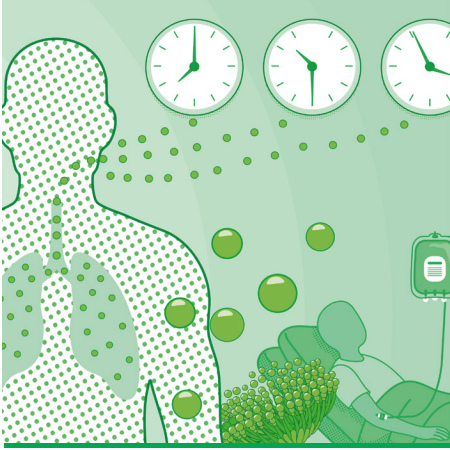




Platelia™ *Aspergillus* Ag Assay Manual Processing Quick Guide

This protocol can be used for Platelia *Aspergillus* Ag testing on Bio-Rad's stand-alone systems:



IPS Incubator



PW 41 Washer

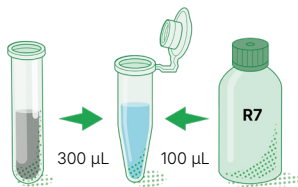
PR 4100 Reader
and Mini-PC

Important: Strict compliance with the technical procedure is essential for the success of this test.
Bring reagents to room temperature (18–30°C) for at least 30 minutes before use.

Step 1. Sample/Control Treatment

- 1 Mix 300 µL control sera and patient samples (sera and/or BAL fluids) with 100 µL R7.

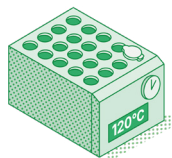
Control sera or patient sample



- 2 Tightly close tubes to prevent opening during heating.

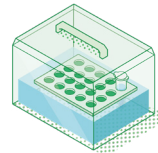
Heat Block

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. Close lid.



< or >

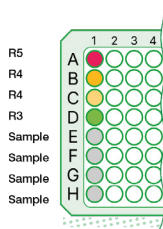
- 3 Centrifuge 10 min at 10,000 g.

Use supernatant in step 6.
It can be stored at 2–8°C for up to 48 hr prior to testing.

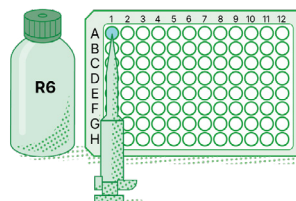


Step 2: EIA Procedure

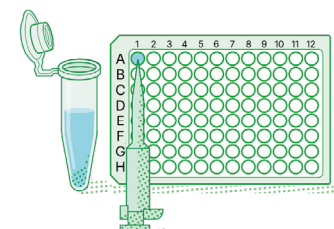
- 4 Refer to this plate map for placement of control sera (R5, R4, R3) and patient samples in step 6.



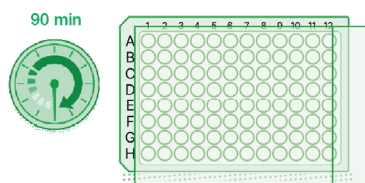
- 5 Dispense 50 µL of ready-to-use conjugate (R6) to each well.



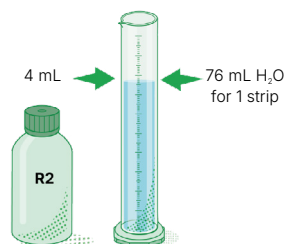
- 6 Add 50 µL supernatant (from treated controls/patient samples) to each well.



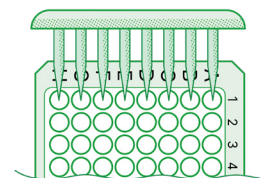
- 7** Apply plate sealer and incubate microplate for 90 ± 5 min at $37 \pm 1^\circ\text{C}$.



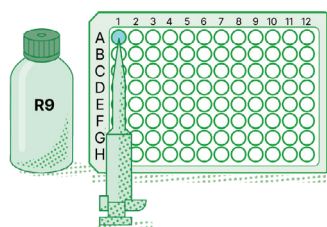
- 8** Prepare the Working Washing Solution: dilute R2 to 1:20 in distilled water.



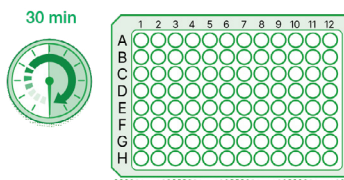
- 9** Use the Microplate Washer to aspirate and wash plate 5 times with Working Washing Solution (800 μL and well-bottom washing is recommended).



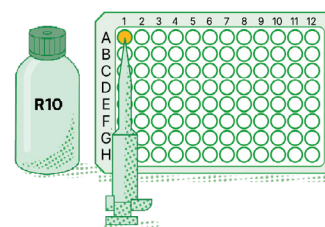
- 10** Quickly add 200 μL of ready-to-use Chromogen TMB Solution (R9) to each well, avoiding exposure to bright light.



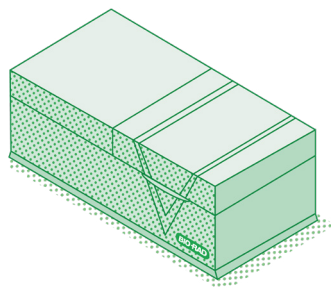
- 11** Incubate microplate **without plate sealer** in the dark at room temperature ($18\text{--}30^\circ\text{C}$) for 30 ± 5 min.



- 12** Add 100 μL of ready-to-use Stopping Solution (R10) to each well. Wipe the microplate bottom carefully.



- 13** Read the OD of wells at 450/620 nm. Microplates must be read within 30 min of adding Stopping Solution.



Ordering Information

Catalog #	Description
12014064	Platelia Aspergillus Ag Assay, 96 tests
62792	Platelia Aspergillus Ag Verification Panel, 3 x 6 tests

Instruments

87530	IPS Microplate Incubator, 1 unit
85508	PW 41 Microplate Washer, 1 unit
91587	PR 4100 Absorbance Microplate Reader, 1 unit
12020451	PR 4100 US Magellan V5.1 Dell 3540 PC, 1 unit

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

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. Platelia is a trademark of Bio-Rad Europe GmbH in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2026 Bio-Rad Laboratories, Inc.

To learn more about the Platelia Aspergillus Ag Assay, contact a Bio-Rad sales representative.

BIO-RAD

**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