



# Platelia Aspergillus Ag Assay Quick Guide Manual Processing

## Bio-Rad Validated Stand-Alone Systems for Manual Processing

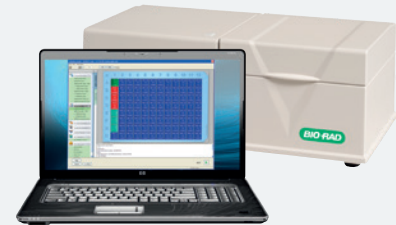
This protocol can be used for Platelia Aspergillus Ag testing on Bio-Rad's stand-alone systems.



IPS Incubator



PW 41 Washer



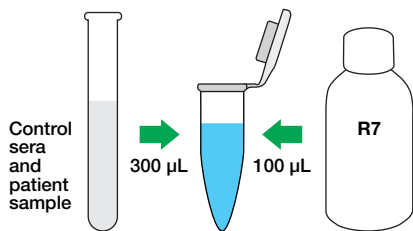
PR 4100 Reader

**Important Note: Strict compliance with the technical procedure is essential for the success of this test.**

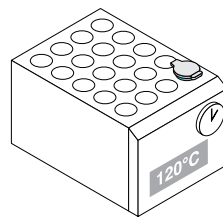
Bring reagents to room temperature (+18–25°C) for at least 30 minutes before use.

### STEP 1: Sample/Control Treatment

- Mix 300  $\mu$ L control sera and patient samples (sera and/or BAL fluids) with 100  $\mu$ L R7



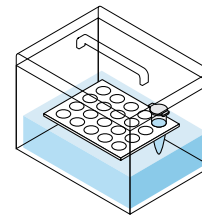
- Tightly close tubes to prevent opening during heating



Heat Block

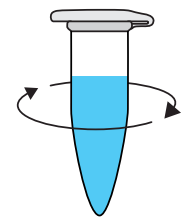
Vortex, then heat tubes for 6 min in a 120°C heat block.

< OR >



Water Bath

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

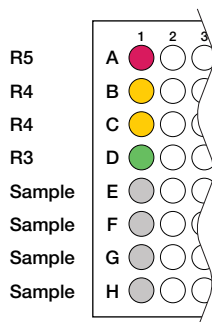


- Centrifuge 10 min at 10,000 g

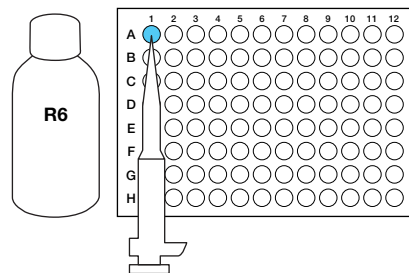
Use supernatant in Step 6. It can be stored at 2–8°C for up to 48h prior to testing.

### STEP 2: EIA Procedure

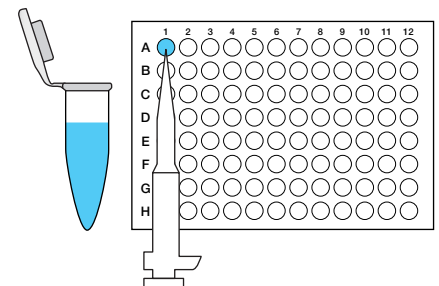
- Refer to this plate map for placement of control sera (R5, R4, R3) and patient samples in Step 6



- Dispense 50  $\mu$ L of ready-to-use conjugate (R6) to each well



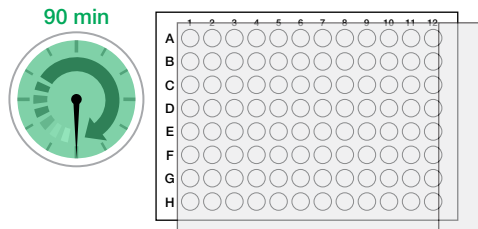
- Add 50  $\mu$ L supernatant (from treated controls/patient samples) to each well



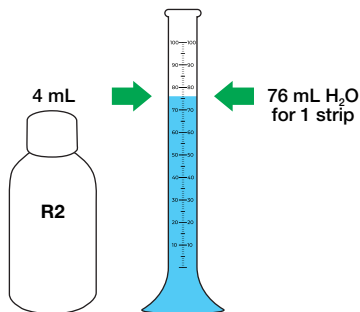
**Important Note: Strict compliance with the technical procedure is essential for the success of this test.**

**STEP 2: EIA Procedure (continued)**

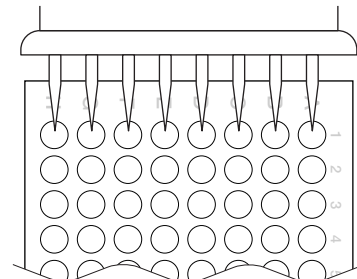
**7** Apply plate sealer and incubate microplate for 90 (±5) min at 37°C (±1°C)



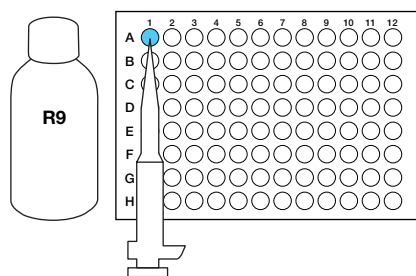
**8** Prepare the Working Washing Solution: dilute R2 to 1:20 in distilled water



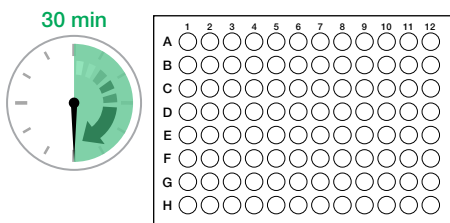
**9** Use the Microplate Washer to aspirate and wash plate 5 times with Working Washing Solution (800 µL and well-bottom washing is recommended)



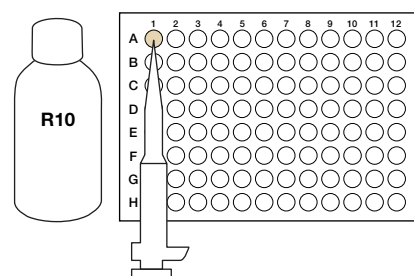
**10** Quickly add 200 µL of ready-to-use Chromogen TMB Solution (R9) to each well, avoiding exposure to bright light



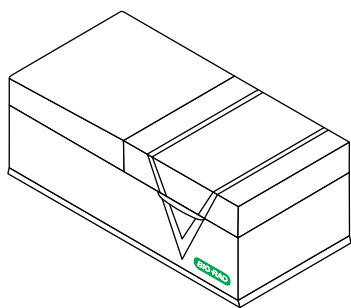
**11** Incubate microplate without plate sealer in the dark at room temperature (18–25°C) for 30 ± 5 min



**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 No.	Description	
12014064	Platelia Aspergillus Ag Assay.....	96 tests
62792	Platelia Aspergillus Ag Verification Panel .....	6-member panel, 3 tests

**Instruments**

87530	IPS Microplate Incubator .....	1 unit
85508	PW 41 Microplate Washer .....	1 unit
91587	PR 4100 Absorbance Microplate Reader.....	1 unit
12020451	PR4100 US Magellan V5.1 Dell 3540 PC.....	1 unit

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc.  
Platelia is a trademark of Bio-Rad Europe GmbH in certain jurisdictions.



**Bio-Rad  
Laboratories, Inc.**

*To learn more about the Platelia Aspergillus Ag Assay, contact a Bio-Rad Sales Representative.*

Clinical  
Diagnostics Group

**Website** www.bio-rad.com/diagnostics **U.S.** 1 800 224 6723 **Australia** +61 (2) 9914 2800 **Austria** +43 (0) 1 877 89 01 9 **Belgium** +32 (0) 3 710 53 00  
**Brazil** +55 11 3065 7550 **Canada** +1 514 334 4372 **China** +86 21 6169 8500 **Czech Republic** +420 241 431 660 **Denmark** +45 44 52 10 00  
**Finland** +358 9 804 22 00 **France** +33 (0) 1 47 95 60 00 **Germany** +49 (0) 89 31884 393 **Greece** +30 210 7774396 **Hong Kong** +85 2 2789 3300  
**Hungary** +36 1 459 6190 **India** +91 124 4029300 **Israel** +972 03 963 6025 **Italy** +39 024 94 86 600 **Japan** +81 3 6361 7070 **Korea** +82 080 007 7373  
**Mexico** +52 (55) 5488 7670 **The Netherlands** +31 (0) 318 540 666 **New Zealand** +64 (9)415 2280 **Norway** +47 23 38 41 30 **Poland** +48 22 331 99 99  
**Portugal** +351 21 47 27 700 **Russia** +7 (495) 721-14-04 **Singapore** +65 6415 3170 **South Africa** +27 11 442 8508 **Spain** +34 91 490 6580  
**Sweden** +46 844 98053 **Switzerland** +41 (0) 61 717 9555 **Taiwan** +886 (2) 2578-7189 **Thailand** (662) 651 8311 **United Kingdom** +44 (0) 1923 471301