



Platelia™ *Aspergillus* Ag Kit

Early detection is key for the diagnosis and treatment of Invasive Aspergillosis

The Galactomannan Test

- Detection of the *Aspergillus* galactomannan antigen, six days before onset of clinical symptoms and 10 days before diagnosis¹
- Results in three hours
- EIA microplate format with ready-to-use reagents and controls

¹ Marr KA, et al. *J Infect Dis* 2004 Aug 1; 190(3):641–649.

FDA-cleared for serum and BAL specimens from adults and pediatric patients

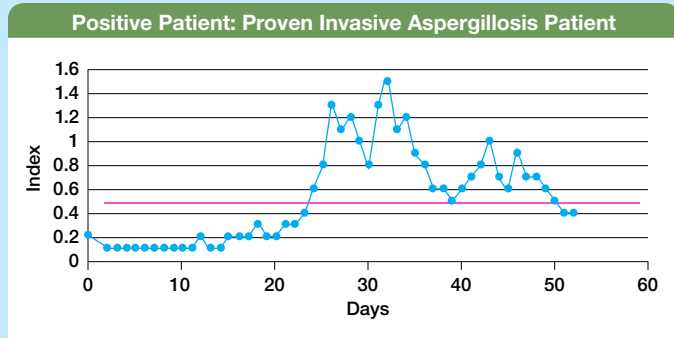


Platelia™ Aspergillus Ag Kit

Testing and Result Interpretation

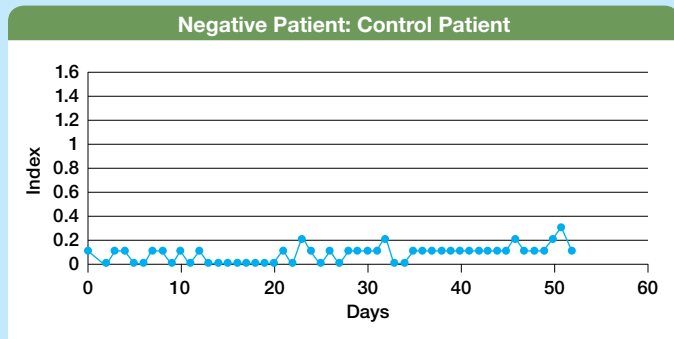
Positive Result = Index ≥ 0.5

- For all positive patient results, re-testing the same Serum/BAL sample (new aliquot) is recommended
- Regular screening (twice weekly) of serum samples of high risk patients is recommended
- Positive results should be considered in conjunction with other diagnostic procedures



Negative Result = Index < 0.5

- Repeat testing is recommended if result is negative but disease is suspected



Convenient

- 96 well microplate with breakable strips
- Ready-to-use colored conjugate

Objective

- Results reported as index

Efficient

- Excellent sensitivity and specificity
- Faster initiation of appropriate treatment

Required Products (not supplied by Bio-Rad)

- Dry heat block or water bath
- Clean work area, away from air vents and open windows
- Centrifuge

Ordering Information

Catalog No. Description

62794	Bio-Rad Platelia™ Aspergillus Ag Kit.....	96 tests
62792	Aspergillus Serum Proficiency Panel.....	6 member panel
25156	Aspergillus Start-Up Package, includes: proficiency panel, microplate assay testing reference manual, material safety data sheet, quick reference guide	1 package



Bio-Rad Laboratories

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

Clinical Diagnostics Group

Website www.bio-rad.com/diagnostics U.S. 1-800-2BIO-RAD Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium 32-9-385-5511 Brazil 5521-3237-9400 Canada 1-514-334-4372 China 86-21-64260808 Czech Republic 420-241-430-532 Denmark +45-4452-1000 Finland 358-9-804-22-00 France 33-1-47-95-60-00 Germany +49-(0)89-318-840 Greece 30-210-7774396 Hong Kong 852-2789-3300 Hungary +36-1-459-6100 India 1-800-180-1224 Israel 972-3-9636050 Italy +39-02-216091 Japan 81-3-6361-7070 Korea 82-2-3473-4460 Mexico +52 (55) 5488-7670 The Netherlands +31-318-540666 New Zealand 64-9-415-2280 Norway 47-23-38-41-30 Poland 48-22-3319999 Portugal 351-21-472-7700 Russia 7-495-721-14-04 Singapore 65-6415-3170 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-(0)20-8328-2000