



BioPlex® 2200 Antiphospholipid Syndrome (APLS) IgG, IgM, IgA Kits

Simultaneous Detection of Autoantibodies for Antiphospholipid Syndrome (APS)

The BioPlex® 2200 offers the highest throughput, fully-automated, random access, multiplex platform for the accurate and timely diagnosis of primary and secondary Antiphospholipid Syndrome (APS).

Bio-Rad's multiplex technology allows consolidated testing of the four first-line serological assays, as recommended in the laboratory criteria for the diagnosis of APS¹—anti-Cardiolipin (aCL) IgG and IgM and anti-Beta-2 Glycoprotein I (a β 2GPI) IgG and IgM. This high level of automation provides the flexibility and improved workflow to combine testing of second-line assays that are frequently detected in specific patient subgroups¹—aCL IgA and a β 2GPI IgA.

Generating a full profile of clinically relevant APS antibodies facilitates identification of patients with multiple APS autoantibody positivity associated with a higher risk of thrombosis¹.

APLS IgG Kit Includes:

- Anti-Cardiolipin IgG
- Anti-Beta-2 Glycoprotein I IgG

APLS IgM Kit Includes:

- Anti-Cardiolipin IgM
- Anti-Beta-2 Glycoprotein I IgM

APLS IgA Kit Includes:

- Anti-Cardiolipin IgA
- Anti-Beta-2 Glycoprotein I IgA

BIO-RAD

BioPlex® 2200 Antiphospholipid Syndrome (APLS) IgG, IgM, IgA Kits

Enhanced Efficiency

- The highest throughput, fully-automated, random access platform available for efficient APS testing
 - Fully-automated, random access processing of up to 200 results per hour provides faster turnaround time by eliminating batched runs
 - One small 5 µL sample generates up to 2 results per kit providing more tests with less sample volume
 - 30-day calibration curve eliminates the need for daily calibration
 - 2 analytes per kit reduces kit inventory
- Multiplexed assays allow for simultaneous detection of multiple antibodies
 - Parallel detection of aCL and aβ2GPI provides physicians with more diagnostic information that aids in the assessment of APS
 - Simultaneous APS antibody measurement facilitates identification of patients with a higher risk of thrombosis through multi-marker positivity
 - Accommodates first-line and second-line APS testing algorithms for more flexible workflow options

Confidence in Patients Results

- Assay cutoffs set at 99th percentile as recommended by laboratory criteria for the diagnosis of APS¹
 - Excellent diagnostic accuracy: high specificity minimizing false-positive results
 - Clinically relevant test results for improved patient management

Excellent Assay Performance

- Excellent precision, total %CV's are less than 10.7% for all 6 assays, providing accurate and reproducible results minimizing repeat testing

¹ Miyakis S. et al. *International Consensus Statement on an Update of the Classification Criteria for Definite Antiphospholipid Syndrome (APS)*. *J Thromb Haemost.* 2006; 4(2): 295-306.

Ordering Information

Catalog No.	Description	
665-1950	BioPlex® 2200 APLS IgG Reagent Pack, 100 samples.....	1 pack
663-1900	BioPlex® 2200 APLS IgG Calibrator Set.....	1 set
663-1930	BioPlex® 2200 APLS IgG Control Set.....	2 sets
665-2050	BioPlex® 2200 APLS IgM Reagent Pack, 100 samples.....	1 pack
663-2000	BioPlex® 2200 APLS IgM Calibrator Set.....	1 set
663-2030	BioPlex® 2200 APLS IgM Control Set.....	2 sets
665-2150	BioPlex® 2200 APLS IgA Reagent Pack, 100 samples.....	1 pack
663-2100	BioPlex® 2200 APLS IgA Calibrator Set.....	1 set
663-2130	BioPlex® 2200 APLS IgA Control Set.....	2 sets



**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. 1-800-224-6723 Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium 32-9-385-5511 Brazil +55 (31)3689-6600 Canada 1-514-334-4372 China 86-21-61698500 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 1800-180-1224 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 (0)26-674-55-05/06 Taiwan 886-2-2578-7189 Thailand 662-651-8311 United Kingdom +44 (0)20-8328-2000