



BioPlex[®] 2200 Celiac IgA and IgG Kits

Comprehensive test kits for confidence in the laboratory assessment of celiac disease

Bio-Rad Laboratories offers state-of-the-art serological tests for the accurate and timely diagnosis of celiac disease. The key diagnostic tests: tissue Transglutaminase (tTG) and deamidated gliadin peptide (DGP) are multiplexed together in two test kits for the simultaneous measurement of IgA or IgG autoantibodies.

The IgA kit includes the unique IgA Verification Bead (AVB) to flag results for samples with deficient IgA levels. This condition is more prevalent in celiac disease and can confound the interpretation of Celiac IgA serology results in other methods.

The Celiac IgA kit includes:

- Anti-tTG IgA
- Anti-DGP IgA
- IgA Verification Bead

The Celiac IgG kit includes:

- Anti-tTG IgG
- Anti-DGP IgG

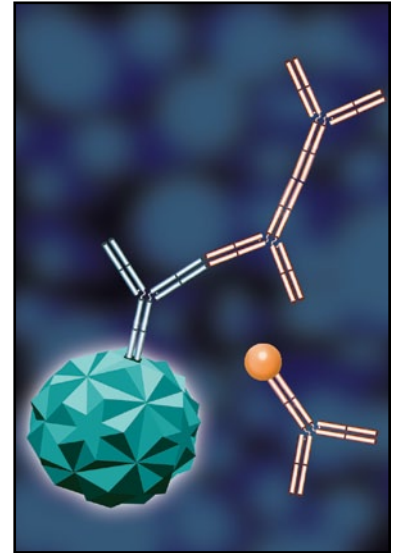
BioPlex® 2200 Celiac IgA and IgG Kits

Confidence in Patient Results

- Unique Celiac IgA Verification Bead (AVB) flags IgA deficient samples for appropriate follow-up testing
- Internal quality control beads (SVB, ISB and AVB) generate a higher level of quality assurance for patient results

Improved Workflow and Random Access

- Random access improves laboratory turnaround time eliminating batch testing
- Multiplexed measurement of Anti-tTG and Anti-DGP
 - Small 5µL sample volume for each kit
 - Reduces reagent kit inventory
- The highest throughput platform available for Celiac testing
 - Up to 200 test results per hour for Celiac testing



AVB facilitates reflex testing to the Celiac IgG kit. Such samples could be falsely interpreted as negative by traditional immunoassays.

Excellent Analytical Performance

- Excellent precision, sensitivity and specificity assures accuracy in patient results and reduces repeat testing

Ordering Information

Catalog No.	Description	
665-2350	BioPlex® 2200 Celiac IgA Reagent Kit	1 pack
663-2300	BioPlex® 2200 Celiac IgA Calibrator Set.....	1 set
663-2330	BioPlex® 2200 Celiac IgA Control Set.....	1 set
665-2250	BioPlex® 2200 Celiac IgG Reagent Kit.....	1 pack
663-2200	BioPlex® 2200 Celiac IgG Calibrator Set	1 set
663-2230	BioPlex® 2200 Celiac IgG Control Set	1 set



**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