

# BioPlex<sup>®</sup> 2200 Autoimmune Panels

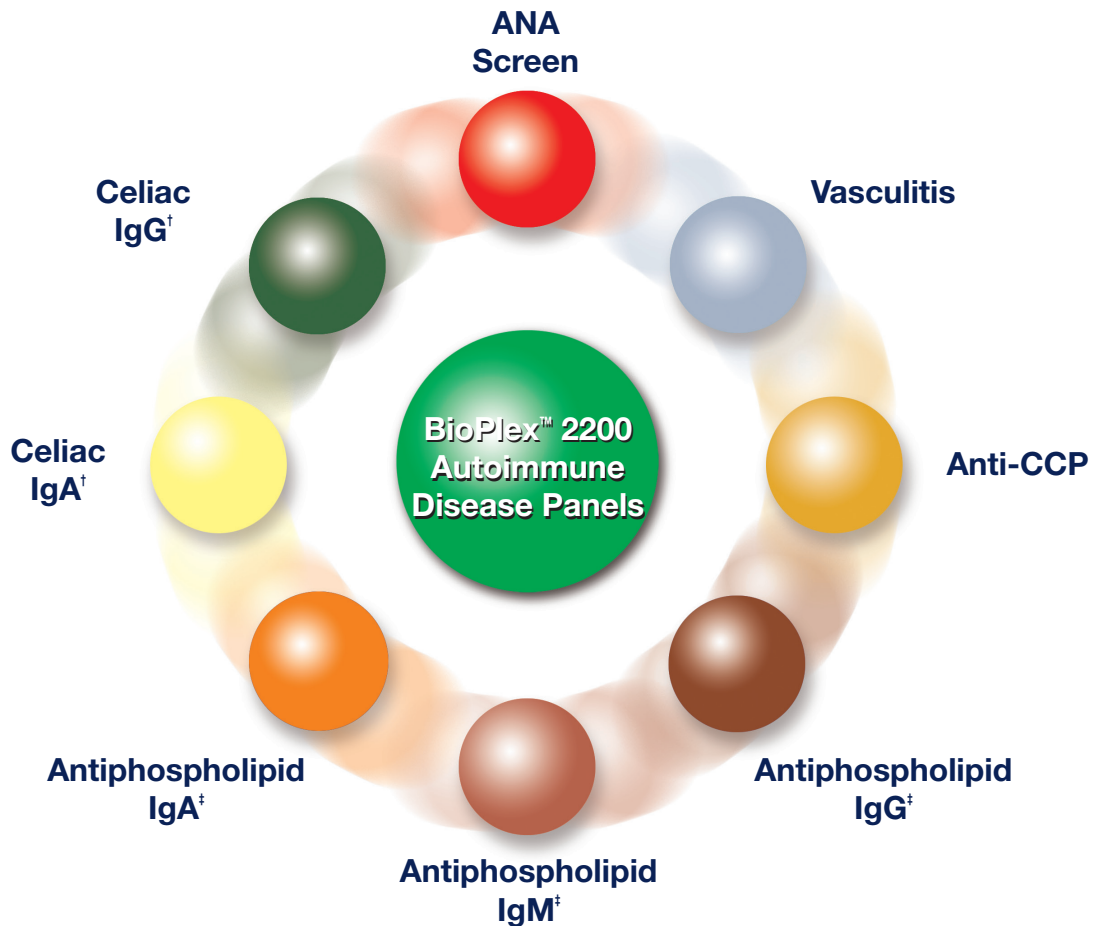
**Bringing a new level of automation and productivity to autoimmune testing with a complete line of disease-specific multiplex panels**

ANA Screen  
Vasculitis  
Anti-CCP  
Antiphospholipid IgG<sup>†</sup>

Antiphospholipid IgM<sup>†</sup>  
Antiphospholipid IgA<sup>‡</sup>  
Celiac IgG<sup>†</sup>  
Celiac IgA<sup>‡</sup>

<sup>†</sup> In development  
<sup>‡</sup> Available outside US only





# BioPlex<sup>®</sup> 2200 Autoimmune Disease Panels

## Leading the way with an expanding multiplexed assay menu

Comprehensive assay menu offering a multiplexed solution for simultaneous detection and identification of the most clinically relevant autoantibodies.

### Offering the benefits of multiplex testing:

- High throughput multiplex testing improves workflow efficiencies and eliminates hands-on steps
- Provides flexibility to screen and confirm specific autoantibodies with one kit using one small patient sample (5  $\mu$ L)
- Consolidates traditional single-analyte tests into one panel
- Time-to-first result is approximately 45 minutes
- Ready-to-use multiplexed calibrators and controls, with curve stability up to 30 days

† In development  
‡ Available outside US only

## Systemic Autoimmune Diseases

### BioPlex® 2200 ANA Screen with MDSS Panel



The BioPlex® 2200 ANA Screen kit will help the clinician to better diagnose systemic autoimmune diseases by detecting only clinically significant autoantibodies.

### BioPlex® 2200 ANA Screen with MDSS

#### Analytes

- ANA Screen\*
- dsDNA (quantitative)
- Chromatin
- Ribosomal Protein
- SS-A (52 and 60)\*\*
- SS-B
- Sm
- Sm/RNP
- RNP (A and 68)\*\*
- Scl-70
- Jo-1
- Centromere B

\*ANA Screen is a composite roll up of all individual results  
\*\* Individual results available outside US

## Systemic Small Vessel Vasculitis

### BioPlex® 2200 Vasculitis Panel



The BioPlex® 2200 Vasculitis kit will help the clinician diagnose some of the most severe autoimmune systemic vasculitic diseases affecting blood vessels in the lungs and kidneys.

### BioPlex® 2200 Vasculitis

#### Analytes

- Anti-MPO
- Anti-PR3
- Anti-GBM

## Rheumatoid Arthritis

### BioPlex® 2200 Anti-CCP



The BioPlex® 2200 Anti-CCP kit aids the clinician in the diagnosis of rheumatoid arthritis by providing important information regarding disease severity and prognosis.

### BioPlex® 2200 Anti-CCP

#### Analytes

- Anti-Cyclic Citrullinated Peptide (CCP) IgG

## Antiphospholipid Syndrome

### BioPlex® 2200 Antiphospholipid Syndrome (APLS) Panels

### BioPlex® 2200 APLS IgG<sup>†</sup>

#### Analytes

- Anti-Cardiolipin IgG
- Anti-Beta 2 Glycoprotein I (β2GPI) IgG

### BioPlex® 2200 APLS IgM<sup>†</sup>

#### Analytes

- Anti-Cardiolipin IgM
- Anti-Beta 2 Glycoprotein I (β2GPI) IgM

### BioPlex® 2200 APLS IgA<sup>†</sup>

#### Analytes

- Anti-Cardiolipin IgA
- Anti-Beta 2 Glycoprotein I (β2GPI) IgA

## Celiac Disease

### BioPlex® 2200 Celiac Panels

### BioPlex® 2200 Celiac IgA<sup>†</sup>

#### Analytes

- Anti-Deamidated Gliadin IgA
- Anti-tissue transglutaminase (tTG) IgA
- IgA Verification Bead - flags IgA deficient samples

### BioPlex® 2200 Celiac IgG<sup>†</sup>

#### Analytes

- Anti-Deamidated Gliadin IgG
- Anti-tissue transglutaminase (tTG) IgG

† In development  
‡ Available outside US only

**Autoimmune Panels**

**Ordering Information**

**Catalog No. Description**

**BioPlex® 2200 ANA Screen with MDSS Panel**

665-1150	BioPlex® 2200 ANA Screen Reagent Pack.....	1 pack
663-1100	BioPlex® 2200 ANA Screen Calibrator Set.....	1 set
663-1130	BioPlex® 2200 ANA Screen Control Set.....	1 set

**BioPlex® 2200 Vasculitis Panel**

665-1850	BioPlex® 2200 Vasculitis Reagent Pack.....	1 pack
663-1800	BioPlex® 2200 Vasculitis Calibrator Set.....	1 set
663-1830	BioPlex® 2200 Vasculitis Control Set.....	1 set

**BioPlex® 2200 Anti-CCP**

665-3250	BioPlex® 2200 Anti-CCP Reagent Pack.....	1 pack
663-3200	BioPlex® 2200 Anti-CCP Calibrator Set.....	1 set
663-3230	BioPlex® 2200 Anti-CCP Control Set.....	1 set

**BioPlex® 2200 Antiphospholipid Syndrome (APLS) Panels**

665-1950	BioPlex® 2200 APLS IgG† Reagent Pack.....	1 pack
663-1900	BioPlex® 2200 APLS IgG† Calibrator Set.....	1 set
663-1930	BioPlex® 2200 APLS IgG† Control Set.....	1 set
665-2050	BioPlex® 2200 APLS IgM† Reagent Pack.....	1 pack
663-2000	BioPlex® 2200 APLS IgM† Calibrator Set.....	1 set
663-2030	BioPlex® 2200 APLS IgM† Control Set.....	1 set
665-2150	BioPlex® 2200 APLS IgA† Reagent Pack.....	1 pack
663-2100	BioPlex® 2200 APLS IgA† Calibrator Set.....	1 set
663-2130	BioPlex® 2200 APLS IgA† Control Set.....	1 set

**BioPlex® 2200 Celiac Panels in Development**

BioPlex® 2200 Celiac IgG†

BioPlex® 2200 Celiac IgA†

† In development  
‡ Available outside US only



**Bio-Rad  
Laboratories**

*For further information, please contact the Bio-Rad office nearest you or visit our website at [www.bio-rad.com/diagnostics](http://www.bio-rad.com/diagnostics)*

*Clinical  
Diagnostics Group*

**Website** [www.bio-rad.com/diagnostics](http://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