

## Efficiency



### High Throughput Capability

Up to 300 results per hour for improved workflow and reduced hands-on time and labor costs.

### The Power of One Panel

Consolidating three traditional single-analyte tests into one test panel significantly reduces turnaround time and reduces the cost per patient result.

Rapid, multi-parameter antibody profiling including STAT capabilities from a single reaction vessel facilitating disease differentiation and treatment.

## Quality



### Antibodies of Clinical Significance

By targeting the most clinically relevant and prevalent antibodies associated with systemic small vessel vasculitis, the BioPlex<sup>™</sup>2200 Vasculitis kit improves clinical specificity, thus reducing the number of false positive results.

### Novel Multiplex Chemistry

Magnetic bead technology enables thorough washing for improved signal to noise ratios and better overall performance compared to homogeneous methods.

## Confidence



### Full Automation

From primary tube to final result, the BioPlex<sup>™</sup>2200 is the only fully-automated, multiplexed Vasculitis testing system featuring:

- Completely automated assay procedure minimizing human errors commonly observed with manual sample and reagent handling
- Standardized assay procedure for improved performance
- Objective results, thus eliminating subjectivity associated with IFA methods

### eFlex<sup>™</sup> Software

The system features a powerful package of user-friendly software tools that provides optimum flexibility.

### Key Features Include:

- “Add-on” of new test orders from previously run patient samples
- Creation of custom Vasculitis panels
- Flexible setup options to create custom result
- Bi-directional LIS connectivity

### Excellent Performance Compared to Traditional EIA and ANCA IFA Methods

The BioPlex<sup>™</sup>2200 Vasculitis kit features the combination of a heterogeneous assay format and a minimum of 150 measurements per analyte per reaction vessel (vs. one photometric EIA measurement) for excellent and consistent lot-to-lot results.

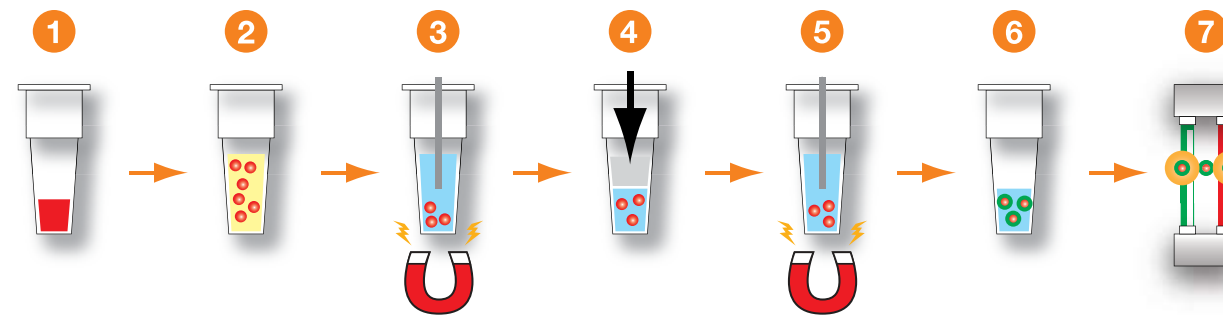
### Internal Quality Controls

For every sample processed, three internal quality controls are also processed, improving confidence in results and as an aid in diagnosis:

- Serum Verification Bead (SVB) confirms presence of serum
- Internal Standard Bead (ISB) standardizes detector performance
- Reagent Blank Bead (RBB) identifies sample problems associated with non-specific binding

## BIOPLEX 2200 SYSTEM

### Vasculitis Assay Procedure



1. Patient sample (5 µL) automatically added to reaction vessel
2. Sample diluent and beads (dilution 1:20) added and incubated at 37°C
3. Wash step
4. Conjugate added and incubated at 37°C
5. Wash step
6. Bead resuspended in wash buffer
7. Flow-based, dual laser detection



### Vasculitis Reagent Pack

Contains all the necessary reagents (sample diluent, beads and conjugate) to process 100 samples and up to 3 results per sample.



### Vasculitis Calibrator Set

Calibrators are ready-to-use and multiplexed. A point-to-point curve fit, using four calibrators, is used to calculate semi-quantitative results.



### Vasculitis Control Set

Quality Controls are ready-to-use and multiplexed. The set includes a negative control as well as a multi-analyte positive control containing antibodies present for analytes within the Reagent Pack.

## Ordering Information

Catalog No.	Description	BioPlex <sup>™</sup> 2200 System and Bulk Reagents
665-1850	Vasculitis Reagent Pack.....	1 pack
663-1800	Vasculitis Calibrator Set.....	1 set
663-1830	Vasculitis Control Set.....	1 set
660-0000	BioPlex <sup>™</sup> 2200.....	1 instrument
660-0800	Sample Racks (30 per box).....	1 box
660-0801	Sample Tray.....	1 tray
660-0817	Instrument Sheath Fluid (2 per box).....	1 box
660-0818	Instrument Wash Buffer.....	1 box
660-2003	Reaction Vessels (1000 per bag).....	1 bag
666-0001	Instrument Detector Calibration Pack....	1 pack
666-0002	Instrument Detector Clean Pack.....	1 pack
666-0003	Instrument Probe Cleaning Solution Bottles (4 per box).....	1 box



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 +49-089-318-840 Greece 30-210-7774395 Hong Kong 852-2789-3300 Hungary +36-1-459-5100 India 91-124-4029300 Israel 972-3-9636050 Italy +39-02-216991 Japan 81-3-6361-7070 Korea 82-2-3473-4460 Mexico 52565200-0520 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 65-6415-3188 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

Printed on recycled paper with soy-based inks

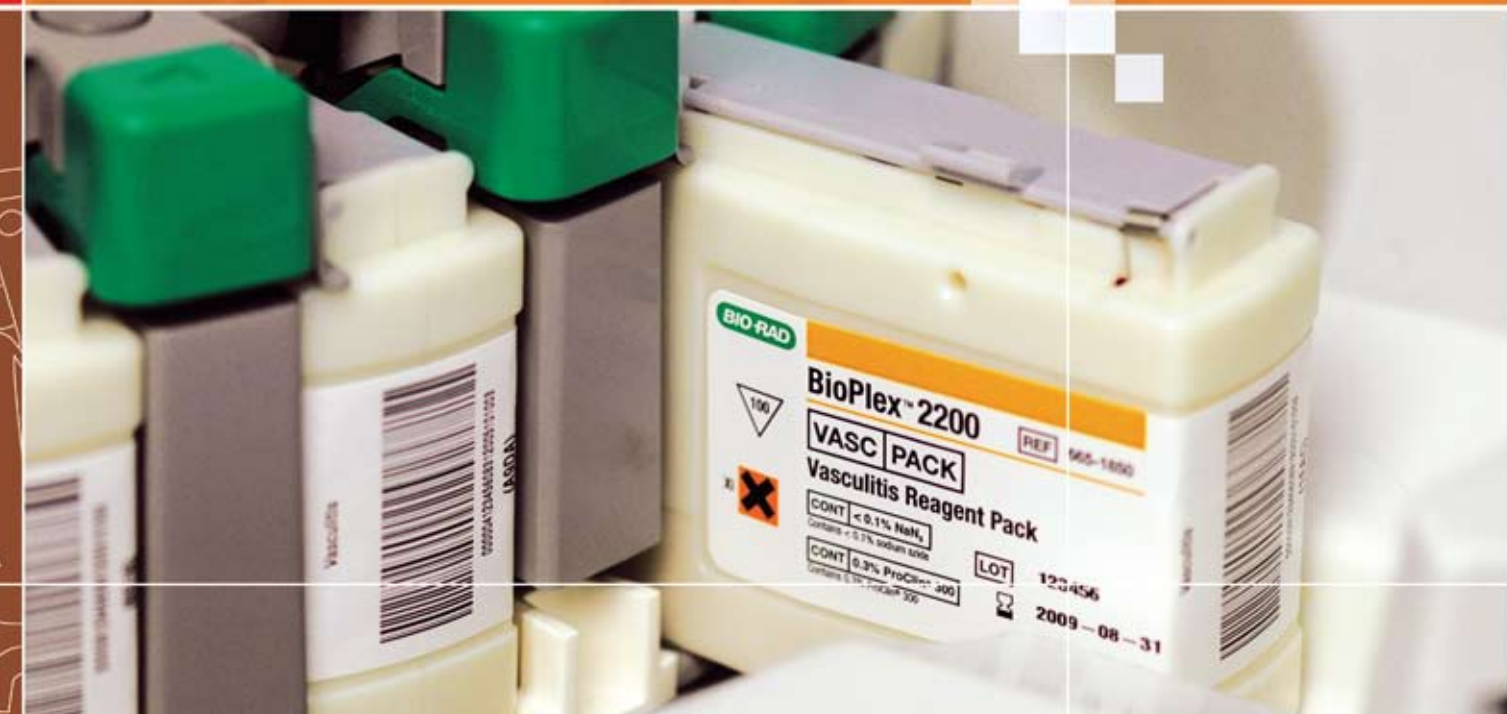
© 2009 Bio-Rad Laboratories, Inc. Printed in USA 09/09 DG09-0758 T-128



# BioPlex<sup>™</sup> 2200 Vasculitis

The first and only fully-automated, random access multiplexed solution for vasculitis testing





Like no other

The BioPlex<sup>™</sup>2200 Vasculitis kit simultaneously detects the three most clinically relevant antibodies.

By detecting the most clinically relevant vasculitis antibodies, the BioPlex<sup>™</sup>2200 Vasculitis kit provides laboratories with more diagnostically useful information for enhanced patient management with highly reproducible results and faster turnaround time.

The BioPlex<sup>™</sup>2200 Vasculitis kit reports results for the following:

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

## Comparison with EIA Methods

### Positive % Agreement

EIA Result	Retrospective Samples Positive for Anti-MPO, Anti-PR3 and Anti-GBM	BioPlex <sup>™</sup> 2200 Vasculitis Anti-MPO Result		BioPlex <sup>™</sup> 2200 Vasculitis Anti-PR3 Result		BioPlex <sup>™</sup> 2200 Vasculitis Anti-GBM Result	
		Positive % Agreement	Overall % Agreement	Positive % Agreement	Overall % Agreement	Positive % Agreement	Overall % Agreement
		93.9% (92/98)	93.0% (93/100)*	100% (79/79)	83.0% (83/100)†	88.9% (16/18)‡	92.6% (25/27)

\* One (1) anti-MPO EIA equivocal result is included in the Overall Agreement.  
 † Ten (10) anti-PR3 EIA equivocal results are included in the Overall Agreement.  
 ‡ Two (2) of the sixteen (16) BioPlex 2200 anti-GBM positive results were weak positive by anti-GBM EIA.  
 § One (1) of the two (2) BioPlex 2200 anti-GBM negative results was a weak positive by anti-GBM EIA

227 retrospective samples positive for anti-MPO (N=100), anti-PR3 (N=100) and anti-GBM (N=27) were tested with the BioPlex<sup>™</sup>2200 Vasculitis kit and the corresponding commercially available microplate EIA methods. The positive % agreement and overall % agreement between the two methods for each analyte are referenced in the table above. The BioPlex<sup>™</sup>2200 Vasculitis kit demonstrated **good positive % agreement** with commercially available EIA assays for each analyte.

### Negative % Agreement

EIA Result	Normal Blood Donors	Unselected Patient Samples*	BioPlex <sup>™</sup> 2200 Vasculitis Anti-MPO Result		BioPlex <sup>™</sup> 2200 Vasculitis Anti-PR3 Result		BioPlex <sup>™</sup> 2200 Vasculitis Anti-GBM Result	
			Negative % Agreement	Overall % Agreement	Negative % Agreement	Overall % Agreement	Negative % Agreement	Overall % Agreement
			100% (293/293)	100% (293/293)	100% (293/293)	100% (293/293)	99.3% (289/291)	98.6% (289/293)
			97.6% (284/291)	96.3% (289/300)	99.0% (292/295)	99.0% (297/300)	99.7% (298/299)	99.3% (298/300)

\*Previously tested with Vasculitis tests

293 samples from normal blood donors and 300 samples from unselected patient samples previously tested with vasculitis tests were tested with the BioPlex<sup>™</sup>2200 Vasculitis kit and the corresponding commercially available microplate EIA methods. The negative % agreement and overall % agreement between the two methods for each analyte are referenced in the table above. The BioPlex<sup>™</sup>2200 Vasculitis kit demonstrated **excellent negative % agreement** with commercially available EIA assays for each analyte.

## Comparison with IFA Method

### Positive % Agreement

pANCA IFA Result	Retrospective Samples Positive for Anti-MPO	BioPlex <sup>™</sup> 2200 Anti-MPO Result		cANCA IFA Result	Retrospective Samples Positive for Anti-PR3	BioPlex <sup>™</sup> 2200 Anti-PR3 Result	
		Positive % Agreement	Overall % Agreement			Positive % Agreement	Overall % Agreement
		93.3% (83/89)	84.0% (84/100)			94.9% (93/98)	93.0% (93/100)

200 retrospective positive samples (100 anti-MPO and 100 anti-PR3) were evaluated by a commercially available ANCA IFA method using ethanol-fixed slides. The positive % agreement and overall % agreement between the two methods are referenced in the adjacent tables. The BioPlex<sup>™</sup>2200 Vasculitis kit demonstrated **good positive % agreement** with the ANCA IFA method.

### BioPlex<sup>™</sup>2200 Vasculitis vs. EIA Methods

Combined p-/c-ANCA IFA Results	Combined BioPlex <sup>™</sup> 2200 Anti-MPO/Anti-PR3 Results		Combined Anti-MPO/Anti-PR3 EIA Results	
	Positive % Agreement	Overall % Agreement	Positive % Agreement	Overall % Agreement
	94.1% (176/187)	88.5% (177/200)	88.2% (165/187)	83.0% (166/200)

For the 200 retrospective positive samples described above, a combined positive agreement of 94.1% was observed between BioPlex<sup>™</sup>2200 Vasculitis anti-MPO/anti-PR3 and ANCA IFA results, compared to a combined positive agreement of 88.2% between anti-MPO/anti-PR3 EIA and ANCA IFA results. Also, a combined overall agreement with ANCA IFA results of 88.5% and 83.0% was observed between BioPlex<sup>™</sup>2200 and EIA results, respectively. Overall, the BioPlex<sup>™</sup>2200 Vasculitis kit demonstrated better concordance with ANCA IFA than the corresponding EIA methods.

## Reproducibility

Positive Control	Sample (N)	Grand Mean AI	Within-Run		Between-Run		Total	
			SD	%CV	SD	%CV	SD	%CV
Anti-MPO	36	2.9	0.198	6.8%	0.000	0.0%	0.223	7.7%
Anti-PR3	36	2.3	0.134	5.8%	0.000	0.0%	0.167	7.3%
Anti-GBM	36	2.8	0.139	4.9%	0.000	0.0%	0.153	5.4%

Following CLSI (formerly NCCLS) EP5-A2 guidelines and ISO/TR 22971:2005, two (2) US testing sites and an internal site performed reproducibility testing using the BioPlex<sup>™</sup>2200 Vasculitis positive control for 3 days with a total of 36 replicates. The overall performance demonstrated **excellent reproducibility** ranging from 0.0% – 7.7%.

The low %CV's on the BioPlex<sup>™</sup>2200 Vasculitis kit can be attributed to:

- Full automation
- Magnetic bead, heterogeneous assay format
- A minimum of 150 measurements per analyte per reaction vessel (up to 3 times more than other multiplex methods)