

BioPlex® 2200 Vasculitis Kit

A key component of the diagnostic picture

Used in conjunction with inflammatory markers, chemistry panels and chest image analysis, autoantibody results are one of the key diagnostic parameters and less invasive laboratory tests for diagnosing some of the most severe autoimmune systemic vasculitis diseases affecting blood vessels in the lungs and kidneys such as:

- Wegener's Granulomatosis
- Churg-Strauss Syndrome
- Microscopic Polyangiitis (MPA)
- Necrotising Glomerulonephritis
- Goodpasture's Syndrome

Consolidated into one test panel

A multiplex flow immunoassay for the semi-quantitative detection in human serum of IgG autoantibodies to:

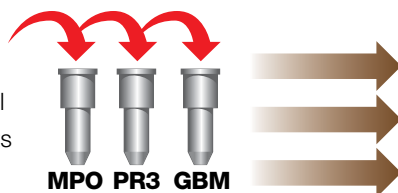
- **MPO** – Myeloperoxidase
- **PR3** – Proteinase 3
- **GBM** – Glomerular Basement Membrane

BIO-RAD

BioPlex® 2200 Vasculitis Kit

Traditional Testing:

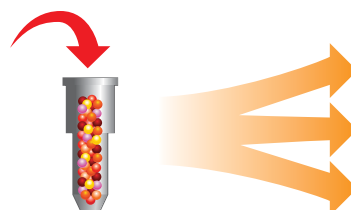
Physician orders 3 Traditional Single-Analyte Vasculitis Tests



3 Reagent equivalents consumed to produce 3 test results

Multiplex Testing:

Physician orders 3 Traditional Single-Analyte Vasculitis Tests



1 Reagent equivalent consumed to produce 3 test results

Efficiency

- Fully-automated, random access, multiplex assay procedure for vasculitis testing from primary tube to final result
- Consolidates 3 traditional single-analyte tests into 1 disease-state panel
- Up to 300 results per hour and up to 3 different reportable test results from 1 small patient sample volume (5 µL)
- Time to first result in approximately 45 minutes
- STAT capability from a single reaction vessel
- Ready-to-use multiplexed calibrators and controls
- Curve stability up to 30 days

Quality

- Excellent negative agreement with commercially available EIA methods resulting in 100%, 100% and 99.3% for anti-MPO, anti-PR3 and anti-GBM, respectively.*
- Excellent positive agreement with ANCA reactivities by commercially available IFA methods resulting in 93.3% and 94.9% for anti-MPO and anti-PR3, respectively.*
- Excellent lot-to-lot reproducibility with %CV ranging from 0.0 – 11.5% for anti-MPO, 0.0 – 18.8% for anti-PR3 and 0.0 – 9.4% for anti-GBM analytes
- Heterogeneous wash format employing novel magnetic bead technology improves assay sensitivity and specificity

* Please refer to Instructions for Use (IFU)

Ordering Information

Catalog No.	Description	
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



**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 (1-800-224-6723) **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