

Bio-Plex Pro Human SARS-CoV-2 N/RBD/S1/S2 4-Plex Panel

High-Performance Multiplex SARS-CoV-2 Serology Assays for Antibody Response Profiling

MAGNETIC SEPARATION ENABLED

IgG antibodies against four SARS-CoV-2 proteins:

- Nucleocapsid
- Receptor binding domain
- Spike 1
- Spike 2

- Acute Phase Response
- Cancer
- Cardiovascular Disease
- Cytokines, Chemokines, and Growth Factors
- Neurology
- Toxicology
- Infectious Disease
- Immunoglobulin Isotyping
- **Serology Assays**
- Signal Transduction



Researchers are interested in measuring antibodies specific to SARS-CoV-2 to determine the exposure rates to the virus in a population, the humoral response to COVID-19 vaccines, and the immune response over time of patients with COVID-19 who have different levels of disease severity. The novelty of symptoms related to COVID-19 inspires a need to understand how long the humoral and cellular adaptive immune response lasts.

New Challenges, New Solution

These assays provide an efficient way to simultaneously profile isotype-specific antibody responses to four SARS-CoV-2 proteins: nucleocapsid, receptor binding domain, spike 1, and spike 2. These qualitative multiplexed immunoassays have been developed and validated to ensure high precision, specificity, and sensitivity.

A negative and a positive control are provided. Sample interpretation guidelines offer median fluorescence intensity (MFI) cutoff values that indicate which samples are positive for antibodies to a specific SARS-CoV-2 protein. The cutoff values provided were based on 282 SARS-CoV-2-negative samples that were collected before December 2015. An isotype-specific positive control with recombinant antibodies specific to the SARS-CoV-2 proteins is included and can be used to compare MFI values across plates.

Assay Features

- Reduced turnaround time
- Reproducible results
- Specific and sensitive for SARS-CoV-2
- Flexible assay configurations
- Antigen-coupled beads and reagents for multispecies studies

Rigorous Assay Validation

- Specificity (cross-reactivity)
- Inter- and intra-assay precision
- Correlation with COVID-19 and healthy serum and plasma samples

For research use only.
Not for use in diagnostic procedures.

Representative assay characteristics.

Parameter	Characteristic
Reactive species	Human
Analytical specificity, % analyte cross-reactivity	<5%
Intra-assay precision, %CV	<10%
Inter-assay precision, %CV	<15%
Clinical specificity* (n = 282)	N = 99%, RBD = 96%, S1 = 99%, S2 = 95%
Clinical sensitivity** (n = 65)	N = 100%, RBD = 100%, S1 = 98%, S2 = 100%
Compatible sample matrices	Serum, plasma

* Clinical specificity was determined by testing 282 SARS-CoV-2–negative samples that were collected earlier than December 2015.

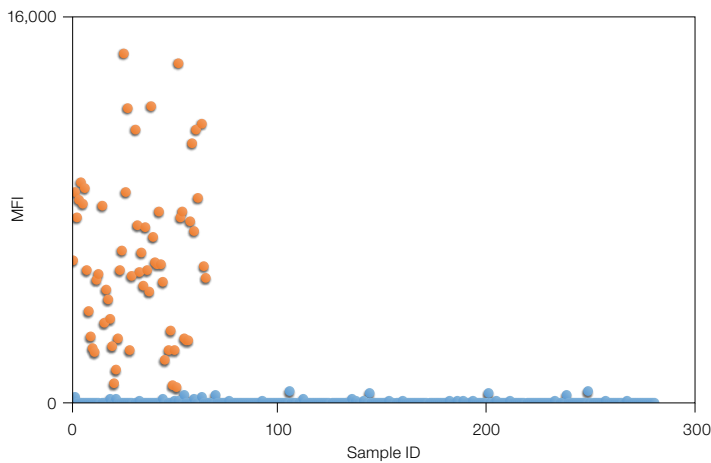
** Clinical sensitivity was determined by running 65 SARS-CoV-2–positive serum and plasma samples that were confirmed to be human IgG anti-SARS-CoV-2 positive.

CV, coefficient of variation; n, number of samples run.

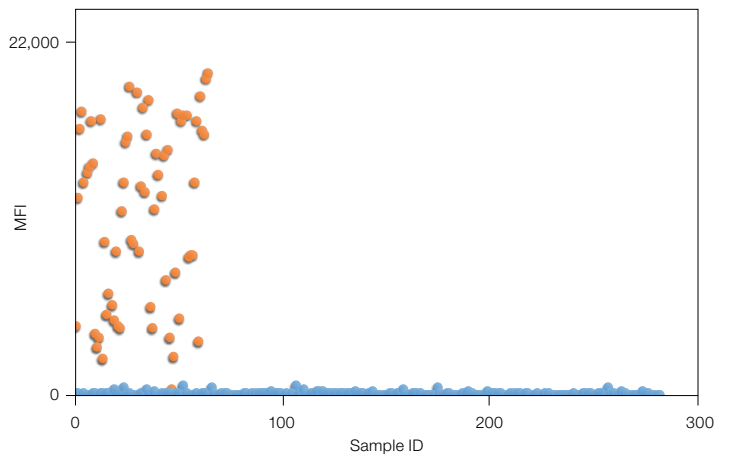
Cross-reactivity results.

Antibody Positive Sera	Number of Samples	Positive	Negative
CMV	3	0	3
EBV	1	0	1
Hepatitis B	1	0	1
Hepatitis C	1	0	1
HIV	1	0	1
HSV	1	0	1
HTLV	1	0	1
Influenza A	2	0	2
Influenza B	2	0	2
RSV	2	0	2
TOXO	1	0	1
WNV	1	0	1

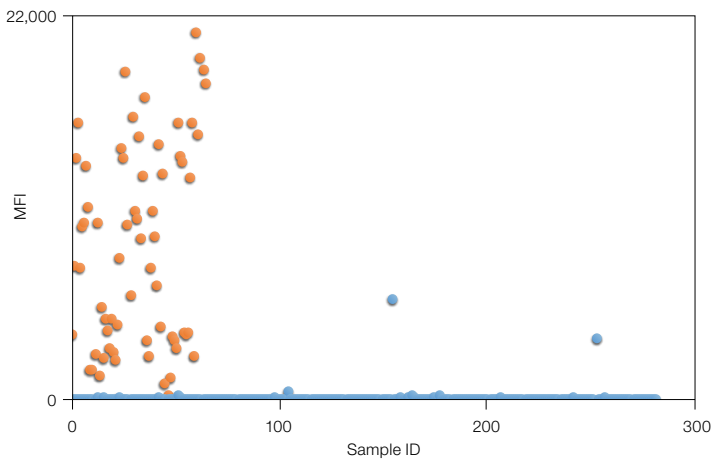
Nucleocapsid



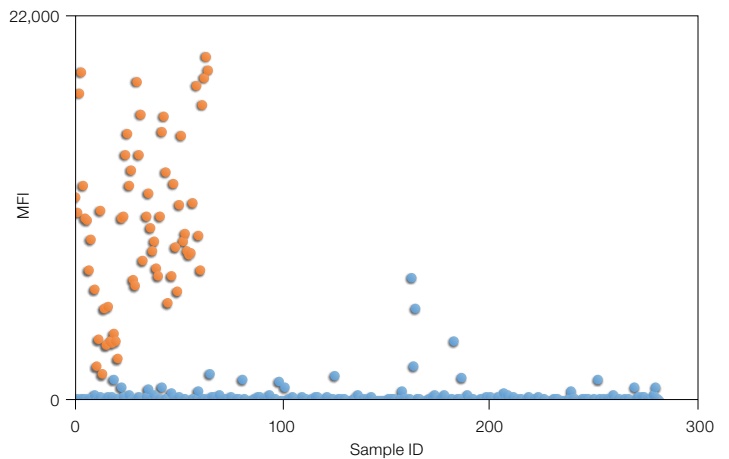
Receptor binding domain



Spike 1



Spike 2



Antibody profile of healthy and diseased samples in COVID-19 infection. The diseased samples (n = 65) were obtained from donors who were SARS-CoV-2 positive (confirmed by PCR testing). The healthy samples (n = 282) were collected from healthy donors prior to December 2015. Both serum and plasma samples were included in the study. Healthy (●); diseased (●).

Ordering Information

Catalog # Description

Multiplex Assays

12014634 **Bio-Plex Pro Human IgG SARS-CoV-2 N/RBD/S1/S2
4-Plex Panel, 1 x 96-well**

Singleplex Items

12014667 **Bio-Plex Pro SARS-CoV-2 N/RBD/S1/S2 4-Plex
Coupled Beads**
12014773 **Bio-Plex Pro SARS-CoV-2 Nucleocapsid Coupled Beads**
12015406 **Bio-Plex Pro SARS-CoV-2 Receptor Binding Domain
Coupled Beads**
12014771 **Bio-Plex Pro SARS-CoV-2 Spike 1 Coupled Beads**
12014772 **Bio-Plex Pro SARS-CoV-2 Spike 2 Coupled Beads**
12014668 **Bio-Plex Pro Human IgG Detection Antibody**
12014774 **Bio-Plex Pro Human IgG SARS-CoV-2 Positive and
Negative Controls**
12014777 **Bio-Plex Pro Human Serology Reagent Kit**

Visit [bio-rad.com/SARS-CoV-2Serology](https://www.bio-rad.com/SARS-CoV-2Serology) for more information.

BIO-RAD and BIO-PLEX are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23
France 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050
Italy 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670
The Netherlands 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23
Russian Federation 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23
Switzerland 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

