

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

Bio-Plex Pro Human Cytokine 48-Plex Screening Panel

Basic FGF, CTACK, eotaxin, G-CSF, GM-CSF, GRO- α , HGF, ICAM-1,* IFN- α 2, IFN- γ , IL-1 α , IL-1 β , IL-1ra, IL-2, IL-2R α , IL-3, IL-4, IL-5, IL-6, IL-7, IL-8, IL-9, IL-10, IL-12 (p40), IL-12 (p70), IL-13, IL-15, IL-16, IL-17A, IL-18, IP-10, LIF, MCP-1 (MCAF), MCP-3, M-CSF, MIF, MIG, MIP-1 α , MIP-1 β , β -NGF, PDGF-BB, RANTES, SCF, SCGF- β , SDF-1 α , TNF- α , TNF- β , TRAIL, VCAM-1,* VEGF-A

* Recommended serum sample dilution: 100-fold. Cannot be multiplexed with analytes in the 48-plex assay.

MAGNETIC SEPARATION ENABLED

- All-in-one 48-plex panel
- Single-level quality control
- Magnetic workflow



High-Performance Multiplex Screening Immunoassays for Research and Drug Discovery

Cytokines and chemokines are extracellular mediators and regulators within a signaling network between cells and are key modulators of inflammation, participating in acute and chronic conditions via a complex network of interactions. This panel integrates a network of biologically relevant cytokines and chemokines in a single assay, enabling you to interrogate 48 biomarkers simultaneously. This tool provides a comprehensive, robust solution for screening 48 biologically relevant analytes in a single well, accelerating breakthrough research and biomarker discovery. All assays in this panel have gone through a vigorous validation process to ensure the data are accurate and reproducible.

Assay Features

- Comprehensive panel coverage
- Broad dynamic range
- Best-in-class performance
- Reduced turnaround time
- Reproducible results
- Streamlined data analysis
- Flexibility

Rigorous Assay Validation

All Bio-Plex Pro Assays undergo rigorous evaluation that includes the following parameters:

- Specificity (cross-reactivity)
- Accuracy (recovery) in key sample matrices
- Inter- and intra-assay precision
- Sensitivity (limit of detection [LOD])
- Assay working range (lower and upper limits of quantification [LLOQ/ULOQ])
- Linearity of dilution
- Parallelism and matrix effect
- Performance characteristics in real samples (Figures 1 and 2)

BIO-RAD

Assay Performance Definitions

The following parameters are indicative of assay performance, as shown in Table 1.

- Assay working range** — the range of concentrations within which the assay is precise and accurate. Boundaries of the assay working range are defined by the LLOQ and ULOQ
- Precision** — the percentage coefficient of variation (%CV) at concentrations within the assay working range
- Accuracy (recovery)** — percentage of the observed concentration relative to the expected concentration of a known amount of analyte within the assay working range
- Sensitivity (LOD)** — the concentration of analyte for which the fluorescence intensity signal is 2 standard deviations above the background signal
- Linearity of dilution** — determines to what extent the dosage response above LLOQ can be diluted and measured accurately within the assay working range (Table 2)
- Parallelism** — determines if a diluent matrix used to construct a standard curve is biologically comparable to the diluent matrix used for sample preparation. This ensures that the recombinant calibrator shares the same binding characteristic to the antibodies as the native molecules (Table 3 and Figure 3)

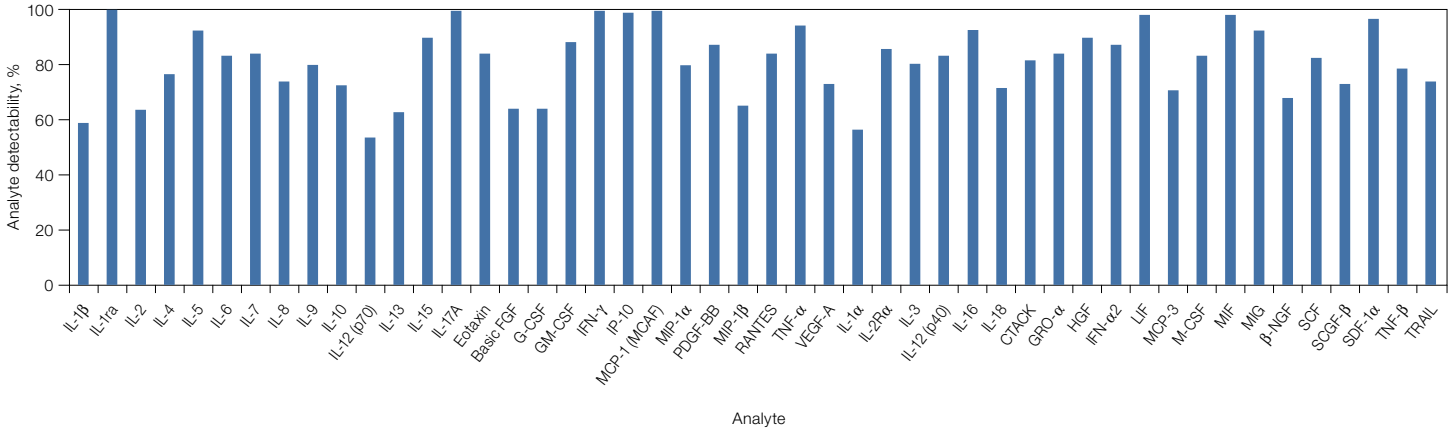


Fig. 1. Analyte detectability. Four hundred and twenty test samples from six laboratories, including culture media, cancer serum, stimulated peripheral blood mononuclear cells, whole blood, and human immunodeficiency virus plasma, were run on the assay. The average analyte detectability for the panel is >80%. Analyte detectability is defined as the percentage of analytes detected within the working assay range for all samples in a study.

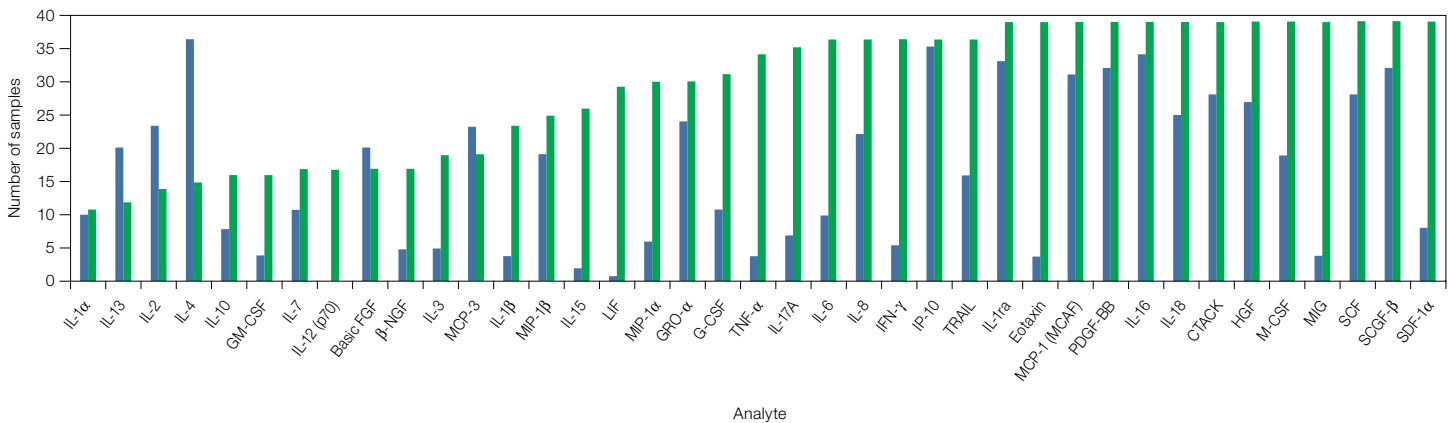


Fig. 2. Competitive comparison. Thirty-nine test samples were run on Vendor R's 39-plex and compared to the same overlapping 39 analytes in the Bio-Plex Pro Human Cytokine 48-Plex Screening Panel. This test configuration generated 1,521 total possible data points (39 samples x 39 analytes). In total, Vendor R's assay detected 637 data points within their working assay range while the 39 analytes from the Bio-Plex Pro Panel detected 1,140 data points. Detected data points divided by total possible data points defines analyte detectability in this comparison. Analyte detectability for Vendor R (■) is 42% while analyte detectability for the Bio-Plex Pro Panel (■) is 75%. Bio-Rad offers 79% improved analyte detectability over Vendor R in the context of these 39 overlapping targets.

Table 1. Representative performance characteristics.

Analyte	Assay Working Range, pg/ml		Assay Sensitivity, pg/ml	Mean Intra-Assay %CV*	Mean Inter-Assay %CV*	Singleplex Bead Region
	LLOQ*	ULOQ*	LOD*			
Basic FGF	3.26	3,341	2.54	3.1	2.4	44
CTACK	2.10	15,656	0.82	2.7	5.2	72
Eotaxin	0.14	2,281	0.05	4.4	1.2	43
G-CSF	6.35	104,106	3.63	3.1	4.0	57
GM-CSF	0.48	7,846	0.19	4.3	2.2	34
GRO- α	21.05	31,255	13.45	2.6	7.9	61
HGF	8.76	143,513	7.09	2.6	2.9	62
ICAM-1	3.84	62,935	1.51	4.4	2.3	12
IFN- α 2	0.95	15,569	0.46	3.3	4.4	20
IFN- γ	1.57	25,665	1.05	3.1	3.6	21
IL-1 α	3.73	61,154	6.65	3.5	4.9	63
IL-1 β	0.29	4,672	0.24	3.6	3.2	39
IL-1ra	6.21	34,949	3.16	4.7	5.1	25
IL-2	1.29	21,178	0.75	1.7	2.5	38
IL-2R α	1.48	24,270	1.65	3.4	4.8	13
IL-3	0.13	2,139	0.13	5.0	3.9	64
IL-4	0.19	3,064	0.09	3.2	1.9	52
IL-5	3.63	59,499	0.86	2.3	2.3	33
IL-6	0.38	6,244	0.34	2.2	3.0	19
IL-7	1.92	31,475	1.22	2.7	3.9	74
IL-8	0.85	13,992	0.36	3.2	2.8	54
IL-9	3.62	31,527	1.08	2.6	7.1	77
IL-10	1.06	17,427	0.69	2.3	3.4	56
IL-12 (p40)	14.68	240,582	6.39	4.5	2.4	28
IL-12 (p70)	1.43	23,425	0.78	3.3	2.9	75
IL-13	0.31	5,157	0.22	3.1	2.7	51
IL-15	12.42	203,426	12.82	2.8	4.1	73
IL-16	1.20	19,639	0.25	2.5	3.0	27
IL-17A	2.44	39,972	1.16	2.4	1.4	76
IL-18	0.66	10,892	0.31	2.9	2.2	42
IP-10	3.41	34,953	1.43	2.8	6.0	48
LIF	3.86	53,806	2.05	2.5	4.7	29
MCP-1 (MCAF)	0.53	8,755	0.44	3.2	3.4	53
MCP-3	0.48	4,899	0.24	4.4	4.2	26
M-CSF	0.75	12,290	0.27	2.4	3.6	67
MIF	2.70	44,168	2.45	3.4	4.7	35
MIG	3.16	32,365	1.39	4.4	4.2	14
MIP-1 α	0.12	1,218	0.06	4.5	4.2	55
MIP-1 β	1.41	1,439	1.41	3.4	2.5	18
β -NGF	0.47	7,655	0.23	2.9	3.9	46
PDGF-BB	7.12	37,133	2.96	3.3	9.7	47
RANTES	16.72	26,467	3.98	3.0	6.7	37
SCF	1.82	29,899	0.99	4.1	2.6	65
SCGF- β	82.11	1,345,200	141.77	2.3	3.8	78
SDF-1 α	7.54	9,381	2.44	2.2	5.4	22
TNF- α	3.33	54,566	1.13	3.5	3.0	36
TNF- β	0.80	13,186	0.38	3.0	4.7	30
TRAIL	1.78	29,188	0.89	3.2	4.5	66
VCAM-1	2.91	47,652	2.05	4.1	3.1	15
VEGF-A	18.01	149,830	10.16	2.8	8.5	45

* The LLOQ, ULOQ, LOD, and inter-assay precision %CV are mean data determined from three independent multiplex assays in a serum-based matrix. Intra-assay %CV is derived from one representative assay. LLOQ and ULOQ are defined as the boundary standard curve points within which the performance specifications of individual standard points were met for a 10% intra-assay CV and recovery range of 70–130%. Data were generated using the magnetic workflow with the Bio-Plex Pro Wash Station.

Table 2. Linearity of dilution.*

Analyte	Linear Dilution Range	Analyte	Linear Dilution Range
Basic FGF	1:9–1:243	IL-13	1–1:2,187
CTACK	1–1:27**	IL-15	1:3–1:81
Eotaxin	1–1:243	IL-16	1–1:243
G-CSF	1–1:243	IL-17A	1–1:243
GM-CSF	1–1:2,187	IL-18	1–1:9
GRO- α	1:27–1:2,187	IP-10	1–1:81
HGF	1–1:81	LIF	1–1:729
IFN- α 2	1:3–1:2,187	MCP-1 (MCAF)	1–1:81
IFN- γ	1–1:2,187	MCP-3	1–1:2,187
IL-1 α	1–1:729	M-CSF	1–1:243
IL-1 β	1–1:2,187	MIF	1:3–1:81
IL-1ra	1:3–1:243	MIG	1–1:81
IL-2	1–1:2,187	MIP-1 α	1:3–1:243
IL-2R α	1–1:81	MIP-1 β	1:3–1:242
IL-3	1–1:729	β -NGF	1–1:729
IL-4	1–1:2,187	PDGF-BB	1:3–1:81
IL-5	1–1:729	RANTES	1–1:9
IL-6	1–1:81	SCF	1–1:243
IL-7	1:9–1:729	SCGF- β	***
IL-8	1–1:243	SDF-1 α	***
IL-9	1:9–1:729	TNF- α	1–1:729
IL-10	1–1:729	TNF- β	1–1:2,187
IL-12 (p40)	1–1:2,187	TRAIL	1–1:243
IL-12 (p70)	1–1:2,187	VEGF-A	1:3–1:243

* Linearity of dilution determines whether the dosage response of the analyte is linear within the working assay range. A 50-plex recombinant standard was spiked into a serum sample which was previously diluted 1:4 with standard diluent HB then serially diluted seven times. Table 2 shows the sample dilution range where the dosage response is linear. The range is defined as the recovery of calculated concentrations within 70–130%.

** Recovery was 69% for CTACK at a 1:9 dilution.

*** Both SCGF- β and SDF-1 α had high endogenous levels in the serum used in this study and were excluded.

Table 3. Parallelism.*

Analyte	%CV
CTACK	21
HGF	9
IL-16	7
IL-18	15
IP-10	12
MCP-1 (MCAF)	13
M-CSF	21
MIF	16
MIG	9
PDGF-BB	22
RANTES	19
SCF	19
SCGF- β	15
TRAIL	11

* Parallelism “is to ascertain that the binding characteristic of the endogenous analyte to the antibodies is the same as for the calibrator” (Andreasson et al. 2015). One human serum sample with high endogenous analytes was serially diluted (1:2). Table 3 shows the %CV of quantitation results from 25% serum samples and the subsequent dilutions.

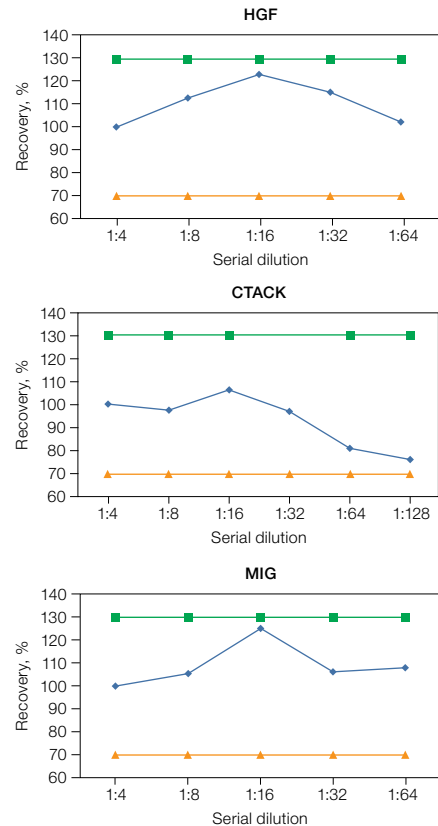


Fig. 3. Parallelism. Representative percentage recovery data from Table 3. Upper limit (■); human serum sample (♦); lower limit (▲).

Ordering Information

Catalog #	Description
12007283	Bio-Plex Pro Human Cytokine 48-Plex Screening Panel , 1 x 96-well, includes coupled magnetic capture beads, premixed detection antibodies, standards, quality control, detection antibody diluent HB, standard diluent HB, sample diluent HB, assay buffer, 10x wash buffer, streptavidin-PE, 96-well flat bottom plate, sealing tape, and instructions, for the detection of basic FGF, CTACK, eotaxin, G-CSF, GM-CSF, GRO- α , HGF, IFN- α 2, IFN- γ , IL-1 α , IL-1 β , IL-1ra, IL-2, IL-2R α , IL-3, IL-4, IL-5, IL-6, IL-7, IL-8, IL-9, IL-10, IL-12 (p40), IL-12 (p70), IL-13, IL-15, IL-16, IL-17A, IL-18, IP-10, LIF, MCP-1 (MCAF), MCP-3, M-CSF, MIF, MIG, MIP-1 α , MIP-1 β , β -NGF, PDGF-BB, RANTES, SCF, SCGF- β , SDF-1 α , TNF- α , TNF- β , TRAIL, VEGF-A

Bio-Plex Pro Human Cytokine Singleplex Assays

171B5016M	Basic FGF Set	171B5012M	IL-13 Set
171B6006M	CTACK Set	171B5013M	IL-15 Set
171B5015M	Eotaxin Set	171B6005M	IL-16 Set
171B5017M	G-CSF Set	171B5014M	IL-17A Set
171B5018M	GM-CSF Set	171B5020M	IP-10 Set
171B6007M	GRO-α Set	171B6011M	LIF Set
171B6008M	HGF Set	171B5021M	MCP-1 (MCAF) Set
171B6009M	ICAM-1 Set*	171B6012M	MCP-3 Set
171B6010M	IFN-α2 Set	171B6013M	M-CSF Set
171B5019M	IFN-γ Set	171B6014M	MIF Set
171B6001M	IL-1α Set	171B6015M	MIG Set
171B5001M	IL-1β Set	171B5022M	MIP-1α Set
171B5002M	IL-1ra Set	171B5023M	MIP-1β Set
171B5003M	IL-2 Set	171B6016M	β-NGF Set
171B6002M	IL-2Rα Set	171B5024M	PDGF-BB Set
171B6003M	IL-3 Set	171B5025M	RANTES Set
171B5004M	IL-4 Set	171B6017M	SCF Set
171B5005M	IL-5 Set	171B6018M	SCGF-β Set
171B5006M	IL-6 Set	171B6019M	SDF-1α Set
171B5007M	IL-7 Set	171B5026M	TNF-α Set
171B5008M	IL-8 Set	171B6020M	TNF-β Set
171B5009M	IL-9 Set	171B6021M	TRAIL Set
171B5010M	IL-10 Set	171B6022M	VCAM-1 Set*
171B6004M	IL-12 (p40) Set	171B5027M	VEGF-A Set
171B5011M	IL-12 (p70) Set		

* Recommended serum sample dilution: 100-fold. Cannot be multiplexed with analytes in the 48-plex assay.

Reagent Kits

171304090M	Bio-Plex Pro Reagent Kit III , 1 x 96-well, includes detection antibody diluent HB, standard diluent HB, sample diluent HB, assay buffer, 10x wash buffer, streptavidin-PE, flat bottom plate, and sealing tape
12005846	Bio-Plex Pro Reagent Kit III , 10 x 96-well, includes detection antibody diluent HB, standard diluent HB, sample diluent HB, assay buffer, 10X wash buffer, streptavidin-PE, flat bottom plate, and sealing tape

Standards

12007919	Bio-Plex Pro Human Cytokine Screening Panel Standards , 1 vial, lyophilized mixture of 50 standard analytes, includes basic FGF, CTACK, eotaxin, G-CSF, GM-CSF, GRO- α , HGF, ICAM-1, IFN- α 2, IFN- γ , IL-1 α , IL-1 β , IL-1ra, IL-2, IL-2R α , IL-3, IL-4, IL-5, IL-6, IL-7, IL-8, IL-9, IL-10, IL-12 (p40), IL-12 (p70), IL-13, IL-15, IL-16, IL-17A, IL-18, IP-10, LIF, MCP-1 (MCAF), MCP-3, M-CSF, MIF, MIG, MIP-1 α , MIP-1 β , β -NGF, PDGF-BB, RANTES, SCF, SCGF- β , SDF-1 α , TNF- α , TNF- β , TRAIL, VCAM-1, VEGF-A
----------	--

Wash Stations and Accessories

30034376	Bio-Plex Pro Wash Station , microplate wash station for magnetic bead-based assays, includes magnetic plate carrier, waste bottle, 2 buffer bottles
171020100	Bio-Plex Handheld Magnetic Washer , includes magnetic washer and adjustment hex tools for use in manual wash steps for all Bio-Plex Magnetic Assays
171025001	Bio-Plex Pro Flat Bottom Plates , pkg of 40 x 96-well plates, for use with Bio-Plex Pro Wash Stations when using magnetic bead-based assays

Catalog #	Description
Software 171001510	Bio-Plex Data Pro Software with Bio-Plex Manager Software , Bio-Plex Data Pro Software (5 seats) for multi-experiment analysis and advanced data visualization and Bio-Plex Manager Software (5 seats) for instrument data evaluation and optimization; CDs and security HASP key included
171001513	Bio-Plex Data Pro Software , 5 seats, for multi-experiment analysis and advanced data visualization
171STND01	Bio-Plex Manager Software , 1 user desktop license, for analysis of Bio-Plex data and generation of protocols; does not operate the instrument

Additional Recommended Bio-Plex Assays for Studying Human Cytokines

171AL001M	Bio-Plex Pro Human Inflammation 37-Plex Panel 1 , for the detection of APRIL/TNFSF13, BAFF/TNFSF13B, sCD30/TNFRSF8, sCD163, chitinase 3-like 1, gp130/sIL-6R β , IFN- α 2, IFN- β , IFN- γ , IL-2, sIL-6R α , IL-8, IL-10, IL-11, IL-12 (p40), IL-12 (p70), IL-19, IL-20, IL-22, IL-26, IL-27 (p28), IL-28A/IFN- λ 2, IL-29/IFN- λ 1, IL-32, IL-34, IL-35, LIGHT/TNFSF14, MMP-1, MMP-2, MMP-3, osteocalcin, osteopontin, pentraxin-3, sTNF-R1, sTNF-R2, TSLP, TWEAK/TNFSF12
171AK99MR2	Bio-Plex Pro Human Chemokine 40-Plex Panel , for the detection of 6CKine/CCL21, BCA-1/CXCL13, CTACK/CCL27, ENA-78/CXCL5, eotaxin/CCL11, eotaxin-2/CCL24, eotaxin-3/CCL26, fractalkine/CX3CL1, GCP-2/CXCL6, GM-CSF, GRO- α /CXCL1, GRO- β /CXCL2, I-309/CCL1, IFN- γ , IL-1 β , IL-2, IL-4, IL-6, IL-8/CXCL8, IL-10, IL-16, IP-10/CXCL10, I-TAC/CXCL11, MCP-1/CCL2, MCP-2/CCL8, MCP-3/CCL7, MCP-4/CCL13, MDC/CCL22, MIF, MIG/CXCL9, MIP-1 α /CCL3, MIP-1 δ /CCL15, MIP-3 α /CCL20, MIP-3 β /CCL19, MPIF-1/CCL23, SCYB16/CXCL16, SDF-1 α + β /CXCL12, TARC/CCL17, TECK/CCL25, TNF- α
171AA001M	Bio-Plex Pro Human Th17 Cytokine 15-Plex Panel , for the detection of IL-1 β , IL-4, IL-6, IL-10, IL-17A, IL-17F, IL-17A/F (singleplex only), IL-21, IL-22, IL-23, IL-25, IL-31, IL-33, IFN- γ , sCD40L, TNF- α
171W4001M	Bio-Plex Pro TGF-β 3-Plex Panel , for the detection of TGF- β 1, TGF- β 2, TGF- β 3

Reference

Andreasson U et al. (2015). A practical guide to immunoassay method validation. *Front Neurol* 6, 179.

Visit bio-rad.com/HCS for more information.

Bio-Plex and Bio-Rad are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions.

The Bio-Plex Suspension Array System includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation.

All trademarks used herein are the property of their respective owner.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00
France 33 01 479 593 00 **Germany** 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300
Israel 972 03 963 6050 **Italy** 39 02 49486600 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666
New Zealand 64 9 415 2280 **Norway** 47 0 233 841 30 **Poland** 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580 **Sweden** 46 08 555 127 00 **Switzerland** 41 0617 17 9555
Taiwan 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 01923 47 1301

