

Ordering Information

Instrument, Software, and Tools

Bio-Plex® System

Bio-Plex system components:

1. Array reader
2. Microplate platform
3. High-throughput fluidics (HTF)
4. Flat-screen monitor
5. PC
6. Reagent kit
7. Cell lysis kit
8. Singleplex or premixed multiplex assays
9. COOH beads
10. Validation kit
11. Maintenance, calibration, and validation (MCV) plate
12. Calibration kit
13. Bio-Plex Pro™ wash station



Catalog # Description

Bio-Plex Systems

- 171-000201* **Bio-Plex 200 System**, 100–240 V, includes array reader, microplate platform, Bio-Plex Manager™ 6.0 software with instrument control license, PC, monitor, calibration kit, validation kit 4.0, maintenance, calibration, and validation (MCV) plate IV, 20 L sheath fluid, spare sample needle
- 171-000205 **Bio-Plex 200 System with HTF**, same as 171-000201 with high-throughput fluidics (HTF)

Software**

- 171-STND01 **Bio-Plex Manager 6.0 Software**, includes 1-user desktop license
- 171-STND05 **Bio-Plex Manager 6.0 Software**, includes 5-user network license
- 171-STND10 **Bio-Plex Manager 6.0 Software**, includes 10-user network license

Software (Security Edition)**

- 171-SCRT00 **Bio-Plex Manager 6.0 Instrument Control License, Security Edition**, upgrades workstation with Bio-Plex Manager 6.0 software, allows compliance with U.S. FDA 21 CFR Part 11 regulations
- 171-SCRT01 **Bio-Plex Manager 6.0 Software, Security Edition**, includes 1-user desktop license, allows compliance with U.S. FDA 21 CFR Part 11 regulations
- 171-SCRT05 **Bio-Plex Manager 6.0 Software, Security Edition**, includes 5-user network license, allows compliance with U.S. FDA 21 CFR Part 11 regulations
- 171-SCRT10 **Bio-Plex Manager 6.0 Software, Security Edition**, includes 10-user network license, allows compliance with U.S. FDA 21 CFR Part 11 regulations

Catalog # Description

Bio-Plex Pro Wash Stations and Accessories

- 30034376 **Bio-Plex Pro Wash Station**, includes magnetic plate carrier, waste bottle, 2 buffer bottles
- 30034377 **Bio-Plex Pro II Wash Station**, includes magnetic plate carrier, vacuum manifold, plate carrier, waste bottle, 2 buffer bottles
- 171-025001*** **Bio-Plex Pro Flat Bottom Plates**, 40 x 96-well plates
- 171-304500 **Bio-Plex Wash Buffer**, 1.5 L

System Tools

- 171-203001 **Bio-Plex Validation Kit 4.0**, includes optics validation, reporter validation, classify validation, and fluidics validation reagents for approximately 50 validation routines using Bio-Plex Manager 4.0 or higher software (standard or Security Edition) and MCV plate IV
- 171-203033 **Bio-Plex MCV Plate IV**, for use with Bio-Plex Manager 4.0 or higher software (standard or Security Edition) and validation kit 4.0
- 171-203050 **Bio-Plex Reservoir**
- 171-203060 **Bio-Plex Calibration Kit**, includes Cal1 and Cal2 calibration beads for approximately 50 daily calibration routines
- 171-000055 **Sheath Fluid**, 1x, 20 L

* Not available in the U.S. and Canada.

** Contact Bio-Rad Technical Support to verify that your current system is compatible with these software packages.

*** Required for washing Bio-Plex Pro assays or other magnetic bead-based assays using the Bio-Plex Pro wash station.

BIO-RAD

Bio-Plex 3D System new



Catalog #	Description
Bio-Plex 3D	Bio-Plex 3D System , includes Bio-Plex 3D suspension array system, xPONENT 4.0 acquisition software, PC, calibration reagents, verification reagents, Bio-Plex Manager 6.0 desktop license
89-20185-00-001*	xPONENT CFR21 Part 11 Software Module , software module to enable CFR21 Part II compliance
89-20187-00-001*	xPONENT Automation Module , software module to enable interfacing with robotics workstations
171-022001*	Swivel Base , swivel mounting to allow rotation of instrument to face robotics workstation and to face forward for maintenance and manual loading
89-20186-00-001*	xPONENT LIS Software Module , software module to enable interfacing with LIS (Laboratory Information Systems) databases
89-20182-00-001*	xPONENT Extra Seat Licenses (3 seats) , 3 additional seats of xPONENT software. Maximum of 15 seats
171-000055	Sheath Fluid , 20 L, 1x bulk fluid in cube container with pour spout, for use with the Bio-Plex 200 and Bio-Plex 3D systems' array reader and high-throughput fluidics (HTF) system

* Orders for these accessories can be placed only at time of Bio-Plex 3D system purchase.

Bio-Plex Pro™ Magnetic Cytokine, Chemokine, and Growth Factor Assays

Bio-Plex Pro cytokine, chemokine, and growth factor assays can be ordered as follows:

Premixed panels — Five human and four mouse premixed assay panels are available in the new all-in-one kit packaging format that includes coupled magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, standard diluent, and sample diluent.

x-Plex assay — Custom premixed assay panels in the new all-in-one kit packaging format can be ordered with any combination of available assays. Two-plate minimum applies if ordering a 6-plex or less. Go to www.bio-rad.com/bio-plex/x-plex/ for details.

Express assay — For faster delivery, customized assay panels in the new all-in-one kit packaging format can be ordered with any combination of available assays up to a 10-plex maximum. Coupled magnetic beads and detection antibodies are not premixed. Go to www.bio-rad.com/bio-plex/x-plex/ for details.

Individual components — Singleplex sets (coupled magnetic beads and detection antibodies), standards, reagent kits (includes diluents), wash buffer, and Bio-Plex Pro flat bottom plates for magnetic bead-based washing are available separately.

For a list of publications using Bio-Plex cytokine assays, request bulletin 5297. For more information, request bulletin 5405.

What to Order



Bio-Plex Pro cytokine, chemokine, and growth factor assays — magnetic bead-based assays built around MagPlex magnetic beads, for the measurement of human and mouse cytokine, chemokine, and growth factor biomarkers. Magnetic beads afford compatibility with Bio-Plex Pro and Bio-Plex Pro II wash stations, providing decreased variability and increased consistency between experiments.

Catalog #	Description
Bio-Plex Pro Premixed Human Cytokine Panels	
M50-000007A	Bio-Plex Pro Human Cytokine 8-Plex Panel , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, standard diluent, sample diluent for the detection of IL-2, IL-4, IL-6, IL-8, IL-10, GM-CSF, IFN- γ , TNF- α
M50-00031YV	Bio-Plex Pro Human Cytokine 17-Plex Panel , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, standard diluent, sample diluent for the detection of IL-1 β , IL-2, IL-4, IL-5, IL-6, IL-7, IL-8, IL-10, IL-12 (p70), IL-13, IL-17, G-CSF, GM-CSF, IFN- γ , MCP-1 (MCAF), MIP-1 β , TNF- α
MF0-005KMII	Bio-Plex Pro Human Cytokine 21-Plex Panel* , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, standard diluent, sample diluent for the detection of IL-1 α , IL-2R α , IL-3, IL-12 (p40), IL-16, IL-18, CTACK, GRO- α , HGF, IFN- α 2, LIF, MCP-3, M-CSF, MIF, MIG, β -NGF, SCF, SCGF- β , SDF-1 α , TNF- β , TRAIL
M50-0KCAF0Y	Bio-Plex Pro Human Cytokine 27-Plex Panel , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, standard diluent, sample diluent for the detection of IL-1 β , IL-1ra, IL-2, IL-4, IL-5, IL-6, IL-7, IL-8, IL-9, IL-10, IL-12 (p70), IL-13, IL-15, IL-17, basic FGF, eotaxin, G-CSF, GM-CSF, IFN- γ , IP-10, MCP-1 (MCAF), MIP-1 α , MIP-1 β , PDGF-BB, RANTES, TNF- α , VEGF
M50-00005L3	Bio-Plex Pro Human Cytokine Th1/Th2 Panel , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, standard diluent, sample diluent for the detection of IL-2, IL-4, IL-5, IL-10, IL-12 (p70), IL-13, GM-CSF, IFN- γ , TNF- α
Bio-Plex Pro Premixed Mouse Cytokine Panels	
M60-000007A	Bio-Plex Pro Mouse Cytokine 8-Plex Panel , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, standard diluent, sample diluent for the detection of IL-1 β , IL-2, IL-4, IL-5, IL-10, GM-CSF, IFN- γ , TNF- α

Catalog #	Description
MD0-00000EL	Bio-Plex Pro Mouse Cytokine 9-Plex Panel , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, standard diluent, sample diluent for the detection of IL-15, IL-18, basic FGF, LIF, M-CSF, MIG, MIP-2, PDGF-BB, VEGF
M60-009RDPD	Bio-Plex Pro Mouse Cytokine 23-Plex Panel , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, standard diluent, sample diluent for the detection of IL-1 α , IL-1 β , IL-2, IL-3, IL-4, IL-5, IL-6, IL-9, IL-10, IL-12 (p40), IL-12 (p70), IL-13, IL-17, eotaxin, G-CSF, GM-CSF, IFN- γ , KC, MCP-1 (MCAF), MIP-1 α , MIP-1 β , RANTES, TNF- α
M60-00003J7	Bio-Plex Pro Mouse Cytokine Th1/Th2 Panel , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, standard diluent, sample diluent for the detection of IL-2, IL-4, IL-5, IL-10, IL-12 (p70), GM-CSF, IFN- γ , TNF- α

* Human ICAM-1 and VCAM-1 require different sample dilutions. It is recommended to order these targets separately or as a 2-plex panel.

Catalog #	Description
Bio-Plex Pro Human Cytokine Singleplex Sets, Group I*	
171-B5001M	Bio-Plex Pro Human Cytokine IL-1 β Set, 1 x 96-well
171-B5002M	Bio-Plex Pro Human Cytokine IL-1ra Set, 1 x 96-well
171-B5003M	Bio-Plex Pro Human Cytokine IL-2 Set, 1 x 96-well
171-B5004M	Bio-Plex Pro Human Cytokine IL-4 Set, 1 x 96-well
171-B5005M	Bio-Plex Pro Human Cytokine IL-5 Set, 1 x 96-well
171-B5006M	Bio-Plex Pro Human Cytokine IL-6 Set, 1 x 96-well
171-B5007M	Bio-Plex Pro Human Cytokine IL-7 Set, 1 x 96-well
171-B5008M	Bio-Plex Pro Human Cytokine IL-8 Set, 1 x 96-well
171-B5009M	Bio-Plex Pro Human Cytokine IL-9 Set, 1 x 96-well
171-B5010M	Bio-Plex Pro Human Cytokine IL-10 Set, 1 x 96-well
171-B5011M	Bio-Plex Pro Human Cytokine IL-12 (p70) Set, 1 x 96-well
171-B5012M	Bio-Plex Pro Human Cytokine IL-13 Set, 1 x 96-well
171-B5013M	Bio-Plex Pro Human Cytokine IL-15 Set, 1 x 96-well
171-B5014M	Bio-Plex Pro Human Cytokine IL-17 Set, 1 x 96-well
171-B5015M	Bio-Plex Pro Human Cytokine Eotaxin Set, 1 x 96-well
171-B5016M	Bio-Plex Pro Human Cytokine Basic FGF Set, 1 x 96-well
171-B5017M	Bio-Plex Pro Human Cytokine G-CSF Set, 1 x 96-well
171-B5018M	Bio-Plex Pro Human Cytokine GM-CSF Set, 1 x 96-well
171-B5019M	Bio-Plex Pro Human Cytokine IFN- γ Set, 1 x 96-well
171-B5020M	Bio-Plex Pro Human Cytokine IP-10 Set, 1 x 96-well
171-B5021M	Bio-Plex Pro Human Cytokine MCP-1 (MCAF) Set, 1 x 96-well
171-B5022M	Bio-Plex Pro Human Cytokine MIP-1 α Set, 1 x 96-well
171-B5023M	Bio-Plex Pro Human Cytokine MIP-1 β Set, 1 x 96-well
171-B5024M	Bio-Plex Pro Human Cytokine PDGF-BB Set, 1 x 96-well
171-B5025M	Bio-Plex Pro Human Cytokine RANTES Set, 1 x 96-well
171-B5026M	Bio-Plex Pro Human Cytokine TNF- α Set, 1 x 96-well
171-B5027M	Bio-Plex Pro Human Cytokine VEGF Set, 1 x 96-well
Bio-Plex Pro Human Cytokine Singleplex Sets, Group II*	
171-B6001M	Bio-Plex Pro Human Cytokine IL-1 α Set, 1 x 96-well
171-B6002M	Bio-Plex Pro Human Cytokine IL-2R α Set, 1 x 96-well
171-B6003M	Bio-Plex Pro Human Cytokine IL-3 Set, 1 x 96-well
171-B6004M	Bio-Plex Pro Human Cytokine IL-12 (p40) Set, 1 x 96-well
171-B6005M	Bio-Plex Pro Human Cytokine IL-16 Set, 1 x 96-well
171-B6006M	Bio-Plex Pro Human Cytokine CTACK Set, 1 x 96-well
171-B6007M	Bio-Plex Pro Human Cytokine GRO- α Set, 1 x 96-well
171-B6008M	Bio-Plex Pro Human Cytokine HGF Set, 1 x 96-well
171-B6009M**	Bio-Plex Pro Human Cytokine ICAM-1 Set, 1 x 96-well
171-B6010M	Bio-Plex Pro Human Cytokine IFN- α 2 Set, 1 x 96-well
171-B6011M	Bio-Plex Pro Human Cytokine LIF Set, 1 x 96-well
171-B6012M	Bio-Plex Pro Human Cytokine MCP-3 Set, 1 x 96-well
171-B6013M	Bio-Plex Pro Human Cytokine M-CSF Set, 1 x 96-well
171-B6014M	Bio-Plex Pro Human Cytokine MIF Set, 1 x 96-well
171-B6015M	Bio-Plex Pro Human Cytokine MIG Set, 1 x 96-well
171-B6016M	Bio-Plex Pro Human Cytokine β -NGF Set, 1 x 96-well
171-B6017M	Bio-Plex Pro Human Cytokine SCF Set, 1 x 96-well
171-B6018M	Bio-Plex Pro Human Cytokine SCGF- β Set, 1 x 96-well
171-B6019M	Bio-Plex Pro Human Cytokine SDF-1 α Set, 1 x 96-well
171-B6020M	Bio-Plex Pro Human Cytokine TNF- β Set, 1 x 96-well
171-B6021M	Bio-Plex Pro Human Cytokine TRAIL Set, 1 x 96-well
171-B6022M**	Bio-Plex Pro Human Cytokine VCAM-1 Set, 1 x 96-well

Catalog #	Description
Bio-Plex Pro Mouse Cytokine Singleplex Sets, Group I*	
171-G5001M	Bio-Plex Pro Mouse Cytokine IL-1 α Set, 1 x 96-well
171-G5002M	Bio-Plex Pro Mouse Cytokine IL-1 β Set, 1 x 96-well
171-G5003M	Bio-Plex Pro Mouse Cytokine IL-2 Set, 1 x 96-well
171-G5004M	Bio-Plex Pro Mouse Cytokine IL-3 Set, 1 x 96-well
171-G5005M	Bio-Plex Pro Mouse Cytokine IL-4 Set, 1 x 96-well
171-G5006M	Bio-Plex Pro Mouse Cytokine IL-5 Set, 1 x 96-well
171-G5007M	Bio-Plex Pro Mouse Cytokine IL-6 Set, 1 x 96-well
171-G5008M	Bio-Plex Pro Mouse Cytokine IL-9 Set, 1 x 96-well
171-G5009M	Bio-Plex Pro Mouse Cytokine IL-10 Set, 1 x 96-well
171-G5010M	Bio-Plex Pro Mouse Cytokine IL-12 (p40) Set, 1 x 96-well
171-G5011M	Bio-Plex Pro Mouse Cytokine IL-12 (p70) Set, 1 x 96-well
171-G5012M	Bio-Plex Pro Mouse Cytokine IL-13 Set, 1 x 96-well
171-G5013M	Bio-Plex Pro Mouse Cytokine IL-17 Set, 1 x 96-well
171-G5014M	Bio-Plex Pro Mouse Cytokine Eotaxin Set, 1 x 96-well
171-G5015M	Bio-Plex Pro Mouse Cytokine G-CSF Set, 1 x 96-well
171-G5016M	Bio-Plex Pro Mouse Cytokine GM-CSF Set, 1 x 96-well
171-G5017M	Bio-Plex Pro Mouse Cytokine IFN- γ Set, 1 x 96-well
171-G5018M	Bio-Plex Pro Mouse Cytokine KC Set, 1 x 96-well
171-G5019M	Bio-Plex Pro Mouse Cytokine MCP-1 (MCAF) Set, 1 x 96-well
171-G5020M	Bio-Plex Pro Mouse Cytokine MIP-1 α Set, 1 x 96-well
171-G5021M	Bio-Plex Pro Mouse Cytokine MIP-1 β Set, 1 x 96-well
171-G5022M	Bio-Plex Pro Mouse Cytokine RANTES Set, 1 x 96-well
171-G5023M	Bio-Plex Pro Mouse Cytokine TNF- α Set, 1 x 96-well
Bio-Plex Pro Mouse Cytokine Singleplex Sets, Group II*	
171-G6001M	Bio-Plex Pro Mouse Cytokine IL-15 Set, 1 x 96-well
171-G6002M	Bio-Plex Pro Mouse Cytokine Basic FGF Set, 1 x 96-well
171-G6003M	Bio-Plex Pro Mouse Cytokine LIF Set, 1 x 96-well
171-G6004M	Bio-Plex Pro Mouse Cytokine M-CSF Set, 1 x 96-well
171-G6005M	Bio-Plex Pro Mouse Cytokine MIG Set, 1 x 96-well
171-G6006M	Bio-Plex Pro Mouse Cytokine MIP-2 Set, 1 x 96-well
171-G6007M	Bio-Plex Pro Mouse Cytokine PDGF-BB Set, 1 x 96-well
171-G6008M	Bio-Plex Pro Mouse Cytokine VEGF Set, 1 x 96-well
Bio-Plex Pro Cytokine Reagent Kits***	
171-304070	Bio-Plex Pro Reagent Kit , 1 x 96-well, includes assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, standard diluent, sample diluent
171-304071	Bio-Plex Pro Reagent Kit , 10 x 96-well, includes assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, standard diluent, sample diluent
Bio-Plex Human Cytokine Standards	
171-D50001	Bio-Plex Human Cytokine Standards Group I , 1-pack of 27 analytes
171-D10501	Bio-Plex Human Cytokine Standards Group I , 50-pack of 27 analytes
171-D60001	Bio-Plex Human Cytokine Standards Group II , 1-pack of 23 analytes
171-D10502	Bio-Plex Human Cytokine Standards Group II , 50-pack of 23 analytes
Bio-Plex Mouse Cytokine Standards	
171-I50001	Bio-Plex Mouse Cytokine Standards Group I , 1-pack of 23 analytes
171-I10501	Bio-Plex Mouse Cytokine Standards Group I , 50-pack of 23 analytes
171-I60001	Bio-Plex Mouse Cytokine Standards Group II , 1-pack of 9 analytes
171-I10502	Bio-Plex Mouse Cytokine Standards Group II , 50-pack of 9 analytes
Bio-Plex Assay Accessories	
171-304500	Bio-Plex Wash Buffer , 1.5 L
171-Q25001	Bio-Plex Pro Plates , 40 pack

* Each singleplex set includes coupled magnetic beads and detection antibodies for the detection of the specified analyte (reagent kits and standards are offered separately).

** Human ICAM-1 and VCAM-1 require different sample dilutions. It is recommended to order these targets separately.

*** Required only with singleplex sets. Reagent kit now contains diluents. Therefore, separate diluent kit is not required.

Note: x-Plex assays or Express assays can be ordered with any combination of analytes within a group, except for ICAM-1 and VCAM-1. For more information, go to www.bio-rad.com/bio-plex/x-plex/.

Available Assays*

Human Group I	8-Plex Panel (M50-000007A)	17-Plex Panel (M50-00031YV)	27-Plex Panel (M50-0KCAF0Y)	Th1/Th2 Panel (M50-00005L3)	Singleplex/x-Plex™ Bead Regions*
IL-1β		●	●		● 39
IL-1ra			●		● 25
IL-2	●	●	●	●	● 38
IL-4	●	●	●	●	● 52
IL-5		●	●	●	● 33
IL-6	●	●	●		● 19
IL-7		●	●		● 74
IL-8	●	●	●		● 54
IL-9			●		● 77
IL-10	●	●	●	●	● 56
IL-12 (p70)		●	●	●	● 75
IL-13		●	●	●	● 51
IL-15			●		● 73
IL-17		●	●		● 76
Basic FGF			●		● 44
Eotaxin			●		● 43
G-CSF		●	●		● 57
GM-CSF	●	●	●	●	● 34
IFN-γ	●	●	●	●	● 21
IP-10			●		● 48
MCP-1 (MCAF)		●	●		● 53
MIP-1α			●		● 55
MIP-1β		●	●		● 18
PDGF-BB			●		● 47
RANTES			●		● 37
TNF-α	●	●	●	●	● 36
VEGF			●		● 45

Human Group II	21-Plex Panel plus ICAM-1 and VCAM-1 (MF0-005KMII)	Singleplex/x-Plex Bead Regions*
IL-1α	●	● 63
IL-2Rα	●	● 13
IL-3	●	● 64
IL-12 (p40)	●	● 28
IL-16	●	● 27
IL-18**	●	● 42
CTACK	●	● 72
GRO-α	●	● 61
HGF	●	● 62
ICAM-1***		● 12
IFN-α2	●	● 20
LIF	●	● 29
MCP-3	●	● 26
M-CSF	●	● 67
MIF	●	● 35
MIG	●	● 14
β-NGF	●	● 46
SCF	●	● 65
SCGF-β	●	● 78
SDF-1α	●	● 22
TNF-β	●	● 30
TRAIL	●	● 66
VCAM-1***		● 15

Mouse Group I	8-Plex Panel (M60-000007A)	23-Plex Panel (M60-009RDPD)	Th1/Th2 Panel (M60-00003J7)	Singleplex/x-Plex Bead Regions*
IL-1α		●		● 53
IL-1β	●	●		● 19
IL-2	●	●	●	● 36
IL-3		●		● 18
IL-4	●	●	●	● 39
IL-5	●	●	●	● 52
IL-6		●		● 38
IL-9		●		● 33
IL-10	●	●	●	● 56
IL-12 (p40)		●		● 76
IL-12 (p70)		●	●	● 78
IL-13		●		● 37
IL-17		●		● 72
Eotaxin		●		● 74
G-CSF		●		● 54
GM-CSF	●	●	●	● 73
IFN-γ	●	●	●	● 34
KC		●		● 57
MCP-1 (MCAF)		●		● 51
MIP-α		●		● 77
MIP-1β		●		● 75
RANTES		●		● 55
TNF-α	●	●	●	● 21

Mouse Group II	9-Plex Panel (MD0-00000EL)	Singleplex/x-Plex Bead Regions*
IL-15	●	● 42
IL-18**	●	● 20
Basic FGF	●	● 25
LIF	●	● 45
M-CSF	●	● 26
MIG	●	● 44
MIP-2	●	● 27
PDGF-BB	●	● 35
VEGF	●	● 47

* You can design an x-Plex assay or Express assay with any combination of analytes within a group, except for ICAM-1 and VCAM-1.

For more information, go to www.bio-rad.com/bio-plex/x-plex/ or contact your local Bio-Rad sales representative.

** IL-18 is not available as a singleplex assay, only as part of an assay panel.

*** This assay requires different sample dilutions.

Note: Red bead regions indicate regions that differ between magnetic and nonmagnetic assay targets.

Bio-Plex Assays

Bio-Plex Rat Cytokine, Chemokine, and Growth Factor Assays

Bio-Plex rat cytokine, chemokine, and growth factor assays can be ordered as a premixed panel, or as an x-Plex assay for any combination of analytes within a group. For more information, go to www.bio-rad.com/bio-plex/x-plex/.

What to Order

1 	2 	3 
Cytokine multiplex assay panel or singleplex assay — contains coupled beads, detection antibody, and lyophilized standards to test specific proteins of interest.	Cytokine reagent kit — includes assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, 96-well filter plate, sealing tape, and instructions required to prepare samples and acquire data.	Diluent kit (optional) — if serum or plasma samples are to be tested, use the species-specific diluent kit for optimum recovery.

Catalog #	Description	Catalog #	Description
Bio-Plex Premixed Rat Cytokine Assay Panel			
171-K11070	Bio-Plex Rat Cytokine 9-Plex A Panel , 1 x 96-well, includes coupled beads, detection antibodies, and standards for the detection of IL-1 α , IL-1 β , IL-2, IL-4, IL-6, IL-10, GM-CSF, IFN- γ , TNF- α	Bio-Plex Reagent Kits	
		171-304000	Bio-Plex Cytokine Reagent Kit , 1 x 96-well
		171-304001	Bio-Plex Cytokine Reagent Kit , 10 x 96-well
		Bio-Plex Rat Serum Diluent Kit (Recommended for serum and plasma samples)	
		171-305008	Bio-Plex Rat Serum Diluent Kit , 1 x 96-well, includes 15 ml sample diluent and 10 ml standard diluent

Bio-Plex Signal Transduction Assays

Available Assays

Assays	Phosphoprotein	Total
Akt Signaling		
Akt (Ser ⁴⁷³)	●	●
GSK-3 α / β (Ser ²¹ /Ser ⁹)	●	
Cell Cycle/Checkpoint Control		
p53 (Ser ¹⁵)	●	●
p53 (Ser ³⁷)	●	●
p53 (Ser ⁴⁶)	●	●
Chromatin Regulation/Acetylation		
Histone H3 (Ser ¹⁰)	●	
Immunology/Inflammation		
I κ B- α (Ser ³² /Ser ³⁶)	●	●
NF- κ B p65 (Ser ⁵³⁶)	●	
STAT2 (Tyr ⁶⁸⁹)	●	
STAT3 (Ser ⁷²⁷)	●	
STAT3 (Tyr ⁷⁰⁵)	●	
STAT6 (Tyr ⁶⁴¹)	●	
Tyk2 (Tyr ¹⁰⁵⁴ /Tyr ¹⁰⁵⁵)	●	
Glucose/Energy Metabolism		
IGF-IR (Tyr ¹¹³¹)	●	
IR- β (Tyr ¹¹⁴⁶)	●	
IRS-1 (Ser ⁶³⁶ /Ser ⁶³⁹)	●	

Assays	Phosphoprotein	Total
MAP Kinase Signaling		
ATF-2 (Thr ⁷¹)	●	●
c-Jun (Ser ⁶³)	●	●
ERK1 (Thr ²⁰² /Tyr ²⁰⁴)	●	
ERK2 (Thr ¹⁸⁵ /Tyr ¹⁸⁷)	●	●
ERK1/2 (Thr ²⁰² /Tyr ²⁰⁴ , Thr ¹⁸⁵ /Tyr ¹⁸⁷)	●	●
HSP27 (Ser ⁷⁸)	●	●
JNK (Thr ¹⁸³ /Tyr ¹⁸⁵)	●	●
MEK1 (Ser ²¹⁷ /Ser ²²¹)	●	●
p38 MAPK (Thr ¹⁸⁰ /Tyr ¹⁸²)	●	●
p90RSK (Thr ³⁵⁹ /Ser ³⁶³)	●	●
Neuroscience		
CREB (Ser ¹³³)	●	●
TrkA (Tyr ⁴⁹⁰)	●	
Translational Control		
p70 S6 kinase (Thr ⁴²¹ /Ser ⁴²⁴)	●	
S6 ribosomal protein (Ser ²³⁵ /Ser ²³⁶)	●	
Tyrosine Kinases		
Bcr-Abl (Tyr ²⁴⁵)	●	●
c-Abl (Tyr ²⁴⁵)	●	
c-Abl (Tyr ⁴¹²)	●	
EGFR (Tyr)	●	
PDGF receptor- β (Tyr ⁷⁵¹)	●	
Src (Tyr ⁴¹⁶)	●	



CST antibodies exclusively developed and validated for Bio-Plex phosphoprotein and total target assays.

Bio-Plex signal transduction assays can be ordered as singleplex assays, premixed panels, or as an x-Plex assay for any combination of analytes within a group. For more information, go to www.bio-rad.com/bio-plex/x-plex/.

What to Order

		
Bio-Plex phosphoprotein assays and Bio-Plex total target assays — includes coupled beads, detection antibodies, and lysates to test for specific proteins of interest in lysate samples.	Bio-Plex cell lysis kit — includes cell lysis buffer, cell wash buffer, factor 1, and factor 2; essential for optimal lysis of your culture or tissue samples.	Bio-Plex phosphoprotein detection reagent kit — includes assay buffer, wash buffer, antibody diluent, resuspension buffer, streptavidin-PE, 96-well filter plate, sealing tape, and instructions required to prepare assays and acquire data.

Catalog #	Description	Catalog #	Description
Bio-Plex Premixed Phosphoprotein Assay Panels		Bio-Plex Total Target Singleplex Assays*	
X70-0000A1B	Bio-Plex Phospho 3-Plex Panel , 1 x 96-well, includes coupled beads, detection antibodies, and lysates for the detection of phosphorylated GSK-3 α/β , p70 S6 kinase, and STAT3	171-V25374**	Bio-Plex Phospho-S6 Ribosomal Protein (Ser²³⁵/Ser²³⁶) Assay , 1 x 96-well
X70-000CLDZ	Bio-Plex Phospho 3-Plex Panel , 1 x 96-well, includes coupled beads, detection antibodies, and lysates for the detection of phosphorylated ATF-2, NF- κ B p65, and p90RSK	171-V23732	Bio-Plex Phospho-STAT2 (Tyr⁶⁸⁹) Assay , 1 x 96-well
X70-00001RD	Bio-Plex Phospho 5-Plex Panel , 1 x 96-well, includes coupled beads, detection antibodies, and lysates for the detection of phosphorylated Akt, ERK1/2, I κ B- α , JNK, and p38 MAPK	171-V26952	Bio-Plex Phospho-STAT3 (Ser⁷²⁷) Assay , 1 x 96-well
		171-V22552	Bio-Plex Phospho-STAT3 (Tyr⁷⁰⁵) Assay , 1 x 96-well
		171-V25772	Bio-Plex Phospho-STAT6 (Tyr⁶⁴¹) Assay , 1 x 96-well
		171-V23973	Bio-Plex Phospho-TrkA (Tyr⁴⁹⁰) Assay , 1 x 96-well
		171-V26533	Bio-Plex Phospho-Tyk2 (Tyr¹⁰⁵⁴/Tyr¹⁰⁵⁵) Assay , 1 x 96-well
Bio-Plex Premixed Total Target Panels		Bio-Plex Reagent Kits	
X90-00001TK	Bio-Plex Total 5-Plex Panel , 1 x 96-well, includes coupled beads, detection antibodies, and lysates for the detection of Akt, ERK1/2, I κ B- α , JNK, and p38 MAPK	171-304004	Bio-Plex Phosphoprotein Detection Reagent Kit , 1 x 96-well
		171-304005	Bio-Plex Phosphoprotein Detection Reagent Kit , 10 x 96-well
Bio-Plex Phosphoprotein Singleplex Assays*		Bio-Plex Sample Preparation Kits	
171-V21075	Bio-Plex Phospho-Akt (Ser⁴⁷³) Assay , 1 x 96-well	171-304011	Bio-Plex Cell Lysis Kit , 1 x 96-well
171-V21620	Bio-Plex Phospho-ATF-2 (Thr⁷¹) Assay , 1 x 96-well	171-304012	Bio-Plex Cell Lysis Kit , 10 x 96-well
171-V27145	Bio-Plex Phospho-Bcr-Abl (Tyr²⁴⁵) Assay , 1 x 96-well		
171-V27545	Bio-Plex Phospho-c-Abl (Tyr²⁴⁵) Assay , 1 x 96-well		
171-V27745	Bio-Plex Phospho-c-Abl (Tyr⁴¹²) Assay , 1 x 96-well		
171-V24356	Bio-Plex Phospho-c-Jun (Ser⁶³) Assay , 1 x 96-well		
171-V26119	Bio-Plex Phospho-CREB (Ser¹³³) Assay , 1 x 96-well		
171-V23120**	Bio-Plex Phospho-EGFR (Tyr) Assay , 1 x 96-well		
171-V21938	Bio-Plex Phospho-ERK1 (Thr²⁰²/Tyr²⁰⁴) Assay , 1 x 96-well		
171-V20438	Bio-Plex Phospho-ERK2 (Thr¹⁸⁵/Tyr¹⁸⁷) Assay , 1 x 96-well		
171-V22238	Bio-Plex Phospho-ERK1/2 (Thr²⁰²/Tyr²⁰⁴, Thr¹⁸⁵/Tyr¹⁸⁷) Assay , 1 x 96-well		
171-V23318	Bio-Plex Phospho-GSK-3α/β (Ser²¹/Ser⁹) Assay , 1 x 96-well		
171-V24777	Bio-Plex Phospho-Histone H3 (Ser¹⁰) Assay , 1 x 96-well		
171-V24551	Bio-Plex Phospho-HSP27 (Ser⁷⁸) Assay , 1 x 96-well		
171-V27343	Bio-Plex Phospho-IGF-IR (Tyr¹¹³¹) Assay , 1 x 96-well		
171-V20758	Bio-Plex Phospho-IκB-α (Ser³²/Ser³⁶) Assay , 1 x 96-well		
171-V27943	Bio-Plex Phospho-IR-β (Tyr¹¹⁴⁶) Assay , 1 x 96-well		
171-V25576	Bio-Plex Phospho-IRS-1 (Ser⁶³⁶/Ser⁶³⁹) Assay , 1 x 96-well		
171-V21034	Bio-Plex Phospho-JNK (Thr¹⁸³/Tyr¹⁸⁵) Assay , 1 x 96-well		
171-V25340	Bio-Plex Phospho-MEK1 (Ser²¹⁷/Ser²²¹) Assay , 1 x 96-well		
171-V24937	Bio-Plex Phospho-NF-κB p65 (Ser⁵³⁶) Assay , 1 x 96-well		
171-V21336	Bio-Plex Phospho-p38 MAPK (Thr¹⁸⁰/Tyr¹⁸²) Assay , 1 x 96-well		
171-V25153	Bio-Plex Phospho-p53 (Ser¹⁵) Assay , 1 x 96-well		
171-V28353	Bio-Plex Phospho-p53 (Ser³⁷) Assay , 1 x 96-well		
171-V28153	Bio-Plex Phospho-p53 (Ser⁴⁶) Assay , 1 x 96-well		
171-V24155	Bio-Plex Phospho-p70 S6 Kinase (Thr⁴²¹/Ser⁴²⁴) Assay , 1 x 96-well		
171-V23535	Bio-Plex Phospho-p90RSK (Thr³⁵⁹/Ser³⁶³) Assay , 1 x 96-well		
171-V25957	Bio-Plex Phospho-PDGF Receptor-β (Tyr⁷⁵¹) Assay , 1 x 96-well		
171-V26741	Bio-Plex Phospho-Src (Tyr⁴¹⁶) Assay , 1 x 96-well		

* For multiplexing capabilities of specific assay combinations, request bulletin 5483, or visit us on the Web at www.bio-rad.com/bio-plex/.

** This assay cannot be multiplexed.

Bio-Plex Pro™ Magnetic Diabetes Assays **new**

Available Assays

Human Assays	2-Plex Panel	10-Plex Panel	Singleplex/Express	Bead Region
Adiponectin*	●		●	64
Adipsin*	●		●	35
C-peptide		●	●	72
Ghrelin (total)		●	●	26
GIP		●	●	14
GLP-1		●	●	27
Glucagon		●	●	15
Insulin		●	●	12
Leptin		●	●	78
PAI-1 (total)		●	●	61
Resistin		●	●	65
Visfatin		●	●	22

Mouse Assays	Singleplex Only	8-Plex Panel	Singleplex/Express	Bead Region
Adiponectin*	●		●	29
Ghrelin (total)		●	●	64
GIP		●	●	46
GLP-1		●	●	62
Glucagon		●	●	63
Insulin		●	●	66
Leptin		●	●	65
PAI-1 (total)		●	●	48
Resistin		●	●	30

* Due to different dilution schemes, adiponectin and adipsin are offered as singleplex assays only.

Bio-Plex Pro diabetes assays can be ordered as follows:

Premixed panels — The new all-in-one kit packaging format includes premixed coupled magnetic beads, detection antibodies (10-plex for human, 8-plex for mouse), standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, standard diluent, and sample diluent.

Express assay — For faster delivery, customized assay panels in the new all-in-one kit packaging format can be ordered with any combination of available assays. Coupled magnetic beads and detection antibodies are not premixed. Go to www.bio-rad.com/bio-plex/x-plex/ for details.

Individual components — Singleplex sets (coupled magnetic beads and detection antibodies), standards, reagent kits, and Bio-Plex Pro flat bottom plates for magnetic bead-based washing are available separately. To run an assay, singleplex sets require standards and reagent kits.

For more information request bulletin 5966 for human diabetes assays or bulletin 5943 for mouse diabetes assays.

What to Order



Bio-Plex Pro diabetes assays — magnetic bead-based assays built around MagPlex magnetic beads, for the measurement of human and mouse diabetes biomarkers. Magnetic beads afford compatibility with Bio-Plex Pro and Bio-Plex Pro II wash stations, providing decreased variability and increased consistency among experiments.

Catalog #	Description
171-A7001M*	Bio-Plex Pro Premixed Human Diabetes Assays Bio-Plex Pro Human Diabetes 10-Plex Panel, 1 x 96-well, includes premixed conjugated magnetic beads and detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, standard diluent, and sample diluent for the detection of C-peptide, ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, resistin, and visfatin
171-A7010M*	Bio-Plex Pro Human Diabetes 10-Plex Panel , 10 x 96-well, includes premixed conjugated magnetic beads and detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plates, sealing tape, standard diluent, and sample diluent for the detection of C-peptide, ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, resistin, and visfatin

Bio-Plex Pro Premixed Human Diabetes Assays

171-A7001M* **Bio-Plex Pro Human Diabetes 10-Plex Panel**, 1 x 96-well, includes premixed conjugated magnetic beads and detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, standard diluent, and sample diluent for the detection of C-peptide, ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, resistin, and visfatin

171-A7010M* **Bio-Plex Pro Human Diabetes 10-Plex Panel**, 10 x 96-well, includes premixed conjugated magnetic beads and detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plates, sealing tape, standard diluent, and sample diluent for the detection of C-peptide, ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, resistin, and visfatin

Catalog #	Description	Catalog #	Description
171-A7002M	Bio-Plex Pro Human Diabetes Adipsin and Adiponectin Assays , 1 x 96-well, includes conjugated magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, and serum-based standard diluent for the detection of adiponectin and adipsin. These are offered as singleplex assays; customers create the 2-plex	171-B7012M	Bio-Plex Pro Human Diabetes Visfatin Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human visfatin, requires reagent kit and standards
171-A7003M	Bio-Plex Pro Human Diabetes Adiponectin Assay , 1 x 96-well, includes conjugated magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, and serum-based standard diluent for the detection of adiponectin	Bio-Plex Pro Mouse Diabetes Singleplex Sets, 1 x 96	
171-A7004M	Bio-Plex Pro Human Diabetes Adipsin Assay , 1 x 96-well, includes conjugated magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, and serum-based standard diluent for the detection of adipsin	171-G7002M	Bio-Plex Pro Mouse Diabetes Ghrelin Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of mouse ghrelin, requires reagent kit and standards
Bio-Plex Pro Premixed Mouse Diabetes Assays		171-G7003M	Bio-Plex Pro Mouse Diabetes GIP Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of mouse GIP, requires reagent kit and standards
171-F7001M*	Bio-Plex Pro Mouse Diabetes 8-Plex Panel , 1 x 96-well, includes premixed conjugated magnetic beads and detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, standard diluent, and sample diluent for the detection of ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, and resistin	171-G7004M	Bio-Plex Pro Mouse Diabetes GLP-1 Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of mouse GLP-1, requires reagent kit and standards
171-F7010M*	Bio-Plex Pro Mouse Diabetes 8-Plex Panel , 10 x 96-well, includes premixed conjugated magnetic beads and detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plates, sealing tape, standard diluent, and sample diluent for the detection of ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, and resistin	171-G7005M*	Bio-Plex Pro Mouse Diabetes Glucagon Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of mouse glucagon, requires reagent kit and standards
171-F7002M	Bio-Plex Pro Mouse Diabetes Adiponectin Assay , 1 x 96-well, includes conjugated magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, and serum-based standard diluent for the detection of adiponectin	171-G7006M	Bio-Plex Pro Mouse Diabetes Insulin Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of mouse insulin, requires reagent kit and standards
Bio-Plex Pro Human Diabetes Singleplex Sets, 1 x 96		171-G7007M	Bio-Plex Pro Mouse Diabetes Leptin Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of mouse leptin, requires reagent kit and standards
171-B7003M	Bio-Plex Pro Human Diabetes C-Peptide Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human C-peptide, requires reagent kit and standards	171-G7008M	Bio-Plex Pro Mouse Diabetes PAI-1 Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of mouse PAI-1, requires reagent kit and standards
171-B7004M	Bio-Plex Pro Human Diabetes Ghrelin Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human ghrelin, requires reagent kit and standards	171-G7009M	Bio-Plex Pro Mouse Diabetes Resistin Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of mouse resistin, requires reagent kit and standards
171-B7005M	Bio-Plex Pro Human Diabetes GIP Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human GIP, requires reagent kit and standards		* Coming soon
171-B7006M	Bio-Plex Pro Human Diabetes GLP-1 Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human GLP-1, requires reagent kit and standards	Bio-Plex Pro Human Diabetes Singleplex Sets, 10 x 96	
171-B7007M*	Bio-Plex Pro Human Diabetes Glucagon Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human glucagon, requires reagent kit and standards	171-B7103M	Bio-Plex Pro Human Diabetes C-Peptide Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human C-peptide, requires reagent kit and standards
171-B7008M	Bio-Plex Pro Human Diabetes Insulin Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human insulin, requires reagent kit and standards	171-B7104M	Bio-Plex Pro Human Diabetes Ghrelin Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human ghrelin, requires reagent kit and standards
171-B7009M	Bio-Plex Pro Human Diabetes Leptin Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human leptin, requires reagent kit and standards	171-B7105M	Bio-Plex Pro Human Diabetes GIP Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human GIP, requires reagent kit and standards
171-B7010M	Bio-Plex Pro Human Diabetes PAI-1 Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human PAI-1, requires reagent kit and standards	171-B7106M	Bio-Plex Pro Human Diabetes GLP-1 Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human GLP-1, requires reagent kit and standards
171-B7011M	Bio-Plex Pro Human Diabetes Resistin Set , 1 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human resistin, requires reagent kit and standards	171-B7107M*	Bio-Plex Pro Human Diabetes Glucagon Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human glucagon, requires reagent kit and standards
		171-B7108M	Bio-Plex Pro Human Diabetes Insulin Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human insulin, requires reagent kit and standards
		171-B7109M	Bio-Plex Pro Human Diabetes Leptin Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human leptin, requires reagent kit and standards
		171-B7110M	Bio-Plex Pro Human Diabetes PAI-1 Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human PAI-1, requires reagent kit and standards

Catalog #	Description	Catalog #	Description
171-B7111M	Bio-Plex Pro Human Diabetes Resistin Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human resistin, requires reagent kit and standards	171-G7108M	Bio-Plex Pro Mouse Diabetes PAI-1 Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of mouse PAI-1, requires reagent kit and standards
171-B7112M	Bio-Plex Pro Human Diabetes Visfatin Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of human visfatin, requires reagent kit and standards	171-G7109M	Bio-Plex Pro Mouse Diabetes Resistin Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of mouse resistin, requires reagent kit and standards
Bio-Plex Pro Mouse Diabetes Singleplex Sets, 10 x 96			
171-G7102M	Bio-Plex Pro Mouse Diabetes Ghrelin Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of mouse ghrelin, requires reagent kit and standards	171-304070	Bio-Plex Pro Reagent Kits** Bio-Plex Pro Reagent Kit , 1 x 96-well, includes assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, standard diluent, and sample diluent
171-G7103M	Bio-Plex Pro Mouse Diabetes GIP Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of mouse GIP, requires reagent kit and standards	171-304071	Bio-Plex Pro Reagent Kit , 10 x 96-well, includes assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plates, sealing tape, standard diluent, and sample diluent
171-G7104M	Bio-Plex Pro Mouse Diabetes GLP-1 Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of mouse GLP-1, requires reagent kit and standards	171-305042	Bio-Plex Pro Diabetes Serum-Based Diluent , 1 x 96-well, 70 ml serum-based diluent for adiponectin and adipsin assays
171-G7105M*	Bio-Plex Pro Mouse Diabetes Glucagon Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of mouse glucagon, requires reagent kit and standards	Bio-Plex Pro Diabetes Standards	
171-G7106M	Bio-Plex Pro Mouse Diabetes Insulin Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of mouse insulin, requires reagent kit and standards	171-D70001	Bio-Plex Pro Human Diabetes Standards , 1-pack of 12 analytes, for adiponectin, adipsin, C-peptide, ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, resistin, and visfatin
171-G7107M	Bio-Plex Pro Mouse Diabetes Leptin Set , 10 x 96-well, includes conjugated magnetic beads and detection antibodies for the detection of mouse leptin, requires reagent kit and standards	171-D70050	Bio-Plex Pro Human Diabetes Standards , 50-pack of 12 analytes, for adiponectin, adipsin, C-peptide, ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, resistin, and visfatin
		171-I70001	Bio-Plex Pro Mouse Diabetes Standards , 1-pack of 9 analytes, for adiponectin, ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, and resistin
		171-I70050	Bio-Plex Pro Mouse Diabetes Standards , 50-pack of 9 analytes, for adiponectin, ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, and resistin

* Coming soon

**Required only with singleplex sets. Same reagent kit can be used for both diabetes and cytokine assays.

Bio-Plex Pro Disease Research Assays

Available Targets

Human Acute Phase	Human Angiogenesis	Human Isotyping	Bio-Plex Precision Pro™ Human Cytokines
α -2-macroglobulin	Angiopoeitin-2	IgG ₁	IL-1 β
CRP	Follistatin	IgG ₂	IL-2
Ferritin	G-CSF	IgG ₃	IL-4
Fibrinogen	HGF	gG ₄	IL-5
Haptoglobin	IL-8	IgM	IL-6
Procalcitonin	Leptin	IgA	IL-10
SAA	PDGF-BB	IgE	IL-12 (p70)
SAP	PECAM-1		IL-13
Tissue plasminogen activator	VEGF		IFN- γ
			TNF- α

What to Order

1 	2 	3 
Bio-Plex Pro assay panel — 1 x 96-well and 10 x 96-well, includes coupled magnetic beads, detection antibodies, standards, and controls.	Bio-Plex Pro reagent kit — includes assay buffer, wash buffer, streptavidin-PE, 96-well filter plate, sealing tape, and panel-specific instructions required to prepare assays and acquire data.	Bio-Plex diluent kit — choose the appropriate diluent kit for the panel for diluting samples, standards, and controls.

Bio-Plex Pro Human Acute Phase Assay Panel

Catalog #	Description
Bio-Plex Pro Human Acute Phase Complete Kits*	
171-A4S07M	Bio-Plex Pro Human Acute Phase 5 + 4-Plex Panel Complete Kit , 1 x 96-well
171-A4C09M	Bio-Plex Pro Human Acute Phase 4-Plex Panel Complete Kit , 1 x 96-well
171-A4C07M	Bio-Plex Pro Human Acute Phase 5-Plex Panel Complete Kit , 1 x 96-well
Bio-Plex Pro Human Acute Phase Assay Kits**	
171-A4007M	Bio-Plex Pro Human Acute Phase 5-Plex Panel , 1 x 96-well, includes coupled beads, detection antibodies, and standards for the detection of ferritin, fibrinogen, procalcitonin, serum amyloid A, tissue plasminogen activator
171-A4008M	Bio-Plex Pro Human Acute Phase 5-Plex Panel , 10 x 96-well
171-A4009M	Bio-Plex Pro Human Acute Phase 4-Plex Panel , 1 x 96-well, includes coupled beads, detection antibodies, and standards for the detection of α -2-macroglobulin, CRP, haptoglobin, serum amyloid P
171-A4010M	Bio-Plex Pro Human Acute Phase 4-Plex Panel , 10 x 96-well
Acute Phase Reagent Kits	
171-304050	Bio-Plex Pro Human Acute Phase Reagent Kit , 1 x 96-well, includes Bio-Plex reagent kit and acute phase instruction manual
171-304051	Bio-Plex Pro Human Acute Phase Reagent Kit , 10 x 96-well
Acute Phase Diluent Kit***	
171-305050	Bio-Plex Pro Human Acute Phase Diluent Kit , 1 x 96-well, includes 15 ml serum-based diluent and 300 ml serum-free diluent

* For simplified ordering, the complete kits include the acute phase panel, acute phase reagent kit, and acute phase diluent kit.

** One acute phase reagent kit is required for use with each assay kit.

*** One acute phase diluent kit is sufficient to perform the 5-plex and the 4-plex assays.

Bio-Plex Pro Human Angiogenesis Assay Panel

Catalog #	Description
Bio-Plex Pro Human Angiogenesis Complete Kit*	
171-A4S11M	Bio-Plex Pro Human Angiogenesis 9-Plex Panel Complete Kit , 1 x 96-well
Bio-Plex Pro Human Angiogenesis Panels	
171-A4011M	Bio-Plex Pro Human Angiogenesis 9-Plex Panel , 1 x 96-well, includes coupled beads, detection antibodies, and standards for the detection of angiopoietin-2, follistatin, G-CSF, HGF, IL-8, leptin, PDGF-BB, PECAM-1, and VEGF
171-A4012M	Bio-Plex Pro Human Angiogenesis 9-Plex Panel , 10 x 96-well
Angiogenesis Reagent Kits	
171-304060	Bio-Plex Pro Angiogenesis Reagent Kit , 1 x 96-well, includes Bio-Plex reagent kit and angiogenesis instruction manual
171-304061	Bio-Plex Pro Angiogenesis Reagent Kit , 10 x 96-well
Angiogenesis Diluent Kit	
171-305000	Bio-Plex Human Serum Diluent Kit , 1 x 96-well, includes 15 ml sample diluent and 10 ml standard diluent

* For simplified ordering, the complete kits include the angiogenesis panel, angiogenesis reagent kit, and human serum diluent kit.

Bio-Plex Pro Human Isotyping Assay Panel

Catalog #	Description
Bio-Plex Pro Human Isotyping Complete Kit*	
171-A3S01M	Bio-Plex Pro Human Isotyping Panel Complete Kit , 1 x 96-well
Bio-Plex Pro Human Isotyping Assay Panels	
171-A3001M	Bio-Plex Pro Human Isotyping 7-Plex Panel , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, and controls for the detection of IgG ₁ , IgG ₂ , IgG ₃ , IgG ₄ , IgA, IgE, IgM
171-A3002M	Bio-Plex Pro Human Isotyping 7-Plex Panel , 10 x 96-well
Isotyping Reagent Kits	
171-304020	Bio-Plex Pro Isotyping Reagent Kit , 1 x 96-well, includes assay buffer, wash buffer, streptavidin-PE, 96-well filter plate, sealing tape, instructions
171-304021	Bio-Plex Pro Isotyping Reagent Kit , 10 x 96-well
Isotyping Diluent Kits	
171-305030	Bio-Plex Pro Isotyping Diluent , 1 x 96-well
171-305031	Bio-Plex Pro Isotyping Diluent , 10 x 96-well

* For simplified ordering, the complete kit includes the isotyping panel, isotyping reagent kit, and isotyping serum diluent.

Bio-Plex® Precision Pro™ Human Cytokine 10-Plex Assay Panel

What to Order



Bio-Plex Precision Pro human cytokine 10-plex panel — highly sensitive, highly reproducible magnetic bead-based assays allow accurate measurement of low levels of human cytokines, while offering low intra- and inter-assay percent coefficient of variation (%CV). Convenient kit format includes assay, reagent, and diluent components all in a single box.

Catalog #	Description
Bio-Plex Precision Pro Human Cytokine Assays	
171-A1001P*	Bio-Plex Precision Pro Human Cytokine 10-Plex Panel , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, controls, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, serum standard diluent, plasma standard diluent, and sample diluent for the detection of IL-1 β , IL-2, IL-4, IL-5, IL-6, IL-10, IL-12 (p70), IL-13, IFN- γ , TNF- α

* This all-in-one assay does not require separate reagent and diluent kits. Not available as x-Plex assays.

Bio-Plex Assay Standards



Bio-Plex assay standards are available in packages of 1 or 50 for all human (groups I and II) and mouse (groups I and II) cytokine assays, and in packages of 2 or 50 for all human acute phase, human angiogenesis, human diabetes, and human isotyping assays.

Available Targets

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

Disease Research				
Human Acute Phase	Human Angiogenesis	Human Diabetes	Mouse Diabetes	Human Isotyping
α -2-macroglobulin	Angiopoeitin-2	Adiponectin	Adiponectin	IgG ₁
CRP	Follistatin	Adipsin	Ghrelin	IgG ₂
Ferritin	G-CSF	C-peptide	GIP	IgG ₃
Fibrinogen	HGF	Ghrelin	GLP-1	IgG ₄
Haptoglobin	IL-8	GIP	Glucagon	IgM
Procalcitonin	Leptin	GLP-1	Insulin	IgA
SAA	PDGF-BB	Glucagon	Leptin	IgE
SAP	PECAM-1	Insulin	PAI-1	
Tissue plasminogen activator	VEGF	Leptin	Resistin	
		PAI-1		
		Resistin		
		Visfatin		

Catalog # Description

Bio-Plex Human Cytokine Standards

171-D50001 **Bio-Plex Human Cytokine Standards Group I**, 1-pack of 27 analytes
 171-D10501 **Bio-Plex Human Cytokine Standards Group I**, 50-pack of 27 analytes
 171-D60001 **Bio-Plex Human Cytokine Standards Group II**, 1-pack of 23 analytes
 171-D10502 **Bio-Plex Human Cytokine Standards Group II**, 50-pack of 23 analytes

Bio-Plex Mouse Cytokine Standards

171-I50001 **Bio-Plex Mouse Cytokine Standards Group I**, 1-pack of 23 analytes
 171-I10501 **Bio-Plex Mouse Cytokine Standards Group I**, 50-pack of 23 analytes
 171-I60001 **Bio-Plex Mouse Cytokine Standards Group II**, 1-pack of 9 analytes
 171-I10502 **Bio-Plex Mouse Cytokine Standards Group II**, 50-pack of 9 analytes

Bio-Plex Disease Research

171-D40002 **Bio-Plex Human Acute Phase Standards**, 2-pack of 9 analytes
 171-D40006 **Bio-Plex Human Acute Phase Standards**, 50-pack of 9 analytes
 171-D40003 **Bio-Plex Human Angiogenesis Standards**, 2-pack of 9 analytes
 171-D40004 **Bio-Plex Human Angiogenesis Standards**, 50-pack of 9 analytes

Catalog # Description

171-D70001 **Bio-Plex Pro Human Diabetes Standards**, 1-pack of 12 analytes
 171-D70050 **Bio-Plex Pro Human Diabetes Standards**, 50-pack of 12 analytes
 171-I70001 **Bio-Plex Pro Mouse Diabetes Standards**, 1-pack of 9 analytes
 171-I70050 **Bio-Plex Pro Mouse Diabetes Standards**, 50-pack of 9 analytes
 171-D30001 **Bio-Plex Human Isotyping Standards**, 2-pack of 7 analytes
 171-D30501 **Bio-Plex Human Isotyping Standards**, 50-pack of 7 analytes

Bio-Plex Coupling Reagents

Bio-Plex Pro™ Magnetic COOH Beads **new**

Bead Region Selection Guide

Bead Types	1 ml Regions Available Off-the-Shelf*
Bio-Plex Pro Magnetic COOH Beads**	26–29, 34–37, 43–46, 52–55, 62–65

* All 1 ml regions not listed above and 4 ml regions 1–150 are available made to order. Go to www.bio-rad.com for which regions are compatible with the Bio-Plex/Luminex 200 system.

** Catalog #MC1-00XX-01, where xx = 2-digit number for bead region (for example, #MC1-0026-01 to order bead region 26).

Bio-Plex COOH Beads

Bead Region Selection Guide

Bead Types	1 ml Regions Available Off-the-Shelf*
Bio-Plex COOH Beads**	11, 16, 18, 20, 24–28, 30, 31, 33, 35, 38, 42, 43–46, 50–53, 56, 66

* All 1 ml regions not listed above and 16 ml regions 1–60 are available made to order. Please contact your local Bio-Rad sales representative for a delivery date.

** Catalog #171-5060xx, where xx = bead region (for example, use #171-506012 to order bead region 12).

Catalog # Description

Beads

See table above **Bio-Plex COOH Beads**, 1 ml

For more information, go to www.bio-rad.com/bio-plex/ or contact your local Bio-Rad sales representative.

Amine Coupling Kit

171-406001 **Bio-Plex Amine Coupling Kit**, 30 reactions, includes bead wash buffer, bead activation buffer, storage buffer, staining buffer coupling reaction tubes

Instruction Manual for Coupling Reagents

Catalog # Description
4110012C **Bio-Plex Pro Magnetic COOH Beads Bio-Plex COOH Beads Amine Coupling Kit Instruction Manual**

The Bio-Plex suspension array system includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation.



CST antibodies exclusively developed and validated for Bio-Plex phosphoprotein and total target assays.

MagPlex and xPONENT are trademarks of Luminex Corporation.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 31 3689 6600 **Canada** 905 364 3435 **China** 86 20 8732 2339 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300 **Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 0508 805 500 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000