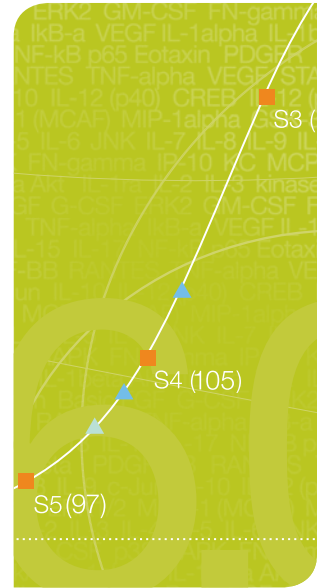
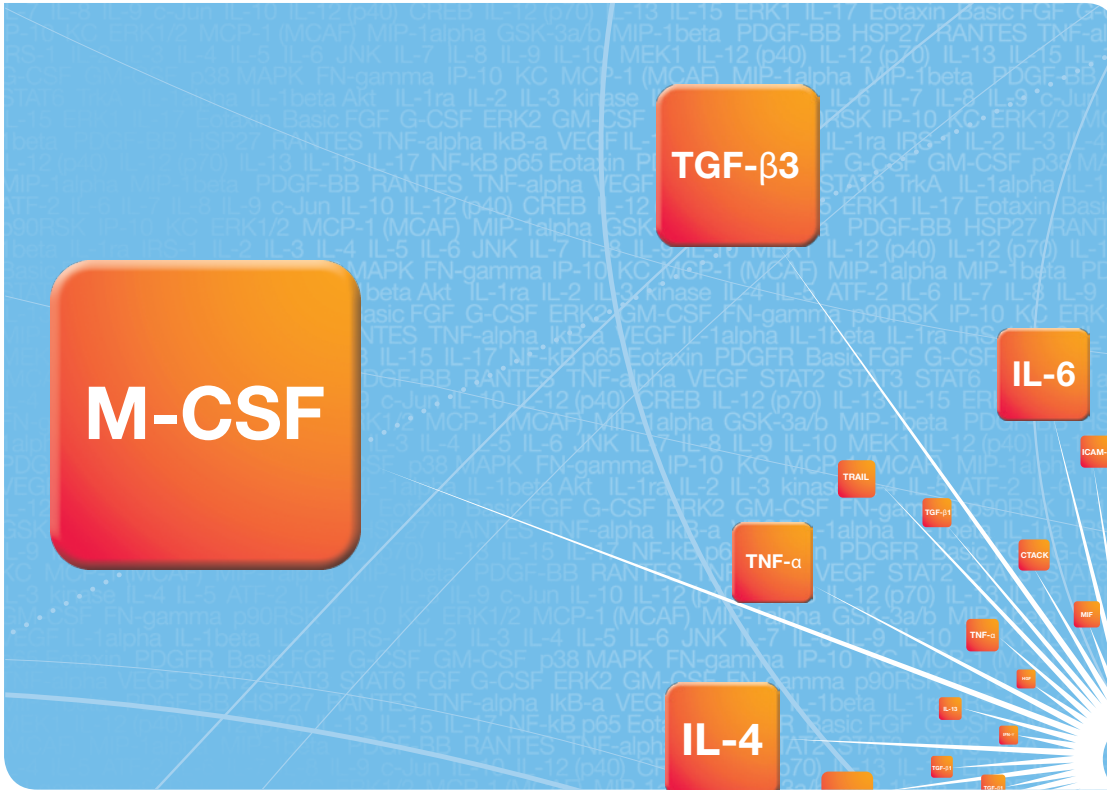


Multiplex Technology



Spring 2013 Ordering Information

Bio-Plex[®] Multiplex System



Table of Contents

Bio-Plex Instruments	2
Bio-Plex 3D Suspension Array System	2
Bio-Plex 200 System	2
System Performance Validation and Calibration Tools	3
Accessories and Replacement Parts	3
Bio-Plex MAGPIX Multiplex Reader	4
Bio-Plex Pro Wash Stations	4
Bio-Plex Software	5
Bio-Plex Software	5
Bio-Plex Assays	6
Bio-Plex Assay Builder for Custom Assays	6
Bio-Plex Pro Cytokine, Chemokine, and Growth Factor Assays	6
Human Cytokines	7
Mouse Cytokines	9
Rat Cytokines	10
TGF- β (for All Species)	11
Cell Signaling Assays	11
Bio-Plex Pro Magnetic Cell Signaling Assays	11
Bio-Plex Nonmagnetic Signal Transduction Assays	13
Kidney Toxicity Assays New	15
Human Kidney Toxicity	15
Canine Kidney Toxicity	15
Rat Kidney Toxicity	15
Disease Research Assays	16
Bio-Plex Pro Human Cancer Biomarker Assays New	16
Bio-Plex Pro Diabetes Assays	17
Human Diabetes	17
Non-Human Primate Diabetes New	18
Mouse Diabetes	18
Rat Diabetes	18
Reagent Kits (for All Species)	18
Bio-Plex Pro Human Acute Phase Assay Panel	19
Bio-Plex Pro Human Angiogenesis Assay Panel	19
Bio-Plex Pro Human Isotyping Assays New	20
Bio-Plex Precision Pro Human Cytokine Assay Panel	20
Beads, Reagents, and Accessories	21
COOH Beads and Related Reagents	21
Bio-Plex Pro COOH Magnetic Beads	21
Bio-Plex COOH Nonmagnetic Beads	21
Bio-Plex Amine Coupling Kit	21
Bio-Plex Streptavidin-PE	21
Accessories	21
Bio-Plex Pro Assay Accessories	21

Bio-Plex Instruments

Bio-Plex® 3D Suspension Array System



Bio-Plex 3D Suspension Array System with Bio-Plex Manager™ Software

Catalog #	Description
Bio-Plex3D	Bio-Plex 3D Suspension Array System , includes Bio-Plex 3D suspension array system, xPONENT acquisition software, PC, calibration and verification reagents, Bio-Plex Manager™ desktop license
8920-18500001*	xPONENT CFR21 Part 11 Software Module , enables U.S. FDA 21 CFR Part 11 compliance
8920-18700001*	xPONENT Automation Module , enables interfacing with robotics workstations
171-022001*	Swivel Base , allows rotation of the instrument away from the robotics workstation for maintenance and sample loading
8920-18600001*	xPONENT LIS Software Module , enables interfacing with LIMS/LIS databases
8920-18200001*	xPONENT Extra Seat Licenses , 3 additional seats of xPONENT software

* All additional accessories may be purchased through Bio-Rad Laboratories, Inc. at the time of system purchase. Post system-purchase accessories must be purchased directly through Luminex Corporation.

For More Information

Web: www.bio-rad.com/bio-plex3D

Bio-Plex 200 System

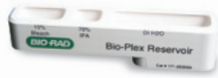


Catalog #	Description
171-000201	Bio-Plex 200 System , 100–240 V, includes array reader, microplate platform, Bio-Plex Manager software with workstation license, PC, monitor, calibration kit, validation kit 4.0, MCV plate IV, Bio-Plex reservoir, 20 L sheath fluid, spare sample needle
171-000205	Bio-Plex 200 System with HTF , same as 171-000201 with high-throughput fluidics (HTF)

For More Information

Web: www.bio-rad.com/bio-plex200

Bio-Plex Instruments



Bio-Plex Reservoir



Bio-Plex MCV Plate IV



Bio-Plex Validation Kit 4.0



Bio-Plex Calibration Kit

Catalog #	Description
-----------	-------------

System Performance Validation and Calibration Tools

171-203001	Bio-Plex Validation Kit 4.0 , includes optics validation, reporter validation, classify validation, and fluidics validation bead sets for approximately 50 validation routines using Bio-Plex Manager 4.0, 4.1.1, 5.0, or 6.0 software (standard or Security Edition) and MCV plate IV
171-203033	Bio-Plex MCV Plate IV , for use with Bio-Plex Manager 4.0, 4.1.1, 5.0, or 6.0 software (standard or Security Edition) and validation kit 4.0
171-203060	Bio-Plex Calibration Kit , includes Cal1 and Cal2 calibration beads for approximately 50 daily calibration routines

For More Information

Web: www.bio-rad.com/bio-plexvalidation

Request or download bulletins: 2879; instruction manual — 4110007

Catalog #	Description
-----------	-------------

Accessories and Replacement Parts

171-203050	Bio-Plex Reservoir
171-000055	Sheath Fluid , 1x, 20 L
171-002001	Communication Cable , 5 ft DB9 cable, connects the microplate platform with the PC serial port
171-002002	Communication Cable , 3 ft CAN BUS cable, connects the array reader with the HTF system
171-002003	Communication Cable , 5 ft USB cable, connects the array reader with a PC USB port
171-002010	Sheath Fluid Bottle , 1 L
171-002012	Sheath Waste Bottle , 1 L
171-002020	Sample Needle , 11.6 cm (4.6")
171-002024	Alignment Guide , aligns the array reader on the microplate platform
171-002026	Needle Adjustment Wrench
171-002030	Protective Shield for Sample Needle
171-002032	Air Intake Filter
171-002033	Syringe Seal
171-002034	Syringe Seal with Cylinder
171-002038	Sheath Fluid Filter with Quick Connect Tubing , removes particles >5 µm in diameter
171-002040	Sheath Cube Filter , 10 µm, removes particles >5 µm in diameter
171-002056	Sheath Cube Filter , 10 µm, includes tubing, removes particles >10 µm in diameter
171-002023	Needle Guide , allows sample needle access to microplates

Bio-Plex Instruments

Bio-Plex® MAGPIX™ Multiplex Reader



Catalog #	Description
171-015001	Bio-Plex MAGPIX Multiplex Reader with Bio-Plex Manager Software , includes Bio-Plex MAGPIX instrument, PC with xPONENT acquisition software installed, Bio-Plex Manager desktop license, calibration kit, verification kit, 2 drive fluid cartridges, 2 waste containers

Kits and Reagents

171-213001	Bio-Plex MAGPIX Calibration Kit , calibration testing kit for Bio-Plex MAGPIX instrument, includes 5 ml calibrator microspheres, twenty-five 8-well strips, CD; 25 applications
171-213002	Bio-Plex MAGPIX Verification Kit , includes MAGPIX verifier, 5 ml microspheres, MAGPIX fluidics 1, MAGPIX fluidics 2, MAGPIX performance verification kit CD, twenty-five 8-well strips; 25 applications
171-213003	Bio-Plex MAGPIX Drive Fluid , pkg of 4, 700 ml drive fluid, for use with Bio-Plex MAGPIX instrument

Accessories

171-012004	Bio-Plex MAGPIX Replacement Waste Fluid Container , pkg of 1, 850 ml container, holds waste fluid, for use with Bio-Plex MAGPIX instrument
171-012005	Bio-Plex MAGPIX Sample Probe Needle , pkg of 1, needle
171-012006	Bio-Plex MAGPIX Sample Probe Height Adjustment Kit , sample probe height adjustment kit
171-012008	Bio-Plex MAGPIX 96-Well Plate Heater Block , pkg of 1, heater block, fits one 96-well plate, for use with Bio-Plex MAGPIX instrument

For More Information

Request or download bulletin: 6005

Bio-Plex Pro™ Wash Stations



Catalog #	Description
300-34376	Bio-Plex Pro Wash Station , includes magnetic plate carrier, waste bottle, 2 buffer bottles
300-34377	Bio-Plex Pro II Wash Station , includes magnetic plate carrier, vacuum manifold plate carrier, waste bottle, 2 buffer bottles
171-020100	Bio-Plex Handheld Magnetic Washer , includes magnetic washer and adjustment hex tools for use in manual wash steps for all Bio-Plex magnetic assays

Accessories

171-025001	Bio-Plex Pro Flat Bottom Plates , 40 x 96-well plates, for washing magnetic beads using the Bio-Plex Pro or Bio-Plex Pro II wash stations
171-304500	Bio-Plex Wash Buffer , 1.5 L
171-304502	Filter Plate , 1 x 96-well filter plate with clear plastic lid, plate holder tray, for washing nonmagnetic beads using the Bio-Plex Pro II wash station, optional vacuum manifold; sealing tape not included

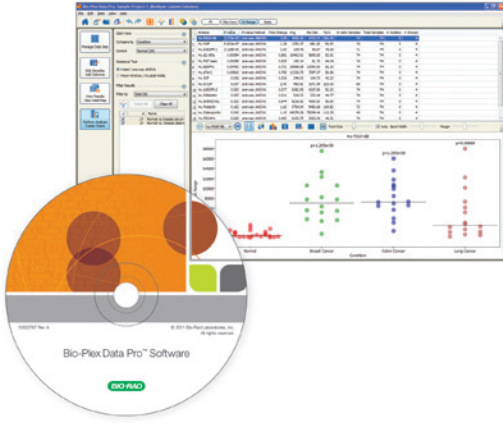
For More Information

Web: www.bio-rad.com/washstation

Request or download bulletin: 5703

Bio-Plex Software

Bio-Plex® Software



Catalog #	Description
171-STND01	Bio-Plex Manager™ Software , includes 1-user desktop license, to analyze Bio-Plex data and generate protocols, does not operate the instrument
171-SCRT00	Bio-Plex Manager Instrument Control License , Security, upgrades workstation with Bio-Plex Manager software, allows compliance with U.S. FDA 21 CFR Part 11 regulations
171-001510	Bio-Plex Data Pro™ Software with Bio-Plex Manager Software , includes 5 Bio-Plex Data Pro software licenses, 5-user desktop licenses for Bio-Plex Manager software
171-001513	Bio-Plex Data Pro Software , includes 5 Bio-Plex Data Pro software licenses

Contact Bio-Rad Technical Support to verify whether your current system is compatible with these software products.

For More Information

Web: www.bio-rad.com/datapro

Request or download bulletin: 5613

Bio-Plex Assays

Bio-Plex® Assay Builder for Custom Assays

Configure a custom assay using our online assay builder.

Design your own assays from our selection of hundreds of analytes in four easy steps. With assay builder you can:

- Get a price quote
- Place your order online
- Look up an existing assay by catalog number

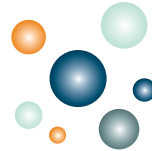
Go to www.bio-rad.com/assaybuilder

x-Plex™ Assay



We Mix

Express Assay



You Mix

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

Available Assays

Analytes	Human	Mouse	Rat
β-NGF	•	—	—
CD40L	—	•	—
CTACK	•	—	—
Eotaxin	•	•	—
EPO	—	—	•
FGF-basic	•	•	—
G-CSF	•	•	•
GM-CSF	•	•	•
GRO/KC	—	•	•
GRO-α	•	—	—
HGF	•	—	—
ICAM-1	•	•	—
IFN-α2	•	—	—
IFN-γ	•	•	•
IL-1α	•	•	•
IL-1β	•	•	•
IL-1ra	•	—	—
IL-2	•	•	•
IL-2Rα	•	—	—
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-17A/F	•	—	—
IL-17e (IL-25)	—	•	—
IL-17F	•	•	—

Available Assays

Analytes	Human	Mouse	Rat
IL-18	•	•	•
IL-21	•	•	—
IL-22	•	•	—
IL-23	•	—	—
IL-23p19	—	•	—
IL-25	•	—	—
IL-27p28	—	•	—
IL-31	•	•	—
IL-33	•	•	—
IP-10	•	—	—
KC	—	•	•
LIF	•	•	—
MCP-1	•	•	•
MCP-3	•	—	—
M-CSF	•	•	•
MIF	•	—	—
MIG	•	•	—
MIP-1α	•	•	•
MIP-1β	•	•	—
MIP-2	—	•	•
MIP-3α	—	•	•
PDGF-BB	•	•	—
RANTES	•	•	•
sCD40L	•	—	—
SCF	•	—	—
SCGF-β	•	—	—
SDF-1α	•	—	—
TGF-β1	•	•	•
TGF-β2	•	•	•
TGF-β3	•	•	•
TNF-α	•	•	•
TNF-β	•	—	—
TRAIL	•	—	—
VCAM-1	•	—	—
VEGF	•	•	•

Bio-Plex Assays



All-in-one Kit

Singleplex Set

Standard

Reagent Kit

Human Cytokines

Catalog #	Description
Premixed All-in-One Kits*	
M50-00007A	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, flat bottom 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-12p70, IL-13, IL-17, G-CSF, GM-CSF, IFN- γ , MCP-1, 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, flat bottom plate, sealing tape, standard diluent, sample diluent for the detection of IL-1 α , IL-2R α , IL-3, IL-12p40, 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-1 α , IL-2, IL-4, IL-5, IL-6, IL-7, IL-8, IL-9, IL-10, IL-12p70, IL-13, IL-15, IL-17, basic FGF, eotaxin, G-CSF, GM-CSF, IFN- γ , IP-10, MCP-1, 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, flat bottom plate, sealing tape, standard diluent, sample diluent, for the detection of IL-2, IL-4, IL-5, IL-10, IL-12p70, IL-13, GM-CSF, IFN- γ , TNF- α
171-AA001M	Bio-Plex Pro Human Th17 Cytokine 15-Plex Panel , 1 x 96-well all-in-one kit, includes premixed coupled magnetic beads, premixed detection antibodies, standards, 2-level quality controls, detection antibody diluent, standard diluent HB, sample diluent HB, assay buffer, wash buffer, streptavidin-PE, 96-well flat bottom plate, sealing tape, and instructions, for the detection of IL-1 β , IL-4, IL-6, IL-10, IL-17A, IL-17F, IL-21, IL-22, IL-23, IL-25, IL-31, IL-33, IFN- γ , sCD40L, TNF- α

* Please contact your local Bio-Rad sales representative for pricing and availability of 10 x 96-well assay kits.

Catalog #	Description	Catalog #	Description
Human Th17 Cytokine Singleplex Sets*, 1 x 96-Well			
171-BA001M	IL-1β Set	171-BA009M	IL-23 Set
171-BA002M	IL-4 Set	171-BA010M	IL-25 Set
171-BA003M	IL-6 Set	171-BA011M	IL-31 Set
171-BA004M	IL-10 Set	171-BA012M	IL-33 Set
171-BA005M	IL-17A Set	171-BA013M	IFN-γ Set
171-BA006M	IL-17F Set	171-BA014M	sCD40L Set
171-BA007M	IL-21 Set	171-BA015M	TNF-α Set
171-BA008M	IL-22 Set	171-BA016M	IL-17A/F Set

* Include coupled magnetic beads and detection antibody. Require Bio-Plex Pro reagent kit II and human Th17 cytokine standards.

Catalog #	Description
Standards	
171-DA0001	Bio-Plex Pro Human Th17 Cytokine 16-plex Standards , pkg of 1 vial, lyophilized mixture of 16 standard analytes
171-DA0501	Bio-Plex Pro Human Th17 Cytokine 16-plex Standards , pkg of 50 lot-matched vials, lyophilized mixture of 16 standard analytes
Reagent Kits*	
171-304055	Bio-Plex Pro Reagent Kit II with Filter Plate , 1 x 96-well, includes detection antibody diluent, standard diluent HB, sample diluent HB, assay buffer, wash buffer, streptavidin-PE, 96-well filter plate, sealing tape, and instructions, for use with vacuum wash methods
171-304055M	Bio-Plex Pro Reagent Kit II with Flat Bottom Plate , 1 x 96-well, includes detection antibody diluent, standard diluent HB, sample diluent HB, assay buffer, wash buffer, streptavidin-PE, 96-well flat bottom plate, sealing tape, and instructions, for use with magnetic wash methods

* Required only with singleplex sets.

Bio-Plex Assays

Catalog #	Description	Catalog #	Description
Human Cytokine Singleplex Sets, Group I*, 1 x 96-Well			
171-B5001M	IL-1β Set	171-B5015M	Eotaxin Set
171-B5002M	IL-1α Set	171-B5016M	Basic FGF Set
171-B5003M	IL-2 Set	171-B5017M	G-CSF Set
171-B5004M	IL-4 Set	171-B5018M	GM-CSF Set
171-B5005M	IL-5 Set	171-B5019M	IFN-γ Set
171-B5006M	IL-6 Set	171-B5020M	IP-10 Set
171-B5007M	IL-7 Set	171-B5021M	MCP-1 Set
171-B5008M	IL-8 Set	171-B5022M	MIP-1α Set
171-B5009M	IL-9 Set	171-B5023M	MIP-1β Set
171-B5010M	IL-10 Set	171-B5024M	PDGF-BB Set
171-B5011M	IL-12p70 Set	171-B5025M	RANTES Set
171-B5012M	IL-13 Set	171-B5026M	TNF-α Set
171-B5013M	IL-15 Set	171-B5027M	VEGF Set
171-B5014M	IL-17A Set		

* Includes coupled magnetic beads and detection antibody. Requires reagent kit and human cytokine standards group I.

Catalog #	Description
Standards	
171-D50001	Bio-Plex Pro Human Cytokine Standards Group I , 1 pack of 27 analytes
171-D10501	Bio-Plex Pro Human Cytokine Standards Group I , 50-pack of 27 analytes

Catalog #	Description	Catalog #	Description
Singleplex Sets, Group II*, 1 x 96-Well			
171-B6001M	IL-1α Set	171-B6012M	MCP-3 Set
171-B6002M	IL-2Rα Set	171-B6013M	M-CSF Set
171-B6003M	IL-3 Set	171-B6014M	MIF Set
171-B6004M	IL-12p40 Set	171-B6015M	MIG Set
171-B6005M	IL-16 Set	171-B6016M	β-NGF Set
171-B6006M	CTACK Set	171-B6017M	SCF Set
171-B6007M	GRO-α Set	171-B6018M	SCGF-β Set
171-B6008M	HGF Set	171-B6019M	SDF-1α Set
171-B6009M	ICAM-1 Set	171-B6020M	TNF-β Set
171-B6010M	IFN-α2 Set	171-B6021M	TRAIL Set
171-B6011M	LIF Set	171-B6022M	VCAM-1 Set

* Includes coupled magnetic beads and detection antibody. Requires reagent kit and human cytokine standards group II.

Catalog #	Description
Standards	
171-D60001	Bio-Plex Pro Human Cytokine Standards Group II , 1 pack of 23 analytes
171-D10502	Bio-Plex Pro Human Cytokine Standards Group II , 50-pack of 23 analytes

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
171-304070M	Bio-Plex Pro Reagent Kit with Flat Bottom Plate , 1 x 96-well, includes flat bottom plate, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, sealing tape, standard diluent, sample diluent, and instructions

Bio-Plex Assays

Mouse Cytokines

Catalog #	Description
Premixed All-in-One Kits*	
M60-00007A	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, flat bottom 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
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, flat bottom 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-12p40, IL-12p70, IL-13, IL-17, eotaxin, G-CSF, GM-CSF, IFN- γ , KC, MCP-1, 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, flat bottom plate, sealing tape, standard diluent, sample diluent for the detection of IL-2, IL-4, IL-5, IL-10, IL-12p70, GM-CSF, IFN- γ , TNF- α
171-FA001M	Bio-Plex Pro Mouse Cytokine Th17 Panel B 8-Plex Group III , 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 CD40L, IL-17F, IL-21, IL-22, IL-23p19, IL-31, IL-33, MIP-3 α
M60-00007NY	Bio-Plex Pro Mouse Cytokine Th17 Panel A 6-Plex Group I , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standard, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, flat bottom plate, sealing tape, standard diluent, sample diluent, and instructions, for the detection of IL-1 β , IL-6, IL-10, IL-17A, IFN- γ , TNF- α
L60-00004C6	Bio-Plex Pro Mouse Cytokine Th1 Panel , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, buffers, diluents, streptavidin-PE, 96-well flat bottom plate, and sealing tape for the detection IL-1 β , IL-2, IL-6, IL-10, IL-12p70, IFN- γ , TNF- α
L60-000UKVT	Bio-Plex Pro Mouse Cytokine Th2 Panel , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, buffers, diluents, streptavidin-PE, 96-well flat bottom plate, and sealing tape for the detection IL-2, IL-4, IL-5, IL-6, IL-9, IL-10, IL-13

* Please contact your local Bio-Rad sales representative for pricing and availability of 10 x 96-well assay kits.

Catalog #	Description	Catalog #	Description
Mouse Cytokine Singleplex Sets, Group I*, 1 x 96-Well			
171-G5001M	IL-1α Set	171-G5013M	IL-17A Set
171-G5002M	IL-1β Set	171-G5014M	Eotaxin Set
171-G5003M	IL-2 Set	171-G5015M	G-CSF Set
171-G5004M	IL-3 Set	171-G5016M	GM-CSF Set
171-G5005M	IL-4 Set	171-G5017M	IFN-γ Set
171-G5006M	IL-5 Set	171-G5018M	KC Set
171-G5007M	IL-6 Set	171-G5019M	MCP-1 Set
171-G5008M	IL-9 Set	171-G5020M	MIP-1α Set
171-G5009M	IL-10 Set	171-G5021M	MIP-1β Set
171-G5010M	IL-12p40 Set	171-G5022M	RANTES Set
171-G5011M	IL-12p70 Set	171-G5023M	TNF-α Set
171-G5012M	IL-13 Set		

* Includes coupled magnetic beads and detection antibody. Requires reagent kit and mouse cytokine standards group I.

Catalog #	Description	Catalog #	Description
Standards			
171-I50001	Bio-Plex Pro Mouse Cytokine Standards Group I , 1 pack of 23 analytes		
171-I10501	Bio-Plex Pro Mouse Cytokine Standards Group I , 50-pack of 23 analytes		
Mouse Cytokine Singleplex Sets, Group II*, 1 x 96-Well			
171-G6001M	IL-15 Set	171-G6005M	MIG Set
171-G6002M	Basic FGF Set	171-G6006M	MIP-2 Set
171-G6003M	LIF Set	171-G6007M	PDGF-BB Set
171-G6004M	M-CSF Set	171-G6008M	VEGF Set

* Includes coupled magnetic beads and detection antibody. Requires reagent kit and mouse cytokine standards group II.

Catalog #	Description
Standards	
171-I60001	Bio-Plex Pro Mouse Cytokine Standards Group II , 1 pack of 9 analytes
171-I10502	Bio-Plex Pro Mouse Cytokine Standards Group II , 50-pack of 9 analytes

Bio-Plex Assays

Catalog #	Description	Catalog #	Description
Mouse Th17 Singleplex Sets, Group III*, 1 x 96-Well			
171-GA001M	CD40L Set	171-GA007M	IL-27p28 Set
171-GA002M	IL-17F Set	171-GA008M	IL-31 Set
171-GA003M	IL-21 Set	171-GA009M	IL-33 Set
171-GA004M	IL-22 Set	171-GA010M	ICAM-1 Set**
171-GA005M	IL-23p19 Set	171-GA011M	MIP-3α Set
171-GA006M	IL-25 Set		

* Includes coupled magnetic beads and detection antibody. Requires reagent kit and mouse cytokine standards group III.

** Requires high dilution reagent kit.

Catalog #	Description
Standards	
171-IA0001	Bio-Plex Pro Mouse Cytokine Standards Group III , pkg of 1 vial, standard for detecting 11 analytes
171-IA0501	Bio-Plex Pro Mouse Cytokine Standards Group III , pkg of 50 lot-matched vials, standard for detecting 11 analytes
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; for vacuum separation methods
171-304070M	Bio-Plex Pro Reagent Kit with Flat Bottom Plate , 1 x 96-well, includes assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, flat bottom plate, sealing tape, standard diluent, sample diluent; for use with magnetic separation methods
171-304080M	Bio-Plex Pro High Dilution Reagent Kit , 1 x 96-well, includes assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, flat bottom plate, filter plate, sealing tape, standard diluent, sample diluent

* Required only with singleplex sets.

Rat Cytokines

Catalog #	Description
Premixed All-in-One Kits*	
171-K1001M	Bio-Plex Pro Rat Cytokine 24-Plex Assay , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, flat bottom plate, sealing tape, standard diluent, sample diluent for the detection of EPO, G-CSF, GM-CSF, GRO/KC, IFN-γ, IL-1α, IL-1β, IL-2, IL-4, IL-5, IL-6, IL-7, IL-10, IL-12p70, IL-13, IL-17A, IL-18, MCP-1, M-CSF, MIP-1α, MIP-3α, RANTES, TNF-α, VEGF
171-K1002M	Bio-Plex Pro Rat Cytokine Th1/Th2 Panel , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standard, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, flat bottom plate, sealing tape, standard diluent, sample diluent and instructions for the detection of IL-α, IL-1β, IL-2, IL-4, IL-5, IL-6, IL-10, IL-12p70, IL-13, GM-CSF, IFN-γ, TNF-α

* Please contact your local Bio-Rad sales representative for pricing and availability of 10 x 96-well assay kits.

Catalog #	Description	Catalog #	Description
Rat Cytokine Singleplex Sets*, 1 x 96-Well			
171-L1002M	EPO Set	171-L1015M	IL-12p40 Set
171-L1003M	G-CSF Set	171-L1016M	IL-12p70 Set
171-L1004M	GM-CSF Set	171-L1017M	IL-13 Set
171-L1005M	GRO/KC Set	171-L1018M	IL-17A Set
171-L1006M	IFN-γ Set	171-L1027M	MCP-1 Set
171-L1007M	IL-1α Set	171-L1020M	M-CSF Set
171-L1008M	IL-1β Set	171-L1021M	MIP-1α Set
171-L1009M	IL-2 Set	171-L1022M	MIP-2 Set
171-L1010M	IL-4 Set	171-L1023M	MIP-3α Set
171-L1011M	IL-5 Set	171-L1024M	RANTES Set
171-L1012M	IL-6 Set	171-L1025M	TNF-α Set
171-L1013M	IL-7 Set	171-L1026M	VEGF Set
171-L1014M	IL-10 Set		

* Includes coupled magnetic beads and detection antibody. Requires reagent kit and rat cytokine standards.

Catalog #	Description
Standards*	
171-NZ0001	Bio-Plex Pro Rat Cytokine Standards , 1 pack of 31 analytes, compatible with rat cytokine and rat diabetes assays
171-NZ0501	Bio-Plex Pro Rat Cytokine Standards , 50-pack of 31 analytes, standard for detecting 31 rat targets. Compatible with rat cytokine and rat diabetes assays

Bio-Plex Assays

Catalog # Description

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
171-304070M	Bio-Plex Pro Reagent Kit with Flat Bottom Plate , 1 x 96-well, includes flat bottom plate, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, sealing tape, standard diluent, sample diluent, and instructions

* Required only with singleplex sets.

TGF- β (for All Species)

Catalog # Description

Premixed All-in-One Kit

171-W4001M	Bio-Plex Pro TGF-β 3-plex Assay , 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 TGF- β 1, TGF- β 2, TGF- β 3
------------	---

Bio-Plex TGF- β Singleplex Sets*, 1 x 96-Well

171-W4001M	TGF-β1 Set
171-V4002M	TGF-β2 Set
171-V4003M	TGF-β3 Set

* Includes coupled magnetic beads and detection antibody. Requires reagent kit and TGF- β standards.

Standards

171-X40001	Bio-Plex Pro TGF-β Standard , 1 pack of 1 TGF- β standard for detecting TGF- β 1, TGF- β 2, and TGF- β 3 analytes
171-X40501	Bio-Plex Pro TGF-β Standard , 1 pack of 50 lot-matched vials, TGF- β standard for detecting TGF- β 1, TGF- β 2, and TGF- β 3 analytes

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, for vacuum separation methods
171-304070M	Bio-Plex Pro Reagent Kit with Flat Bottom Plate , 1 x 96-well, includes assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, flat bottom plate, sealing tape, standard diluent, sample diluent, for use with magnetic separation methods

Note: x-Plex™ assay panels can be ordered with any combination of available assays. For more information, go to www.bio-rad.com/assaybuilder.

For More Information

Web: www.bio-rad.com/CCGFassays

Request or download bulletins: 5828, 6054, 6100, and 6110

Cell Signaling Assays




Bio-Plex Pro Magnetic Cell Signaling Assays

Available Magnetic Assays

Assays	Phosphoprotein	Total
Akt (Ser ⁴⁷³)	•	•
Akt (Thr ³⁰⁸)	•	•
c-Jun (Ser ⁶³)	•	•
EGFR (Tyr ¹⁰⁶⁸)	•	—
EGFR (Tyr ¹¹⁷³)	•	—
Erk1/2 (Thr ²⁰² /Tyr ²⁰⁴ , Thr ¹⁸⁵ /Tyr ¹⁸⁷)	•	•
GSK-3 α / β (Ser ²¹ /Ser ⁹)	•	•
HER-2 (Tyr ¹²⁴⁸)	•	•
IGF-IR (Tyr ¹¹³¹)	•	—
I κ B- α (Ser ³² /Ser ³⁶)	•	•
JNK (Thr ¹⁸³ /Tyr ¹⁸⁵)	•	•

Assays	Phosphoprotein	Total
MEK1 (Ser ²¹⁷ /Ser ²²¹)	•	•
NF- κ B p65 (Ser ⁵³⁶)	•	—
p38 MAPK (Thr ¹⁸⁰ /Tyr ¹⁸²)	•	•
p70 S6 Kinase (Thr ⁴²¹ /Ser ⁴²⁴)	•	•
p70 S6 Kinase (Thr ³⁸⁹)	•	•
PDGFR- α (Tyr ⁷⁵⁴)	•	—
PDGFR- β (Tyr ⁷⁵¹)	•	—
Smad2 (Ser ⁴⁶⁵ /Ser ⁴⁶⁷)	•	•
Stat1 (Tyr ⁷⁰¹)	•	—
Stat3 (Ser ⁷²⁷)	•	—
Stat3 (Tyr ⁷⁰⁵)	•	—
VEGFR-2 (Tyr ¹¹⁷⁵)	•	—

Bio-Plex Assays

		
<p>Bio-Plex Pro cell signaling singleplex sets* — include coupled beads and detection antibodies. Users can mix compatible sets to create multiplex assays.</p>	<p>Bio-Plex Pro cell signaling reagent kit** — includes cell lysis buffer, cell lysis factor QG, cell wash buffer, bead resuspension buffer, detection antibody diluent, wash buffer, streptavidin-PE, 96-well flat bottom plate, sealing tape, and instructions.</p>	<p>Bio-Plex Pro cell lysate controls*** — refer to the lysate controls selection guide to identify the appropriate lysate for your assay.</p>

* Required for detecting your phosphoprotein or total protein of interest.

** Required for preparing lysates and running the assay.

*** Recommended for use as qualitative phosphoprotein and total protein assay controls.

Catalog #	Description	Catalog #	Description
Phosphoprotein Singleplex Sets*, 1 x 96-Well			
171-V50001M	Akt (Ser⁴⁷³)	171-V50013M	NF-κB p65 (Ser⁵³⁶)
171-V50002M	Akt (Thr³⁰⁸)	171-V50014M	p38 MAPK (Thr¹⁸⁰/Tyr¹⁸²)
171-V50003M	c-Jun (Ser⁶³)	171-V50015M	p70 S6 Kinase (Thr⁴²¹/Ser⁴²⁴)
171-V50004M	EGFR (Tyr¹⁰⁶⁸)	171-V50016M	p70 S6 Kinase (Thr³⁸⁹)
171-V50005M	EGFR (Tyr¹¹⁷³)	171-V50017M	PDGFR-α (Tyr⁷⁵⁴)
171-V50006M	Erk1/2 (Thr²⁰²/Tyr²⁰⁴, Thr¹⁸⁵/Tyr¹⁸⁷)	171-V50018M	PDGFR-β (Tyr⁷⁵¹)
171-V50007M	GSK-3α/β (Ser²¹/Ser⁹)	171-V50019M	Smad2 (Ser⁴⁶⁵/Ser⁴⁶⁷)
171-V50008M	HER-2 (Tyr¹²⁴⁸)	171-V50020M	Stat1 (Tyr⁷⁰¹)
171-V50009M	IGF-IR (Tyr¹¹³¹)	171-V50021M	Stat3 (Ser⁷²⁷)
171-V50010M	IκB-α (Ser³²/Ser³⁶)	171-V50022M	Stat3 (Tyr⁷⁰⁵)
171-V50011M	JNK (Thr¹⁸³/Tyr¹⁸⁵)	171-V50023M	VEGFR-2 (Tyr¹¹⁷⁵)
171-V50012M	MEK1 (Ser²¹⁷/Ser²²¹)		

* Includes coupled magnetic beads and detection antibodies for detecting phosphorylated protein. Requires cell signaling reagent kit and lysate control.

Catalog #	Description	Catalog #	Description
Total Target Singleplex Sets			
171-V60001M	Akt	171-V60007M	JNK
171-V60002M	c-Jun	171-V60008M	MEK1
171-V60003M	Erk1/2	171-V60009M	p38 MAPK
171-V60004M	GSK-3β	171-V60010M	p70 S6 Kinase
171-V60005M	HER-2	171-V60011M	Smad2
171-V60006M	IκB-α		
Catalog #	Description	Catalog #	Description

Lysate Controls*

171-YZ0001	EGF-Treated HEK-293 Lysate	171-YZ0008	TNF-α-Treated HeLa Lysate
171-YZ0002	EGF-Treated HeLa Lysate	171-YZ0009	UV-Treated HEK-293 Lysate
171-YZ0003	EGF-Treated SK-Br3 Lysate	171-YZ0010	VEGF-Treated HUVEC Lysate
171-YZ0004	IFN-α-Treated HeLa Lysate	171-YZT001	Untreated HEK-293
171-YZ0005	IGF-1-Treated HEK-293 Lysate	171-YZT002	Untreated HeLa Lysate
171-YZ0006	NGFβ-Treated PC12 Lysate	171-YZB001	Phosphatase HeLa Lysate**
171-YZ0007	PDGF-Treated NIH3T3 Lysate		

* All controls are shipped as one vial of lyophilized cell lysate.

** Negative control.

Catalog #	Description
Reagents	
171-304006M	Bio-Plex Pro Cell Signaling Reagent Kit , 1 x 96-well, includes cell lysis buffer, cell lysis factor QG, cell wash buffer, bead resuspension buffer, detection antibody diluent, wash buffer, streptavidin-PE, 96-well flat bottom plate, sealing tape, and instructions
171-304515	Bio-Plex Pro Signaling Wash Buffer , 330 ml, for 1 x 96-well assay, for use with Bio-Plex Pro cell signaling assays only, compatible with both magnetic and vacuum separation methods

For More Information

Request or download bulletin: 6285

Bio-Plex Assays

Bio-Rad Cell Lysate Controls* — Selection Guide for Use with Bio-Plex Pro Magnetic Cell Signaling Assays

Phosphoprotein	Lysate Control	Catalog #	Phosphoprotein	Lysate Control	Catalog #
Akt (Ser ⁴⁷³)			PDGFR- α (Tyr ⁷⁵⁴)	PDGF-treated NIH3T3	171-YZ0007
Akt (Thr ³⁰⁸)			PDGFR- β (Tyr ⁷⁵¹)		
Erk1/2 (Thr ²⁰² /Tyr ²⁰⁴ , Thr ¹⁸⁵ /Tyr ¹⁸⁷)	EGF-treated HEK-293	171-YZ0001	Stat1 (Tyr ⁷⁰¹)		
GSK-3 α / β (Ser ²¹ /Ser ⁹)			Stat3 (Ser ⁷²⁷)	IFN- α -treated HeLa	171-YZ0004
MEK1 (Ser ²¹⁷ /Ser ²²¹)			Stat3 (Tyr ⁷⁰⁵)		
c-Jun (Ser ⁶³)			VEGFR-2 (Tyr ¹¹⁷⁵)	VEGF-treated HUVEC	171-YZ0010
JNK (Thr ¹⁸³ /Tyr ¹⁸⁵)			Negative control for all phosphoprotein assays	Phosphatase-treated HeLa	171-YZB001
p38 MAPK (Thr ¹⁸⁰ /Tyr ¹⁸²)	UV-treated HEK-293	171-YZ0009	Total Target	Lysate Control	Catalog #
EGFR (Tyr ¹⁰⁶⁸)			All total assays except HER-2 and c-Jun	Untreated HeLa	171-YZT002
EGFR (Tyr ¹¹⁷³)	EGF-treated HeLa	171-YZ0002	Total c-Jun	Untreated HEK-293	171-YZT001
HER-2 (Tyr ¹²⁴⁸)	EGF-treated SK-Br3	171-YZ0003	Total HER-2	EGF-treated SK-Br3	171-YZ0003
IGF-IR (Tyr ¹¹³¹)	IGF-1-treated HEK-293	171-YZ0005	Negative control for all total target assays	Detection antibody diluent**	
I κ B- α (Ser ³² /Ser ³⁶)					
NF- κ B p65 (Ser ⁵³⁶)	TNF- α -treated HeLa	171-YZ0008			
Smad2 (Ser ⁴⁶⁵ /Ser ⁴⁶⁷)					
p70 S6 Kinase (Thr ⁴²¹ /Ser ⁴²⁴)					
p70 S6 Kinase (Thr ³⁸⁹)	NGF β -treated PC-12	171-YZ0006			

* All controls are shipped as one vial of lyophilized cell lysate.

** Included with the Bio-Plex Pro cell signaling reagent kit (171-304006M).

Bio-Plex Nonmagnetic Signal Transduction Assays

Available Nonmagnetic 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)	•	•
ERK1/2 (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)	•	—

Bio-Plex Assays

		
<p>Bio-Plex phosphoprotein and total target assays — include coupled beads, detection antibodies, and lysates to test for specific proteins of interest in lysate samples.</p>	<p>Bio-Plex cell lysis kit* — includes cell lysis buffer, cell wash buffer, factor 1, and factor 2.</p>	<p>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.</p>

* Essential for optimal lysis of your cell culture or tissue samples.

** Required to prepare assays and acquire data.

Catalog #	Description	Catalog #	Description
Phosphoprotein Singleplex Assays*			
171-V21075	Akt (Ser ⁴⁷³)	171-V25340	MEK1 (Ser ²¹⁷ /Ser ²²¹)
171-V21620	ATF-2 (Thr ⁷¹)	171-V24937	NF-κB p65 (Ser ⁵³⁶)
171-V27145	Bcr-Abl (Tyr ²⁴⁵)	171-V21336	p38 MAPK (Thr ¹⁸⁰ /Tyr ¹⁸²)
171-V27545	c-Abl (Tyr ²⁴⁵)	171-V25153	p53 (Ser ¹⁵)
171-V27745	c-Abl (Tyr ⁴¹²)	171-V28353	p53 (Ser ³⁷)
171-V24356	c-Jun (Ser ⁶³)	171-V28153	p53 (Ser ⁴⁶)
171-V26119	CREB (Ser ¹³³)	171-V24155	p70 S6 Kinase (Thr ⁴²¹ /Ser ⁴²⁴)
171-V23120**	EGFR (Tyr)	171-V23535	p90RSK (Thr ³⁵⁹ /Ser ³⁶³)
171-V22238	ERK1/2 (Thr ²⁰² /Tyr ²⁰⁴ , Thr ¹⁸⁵ /Tyr ¹⁸⁷)	171-V25957	PDGF Receptor-β (Tyr ⁵¹)
171-V23318	GSK-3α/β (Ser ²¹ /Ser ⁹)	171-V25374**	S6 Ribosomal Protein (Ser ²³⁵ /Ser ²³⁶)
171-V24777	Histone H3 (Ser ¹⁰)	171-V26741	Src (Tyr ⁴¹⁶)
171-V24551	HSP27 (Ser ⁷⁸)	171-V23732	STAT2 (Tyr ⁶⁸⁹)
171-V27343	IGF-IR (Tyr ¹¹³¹)	171-V26952	STAT3 (Ser ⁷²⁷)
171-V20758	IκB-α (Ser ³² /Ser ³⁶)	171-V22552	STAT3 (Tyr ⁷⁰⁵)
171-V27943	IR-β (Tyr ¹¹⁴⁶)	171-V25772	STAT6 (Tyr ⁶⁴¹)
171-V25576	IRS-1 (Ser ⁶³⁶ /Ser ⁶³⁹)	171-V23973	TrkA (Tyr ⁴⁹⁰)
171-V21034	JNK (Thr ¹⁸³ /Tyr ¹⁸⁵)	171-V26533	Tyk2 (Tyr ¹⁰⁵⁴ /Tyr ¹⁰⁵⁵)

* Requires cell lysis kit and phosphoprotein detection reagent kit.

** This assay cannot be multiplexed.

Catalog #	Description	Catalog #	Description
Total Target Singleplex Assays*			
171-V31075	Akt	171-V30758	IκB-α
171-V31620	ATF-2	171-V31034	JNK
171-V37145	Bcr-Abl	171-V35340	MEK1
171-V34356	c-Jun	171-V31336	p38 MAPK
171-V32238	ERK1/2	171-V35153	p53
171-V34551	HSP27	171-V33535	p90RSK

* Requires cell lysis kit and phosphoprotein detection reagent kit.

Catalog #	Description
171-304004	Bio-Plex Phosphoprotein Detection Reagent Kit , 1 x 96-well, includes assay buffer, wash buffer, antibody diluent, resuspension buffer, streptavidin-PE, filter plate, sealing tape
171-304005	Bio-Plex Phosphoprotein Detection Reagent Kit , 10 x 96-well, includes assay buffer, wash buffer, antibody diluent, resuspension buffer, streptavidin-PE, filter plate, sealing tape

Sample Preparation Kits

171-304011	Bio-Plex Cell Lysis Kit , 1 x 96-well, includes cell lysis and wash buffers, factor 1 and factor 2
171-304012	Bio-Plex Cell Lysis Kit , 10 x 96-well, includes cell lysis and wash buffers, factor 1 and factor 2

Note: Customized Bio-Plex assays can be ordered with any combination of available analytes. Request bulletin 5483 for multiplexing capabilities of specific assay combinations. For more information, go to www.bio-rad.com/bio-plex/assaybuilder.

For More Information

Web: www.bio-rad.com/phosphoprotein

Request or download bulletins: 5405 and 5483

Bio-Plex Assays

Kidney Toxicity Assays New

Available Bio-Plex Pro RBM Kidney Toxicity Kits

Human Kidney Toxicity Panel 1	Human Kidney Toxicity Panel 2
Calbindin	Albumin
Clusterin	B2M
GST- π	Cystatin C
IL-18	NGAL
KIM-1	Osteopontin
MCP-1	TFF3



MYRIAD RBM
Manufactured by Myriad RBM®

Rat Kidney Toxicity Albumin Kit	Rat Kidney Toxicity Panel 1	Rat Kidney Toxicity Panel 2
Albumin	Clusterin	B2M
	IL-18	Calbindin
	KIM-1	Cystatin C
	MCP-1	NGAL
	Osteopontin	
Canine Kidney Toxicity Albumin Kit	Canine Kidney Toxicity Panel 1	
Albumin	Clusterin	
	KIM-1	
	MCP-1	
	NGAL	

Human Kidney Toxicity

Catalog #	Description
Premixed All-in-One Kit	
171-ATR1CK	Bio-Plex Pro RBM Human Kidney Toxicity Panel 1 , 1 x 96-well, includes premixed magnetic capture beads and detection antibodies, standards, 2-level controls, standard diluent, buffers (blocking, sample dilution, and 10x assay), 10x streptavidin-PE, flat bottom plate/seals, and instructions for the detection of the following analytes: calbindin, clusterin, GST- π , IL-18, KIM-1, MCP-1
171-ATR2CK	Bio-Plex Pro RBM Human Kidney Toxicity Panel 2 , 1 x 96-well, includes premixed magnetic capture beads and detection antibodies, standards, 2-level controls, standard diluent, buffers (blocking, sample dilution, and 10x assay), 10x streptavidin-PE, flat bottom plate/seals, and instructions for the detection of the following analytes: albumin, B2M, cystatin C, NGAL, osteopontin, TFF3

Rat Kidney Toxicity

Catalog #	Description
Premixed All-in-One Kit	
171-KTR1CK	Bio-Plex Pro RBM Rat Kidney Toxicity Panel 1 , 1 x 96-well, includes premixed magnetic capture beads and detection antibodies, standards, 2-level controls, standard diluent, buffers (blocking, sample dilution, and 10x assay), 10x streptavidin-PE, flat bottom plate/seals, and instructions for the detection of the following analytes: clusterin, IL-18, KIM-1, MCP-1, osteopontin
171-KTR2CK	Bio-Plex Pro RBM Rat Kidney Toxicity Panel 2 , 1 x 96-well, includes premixed magnetic capture beads and detection antibodies, standards, 2-level controls, standard diluent, buffers (blocking, sample dilution, and 10x assay), 10x streptavidin-PE, flat bottom plate/seals, and instructions for the detection of the following analytes: B2M, calbindin, cystatin C, NGAL
171-KTR3CK	Bio-Plex Pro RBM Rat Kidney Toxicity Albumin Kit , 1 x 96-well, includes magnetic capture beads, detection antibodies, standards, controls, diluent, buffers, streptavidin-PE, flat bottom plate, and plate seals, for the detection of albumin

Canine Kidney Toxicity

Catalog #	Description
Premixed All-in-One Kit	
171-QTR1CK	Bio-Plex Pro RBM Canine Kidney Toxicity Panel 1 , 1 x 96-well, includes premixed magnetic capture beads and detection antibodies, standards, 2-level controls, standard diluent, buffers (blocking, sample dilution, and 10x assay), 10x streptavidin-PE, flat bottom plate/seals, and instructions for the detection of the following analytes: clusterin, KIM-1, MCP-1, NGAL
171-QTR2CK	Bio-Plex Pro RBM Canine Kidney Toxicity Albumin Kit , 1 x 96-well, includes premixed magnetic capture beads and detection antibodies, standards, 2-level controls, standard diluent, buffers (blocking, sample dilution, and 10x assay), 10x streptavidin-PE, flat bottom plate/seals, and instructions for the detection of albumin

Bio-Plex Assays

Disease Research Assays



Premixed All-In-One Kit

Singleplex Set

Standard

Reagent Kit

Diluent Kit

Bio-Plex Pro Human Cancer Biomarker Assays

Available Bio-Plex Pro Human Cancer Biomarker Panel 1 Assays — receptors, cytokines, chemokines, growth factors, and hormones

FGF-basic	PDGF-AB/BB	sHER-2/neu
Follistatin	PECAM-1	sIL-6R α
G-CSF	Prolactin	sTIE-2
HGF	SCF	sVEGFR-1
Leptin	sEGFR	sVEGFR-2
Osteopontin		

New

Available Bio-Plex Pro Human Cancer Biomarker Panel 2 Assays — ligands for the receptors in panel 1, cytokines, chemokines, and growth factors

Angiopoietin-2	IGFBP-1	TGF- α
sCD40L	IL-6	TNF- α
EGF	IL-8	uPA
Endoglin	IL-18	VEGF-A
sFASL	PAI-1	VEGF-C
HB-EGF	PLGF	VEGF-D

Catalog # Description

Premixed All-in-One Kit

171-AC500M	Bio-Plex Pro Human Cancer Biomarker Panel 1 16-Plex , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standard, 2-level controls, detection antibody diluent, standard diluent HB, sample diluent HB, assay buffer, wash buffer, streptavidin-PE, flat bottom plate, sealing tape, and instructions for the detection of sEGFR, FGF-Basic, follistatin, G-CSF, HGF, sHER-2/neu, sIL-6R α , leptin, osteopontin, PECAM-1, PDGF-AB/BB, prolactin, SCF, sTIE-2, sVEGFR-1, and sVEGFR-2
171-AC600M	Bio-Plex Pro Human Cancer Biomarker Panel 2, 18-Plex , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standard, 2-level controls, detection antibody diluent, standard diluent HB, sample diluent HB, assay buffer, wash buffer, streptavidin-PE, flat bottom plate, sealing tape, and instructions for the detection of angiopoietin-2, sCD40L, EGF, endoglin, sFASL, HB-EGF, IGFBP-1, IL-6, IL-8, IL-18, PAI-1, PLGF, TGF- α , TNF- α , uPA, VEGF-A, VEGF-C, VEGF-D

Catalog # Description

Human Cancer Biomarker Panel 1 Singleplex Sets*, 1 x 96-Well**

171-BC501M	sEGFR Set	171-BC509M	Osteopontin Set
171-BC502M	FGF-basic Set	171-BC510M	PECAM-1 Set
171-BC503M	Follistatin Set	171-BC511M	PDGF-AB/BB Set
171-BC504M	G-CSF Set	171-BC512M	Prolactin Set
171-BC505M	HGF Set	171-BC513M	SCF Set
171-BC506M	sHER-2/neu Set	171-BC514M	sTIE-2 Set
171-BC507M	sIL-6Rα Set	171-BC515M	sVEGFR-1 Set
171-BC508M	Leptin Set	171-BC516M	sVEGFR-2 Set

* Requires reagent kit II and human cancer biomarker panel 1 standards.

Catalog # Description

Bio-Plex Pro Human Cancer Biomarker Panel 2 Singleplex Sets**

1 x 96-well, include coupled magnetic beads and detection antibodies. Require Bio-Plex Pro reagent kit II and panel 2 standards (171-DC6000).

171-BC601M	Angiopoietin-2 Set	171-BC611M	PAI-1 Set
171-BC602M	sCD40L Set	171-BC612M	PLGF Set
171-BC603M	EGF Set	171-BC613M	TGF-α Set
171-BC604M	Endoglin Set	171-BC614M	TNF-α Set
171-BC605M	sFASL Set	171-BC615M	uPA Set
171-BC606M	HB-EGF Set	171-BC616M	VEGF-A Set
171-BC607M	IGFBP-1 Set	171-BC617M	VEGF-C Set
171-BC608M	IL-6 Set	171-BC618M	VEGF-D Set

Bio-Plex Assays

171-BC609M **IL-8 Set**

** It is not recommended to mix assays between panel 1 and panel 2.

Catalog # Description

Standards

171-DC5000 **Bio-Plex Pro Human Cancer Biomarker Panel 1 16-Plex Standards**, pkg of 1 vial, lyophilized mixture of 16 standard analytes
 171-DC5001 **Bio-Plex Pro Human Cancer Biomarker Panel 1 16-Plex Standards**, pkg of 50 lot-matched vials, lyophilized mixture of 16 standard analytes
 171-DC6000 **Bio-Plex Pro Human Cancer Biomarker Panel 2, 18-plex Standards**, pkg of 1 vial, lyophilized mixture of 18 standard analytes
 171-DC6001 **Bio-Plex Pro Human Cancer Biomarker Panel 2, 18-plex Standards**, pkg of 50 lot-matched vials, lyophilized mixture of 18 standard analytes

Reagent Kits

171-304055 **Bio-Plex Pro Reagent Kit II with Filter Plate**, 1 x 96-well, includes detection antibody diluent, standard diluent HB, sample diluent HB, assay buffer, wash buffer, streptavidin-PE, 96-well filter plate, sealing tape, and instructions for use with vacuum separation methods
 171304055M **Bio-Plex Pro Reagent Kit II with Flat Bottom Plate**, 1 x 96-well, includes detection antibody diluent, standard diluent HB, sample diluent HB, assay buffer, wash buffer, streptavidin-PE, 96-well flat bottom plate, sealing tape, and instructions for use with magnetic separation methods

For More Information

Bio-Plex Pro Diabetes Assays



Premixed All-In-One Kit

New Available Bio-Plex Pro Diabetes Assays

Human	Non-Human Primate	Mouse	Rat
Adiponectin	Adiponectin	Adiponectin	Ghrelin
Adipsin	Adipsin	Ghrelin	GLP-1
C-peptide	C-peptide	GIP	Glucagon
Ghrelin	Ghrelin	GLP-1	Leptin
GIP	GIP	Glucagon	PAI-1
GLP-1	GLP-1	Insulin	
Glucagon	Glucagon	Leptin	
Insulin	Insulin	PAI-1	
Leptin	Leptin	Resistin	
PAI-1	PAI-1		
Resistin	Resistin		
Visfatin	Visfatin		

Web: www.bio-rad.com/bio-plexprocancer

Request or download bulletins: 6156 and 6159

Human Diabetes

Catalog # Description

Premixed All-in-One Kits

171-A7001M **Bio-Plex Pro Human Diabetes 10-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 C-peptide, ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, resistin, visfatin
 171-A7002M **Bio-Plex Pro Human Diabetes Adipsin and Adiponectin Assays**, 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, serum-based diluent for the detection of adipsin and adiponectin
 171-A7003M **Bio-Plex Pro Human Diabetes Adiponectin Assay**, 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, serum-based diluent for the detection of adiponectin
 171-A7004M **Bio-Plex Pro Human Diabetes Adipsin Assay**, 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape, serum-based diluent for the detection of adipsin

Catalog # Description

Catalog # Description

Human Diabetes Singleplex Sets*, 1 x 96-Well

171-B7003M C-Peptide Set	171-B7008M Insulin Set
171-B7004M Ghrelin Set	171-B7009M Leptin Set
171-B7005M GIP Set	171-B7010M PAI-1 Set
171-B7006M GLP-1 Set	171-B7011M Resistin Set
171-B7007M Glucagon Set	171-B7012M Visfatin Set

* Requires reagent kit and human diabetes standards.

Catalog # Description

Standards

171-D70001 **Bio-Plex Pro Human Diabetes Standards**, pkg of 1 vial, lyophilized mixture of 12 standard analytes, including adiponectin, adipsin, C-peptide, ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, resistin, and visfatin
 171-D70050 **Bio-Plex Pro Human Diabetes Standards**, pkg of 50 lot-matched vials, lyophilized mixture of 12 standard analytes, including adiponectin, adipsin, C-peptide, ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, resistin, and visfatin

Bio-Plex Assays

Non-Human Primate Diabetes New

Catalog #	Description
-----------	-------------

Premixed All-in-One Kits

171-W7001M	Bio-Plex Pro Non-Human Primate Diabetes 11-Plex Panel , 1 x 96-well, includes premixed coupled magnetic beads, premixed detection antibodies, standard, detection antibody diluent, standard diluent, sample diluent, assay buffer, wash buffer, streptavidin-PE, a 96-well flat bottom plate, sealing tape, and quick guide for the detection of adipsin, C-peptide, ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, resistin, and visfatin
171-W7002M	Bio-Plex Pro Non-Human Primate Diabetes Adiponectin Kit , 1 x 96-well, includes premixed coupled magnetic beads, detection antibodies, standard, reagents, flat bottom plate, and sealing tape for the detection of adiponectin

Catalog #	Description	Catalog #	Description
-----------	-------------	-----------	-------------

Bio-Plex Pro Non-Human Primate Diabetes Singleplex Sets

171-W7003M	Adipsin	171-W7009M	Insulin
171-W7004M	C-Peptide	171-W7010M	Leptin
171-W7005M	Ghrelin	171-W7011M	PAI-1
171-W7006M	GIP	171-W7012M	Resistin
171-W7007M	GLP-1	171-W7013M	Visfatin
171-W7008M	Glucagon		

Bio-Plex Pro Non-Human Primate Diabetes Standards

Standards are for the detection of adiponectin, adipsin, C-peptide, ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, resistin, visfatin

171-D70001	Bio-Plex Pro Diabetes Standards , Pkg of 1 vial, lyophilized mixture of 11 analytes for use with Bio-Plex Pro non-human primate diabetes assays
------------	--

Mouse Diabetes

Catalog #	Description
-----------	-------------

Premixed All-in-One Kits

171-F7001M	Bio-Plex Pro Mouse Diabetes 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, flat bottom plate, sealing tape, standard diluent, sample diluent for the detection of ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, resistin
171-F7002M	Bio-Plex Pro Mouse Diabetes Adiponectin Assay , 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, flat bottom plate, sealing tape, serum-based diluent for the detection of adiponectin

Catalog #	Description	Catalog #	Description
-----------	-------------	-----------	-------------

Mouse Diabetes Singleplex Sets*, 1 x 96-Well

171-G7002M	Ghrelin Set	171-G7006M	Insulin Set
171-G7003M	GIP Set	171-G7007M	Leptin Set
171-G7004M	GLP-1 Set	171-G7008M	PAI-1 Set
171-G7005M	Glucagon Set	171-G7009M	Resistin Set

* Requires reagent kit and mouse diabetes standards. Due to the different dilution scheme, adiponectin is offered in an all-in-one format only.

Catalog #	Description
-----------	-------------

Standards

171-I70001	Bio-Plex Pro Mouse Diabetes Standards , pkg of 1 vial, lyophilized mixture of 9 standard analytes, including adiponectin, ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, and resistin
171-I70050	Bio-Plex Pro Mouse Diabetes Standards , pkg of 50 lot-matched vials, lyophilized mixture of 9 standard analytes, including adiponectin, ghrelin, GIP, GLP-1, glucagon, insulin, leptin, PAI-1, and resistin

Rat Diabetes

Catalog #	Description
-----------	-------------

Rat Diabetes Singleplex Sets*, 1 x 96-Well

171-L7001M	Ghrelin Set
171-L7003M	GLP-1 Set
171-L7004M	Glucagon Set
171-L7006M	Leptin Set
171-L7007M	PAI-1 Set

* Requires reagent kit and rat standards.

Standards

171-NZ0001	Bio-Plex Pro Rat Standards , pkg of 1 vial, lyophilized mixture of 31 standard analytes. Compatible with rat cytokine and rat diabetes assays
171-NZ0501	Bio-Plex Pro Rat Standards , pkg of 50 lot-matched vials, lyophilized mixture of 31 standard analytes. Compatible with rat cytokine and rat diabetes assays

Reagent Kits* (for All Species)

Catalog #	Description
-----------	-------------

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
171-304070M	Bio-Plex Pro Reagent Kit with Flat Bottom Plate , 1 x 96-well, includes flat bottom plate, assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, sealing tape, standard diluent, sample diluent, and instructions

* Same reagent kit can be used for both diabetes and cytokine assays.

For More Information

Web: www.bio-rad.com/bio-plexprodiabetes

Request or download bulletins: 6342 and 6119

Bio-Plex Assays

Bio-Plex Pro Human Acute Phase Assay Panel

Available Bio-Plex Pro Human Acute Phase Assays

α -2-macroglobulin	Procalcitonin
CRP	SAA
Ferritin	SAP
Fibrinogen	Tissue plasminogen activator
Haptoglobin	

Catalog # Description

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

Assay Kits**

171-A4007M	Bio-Plex Pro Human Acute Phase 5-Plex Panel , 1 x 96-well, includes premixed coupled magnetic beads, detection antibodies, standards, and controls for the detection of ferritin, fibrinogen, procalcitonin, serum amyloid A, tissue plasminogen activator, requires acute phase reagent kit and acute phase diluent kit
171-A4009M	Bio-Plex Pro Human Acute Phase 4-Plex Panel , 1 x 96-well, includes premixed coupled magnetic beads, detection antibodies, standards, and controls for the detection of α -2-macroglobulin, CRP, haptoglobin, serum amyloid P, requires acute phase reagent kit and acute phase diluent kit

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

** Requires acute phase reagent kit and acute phase diluent kit.

Standards

171-D40002	Bio-Plex Pro Human Acute Phase Standards , pkg of 2 vials, lyophilized mixture of 9 standard analytes
171-D40006	Bio-Plex Pro Human Acute Phase Standards , pkg of 50 lot-matched vials, lyophilized mixture of 9 standard analytes

Reagent Kit

171-304050	Bio-Plex Pro Human Acute Phase Reagent Kit , 1 x 96-well, includes assay buffer, wash buffer, detection antibody diluent, streptavidin-PE, filter plate, sealing tape
------------	--

Diluent Kit*

171-305050	Bio-Plex Pro Human Acute Phase Diluent Kit , 1 x 96-well, includes 10 ml serum-based diluent and 300 ml serum-free diluent
------------	---

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

For More Information

Web: www.bio-rad.com/bio-plexproacutephase

Request or download bulletin: 5650

Bio-Plex Pro Human Angiogenesis Assay Panel

Available Bio-Plex Pro Human Angiogenesis Assay Panel

Angiopoietin-2	HGF	PDGF-BB
Follistatin	IL-8	PECAM-1
G-CSF	Leptin	VEGF

Catalog # Description

Complete Kit*

171-A4S11M	Bio-Plex Pro Human Angiogenesis 9-Plex Panel Complete Kit , 1 x 96-well
------------	--

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

Assay Kit

171-A4011M	Bio-Plex Pro Human Angiogenesis 9-Plex Panel , 1 x 96-well, includes premixed coupled magnetic beads, detection antibodies, standards, and controls for the detection of angiopoietin-2, follistatin, G-CSF, HGF, IL-8, leptin, PDGF-BB, PECAM-1, and VEGF; requires angiogenesis reagent kit and human serum diluent kit
------------	--

Standards

171-D40003	Bio-Plex Pro Human Angiogenesis Standards , pkg of 2 vials, lyophilized mixture of 9 standard analytes
171-D40004	Bio-Plex Pro Human Angiogenesis Standards , pkg of 50 lot-matched vials, lyophilized mixture of 9 standard analytes

Reagent Kit

171-304060	Bio-Plex Pro Angiogenesis Reagent Kit , 1 x 96-well
------------	--

Diluent Kit

171-305000	Bio-Plex Human Serum Diluent Kit , 1 x 96-well, includes 15 ml sample diluent, 10 ml standard diluent
------------	--

For More Information

Web: www.bio-rad.com/bio-plexproangiogenesis

Request or download bulletins: 5653

Bio-Plex Assays

Bio-Plex Pro Human Isotyping Assays

New

Available Bio-Plex Pro Human Isotyping Assay Panel

IgG ₁	IgG ₄	IgM
IgG ₂	IgA	IgE
IgG ₃		IgG Total



Premixed All-In-One Kit

Catalog #	Description
-----------	-------------

Bio-Plex Pro Human Isotyping Panel Premixed All-In-One Multiplex Kit

171-A3100M **Bio-Plex Pro Human Isotyping Panel, 6-plex**, 1 x 96-well, includes coupled magnetic beads, detection antibodies, standards, 1-level quality controls, detection antibody diluent, isotyping diluent (for use with samples, standards, and controls), assay buffer, wash buffer, streptavidin-PE, 96-well flat bottom plate, sealing tape, assay quick guide, and product data sheet; for the detection of IgG₁, IgG₂, IgG₃, IgG₄, IgA, and IgM.

Bio-Plex Pro Human Isotyping All-In-One Singleplex Kits

Singleplex kits are specific for each assay indicated below and include coupled magnetic beads, detection antibodies, standards, 1-level quality controls, detection antibody diluent, isotyping diluent (for use with samples, standards, and controls), assay buffer, wash buffer, streptavidin-PE, 96-well flat bottom plate, sealing tape, assay quick guide, and product data sheet.

171-A3101M	IgA Kit , 1 x 96-well
171-A3104M	IgM Kit , 1 x 96-well
171-A3102M	IgE Kit , 1 x 96-well
171-A3103M	IgG Total Kit , 1 x 96-well

Wash Stations and Accessories

300-34376	Bio-Plex Pro Wash Station , includes magnetic plate carrier, waste bottle, 2 buffer bottles
300-34377	Bio-Plex Pro II Wash Station , includes magnetic plate carrier, vacuum manifold plate carrier, waste bottle, 2 buffer bottles
171-025001	Bio-Plex Flat Bottom Plates , 40 x 96-well plates
171-304500	Bio-Plex Wash Buffer , 1.5 L
171-020100	Bio-Plex Handheld Magnetic Washer , includes magnetic washer and adjustment hex tools for use in manual wash steps for all Bio-Plex magnetic assays
171-304502	Filter Plate , pkg of 1, 96-well plate, with clear plastic lid and tray, for Bio-Plex assays using the vacuum wash method, sealing tape not included

Software

171-001510	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
171-001513	Bio-Plex Data Pro Software , (5 seats), for multi-experiment analysis and advanced data visualization
171-001523	Bio-Plex Data Pro Plus Software , contains all the features of Bio-Plex Data Pro Software with added visualization, sharing, and analysis functionality
171-STND01	Bio-Plex Manager Software , includes 1 user desktop license, to analyze Bio-Plex data and generate protocols, does not operate the instrument

For More Information

Web: www.bio-rad.com/bio-plexproisotyping

Request or download bulletin: 6344

Bio-Plex® Precision Pro™ Human Cytokine Assay Panel



Premixed All-In-One Kit

Available Bio-Plex Precision Pro Human Cytokine Assay Panel

IL-1 β	IL-6	IL-13
IL-2	IL-10	IFN- γ
IL-4	IL-12p70	TNF- α
IL-5		

Catalog #	Description
-----------	-------------

171-A1001P **Bio-Plex Precision Pro Human Cytokine 10-Plex Panel**, 1 x 96-well, includes premixed 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-12p70, IL-13, IFN- γ , TNF- α

For More Information

Web: www.bio-rad.com/bio-plexprecisionpro

Request or download bulletin: 5629

Beads, Reagents and Accessories

COOH Beads and Related Reagents

Bio-Plex Pro™ COOH Magnetic Beads



Catalog #	Description
MC100xx-01*	Bio-Plex Pro Magnetic COOH Beads , 1 ml, 1.25×10^7 beads/ml
MC100xx-04*	Bio-Plex Pro Magnetic COOH Beads , 4 ml, 1.25×10^7 beads/ml

* xx = bead region. Bead regions available, 26–29, 34–37, 43–46, 52–55, 62–65.

For More Information

Web: www.bio-rad.com/COOH

Bio-Plex® COOH Nonmagnetic Beads



Catalog #	Description
171-5060xx*	Bio-Plex COOH Beads , 1 ml, 1.25×10^7 beads/ml
171-6060xx*	Bio-Plex COOH Beads , 16 ml, 1.25×10^7 beads/ml

* xx = bead region. Bead regions available, 11, 16, 18, 20, 24–28, 30, 31, 33, 35, 38, 42–46, 50–53, 56, 66.

For More Information

Web: www.bio-rad.com/COOH

Bio-Plex Amine Coupling Kit



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

Bio-Plex Streptavidin-PE*



Catalog #	Description
Bio-Plex Streptavidin-PE*	
171-304501	Bio-Plex Streptavidin-PE , streptavidin-phycoerythrin for 10 x 96-well reactions

* Streptavidin-PE is a component of all reagent kits and is available separately.

For More Information

Web: www.bio-rad.com/COOH

Accessories

Bio-Plex Pro Assay Accessories

Catalog #	Description
171-304500	Bio-Plex Wash Buffer , 1.5 L
171-025001*	Bio-Plex Pro Flat Bottom Plates , 40 x 96-well plates
171-304502	Filter Plate , 1 x 96-well filter plate with clear plastic lid and tray. Sealing tape not included

* Required for washing Bio-Plex Pro assays or other magnetic bead-based assays using Bio-Plex Pro and Bio-Plex Pro II wash stations.

xPONENT is a trademark of the Luminex Corporation.

The Bio-Plex suspension array system includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the 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 11 5044 5699
Canada 905 364 3435 **China** 86 21 6169 8500 **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 9532 220 **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 64 9 415 2280 **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** 026 674 55 05
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000