

Bio-Plex™ Human Cytokine
Singleplex Assay
(1 x 96-well)
Product Insert

Requires manual and reagents in
Bio-Plex Cytokine Reagent Kit (1x96)
Catalog # 171-304000

For research use only.
Not for diagnostic procedures.

BIO-RAD

- ☐ This kit includes 2 vials of Human Cytokine Standard containing the following premixed cytokines:

IL-1 β	IL-8	Eotaxin	MIP-1 α
IL-1ra	IL-9	FGF basic	MIP-1 β
IL-2	IL-10	G-CSF	PDGF bb
IL-4	IL-12 (p70)	GM-CSF	RANTES
IL-5	IL-13	IFN- γ	TNF- α
IL-6	IL-15	IP-10	VEGF
IL-7	IL-17	MCP-1 (MCAF)	

Note: To increase 'plexes', refer to the Cytokine Assay Instruction Manual.

 Singleplex Assay Product list

Human Cytokine	Bead Region	Catalog #
IL-1 β	32	171-B12832
IL-1ra	25	x-plex only
IL-2	38	171-B10138
IL-4	52	171-B10452
IL-5	33	171-B12533
IL-6	19	171-B10719
IL-7	74	171-B13174
IL-8	54	171-B11054
IL-9	77	x-plex only
IL-10	56	171-B11356
IL-12 (p70)	75	171-B13475
IL-13	51	171-B13751
IL-15	73	x-plex only
IL-17	76	171-B14076

Human Cytokine	Bead Region	Catalog #
Eotaxin	43	x-plex only
FGF basic	44	x-plex only
G-CSF	57	171-B14357
GM-CSF	34	171-B11634
IFN- γ	21	171-B11921
IP-10	41	x-plex only
MCP-1 (MCAF)	53	171-B14653
MIP-1 α	55	x-plex only
MIP-1 β	18	171-B14918
PDGF bb	47	x-plex only
RANTES	37	x-plex only
TNF- α	36	171-B12236
VEGF	45	x-plex only

For x-plex assays visit www.bio-rad.com/bio-plex/x-plex

The concentration of the singleplex detection antibody is listed on the vial.

Product Specifications:

Storage 2 to 8°C

Stability 6 to 12 months

Instructions Please refer to the Cytokine Assay Instruction Manual in Catalog #171-304000

For Technical Specifications refer to bulletin 3157.

Sample Matrix	Diluent Kit Catalog #
Serum	171-305000

For technical service, call your local Bio-Rad office, or in the US, call 1-800-4BIORAD (1-800-424-6723)

4110005

Rev G