

Bio-Plex™ Human Cytokine
17-Plex Panel
(1 x 96-well)
Product Insert

Requires the 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.

- The Human Cytokine 17-Plex Panel contains the following beads and detection antibodies:

Target	Bead Region
IL-1 β	32
IL-2	38
IL-4	52
IL-5	33
IL-6	19
IL-7	74
IL-8	54
IL-10	56
IL-12 (p70)	75
IL-13	51

Target	Bead Region
IL-17	76
G-CSF	57
GM-CSF	34
IFN- γ	21
MCP-1 (MCAF)	53
MIP-1 β	18
TNF- α	36

The concentration of Human Cytokine 17-Plex Panel 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)

4110059

Rev D