

Bio-Plex™ Mouse Cytokine
18-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 Mouse Cytokine Standard (25 ng/cytokine), containing the following 18 premixed cytokines:

IL-1 α	IL-12 (p70)
IL-1 β	IL-17
IL-2	G-CSF
IL-3	GM-CSF
IL-4	IFN- γ
IL-5	KC
IL-6	MIP-1 α
IL-10	RANTES
IL-12 (p40)	TNF- α

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

- The Mouse Cytokine 18-Plex Panel contains the following beads and detection antibodies:

Target	Bead Region
IL-1 α	53
IL-1 β	19
IL-2	36
IL-3	18
IL-4	32
IL-5	52
IL-6	38
IL-10	56
IL-12 (p40)	76

Target	Bead Region
IL-12 (p70)	58
IL-17	72
G-CSF	54
GM-CSF	73
IFN- γ	34
KC	57
MIP-1 α	77
RANTES	55
TNF- α	21

The concentration of Mouse Cytokine 18-Plex Panel detection antibody is 25x. Please check the concentration listed on the vial before mixing with other assay kits.

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

Sensitivity <10 pg/ml
(Based on detectable signal higher than 2 SD above background)

Precision
Intra- and Interassay CV<10%

Assay range 1–32,000 pg/ml
(Based on 4 or 5 PL curve fit)

% Recovery

Serum	80–120%
Plasma	70–130%
Cell Culture	90–110%

Cross-reactivity Negligible

Sample Matrix	Diluent Kit Catalog #
Serum	171-305004

For Technical Service
Call Your Local Bio-Rad Office or
In the U.S. call 1-800-424-6723