

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

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

For use with Bio-Plex reagent kit A
Catalog #171-304000

Serum samples also require
Bio-Plex diluent kit B
Catalog #171-305004

For research use only.
Not for diagnostic procedures.

4110010 Rev B 0101



Mouse Cytokine	Bead Region	Catalog #
IL-1 β	19	171-G12819
IL-2	36	171-G10136
IL-4	32	171-G10432
IL-5	52	171-G12552
IL-10	56	171-G11356
GM-CSF	73	171-G11673
IFN- γ	34	171-G11934
TNF- α	21	171-G12221
8-Plex B	all above premixed	171-F11080

Storage 2–8°C
Stability 6 months
Instructions See Bio-Plex reagent kit manual

Limit of detection <10 pg/ml
(based on detectable signal higher
than 2 SD above background)

Intra-assay CV <10%

Inter-assay CV <10%

Assay range 10–32,000 pg/ml
(based on 4 PL curve fit)

% Recovery 80–120%

Cross-reactivity <10%

Matrices Serum and tissue
culture media

Multiplex
compatibility Mouse IL-1 β , IL-2,
IL-4, IL-5, IL-10,
TNF- α , IFN- γ , GM-CSF