

Bio-Plex™ Rat
Cytokine 9-Plex A Panel
(1 x 96-Well)
Product Insert

Requires the use of
Bio-Plex Cytokine Reagent Kit
Catalog #171-304000

For research use only.
Not for diagnostic procedures.

BIO-RAD

Rat Cytokine	Bead Region
IL-1 α	21
IL-1 β	35
IL-2	18
IL-4	32
IL-6	55
IL-10	20
GM-CSF	37
IFN- γ	34
TNF- α	72

Product Specifications:

Storage

2 to 8°C

Stability	6 to 12 months
Instructions	See Bio-Plex reagent kit manual
Sensitivity (based on detectable signal higher than 2 SD above background)	<10 pg/ml
Precision	Intra- and Interassay CV<10%
Assay range (based on 4 or 5 PL curve fit)	1–32,000 pg/ml
% Recovery	
Serum	80–120%
Plasma	70–130%
Cell culture	90–110%
Cross-reactivity	Negligible

Sample Matrix	Diluent Kit Catalog #
Serum	171-305008

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