

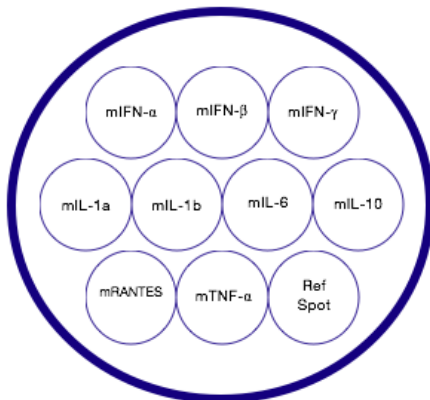
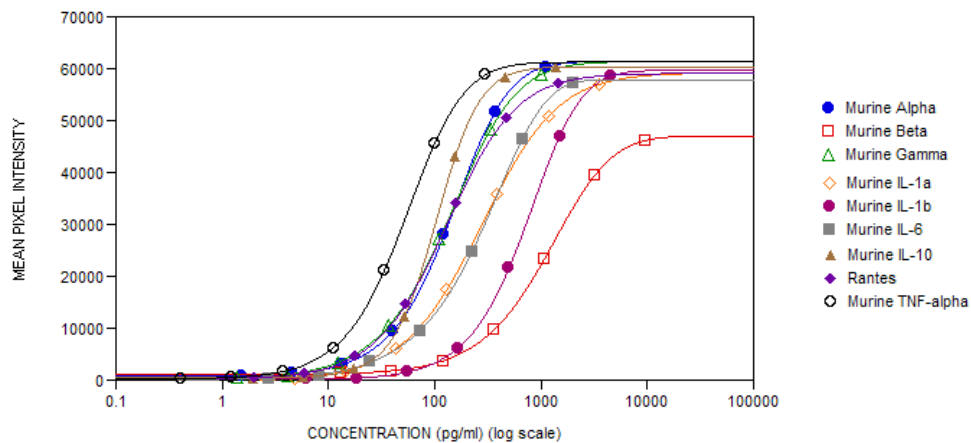


VeriPlex Mouse Cytokine 9-Plex ELISA

Catalog No. 52500

- Simultaneous quantification of Type I and Type II Mouse IFNs and cytokines:
 - IFN- α , IFN- β , IFN- γ , IL-1 α , IL-1 β , IL-6, IL-10, RANTES and TNF- α
- Sample volume of only 50 μ l needed to measure all 9 analytes in a single assay
- Compatible with sera, plasma, tissue culture media (TCM) and buffers
- No cross-reactivity detected with Human IFN- α 2a, IFN- β 1a, IFN- γ , IFN- λ 1, 2 or 3; Rat IFN- α , IFN- β

Figure 1. Representative Analyte Standard Curves and Spot Location in each well



The VeriPlex Mouse Cytokine 9-Plex ELISA is a multiplex immunoassay with the functional format of a sandwich ELISA with a chemiluminescent output. It has been developed to simultaneously detect key cytokines released upstream and downstream of interferon signaling using the Q-Plex™ array spotting technology in which capture antibodies to the specific analytes are spotted in a single well in a defined array. Each *spot* represents a distinct capture antibody population. It is compatible with multiple matrices including tissue culture media, mouse serum, mouse plasma and buffers.

This kit provides a cost-effective option to profile the biological complexity of interferon-related events by simultaneously quantifying several cytokines within a single sample. The ability to assess multiple analytes in one assay means less cost per data point and the ability to generate unique cytokine profiles or “fingerprints” among different patient samples or disease states.

Figure 2. Lower Limit of Detection (LLOD) and Lower Limit of Quantitation (LLOQ)

Analyte	Curve Range (pg/ml)	LLOD (pg/ml)	LLOQ (pg/ml)
Mouse IFN- α	1100 to 1.5	0.5	2.1
Mouse IFN- β	9650 to 13.2	6.5	19.9
Mouse IFN- γ	1000 to 1.4	0.6	1.4
Mouse IL-1 α	3500 to 4.8	1.0	4.8
Mouse IL-1 β	4500 to 6.2	11.0	20.4
Mouse IL-6	2000 to 2.7	1.0	3.7
Mouse IL-10	1400 to 1.9	1.4	4.7
Mouse RANTES	1450 to 2.0	0.8	3.4
Mouse TNF- α	300 to 0.4	0.4	0.5

*LLOD and LLOQ values may vary from lot to lot.



PBL recommends using the Quansys Q-View™ imager in conjunction with the latest version of Q-View™ software for optimal image analysis.

Product Information*

Catalog No.	52500-1
Description	VeriPlex Mouse Cytokine 9-Plex ELISA
Size	1 x 96-well plate
Compatibility	Serum, Plasma, Cell Culture Supernatant & Buffers
Assay Range	2.86 log with LLOD & LLOQ in the low pg/ml range

* Powered by Quansys Biosciences Q-Plex™ Technology **QUANSYS** BIOSCIENCES

Q-Plex, Q-View, and Quansys trademark under Quansys Biosciences, Utah, USA.