

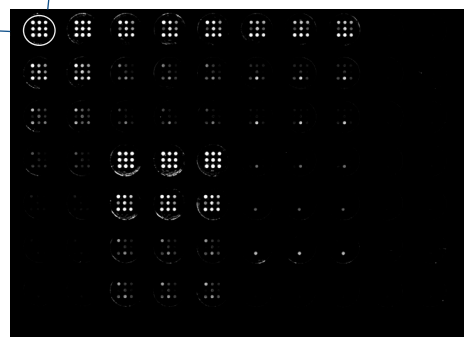
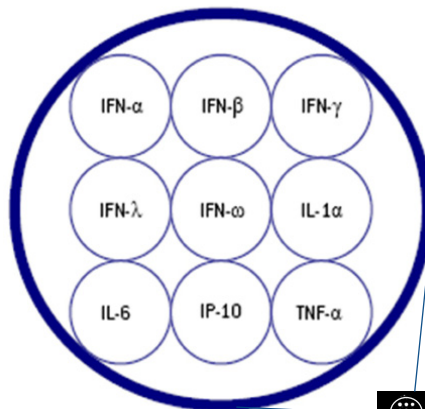


VeriPlex

Human Interferon 9-Plex ELISA

Catalog No. 51500

- Simultaneous measurement of various Type I, II and III Human IFNs and related pro-inflammatory cytokines:
 - IFN- α , IFN- β , IFN- γ , IFN- λ , IFN- ω , IL-1 α , IL-6, IP-10 and TNF- α
- Quantify all 9 analytes in a single well using a sample volume of only 50 μ l
- Compatible with autoimmune sera, plasma, tissue culture media, and buffers
- Specificity:
 - Detects multiple subtypes of IFN- α and all 3 subtypes of IFN- λ
 - Minimal cross-reactivity detected with Mouse IFN- α ; No cross-reactivity detected with Mouse IFN- β , IFN- γ , and IFN- λ 3
 - Low cross-reactivity detected with Rhesus and Cynomolgus monkey IFN- α ; Detects Cynomolgus monkey IFN- β



The VeriPlex Human Interferon 9-Plex ELISA has been developed to simultaneously detect Type I, II and III IFNs and related pro-inflammatory cytokines. This immunoassay employs an established ELISA-based multiplex technology in which 9 distinct capture antibodies have been adsorbed to each well of a 96-well plate in a defined array. Each spot represents a different distinct capture antibody population.

This ELISA kit provides an easy and cost-effective way to profile the biological complexity of interferon-related events by simultaneously quantifying the amount of human IFN- α , IFN- β , IFN- ω , IFN- γ , and IFN- λ in a single sample. The ability to assess multiple interferons 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, even autoimmune diseases.

Figure 1. Lower Limit of Detection (LLOD) and Lower Limit of Quantitation (LLOQ)

Analyte	Curve Range (pg/ml)	LLOD (pg/ml)	LLOQ (pg/ml)
Human IFN- α	3500 to 4.8	2.92	8.38
Human IFN- β	19,000 to 26.06	36.8	82.8
Human IFN- γ	5400 to 7.41	25.3	58.8
Human IFN- λ 1/2	8000 to 10.97	103.0	186.1
Human IFN- ω	4050 to 5.56	14.5	34.9
Human IL-1 α	1100 to 1.51	5.7	15.5
Human IL-6	20000 to 27.43	53.5	131.5
Human IP-10	1500 to 2.06	3.2	9.4
Human TNF- α	3000 to 4.12	7.5	17.6

*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.	51500-1
Description	VeriPlex Human Interferon 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.