

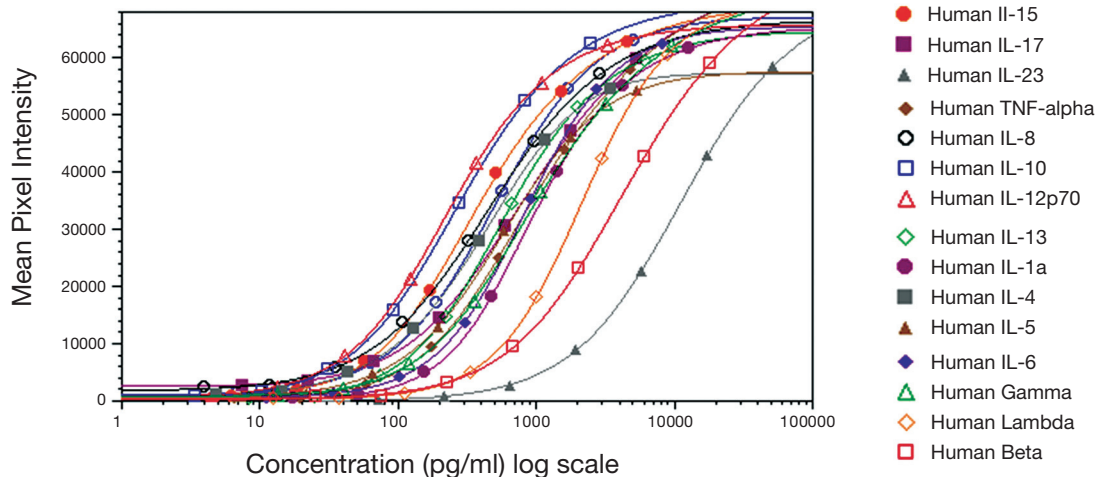


VeriPlex Human Cytokine 16-Plex ELISA

Catalog No. 51510

- Enables simultaneous quantification of Type I, II & III IFNs and other key innate and adaptive immunity cytokines in a single sample:
 - IFN- α , - β , - γ , - λ ; IL-1 α , -4, -5, -6, -8, -10, -12p70, -13, -15, -17, -23; TNF- α
- Utilizes easy to use 96-well plate-based ELISA technology and only requires a sample volume of 50 μ l per well
- Provides analyte sensitivity to as low as 5 pg/ml
- Achieves reproducible results with < 10% inter and intra-assay CVs in various sample matrices, including autoimmune sera and plasma

Figure 1. Representative standard curves



The Human Cytokine 16-Plex ELISA has been developed to simultaneously quantify 16 human cytokines in a single sample in low volumes of serum, plasma, and tissue culture media. This immunoassay employs an ELISA-based multiplex technology in which 16 distinct capture antibodies are spotted in each well of a 96-well plate in a defined 4x4 array. This ELISA kit provides unique chemiluminescent results on each analyte within a sample and allows for a global understanding of ongoing immune responses in diseases ranging from cancer to autoimmunity.

Product Information*

Catalog No.	51510-1
Description	VeriPlex Human Cytokine 16-Plex ELISA
Size	1 x 96-well plate
Compatibility	Serum, Plasma, Tissue Culture Media & Buffers
Assay Range	2.86 log with LLOD & LLOQ in the low pg/ml range

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

Analyte	LLOD (pg/ml)	LLOQ (pg/ml)	Assay Range (pg/ml)
IFN- α	2.41	7.00	5100 - 7.00
IFN- β	19.73	28.78	18300 - 25.10
IFN- γ	5.65	13.31	9700 - 13.31
IFN- λ	36.52	77.17	9000 - 12.35
IL-1 α	14.75	17.42	12700 - 17.42
IL-4	1.96	5.26	3500 - 4.80
IL-5	0.63	7.27	5300 - 7.27
IL-6	11.39	13.77	8300 - 11.39
IL-8	16.35	37.18	2900 - 3.98
IL-10	2.45	4.62	2500 - 3.43
IL-12p70	0.07	4.53	3300 - 4.53
IL-13	7.98	14.90	6000 - 8.23
IL-15	2.91	6.31	4600 - 6.31
IL-17	24.17	63.24	5400 - 7.41
IL-23	96.93	171.05	52000 - 71.33
TNF- α	4.68	9.79	4800 - 6.58

Figure 3. Human Cytokine 16-Plex ELISA IFN Alpha Subtypes Cross-Reactivity

Analyte	% Cross-Reactivity
Alpha J1	187.86%
Alpha 4a	159.58%
Alpha C	149.08%
Alpha 2b	133.04%
Alpha A 2a	114.67%
Alpha D (Alpha 1 [Ala114])	69.57%
Alpha D (Alpha 1 [Val114])	49.33%
Alpha K	37.58%
Alpha WA	19.69%
Alpha 4b	19.02%
Alpha G	2.80%
Alpha I	2.04%
Alpha H2	1.29%
Alpha B2	1.28%
Alpha F	0.00%

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



* Powered by Quansys Biosciences
Q-Plex™ Technology

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