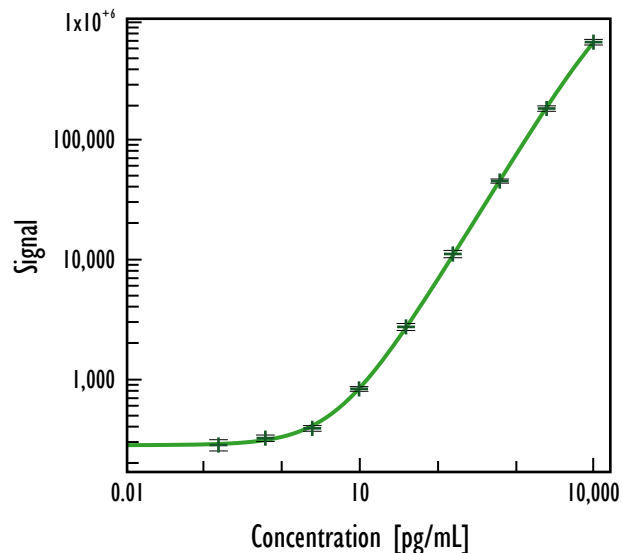


MULTI-ARRAY[®] Human IL-17 Tissue Culture Assay

Detection of IL-17 in Tissue Culture Samples

Typical Standard Curve Data (TCM)



Human IL-17		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	289	9.6
0.15	283	10.8
0.61	323	6.1
2.4	390	5.4
9.8	832	4.5
39	2,747	6.7
156	11,182	7.1
625	45,497	4.0
2,500	183,875	5.2
10,000	659,718	5.3

Typical Detection Parameters

Protocol	LLOD (pg/mL)
Tissue Culture	0.4

LLOD = Lower Limit of Detection is 2.5 stdev above the background

Note:

The MSD IL-17 assay detects IL-17A isoform and not the IL-17B, IL-17C, IL-17E and IL17F isoforms.

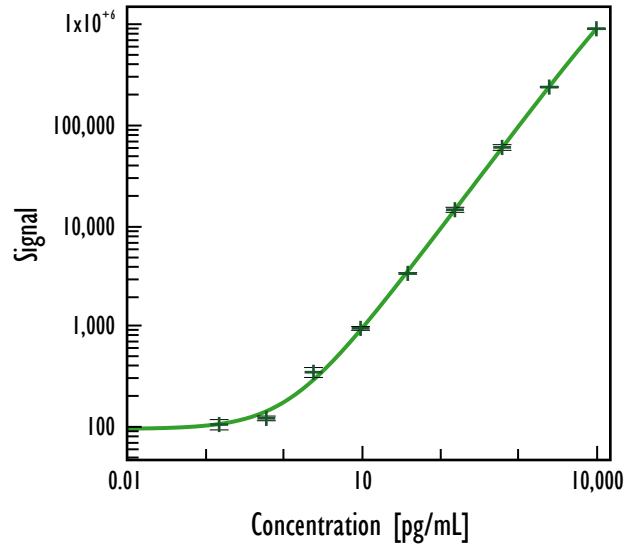
The IL-17A assay can be multiplexed with most other MSD cytokine assays.

Kit Model	Kit Size	Catalog Number
Tissue Culture	1 plate	K15IATB-1
Tissue Culture	5 plates	K15IATB-2
Tissue Culture	25 plates	K15IATB-4
Ultra-Sensitive	1 plate	K15IATC-1
Ultra-Sensitive	5 plates	K15IATC-2
Ultra-Sensitive	25 plates	K15IATC-4
Base	20 plates	K15IATA-4
Base	250 plates	K15IATA-5

MULTI-ARRAY[®] Human IL-17 Ultra-Sensitive Assay

Detection of IL-17 in Human Serum and Plasma Samples

Serum/Plasma Assay



Human IL-17		
Concentration (pg/mL)	Signal	
	Mean	%CV
0	108	29.2
0.15	106	11.8
0.61	122	5.4
2.4	351	11.3
9.8	958	4.3
39	3,381	1.8
156	14,518	5.8
625	60,521	6.4
2,500	241,099	1.3
10,000	916,742	1.2

Typical Detection Parameters

Protocol	LLOD (pg/mL)
Serum/Plasma	0.2

LLOD = Lower Limit of Detection
is 2.5 stdev above the background

Spike Recoveries

Sample Type	Average % Recovery
Serum	94
EDTA Plasma	85
Heparin Plasma	97

$$\% \text{ recovery} = \frac{\text{measured value} * 100}{\text{expected value}}$$

Spike recoveries were determined in each matrix over a range of spike levels.

Endogenous Levels

Sample Type	Samples Below LLOD
Serum	8

Eight human serum samples were measured in the IL-17 assay.

Note:

The MSD IL-17 assay detects IL-17A isoform and not the IL-17B, IL-17C, IL-17E and IL17F isoforms.
The IL-17A assay can be multiplexed with most other MSD cytokine assays.