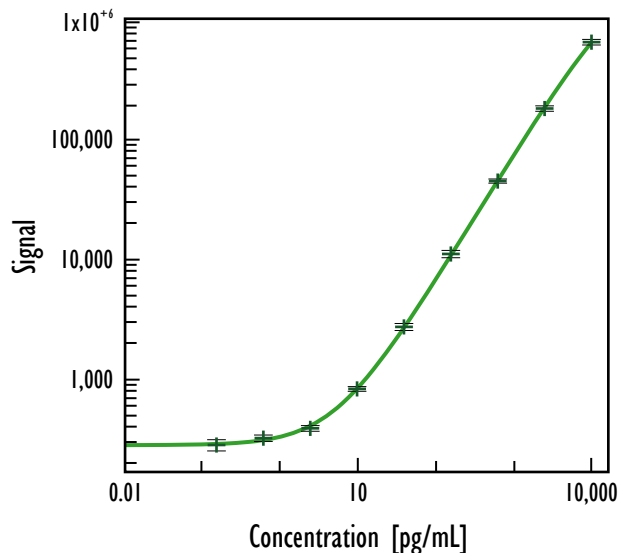


# MULTI-ARRAY<sup>®</sup> Human IL-17 Tissue Culture Assay

Detection of IL-17 in Tissue Culture Samples

## Typical Standard Curve Data (TCM)



Human IL-7		
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.

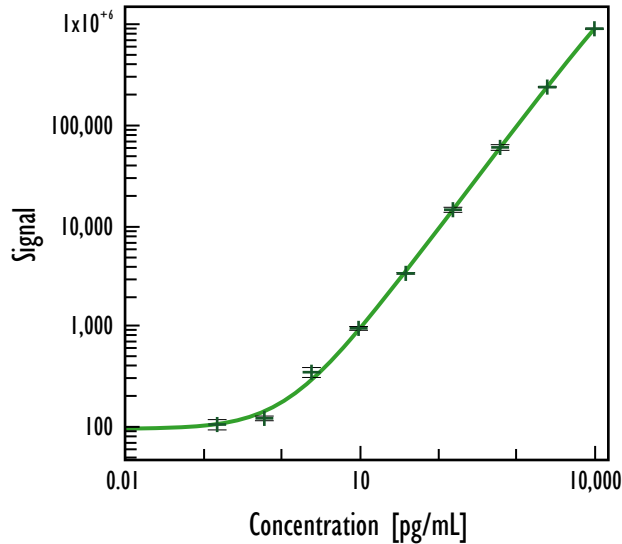
The IL-17A assay can be multiplexed with most other MSD cytokine assays.

Kit Model	Kit Size	SI2400	SI6000
Tissue Culture	1 plate	K151ATB-1	K111ATB-1
Tissue Culture	5 plates	K151ATB-2	K111ATB-2
Tissue Culture	25 plates	K151ATB-4	K111ATB-4
Ultra-Sensitive	1 plate	K151ATC-1	K111ATC-1
Ultra-Sensitive	5 plates	K151ATC-2	K111ATC-2
Ultra-Sensitive	25 plates	K151ATC-4	K111ATC-4
Base	20 plates	K151ATA-4	K111ATA-4
Base	250 plates	K151ATA-5	K111ATA-5

# MULTI-ARRAY<sup>®</sup> Human IL-17 Ultra-Sensitive Assay

## Detection of IL-17 in Human Serum and Plasma Samples

### Serum/Plasma Assay



Human IL-7		
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.