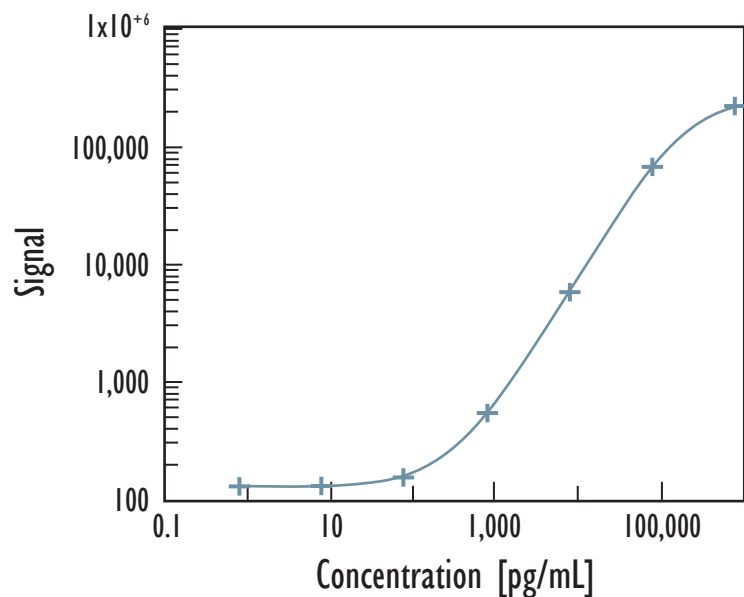
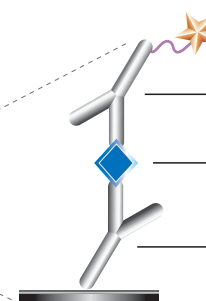
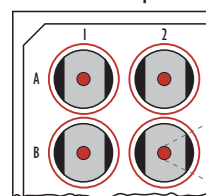


MULTI-ARRAY[®] Human TGF- β 1 Assay

Detection of Transforming Growth Factor β 1 (TGF- β 1) in Human Serum Samples



MSD MULTI-ARRAY[®]
96-Well Small Spot Plate



anti-TGF- β 1 SULFO-TAG[™] antibody

TGF- β 1

anti-TGF- β 1 capture antibody

Concentration (pg/mL)	Signal
0	138
0.1	129
1	135
10	157
100	565
1,000	5,853
10,000	67,476
100,000	218,841

TGF- β 1 LLOD	17 (pg/mL)
---------------------	------------

LLOD (Lower Limit of Detection) is defined as 2.5x stdev above the background

Kit Size	SI2400	SI6000
1 plate	K151IUC-1	K111IUC-1
5 plates	K151IUC-2	K111IUC-2
20 plates	K151IUC-3	K111IUC-3
20 plates (Base)	K151IUA-3	K111IUA-3

Above is representative calibration curve data

Serum and platelet —poor EDTA samples should be activated (acid treatment followed by neutralization) and then diluted 1:4 prior to use

MULTI-ARRAY[®] Human TGF- β 1 Assay

Detection of Transforming Growth Factor β 1 (TGF- β 1) in Human Serum Samples

Dilutional Linearity

- Serum samples from 5-6 individuals each were diluted with the appropriate calibrator diluents. Recoveries were measured at three dilution levels – the average recovery at each dilution level is tabulated.

$$\% \text{ recovery} = \frac{(\text{measured value} * \text{dilution factor} * 100)}{\text{predicted value}}$$

	Average Recovery at 3/4 Dilution Factor	Average Recovery at 1/2 Dilution Factor	Average Recovery at 1/4 Dilution Factor
TGF- β 1	99%	107%	133%

Spike Recovery

- Serum samples from 16 different individuals were spiked with known amounts of TGF- β 1 at defined levels (2-20 fold above endogenous) and measured. The average spike recovery is tabulated.

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

	Serum
TGF- β 1	92%

Endogenous Levels in Human Samples

- 14 normal human serum samples were measured.

	N	Mean	Median	Range
TGF- β 1 (pg/mL)	14	3,132	2,035	<1,307 - 13,403