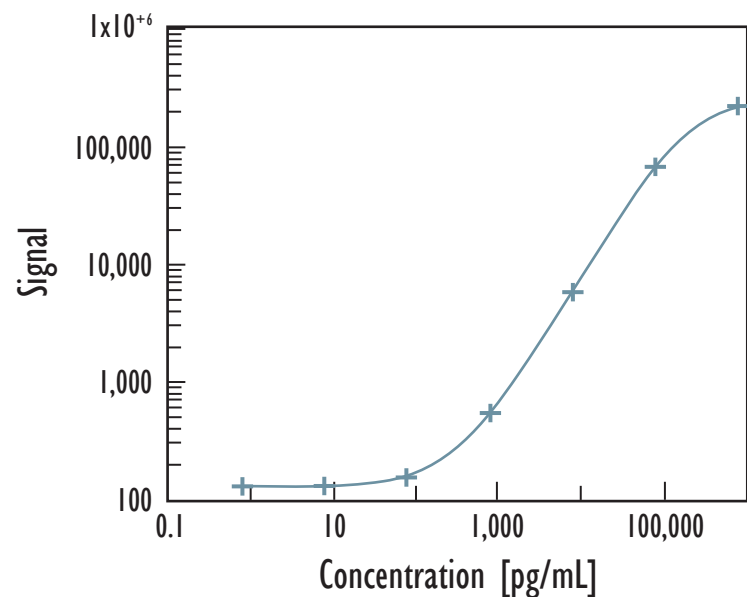
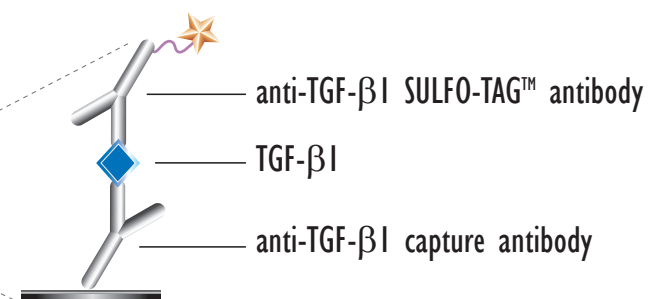
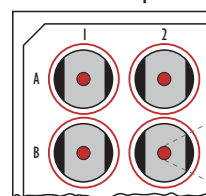


MULTI-ARRAY[®] Human TGF- β 1 Assay

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



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



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

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

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

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

Kit Size	Catalog Number
1 plate	K1511UC-1
5 plates	K1511UC-2
20 plates	K1511UC-3
20 plates (Base)	K1511UA-3

MULTI-ARRAY[®] Human TGF-β1 Assay

Detection of Transforming Growth Factor β1 (TGF-β1) 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-β1	99%	107%	133%

Spike Recovery

- Serum samples from 16 different individuals were spiked with known amounts of TGF-β1 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-β1	92%

Endogenous Levels in Human Samples

- 14 normal human serum samples were measured.

	N	Mean	Median	Range
TGF-β1 (pg/mL)	14	3,132	2,035	<1,307 - 13,403