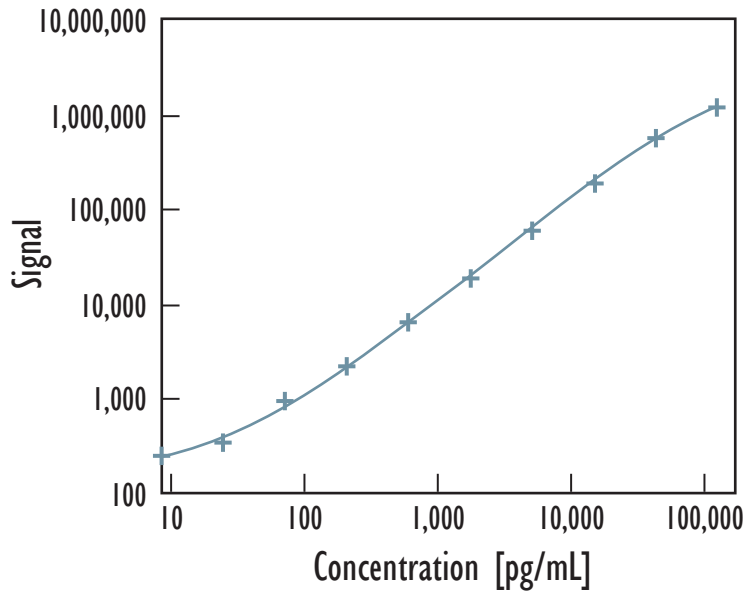


MULTI-ARRAY[®] Human c-Kit Assay

Detection of Kit Ligand Receptor (c-Kit) in Human Serum and Plasma Samples

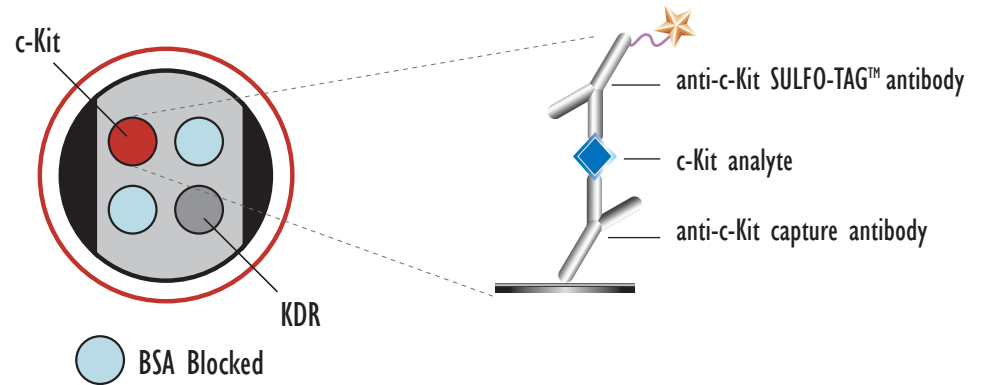


Concentration (pg/mL)	Average	%CV
0	205	6
8	279	6
23	385	5
69	1,047	7
210	2,381	7
620	6,848	6
1,900	20,203	3
5,600	65,068	1
17,000	206,719	3
50,000	614,835	4
150,000	1,276,402	4

Above is representative calibration curve data

50X dilution of samples is recommended

MSD MULTI-SPOT[®]
96-Well 4-Spot Plate



c-Kit LLOD	5 (pg/mL)
------------	-----------

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

Kit Size	SI2400	SI6000
1 plate	KI51BPC-1	KI11BPC-1
5 plates	KI51BPC-2	KI11BPC-2
20 plates	KI51BPC-3	KI11BPC-3
20 plates (Base)	KI51BPA-3	KI11BPA-3

MULTI-ARRAY[®] Human c-Kit Assay

Detection of Kit Ligand Receptor (c-Kit) in Human Serum and Plasma Samples

Dilutional Linearity

- Measured analyte levels in samples diluted in assay diluent

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

- Values presented are from pooled samples
- Recommend 1:50 dilution for samples; dilution factors indicated are in addition to 1:50 dilution

	Dilution Factor	c-Kit
Serum Pool 1	1:2	112
	1:4	102
	1:8	104
Serum Pool 2	1:2	120
	1:4	111
	1:8	108

Endogenous Levels in Human Samples

- Normal pooled samples
- Detected level is above LLOQ
- Average CVs for measured samples was less than 8%
- Samples diluted 1:50 for use in assay

	c-Kit (ng/mL)
Serum	165
EDTA Plasma	140

Spike Recovery

- Measured analyte spiked into human samples
- Samples were diluted 50X and then measured

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

	Spike Level (ng/mL)	c-Kit
Plasma 1	156	94
	625	98
	2,500	104
Plasma 2	156	84
	625	82
	2,500	103
Plasma 3	156	99
	625	103
	2,500	113
Plasma 4	156	95
	625	100
	2,500	110
Plasma 5	156	120
	625	113
	2,500	112