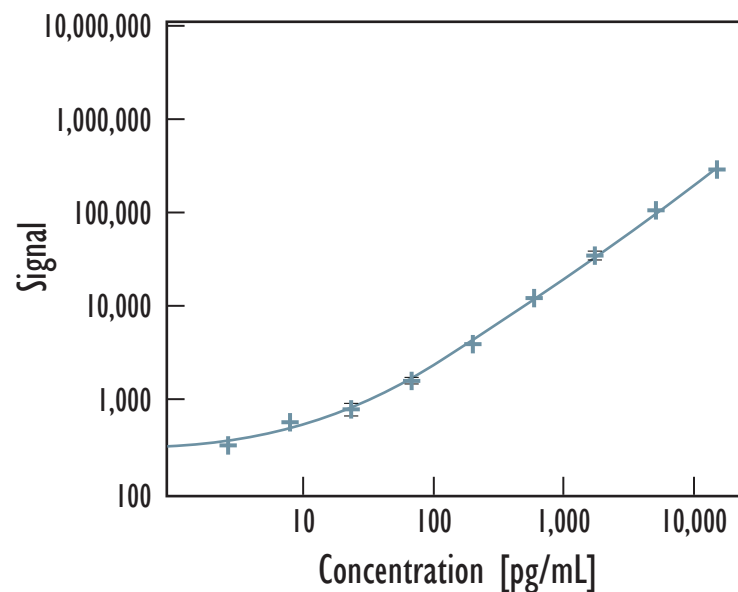


MULTI-ARRAY[®] Human KDR Assay

Detection of Kinase Domain Insert (KDR) in Human Serum and Plasma Samples

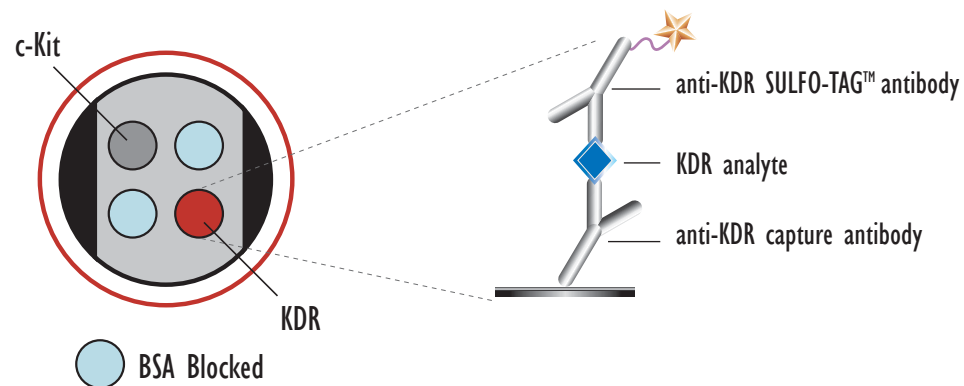


Concentration (pg/mL)	Average	%CV
0	276	6
0.8	305	1
2.3	321	5
6.9	541	1
21	736	15
62	1,480	7
190	3,624	5
560	10,974	2
1,700	30,472	10
5,000	91,732	2
15,000	248,936	1

Above is representative calibration curve data

50X dilution of samples is recommended

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



KDR LLOD	2 (pg/mL)
----------	-----------

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

Kit Size	SI2400	SI6000
1 plate	KI51BOC-1	KI11BOC-1
5 plates	KI51BOC-2	KI11BOC-2
20 plates	KI51BOC-3	KI11BOC-3
20 plates (Base)	KI51BOA-3	KI11BOA-3

MULTI-ARRAY[®] Human KDR Assay

Detection of Kinase Domain Insert (KDR) 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	KDR
Serum Pool 1	1:2	103
	1:4	97
	1:8	106
Serum Pool 2	1:2	111
	1:4	107
	1:8	101

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

	KDR (ng/mL)
Serum	31
EDTA Plasma	28

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)	KDR
Plasma 1	156	93
	625	96
	2,500	106
Plasma 2	156	82
	625	78
	2,500	105
Plasma 3	156	105
	625	100
	2,500	119
Plasma 4	156	96
	625	97
	2,500	114
Plasma 5	156	93
	625	97
	2,500	108