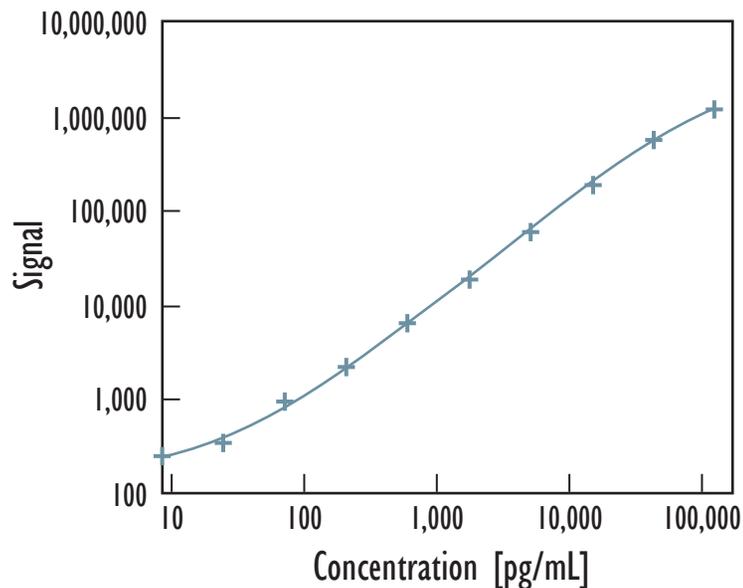


MULTI-ARRAY[®] Human c-Kit Assay

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

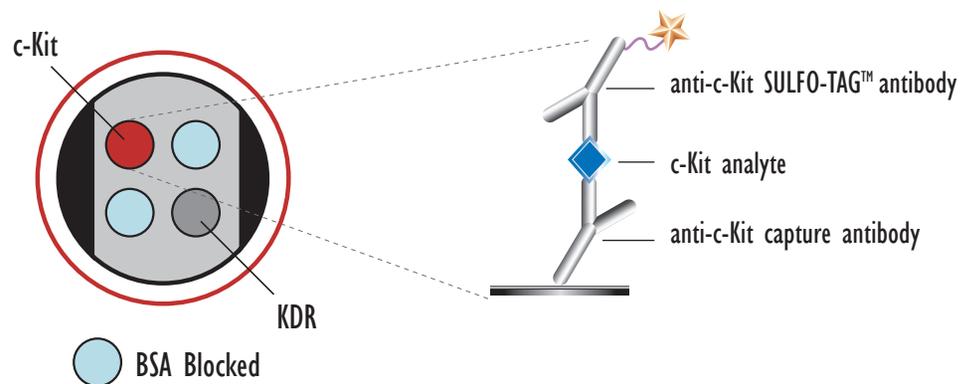


| Concentration (pg/mL) | Average | %CV |
|-----------------------|-----------|-----|
| 0 | 205 | 6 |
| 8 | 279 | 6 |
| 23 | 385 | 5 |
| 69 | 1,047 | 7 |
| 210 | 2,381 | 7 |
| 620 | 6,848 | 6 |
| 1,900 | 20,203 | 3 |
| 5,600 | 65,068 | 1 |
| 17,000 | 206,719 | 3 |
| 50,000 | 614,835 | 4 |
| 150,000 | 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 | Catalog Number |
|------------------|----------------|
| 1 plate | KI51BPC-1 |
| 5 plates | KI51BPC-2 |
| 20 plates | KI51BPC-3 |
| 20 plates (Base) | KI51BPA-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 |