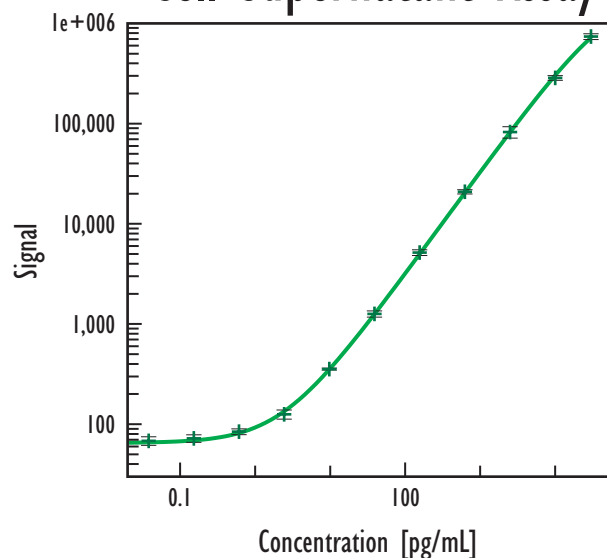


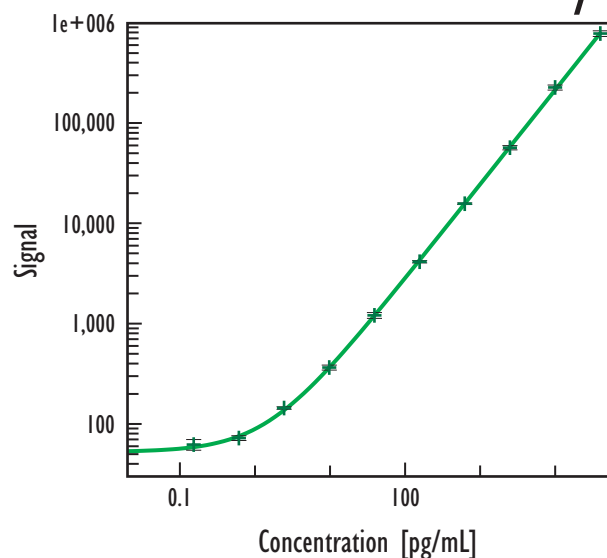
Human IL-6 Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



| Human IL-6 | | |
|-----------------------|---------|------|
| Concentration (pg/mL) | Signal | |
| | Mean | %CV |
| 0 | 59 | 9.0 |
| 0.010 | 68 | 7.1 |
| 0.038 | 68 | 9.8 |
| 0.15 | 72 | 8.6 |
| 0.61 | 84 | 6.6 |
| 2.4 | 126 | 10.5 |
| 9.8 | 355 | 2.2 |
| 39 | 1,262 | 7.1 |
| 156 | 5,165 | 6.5 |
| 625 | 20,875 | 4.8 |
| 2,500 | 82,404 | 13.0 |
| 10,000 | 285,888 | 5.4 |
| 40,000 | 738,908 | 6.6 |

Serum / Plasma Assay



| Human IL-6 | | |
|-----------------------|---------|------|
| Concentration (pg/mL) | Signal | |
| | Mean | %CV |
| 0 | 51 | 11.0 |
| 0.15 | 62 | 12.3 |
| 0.61 | 72 | 5.5 |
| 2.4 | 144 | 2.9 |
| 9.8 | 365 | 5.7 |
| 39 | 1,209 | 6.6 |
| 156 | 4,140 | 2.7 |
| 625 | 15,700 | 1.7 |
| 2,500 | 56,809 | 4.9 |
| 10,000 | 226,443 | 5.8 |
| 40,000 | 781,884 | 6.5 |

Detection Limits

| Sample Type | Detection Limit (pg/mL) |
|------------------|-------------------------|
| Cell Supernatant | 0.9 |
| Serum | 0.7 |
| Plasma | 1.6 |

Detection Limits were determined across multiple runs using 2.5 standard deviations above the background.

Recoveries

| Sample Type | Average % Recovery |
|------------------|--------------------|
| Cell Supernatant | 108 |
| Serum | 86 |
| EDTA Plasma | 102 |
| Heparin Plasma | 99 |

Spike recoveries were determined in each matrix over a range of spike levels from 9.8 to 313 pg/mL. Each spike was tested in ≥ 3 replicates.

Endogenous Levels

| Sample Type (# of unique samples) | Samples Below Detectable Range | Range |
|-----------------------------------|--------------------------------|------------|
| Serum (n=9) | 0 | 1.2 - 15.3 |
| EDTA Plasma (n=5) | 1 | n/d - 4.7 |
| Heparin Plasma (n=5) | 0 | 1.9 - 3.9 |

Endogenous cytokine levels (pg/mL) were determined for different sample types across multiple samples. (n/d indicates cytokine level below the detection limit)

Representative data shown above is for demonstration purposes only. Individual results may vary depending upon samples tested.