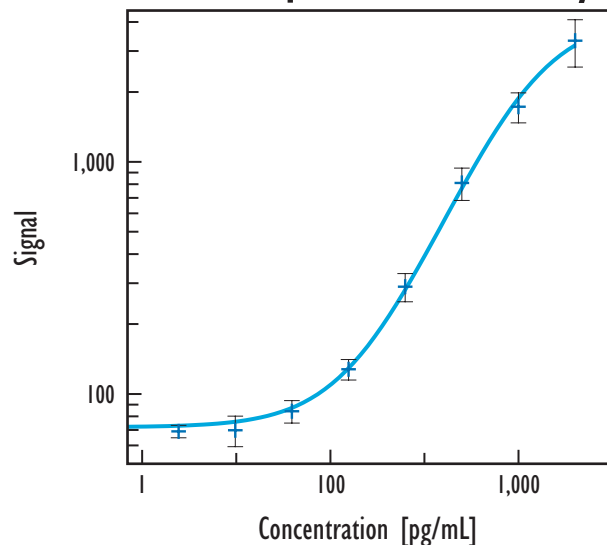


Mouse IL-12p70 Cytokine Assay (MULTI-ARRAY™ 96-Well Small Spot Plate)

Cell Supernatant Assay



| Mouse IL-12p70 | | |
|-----------------------|--------|------|
| Concentration (pg/mL) | Signal | |
| | Mean | %CV |
| 0 | 66 | 10.8 |
| 0.010 | 79 | 7.5 |
| 0.038 | 82 | 10.7 |
| 0.15 | 75 | 6.6 |
| 0.61 | 75 | 3.9 |
| 2.4 | 69 | 6.1 |
| 9.8 | 70 | 15.0 |
| 39 | 84 | 11.1 |
| 156 | 128 | 10.2 |
| 625 | 290 | 13.9 |
| 2,500 | 756 | 11.0 |
| 10,000 | 1,629 | 12.3 |
| 40,000 | 2,958 | 8.8 |

Detection Limits

| Sample Type | Detection Limit (pg/mL) |
|------------------|-------------------------|
| Cell Supernatant | 65 |
| Serum / Plasma | 346 |

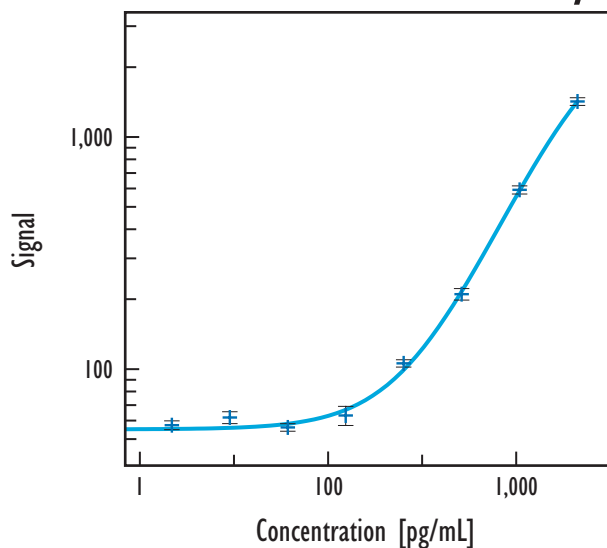
Detection Limits were determined across multiple runs using 2.5 standard deviations above the background.

Recoveries

| Sample Type | Average % Recovery |
|------------------|--------------------|
| Cell Supernatant | 103 |
| Serum | 88 |
| EDTA Plasma | 87 |

Spike recoveries were determined in each matrix over a range of spike levels. Each spike was tested in triplicate.

Serum / Plasma Assay



| Mouse IL-12p70 | | |
|-----------------------|--------|-----|
| Concentration (pg/mL) | Signal | |
| | Mean | %CV |
| 0 | 50 | 7.4 |
| 0.15 | 61 | 4.8 |
| 0.61 | 52 | 7.4 |
| 2.4 | 57 | 4.4 |
| 9.8 | 61 | 5.9 |
| 39 | 56 | 3.8 |
| 156 | 63 | 9.5 |
| 625 | 105 | 3.6 |
| 2,500 | 209 | 5.7 |
| 10,000 | 590 | 4.0 |
| 40,000 | 1,421 | 3.8 |

Endogenous Levels

| Sample Type (# of unique samples) | Samples Below Detectable Range | Range |
|-----------------------------------|--------------------------------|------------|
| Serum (n=8) | 0 | 471 - 1003 |
| Heparin Plasma (n=8) | 5 | n/d - 683 |

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.