

| | |
|-----------------|-------------|
| Product Number: | 85-0329 |
| Lot Number: | 319713 |
| Expiration: | 10-May-2023 |
| Platform(s): | SP-X |

| Component | Part Number | Lot Number |
|--|-------------|------------|
| CorPlex Cytokine Plate | 97-0329 | 319713P |
| Biotinylated Antibody Rgt | 98-0329 | 319713B |
| Sample Diluent 21 | 1863329 | 319579 |
| IL-12p70 Calibrator | 18910034 | 319156 |
| IL-1 β Calibrator | 18910040 | 319564 |
| IL-4 Calibrator | 1858452 | 319458 |
| IL-5 Calibrator | 1858462 | 319223 |
| IFN γ Calibrator | 1858260 | 318898 |
| IL-6 Calibrator | 18910044 | 319484 |
| IL-8 Calibrator | 1858454 | 319436 |
| IL-22 Calibrator | 18910036 | 319459 |
| TNF α Calibrator | 1880012 | 319460 |
| IL-10 Calibrator | 18910043 | 319215 |
| Wash Buffer (25X) ¹ | 1860055 | N/A |
| Streptavidin-HRP Reagent ¹ | 1858263 | N/A |
| Super Signal [®] Luminol* ¹ | 1858089 | N/A |
| Super Signal [®] Peroxide* ¹ | 1858090 | N/A |



Data below represents results generated on the SP-X Imaging and Analysis System

¹ Reagents are not Kit Lot Specific

*SuperSignal[®] technology is protected by US Patent #6,432,662 and is a registered trademark of ThermoFisher Scientific.

All reagents are specific for use on the SP-X platform - refer to the CorPlex Cytokine Kit Instructions (85A-0329).

Release Materials

IL-12p70

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|---------|--------|----------------|--------------------|
| Panel 1 | 318935 | 167.6 | 133.1 - 204.3 |
| Panel 2 | 318935 | 51.74 | 37.09 - 66.72 |
| Panel 3 | 318935 | 17.29 | 13.48 - 21.97 |
| Panel 4 | 318935 | 5.157 | 3.812 - 6.073 |
| Panel 5 | 318935 | 1.613 | 1.188 - 2.091 |

IL-1 β

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|---------|--------|----------------|--------------------|
| Panel 1 | 318935 | 27.48 | 12.49 - 32.68 |
| Panel 2 | 318935 | 8.150 | 4.122 - 9.719 |
| Panel 3 | 318935 | 2.790 | 1.349 - 3.475 |
| Panel 4 | 318935 | 0.830 | 0.418 - 0.968 |
| Panel 5 | 318935 | 0.282 | 0.133 - 0.345 |

IL-4

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|---------|--------|----------------|--------------------|
| Panel 1 | 318935 | 69.35 | 51.07 - 77.41 |
| Panel 2 | 318935 | 20.91 | 14.17 - 24.05 |
| Panel 3 | 318935 | 7.526 | 5.217 - 8.887 |
| Panel 4 | 318935 | 2.294 | 1.455 - 2.760 |
| Panel 5 | 318935 | 0.882 | 0.418 - 1.250 |

IL-5

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|---------|--------|----------------|--------------------|
| Panel 1 | 318935 | 60.47 | 43.59 - 101.9 |
| Panel 2 | 318935 | 22.03 | 13.43 - 32.50 |
| Panel 3 | 318935 | 6.604 | 4.948 - 11.09 |
| Panel 4 | 318935 | 2.196 | 1.317 - 3.402 |
| Panel 5 | 318935 | 0.778 | 0.473 - 1.224 |

Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Release Materials

IFN γ

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|---------|--------|----------------|--------------------|
| Panel 1 | 318935 | 9.877 | 8.190 - 14.53 |
| Panel 2 | 318935 | 3.106 | 2.393 - 4.771 |
| Panel 3 | 318935 | 1.001 | 0.855 - 1.720 |
| Panel 4 | 318935 | 0.330 | 0.277 - 0.494 |
| Panel 5 | 318935 | 0.141 | 0.114 - 0.221 |

IL-6

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|---------|--------|----------------|--------------------|
| Panel 1 | 318935 | 172.8 | 142.9 - 214.3 |
| Panel 2 | 318935 | 77.05 | 62.87 - 99.85 |
| Panel 3 | 318935 | 15.76 | 13.37 - 20.05 |
| Panel 4 | 318935 | 7.599 | 5.951 - 8.927 |
| Panel 5 | 318935 | 2.452 | 1.992 - 3.084 |

IL-8

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|---------|--------|----------------|--------------------|
| Panel 1 | 318935 | 460.7 | 365.0 - 604.4 |
| Panel 2 | 318935 | 129.6 | 92.02 - 191.0 |
| Panel 3 | 318935 | 48.59 | 32.36 - 67.44 |
| Panel 4 | 318935 | 13.55 | 9.991 - 18.54 |
| Panel 5 | 318935 | 4.693 | 3.355 - 6.599 |

IL-22

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|---------|--------|----------------|--------------------|
| Panel 1 | 318935 | 69.95 | 50.55 - 160.4 |
| Panel 2 | 318935 | 13.08 | 7.411 - 28.41 |
| Panel 3 | 318935 | 3.989 | 2.237 - 9.205 |
| Panel 4 | 318935 | 1.312 | 0.738 - 2.619 |
| Panel 5 | 318935 | 0.386 | 0.226 - 0.888 |

TNF α

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|---------|--------|----------------|--------------------|
| Panel 1 | 318935 | 304.4 | 250.8 - 453.3 |
| Panel 2 | 318935 | 94.37 | 68.46 - 156.1 |
| Panel 3 | 318935 | 33.15 | 25.97 - 51.08 |
| Panel 4 | 318935 | 10.32 | 7.455 - 14.81 |
| Panel 5 | 318935 | 3.364 | 2.339 - 4.887 |

IL-10

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|---------|--------|----------------|--------------------|
| Panel 1 | 318935 | 104.4 | 76.56 - 126.0 |
| Panel 2 | 318935 | 31.73 | 21.73 - 39.58 |
| Panel 3 | 318935 | 10.34 | 7.682 - 13.72 |
| Panel 4 | 318935 | 3.267 | 2.408 - 3.832 |
| Panel 5 | 318935 | 1.220 | 0.837 - 1.442 |

Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Review/Approval

Shivani Goel

Quality Control
Manager

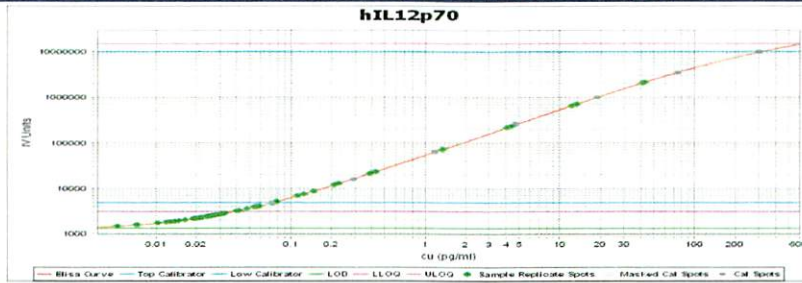
Shivani Goel 12 May 2022

Name

Title

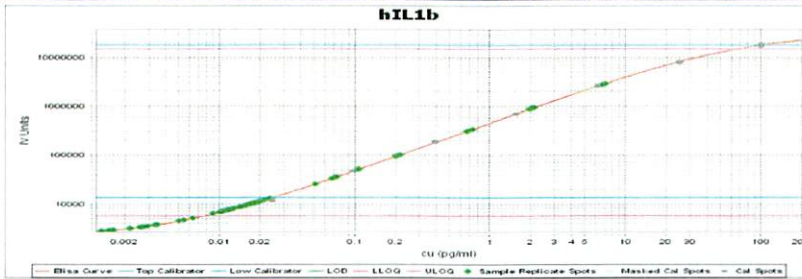
Signature/Date

IL-12p70 Calibration Curve



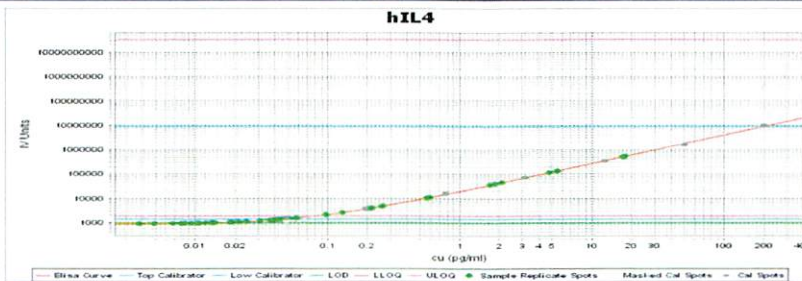
| Calibrator Levels (pg/ml) | |
|---------------------------|-------|
| 1 | 300 |
| 2 | 75 |
| 3 | 18.75 |
| 4 | 4.688 |
| 5 | 1.172 |
| 6 | 0.293 |
| 7 | 0.073 |
| 8 | 0.000 |

IL-1β Calibration Curve



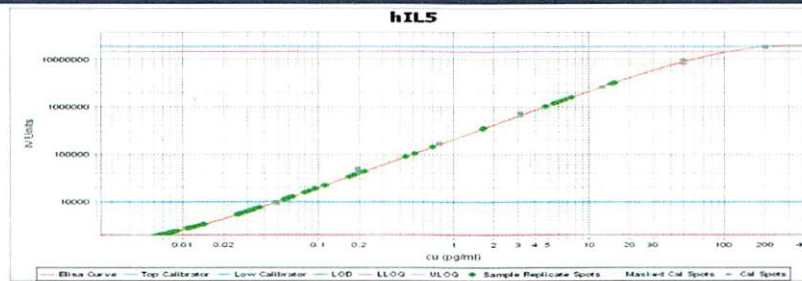
| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| 1 | 100 |
| 2 | 25 |
| 3 | 6.25 |
| 4 | 1.563 |
| 5 | 0.391 |
| 6 | 0.098 |
| 7 | 0.024 |
| 8 | 0.000 |

IL-4 Calibration Curve



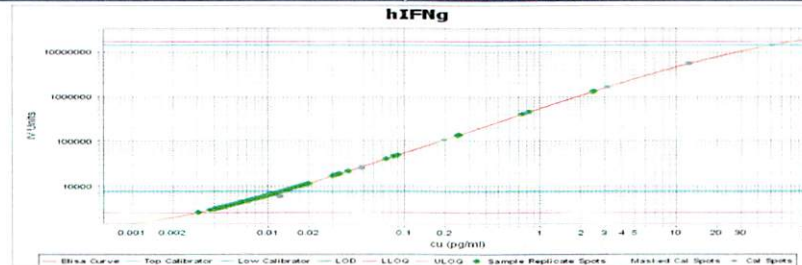
| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| 1 | 200 |
| 2 | 50 |
| 3 | 12.50 |
| 4 | 3.125 |
| 5 | 0.781 |
| 6 | 0.195 |
| 7 | 0.049 |
| 8 | 0.000 |

IL-5 Calibration Curve



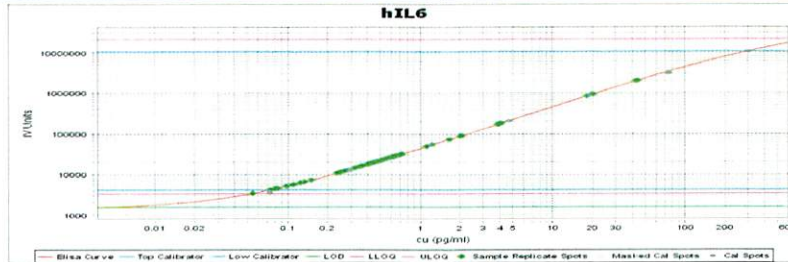
| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| 1 | 200 |
| 2 | 50 |
| 3 | 12.50 |
| 4 | 3.125 |
| 5 | 0.781 |
| 6 | 0.195 |
| 7 | 0.049 |
| 8 | 0.000 |

IFNγ Calibration Curve



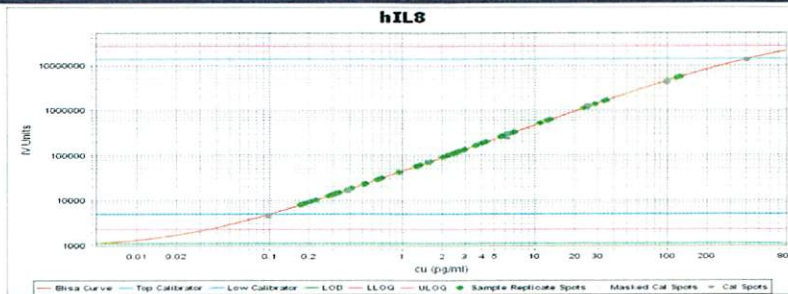
| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| 1 | 50 |
| 2 | 12.5 |
| 3 | 3.13 |
| 4 | 0.781 |
| 5 | 0.195 |
| 6 | 0.049 |
| 7 | 0.012 |
| 8 | 0.000 |

IL-6 Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| 1 | 300 |
| 2 | 75 |
| 3 | 18.75 |
| 4 | 4.688 |
| 5 | 1.172 |
| 6 | 0.293 |
| 7 | 0.073 |
| 8 | 0.000 |

IL-8 Calibration Curve



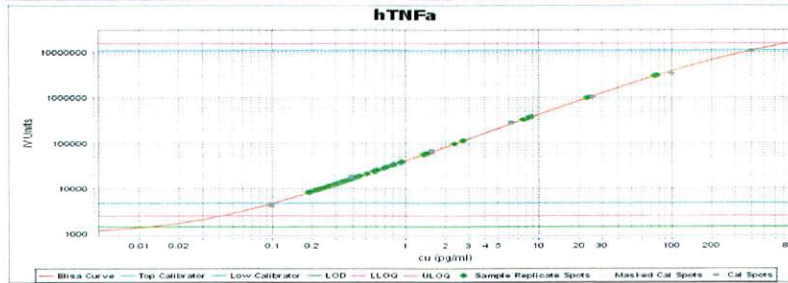
| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| 1 | 400 |
| 2 | 100 |
| 3 | 25.00 |
| 4 | 6.250 |
| 5 | 1.563 |
| 6 | 0.391 |
| 7 | 0.098 |
| 8 | 0.000 |

IL-22 Calibration Curve



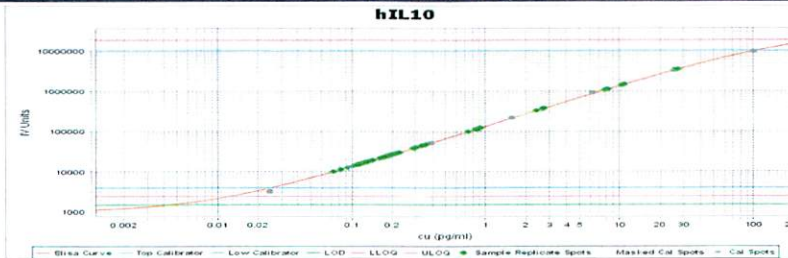
| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| 1 | 100 |
| 2 | 25 |
| 3 | 6.25 |
| 4 | 1.563 |
| 5 | 0.391 |
| 6 | 0.098 |
| 7 | 0.024 |
| 8 | 0.000 |

TNFα Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| 1 | 400 |
| 2 | 100 |
| 3 | 25.00 |
| 4 | 6.250 |
| 5 | 1.563 |
| 6 | 0.391 |
| 7 | 0.098 |
| 8 | 0.000 |

IL-10 Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| 1 | 100 |
| 2 | 25 |
| 3 | 6.25 |
| 4 | 1.563 |
| 5 | 0.391 |
| 6 | 0.098 |
| 7 | 0.024 |
| 8 | 0.000 |