

Product Number:	85-0329
Lot Number:	319694
Expiration:	10-May-2023
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	319694P
Biotinylated Antibody Rgt	98-0329	319694B
Sample Diluent 21	1863329	319579
IL-12p70 Calibrator	18910034	319156
IL-1 $\beta$ Calibrator	18910040	319564
IL-4 Calibrator	1858452	319458
IL-5 Calibrator	1858462	319223
IFN $\gamma$ Calibrator	1858260	318886
IL-6 Calibrator	18910044	319484
IL-8 Calibrator	1858454	319436
IL-22 Calibrator	18910036	319176
TNF $\alpha$ Calibrator	1880012	319460
IL-10 Calibrator	18910043	319215
Wash Buffer (25X) <sup>1</sup>	1860055	N/A
Streptavidin-HRP Reagent <sup>1</sup>	1858263	N/A
Super Signal <sup>®</sup> Luminol* <sup>1</sup>	1858089	N/A
Super Signal <sup>®</sup> Peroxide* <sup>1</sup>	1858090	N/A



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

<sup>1</sup> Reagents are not Kit Lot Specific

\*SuperSignal<sup>®</sup> 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	141.8	133.1 - 204.3
Panel 2	318935	45.49	37.09 - 66.72
Panel 3	318935	16.05	13.48 - 21.97
Panel 4	318935	4.850	3.812 - 6.073
Panel 5	318935	1.650	1.188 - 2.091

**IL-1 $\beta$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	22.95	12.49 - 32.68
Panel 2	318935	6.944	4.122 - 9.719
Panel 3	318935	2.519	1.349 - 3.475
Panel 4	318935	0.753	0.418 - 0.968
Panel 5	318935	0.288	0.133 - 0.345

**IL-4**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	53.42	51.07 - 77.41
Panel 2	318935	16.84	14.17 - 24.05
Panel 3	318935	6.440	5.217 - 8.887
Panel 4	318935	2.058	1.455 - 2.760
Panel 5	318935	0.827	0.418 - 1.250

**IL-5**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	55.27	43.59 - 101.9
Panel 2	318935	18.53	13.43 - 32.50
Panel 3	318935	6.447	4.948 - 11.09
Panel 4	318935	2.240	1.317 - 3.402
Panel 5	318935	0.759	0.473 - 1.224

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

**Release Materials**

**IFN $\gamma$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	10.79	8.190 - 14.53
Panel 2	318935	3.535	2.393 - 4.771
Panel 3	318935	1.248	0.855 - 1.720
Panel 4	318935	0.410	0.277 - 0.494
Panel 5	318935	0.185	0.114 - 0.221

**IL-6**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	176.9	142.9 - 214.3
Panel 2	318935	83.11	62.87 - 99.85
Panel 3	318935	17.03	13.37 - 20.05
Panel 4	318935	8.338	5.951 - 8.927
Panel 5	318935	2.923	1.992 - 3.084

**IL-8**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	421.4	365.0 - 604.4
Panel 2	318935	124.0	92.02 - 191.0
Panel 3	318935	45.68	32.36 - 67.44
Panel 4	318935	15.07	9.991 - 18.54
Panel 5	318935	5.260	3.355 - 6.599

**IL-22**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	95.28	50.55 - 160.4
Panel 2	318935	15.98	7.411 - 28.41
Panel 3	318935	5.473	2.237 - 9.205
Panel 4	318935	1.765	0.738 - 2.619
Panel 5	318935	0.652	0.226 - 0.888

**TNF $\alpha$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	254.9	250.8 - 453.3
Panel 2	318935	82.41	68.46 - 156.1
Panel 3	318935	30.03	25.97 - 51.08
Panel 4	318935	9.639	7.455 - 14.81
Panel 5	318935	3.318	2.339 - 4.887

**IL-10**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	90.47	76.56 - 126.0
Panel 2	318935	27.27	21.73 - 39.58
Panel 3	318935	9.788	7.682 - 13.72
Panel 4	318935	3.127	2.408 - 3.832
Panel 5	318935	1.188	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* 09 May 2022

Name

Title

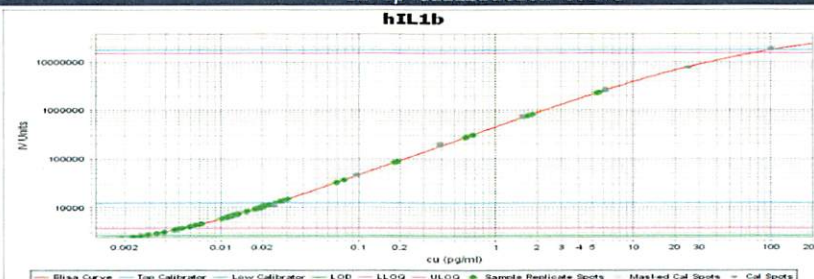
Signature/Date

**IL-12p70 Calibration Curve**



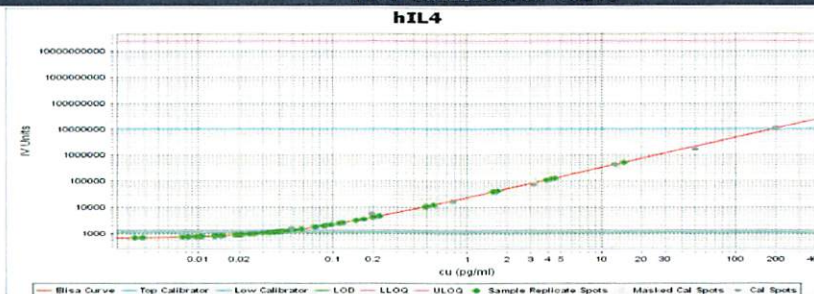
Calibrator Levels (pg/ml)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

**IL-1β Calibration Curve**



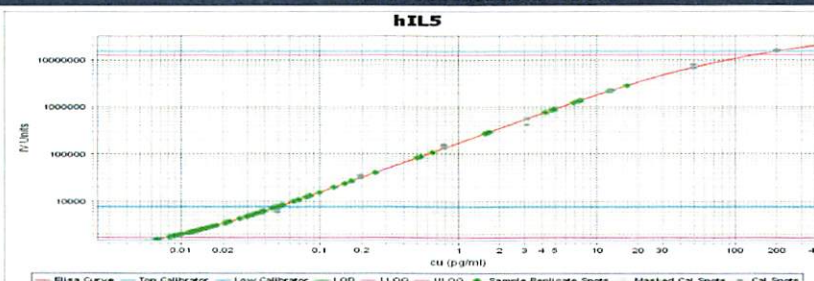
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

**IL-4 Calibration Curve**



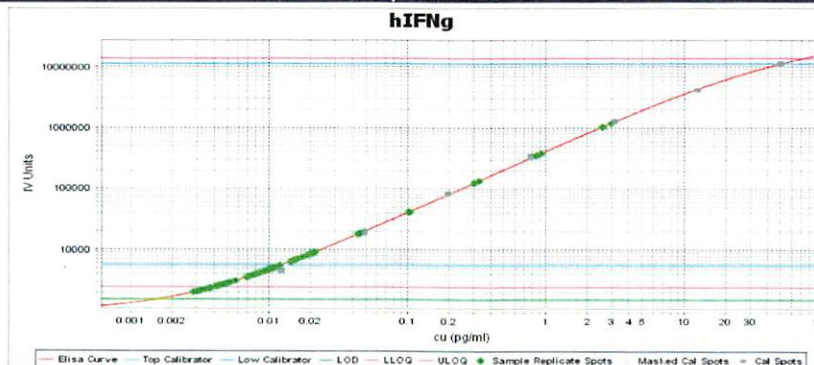
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

**IL-5 Calibration Curve**



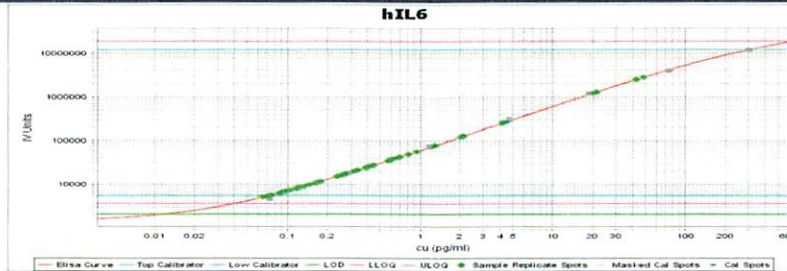
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

**IFNγ Calibration Curve**



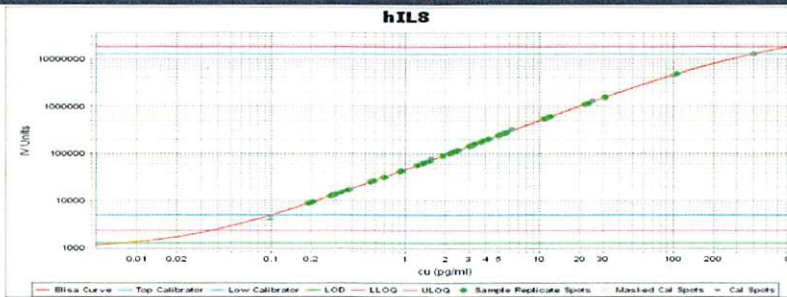
Calibrator Levels (pg/mL)	
1	50
2	12.5
3	3.13
4	0.781
5	0.195
6	0.049
7	0.012
8	0.000

**IL-6 Calibration Curve**



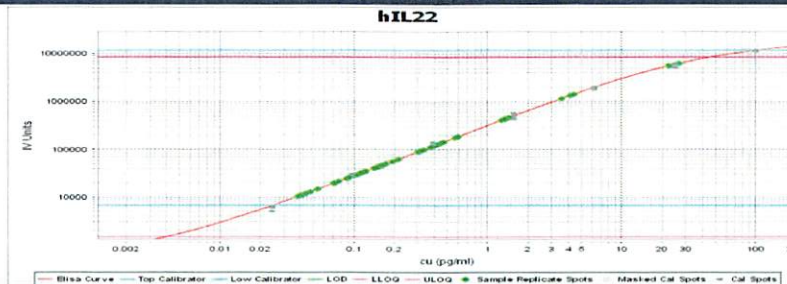
Calibrator Levels (pg/mL)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

**IL-8 Calibration Curve**



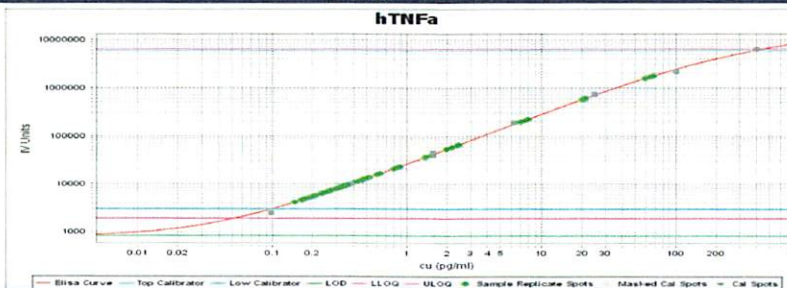
Calibrator Levels (pg/mL)	
1	400
2	100
3	25.00
4	6.250
5	1.563
6	0.391
7	0.098
8	0.000

**IL-22 Calibration Curve**



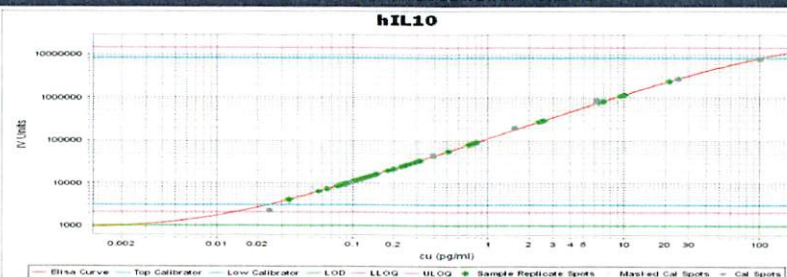
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

**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