

Product Number:	85-0329
Lot Number:	319275
Expiration:	27-Nov-2021
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	319275P
Biotinylated Antibody Rgt	98-0329	319275B
Sample Diluent 21	1863329	319289
IL-12p70 Calibrator	18910034	318756
IL-1 $\beta$ Calibrator	18910040	318728
IL-4 Calibrator	1858452	318580
IL-5 Calibrator	1858462	318837
IFN $\gamma$ Calibrator	1858260	318886
IL-6 Calibrator	18910044	318724
IL-8 Calibrator	1858454	318500
IL-22 Calibrator	18910036	318813
TNF $\alpha$ Calibrator	1880012	318883
IL-10 Calibrator	18910043	318755
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	193.3	140.6 - 239.0
Panel 2	318935	58.37	44.36 - 73.70
Panel 3	318935	21.22	15.74 - 23.61
Panel 4	318935	5.645	4.400 - 6.601
Panel 5	318935	1.562	1.243 - 2.397

**IL-1 $\beta$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	22.04	16.60 - 33.00
Panel 2	318935	7.084	5.060 - 10.28
Panel 3	318935	2.383	1.780 - 3.430
Panel 4	318935	0.803	0.496 - 1.030
Panel 5	318935	0.233	0.151 - 0.367

**IL-4**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	71.07	53.27 - 82.46
Panel 2	318935	21.30	16.02 - 24.04
Panel 3	318935	8.046	6.011 - 9.016
Panel 4	318935	2.381	1.729 - 2.776
Panel 5	318935	0.948	0.434 - 1.312

**IL-5**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	87.51	54.07 - 121.1
Panel 2	318935	26.82	18.39 - 34.26
Panel 3	318935	9.31	6.909 - 11.42
Panel 4	318935	2.688	2.007 - 3.306
Panel 5	318935	0.935	0.615 - 1.271

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	12.48	9.152 - 15.15
Panel 2	318935	3.890	2.766 - 5.014
Panel 3	318935	1.385	1.021 - 1.670
Panel 4	318935	0.435	0.320 - 0.515
Panel 5	318935	0.173	0.142 - 0.223

**IL-6**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	191.2	140.8 - 249.6
Panel 2	318935	86.01	68.32 - 107.0
Panel 3	318935	17.34	14.11 - 21.96
Panel 4	318935	8.078	6.379 - 9.569
Panel 5	318935	2.511	1.951 - 3.380

**IL-8**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	490.1	287.6 - 812.3
Panel 2	318935	147.5	99.01 - 213.4
Panel 3	318935	50.64	33.87 - 79.01
Panel 4	318935	15.26	10.72 - 20.44
Panel 5	318935	5.042	3.706 - 7.504

**IL-22**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	131.2	73.64 - 264.5
Panel 2	318935	23.07	13.35 - 41.63
Panel 3	318935	7.407	6.048 - 11.67
Panel 4	318935	2.148	1.025 - 3.238
Panel 5	318935	0.665	0.478 - 1.052

**TNF $\alpha$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	403.2	213.6 - 599.8
Panel 2	318935	129.8	81.39 - 179.7
Panel 3	318935	46.07	30.07 - 60.48
Panel 4	318935	12.29	9.413 - 16.87
Panel 5	318935	3.644	2.187 - 6.149

**IL-10**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	104.8	86.85 - 130.3
Panel 2	318935	32.00	25.35 - 40.54
Panel 3	318935	11.01	9.246 - 13.87
Panel 4	318935	3.122	2.649 - 4.032
Panel 5	318935	1.042	0.890 - 1.546

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 26 May 2021

Name

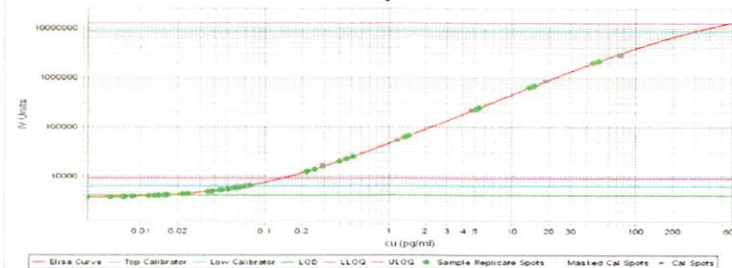
Title

Signature/Date



**IL-12p70 Calibration Curve**

**hIL12p70**

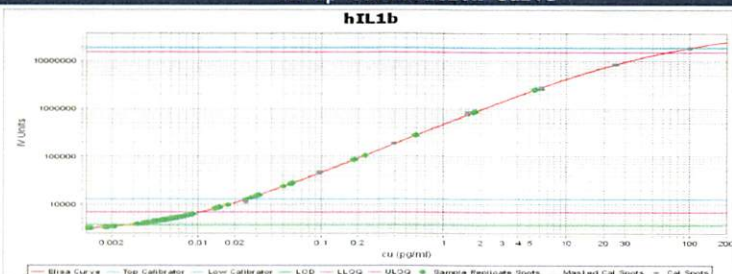


**Calibrator Levels (pg/ml)**

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**

**hIL1b**

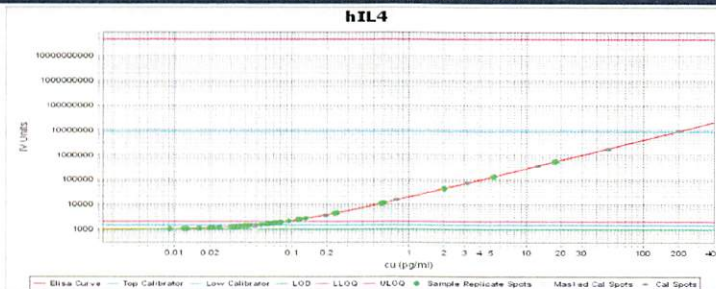


**Calibrator Levels (pg/mL)**

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**

**hIL4**

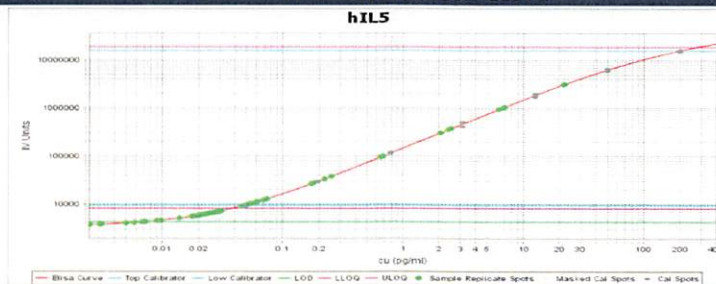


**Calibrator Levels (pg/mL)**

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**

**hIL5**

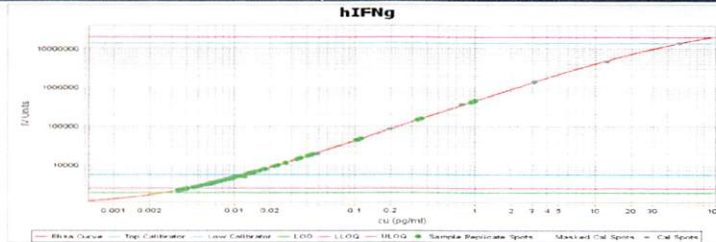


**Calibrator Levels (pg/mL)**

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**

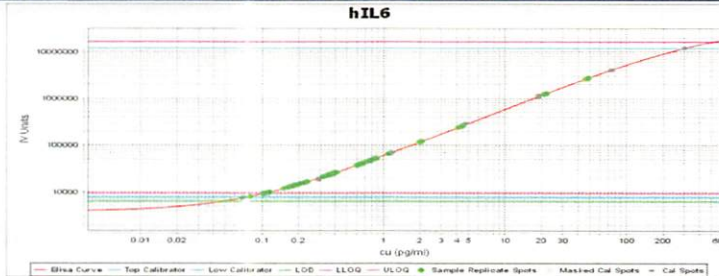
**hIFNγ**



**Calibrator Levels (pg/mL)**

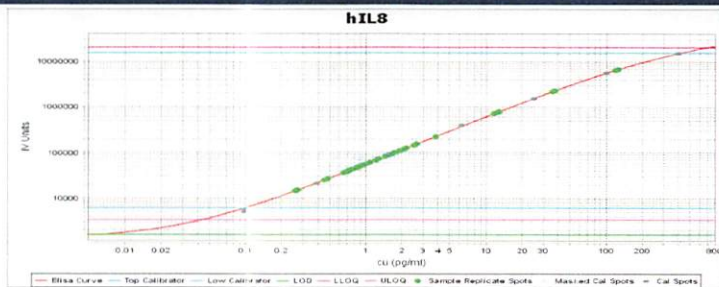
1	50
2	13
3	3.13
4	0.781
5	0.195
6	0.049
7	0.012
8	0.000

**IL-6 Calibration Curve**



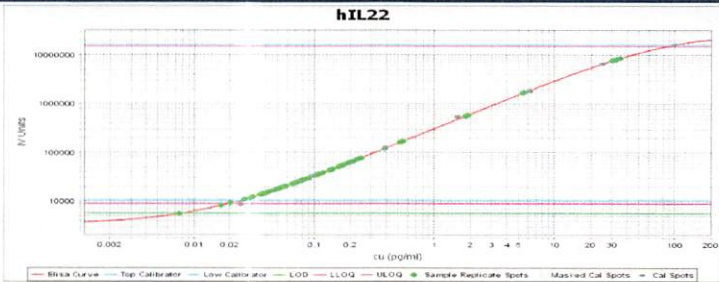
Calibrator Levels (pg/mL)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

**IL-8 Calibration Curve**



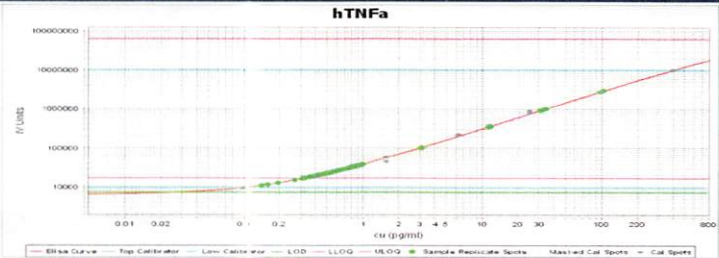
Calibrator Levels (pg/mL)	
1	400
2	100
3	25.00
4	6.250
5	1.563
6	0.391
7	0.098
8	0.000

**IL-22 Calibration Curve**



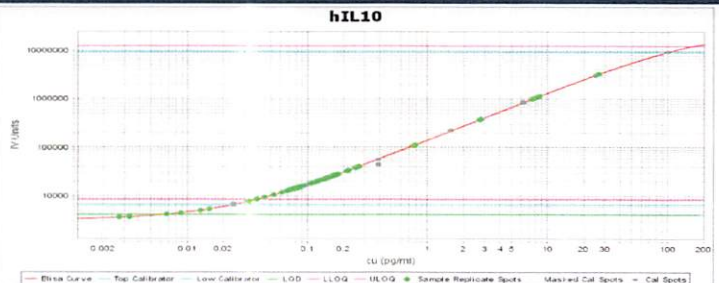
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

**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