

Product Number:	85-0329
Lot Number:	319497
Expiration:	3-Feb-2024
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	319352P
Biotinylated Antibody Rgt	98-0329	319352B
Sample Diluent 21	1863329	319172
IL-12p70 Calibrator	18910034	318756
IL-1 $\beta$ Calibrator	18910040	319165
IL-4 Calibrator	1858452	319016
IL-5 Calibrator	1858462	318837
IFN $\gamma$ Calibrator	1858260	318898
IL-6 Calibrator	18910044	318724
IL-8 Calibrator	1858454	319011
IL-22 Calibrator	18910036	318813
TNF $\alpha$ Calibrator	1880012	318883
IL-10 Calibrator	18910043	318755
Wash Buffer (12.5X) <sup>1</sup>	1863332	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	158.7	131.8 - 233.4
Panel 2	318935	51.89	39.35 - 72.53
Panel 3	318935	15.87	14.88 - 23.02
Panel 4	318935	4.710	4.159 - 6.457
Panel 5	318935	1.653	1.157 - 2.321

**IL-1 $\beta$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	18.24	12.01 - 32.00
Panel 2	318935	6.083	3.739 - 9.953
Panel 3	318935	2.021	1.270 - 3.351
Panel 4	318935	0.605	0.387 - 0.986
Panel 5	318935	0.199	0.111 - 0.348

**IL-4**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	65.29	51.07 - 80.47
Panel 2	318935	20.62	14.40 - 24.06
Panel 3	318935	7.524	5.421 - 8.931
Panel 4	318935	2.319	1.585 - 2.879
Panel 5	318935	0.803	0.478 - 1.311

**IL-5**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	85.52	47.13 - 109.0
Panel 2	318935	29.15	15.94 - 32.60
Panel 3	318935	10.13	5.192 - 11.65
Panel 4	318935	3.272	1.649 - 3.323
Panel 5	318935	1.053	0.535 - 1.226

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	10.90	7.193 - 14.98
Panel 2	318935	3.784	1.985 - 4.964
Panel 3	318935	1.283	0.714 - 1.703
Panel 4	318935	0.403	0.249 - 0.513
Panel 5	318935	0.176	0.103 - 0.227

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	172.5	122.5 - 249.4
Panel 2	318935	81.42	57.98 - 109.1
Panel 3	318935	16.44	12.14 - 22.22
Panel 4	318935	7.554	5.891 - 9.589
Panel 5	318935	2.271	1.746 - 3.433

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	561.1	285.9 - 755.9
Panel 2	318935	166.1	89.15 - 204.6
Panel 3	318935	62.99	31.29 - 74.91
Panel 4	318935	17.72	10.10 - 19.60
Panel 5	318935	6.063	3.241 - 7.262

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	110.9	61.61 - 225.7
Panel 2	318935	20.68	11.69 - 35.93
Panel 3	318935	6.315	4.089 - 11.05
Panel 4	318935	1.871	1.243 - 2.815
Panel 5	318935	0.582	0.443 - 0.957

TNF $\alpha$

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	337.7	189.4 - 566.2
Panel 2	318935	115.6	66.23 - 174.3
Panel 3	318935	40.93	25.41 - 58.84
Panel 4	318935	11.86	7.756 - 16.68
Panel 5	318935	3.667	2.063 - 5.792

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	101.5	75.66 - 129.7
Panel 2	318935	32.18	21.68 - 40.39
Panel 3	318935	11.62	7.652 - 14.16
Panel 4	318935	3.398	2.387 - 4.063
Panel 5	318935	1.130	0.777 - 1.557

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

Review/Approval

Robert Somogyi

Quality Control  
Manager

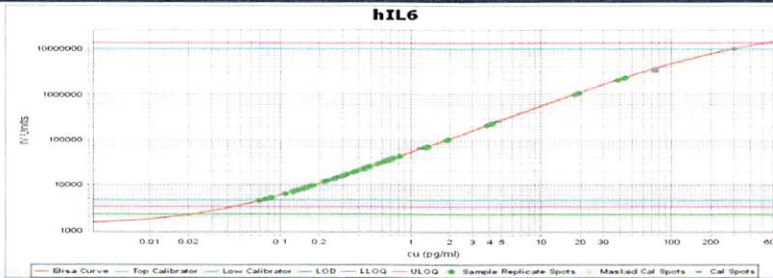
Name

Title

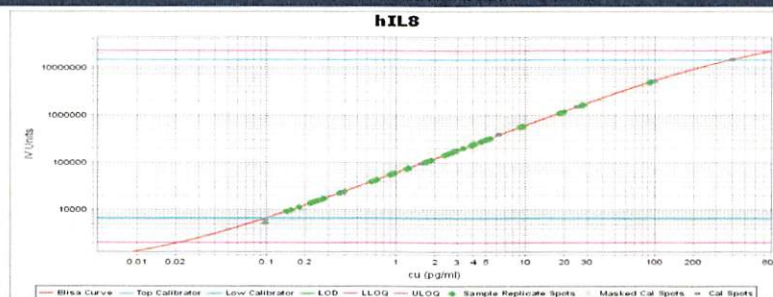
Signature/Date

**IL-6 Calibration Curve**



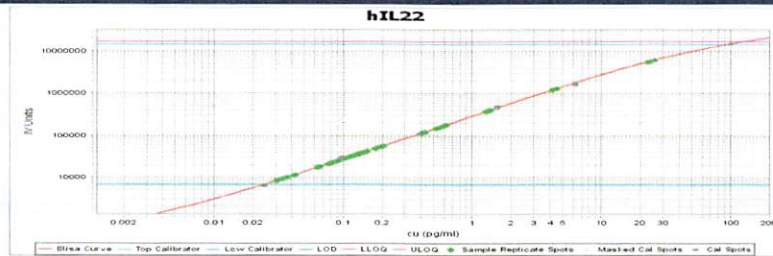
Calibrator Levels (pg/mL)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

**IL-8 Calibration Curve**



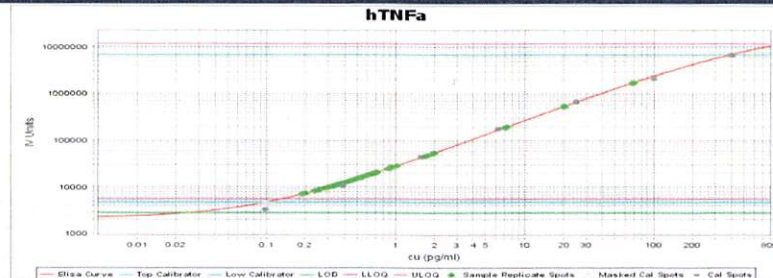
Calibrator Levels (pg/mL)	
1	400
2	100
3	25.00
4	6.250
5	1.563
6	0.391
7	0.098
8	0.000

**IL-22 Calibration Curve**



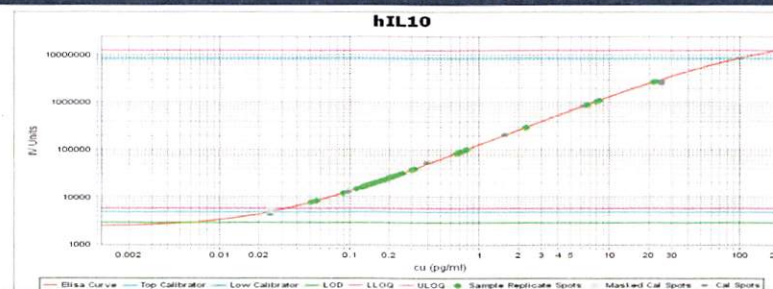
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

**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

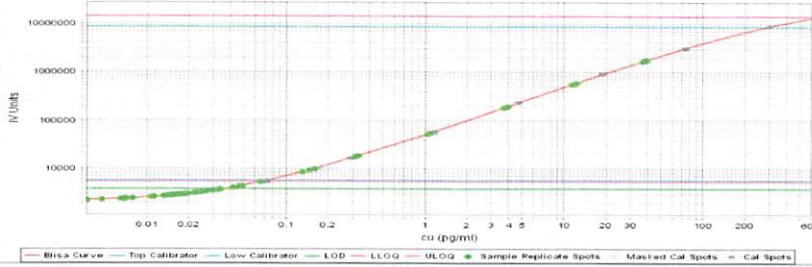
**IL-10 Calibration Curve**



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

**IL-12p70 Calibration Curve**

**hIL12p70**

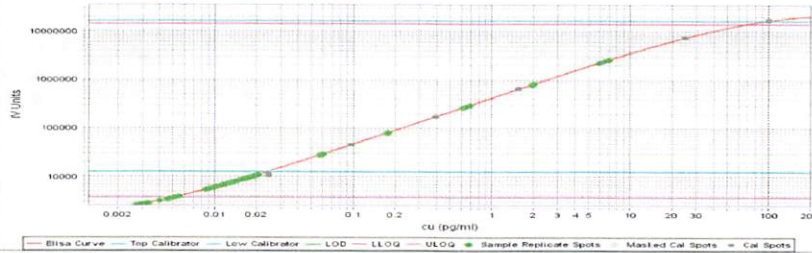


**Calibrator Levels  
(pg/ml)**

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**

**hIL1β**

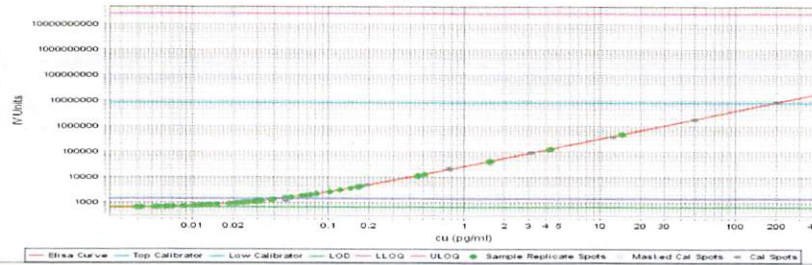


**Calibrator Levels  
(pg/mL)**

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**

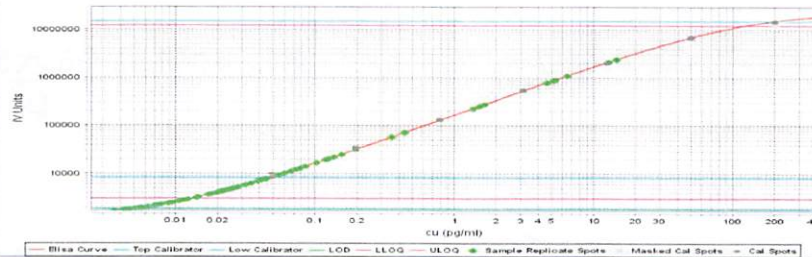


**Calibrator Levels  
(pg/mL)**

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**

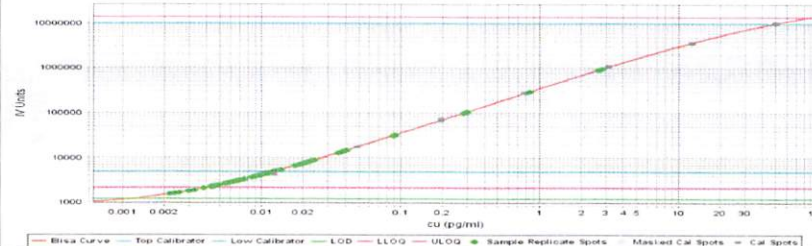


**Calibrator Levels  
(pg/mL)**

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	12.5
3	3.13
4	0.781
5	0.195
6	0.049
7	0.012
8	0.000