

Product Number:	85-0329
Lot Number:	318783
Expiration:	25-Oct-2021
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	318182P
Biotinylated Antibody Rgt	98-0329	318782B
Sample Diluent 21 <sup>1</sup>	1863329	N/A
IL-12p70 Calibrator	18910034	318629
IL-1β Calibrator	18910040	318673
IL-4 Calibrator	1858452	318480
IL-5 Calibrator	1858462	318465
IFNγ Calibrator	1858260	318631
IL-6 Calibrator	18910044	318724
IL-8 Calibrator	1858454	318500
IL-22 Calibrator	18910036	318408
TNFα Calibrator	1880012	318436
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	318641	168.9	148.9 - 240.7
Panel 2	318641	63.79	51.04 - 81.68
Panel 3	318641	22.03	17.95 - 26.93
Panel 4	318641	7.242	5.525 - 9.670
Panel 5	318641	2.234	1.672 - 3.404

**IL-1β**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	23.82	19.65 - 29.48
Panel 2	318641	8.389	6.475 - 10.24
Panel 3	318641	2.914	2.123 - 3.489
Panel 4	318641	1.012	0.238 - 1.776
Panel 5	318641	0.351	0.223 - 0.460

**IL-4**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	83.21	66.79 - 105.1
Panel 2	318641	26.60	20.41 - 35.27
Panel 3	318641	9.984	6.722 - 12.50
Panel 4	318641	3.374	2.180 - 4.624
Panel 5	318641	1.292	0.815 - 1.597

**IL-5**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	64.73	50.55 - 76.98
Panel 2	318641	22.15	16.93 - 25.93
Panel 3	318641	7.992	5.920 - 9.298
Panel 4	318641	2.607	1.943 - 3.293
Panel 5	318641	0.943	0.700 - 1.238

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	318641	12.43	8.541	- 12.86
Panel 2	318641	4.415	2.940	- 4.681
Panel 3	318641	1.652	0.994	- 1.686
Panel 4	318641	0.607	0.348	- 0.647
Panel 5	318641	0.236	0.138	- 0.270

**IL-6**

	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Panel 1	318641	121.3	97.18	- 184.2
Panel 2	318641	3.809	2.780	- 7.433
Panel 3	318641	1.236	0.928	- 2.564
Panel 4	318641	0.411	0.330	- 1.033

**IL-8**

	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Panel 1	318641	470.6	372.9	- 559.4
Panel 2	318641	158.0	109.4	- 188.6
Panel 3	318641	51.83	36.67	- 64.61
Panel 4	318641	20.03	11.28	- 24.27
Panel 5	318641	6.205	3.867	- 7.799

**IL-22**

	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Panel 1	318641	20.40	14.24	- 33.29
Panel 2	318641	6.481	4.355	- 9.706
Panel 3	318641	2.286	1.528	- 3.283
Panel 4	318641	0.744	0.462	- 1.223
Panel 5	318641	0.269	0.156	- 0.428

**TNF $\alpha$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Panel 1	318641	71.36	59.79	- 100.9
Panel 2	318641	24.16	20.44	- 35.53
Panel 3	318641	8.588	6.956	- 12.29
Panel 4	318641	3.007	2.303	- 4.312
Panel 5	318641	0.926	0.547	- 1.821

**IL-10**

	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Panel 1	318641	66.48	48.59	- 73.50
Panel 2	318641	23.44	15.68	- 27.33
Panel 3	318641	8.371	5.500	- 9.707
Panel 4	318641	3.012	2.011	- 3.603
Panel 5	318641	1.040	0.706	- 1.339

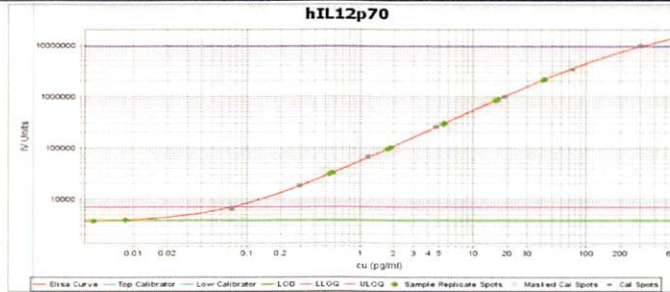
Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Review/Approval

Shivani Goel  
Name Title  
Manager, QC

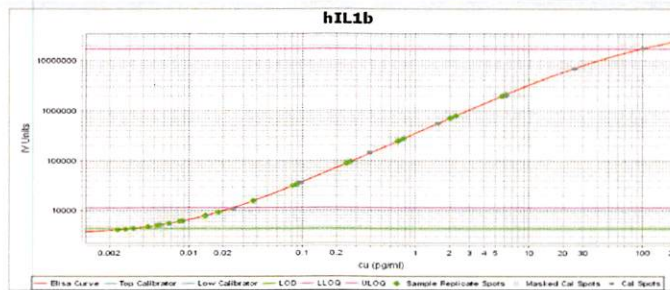
Shivani Goel 2/10/2020  
Signature/Date

**IL-12p70 Calibration Curve**



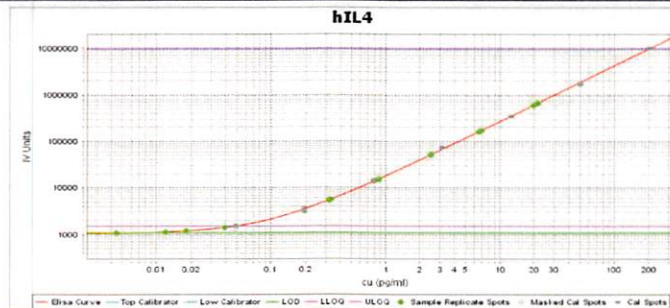
Calibrator Levels (pg/ml)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

**IL-1β Calibration Curve**



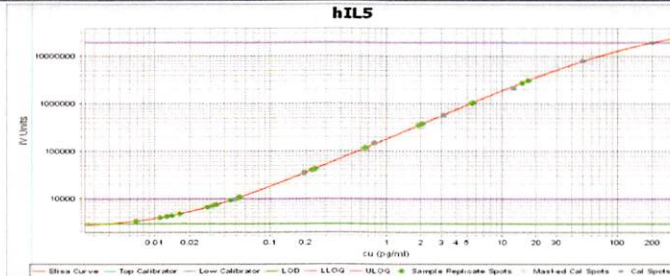
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

**IL-4 Calibration Curve**



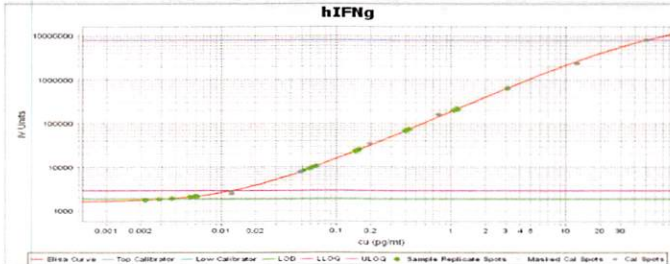
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

**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

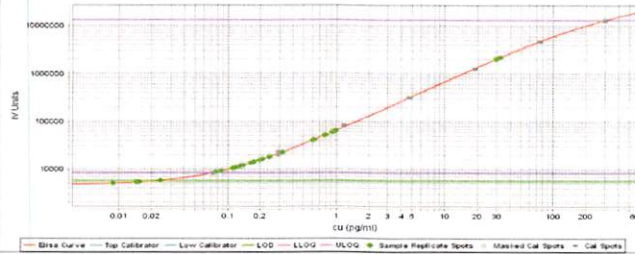
**IFNγ Calibration Curve**



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**

**hIL6**

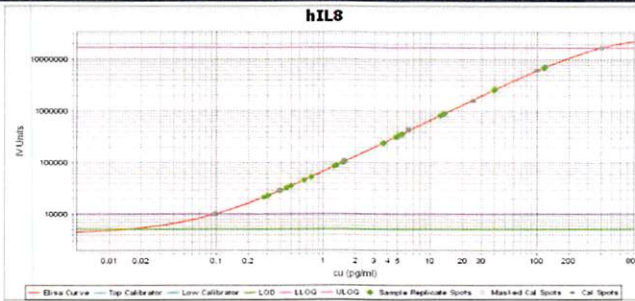


**Calibrator Levels (pg/mL)**

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**

**hIL8**

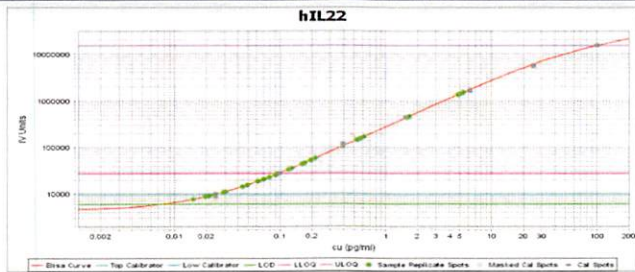


**Calibrator Levels (pg/mL)**

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**

**hIL22**

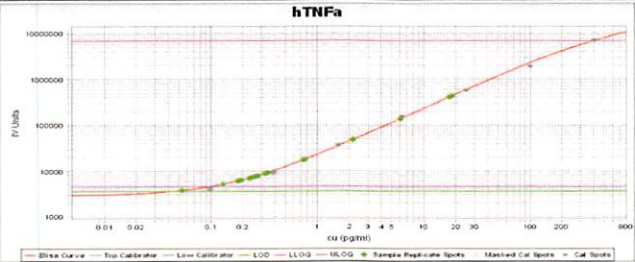


**Calibrator Levels (pg/mL)**

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**

**hTNFα**

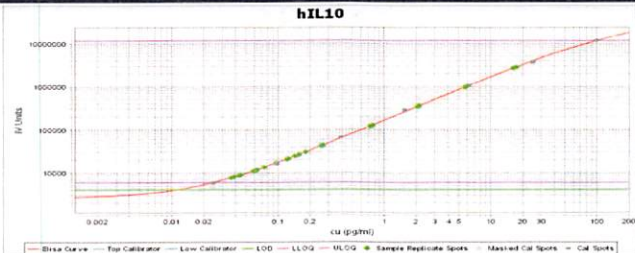


**Calibrator Levels (pg/mL)**

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**

**hIL10**



**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