

Product Number:	85-0329
Lot Number:	318839
Expiration:	27-Nov-2021
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	318839P
Biotinylated Antibody Rgt	98-0329	318839B
Sample Diluent 2l <sup>1</sup>	1863329	N/A
IL-12p70 Calibrator	18910034	318629
IL-1β Calibrator	18910040	318728
IL-4 Calibrator	1858452	318480
IL-5 Calibrator	1858462	318465
IFNγ Calibrator	1858260	318538
IL-6 Calibrator	18910044	318583
IL-8 Calibrator	1858454	318342
IL-22 Calibrator	18910036	318408
TNFα Calibrator	1880012	318807
IL-10 Calibrator	18910043	318630
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

<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).



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

**Release Materials**

**IL-12p70**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	164.3	148.5 - 237.3
Panel 2	318641	53.70	50.80 - 80.57
Panel 3	318641	20.22	17.83 - 26.74
Panel 4	318641	8.087	5.565 - 9.512
Panel 5	318641	2.728	1.710 - 3.318

**IL-1β**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	19.17	17.97 - 29.91
Panel 2	318641	6.423	5.806 - 10.49
Panel 3	318641	2.331	1.923 - 3.568
Panel 4	318641	0.879	0.600 - 1.285
Panel 5	318641	0.285	0.211 - 0.459

**IL-4**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	83.78	65.00 - 104.1
Panel 2	318641	25.64	20.30 - 34.72
Panel 3	318641	9.441	6.646 - 12.28
Panel 4	318641	3.709	2.094 - 4.553
Panel 5	318641	1.287	0.814 - 1.563

**IL-5**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	64.81	50.70 - 76.78
Panel 2	318641	21.45	17.10 - 25.74
Panel 3	318641	7.594	5.870 - 9.215
Panel 4	318641	2.925	1.866 - 3.322
Panel 5	318641	1.028	0.662 - 1.245

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	10.55	8.584 - 12.88
Panel 2	318641	3.590	2.915 - 4.650
Panel 3	318641	1.247	1.005 - 1.662
Panel 4	318641	0.534	0.348 - 0.641
Panel 5	318641	0.203	0.143 - 0.264

**IL-6**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	126.9	96.92 - 180.1
Panel 2	318641	4.189	2.328 - 7.392
Panel 3	318641	1.474	0.715 - 2.582
Panel 4	318641	0.637	0.223 - 1.047

**IL-8**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	484.0	367.9 - 561.8
Panel 2	318641	163.3	111.2 - 187.0
Panel 3	318641	54.85	37.26 - 64.06
Panel 4	318641	22.35	11.52 - 23.94
Panel 5	318641	6.761	3.937 - 7.652

**IL-22**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	20.66	13.41 - 32.67
Panel 2	318641	5.673	3.573 - 9.829
Panel 3	318641	1.983	1.262 - 3.329
Panel 4	318641	0.776	0.353 - 1.238
Panel 5	318641	0.242	0.123 - 0.430

**TNF $\alpha$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	60.69	54.09 - 101.5
Panel 2	318641	20.75	18.91 - 35.47
Panel 3	318641	6.664	6.132 - 12.45
Panel 4	318641	2.636	2.078 - 4.324
Panel 5	318641	0.876	0.540 - 1.765

**IL-10**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	70.69	48.85 - 74.86
Panel 2	318641	23.92	16.26 - 27.13
Panel 3	318641	8.455	5.725 - 9.560
Panel 4	318641	3.412	2.075 - 3.556
Panel 5	318641	1.152	0.738 - 1.323

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

Review/Approval

Patrizia Stadler

Associate  
Director of QC

*P. Stadler*

16 OCT 2020

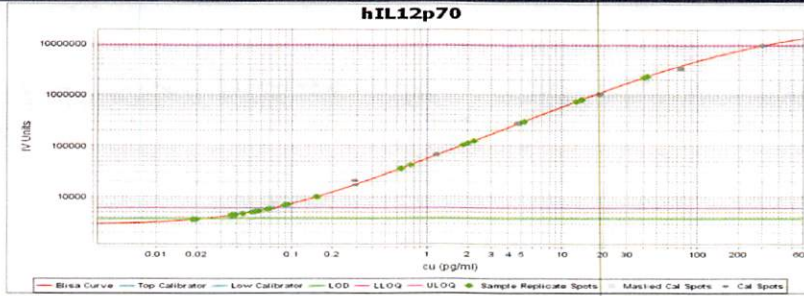
Name

Title

Signature/Date

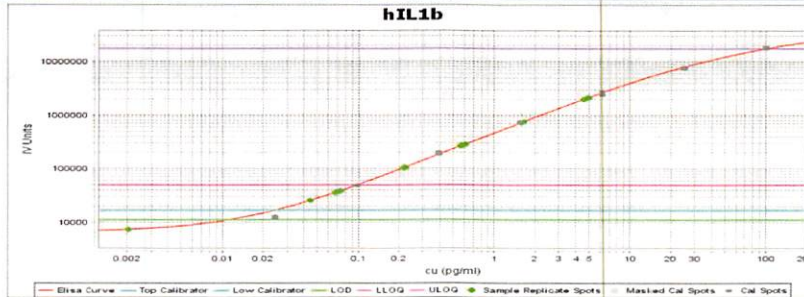


**IL-12p70 Calibration Curve**



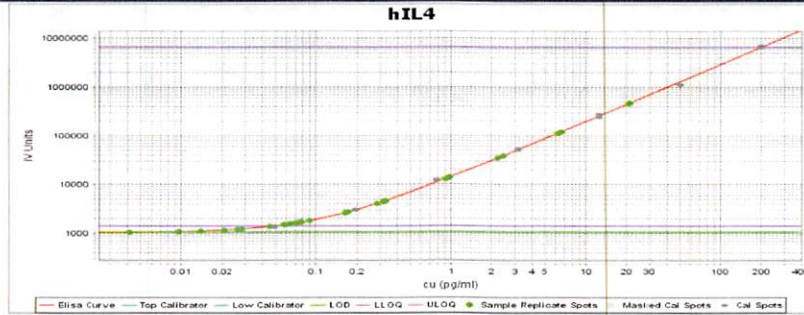
Calibrator Levels (pg/ml)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

**IL-1β Calibration Curve**



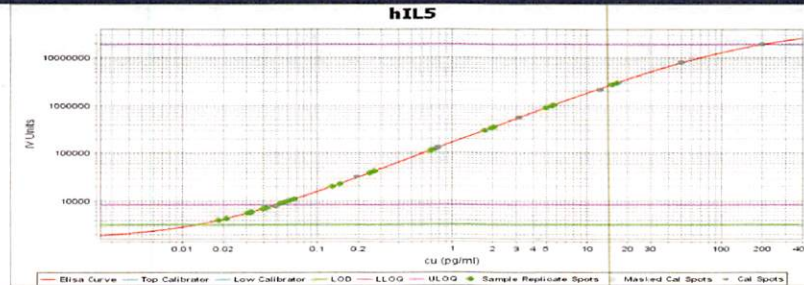
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

**IL-4 Calibration Curve**



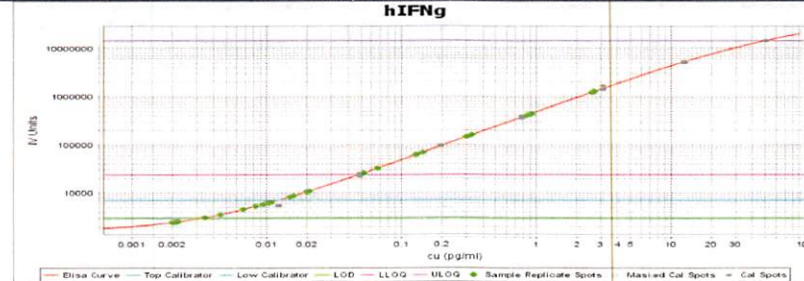
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

**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

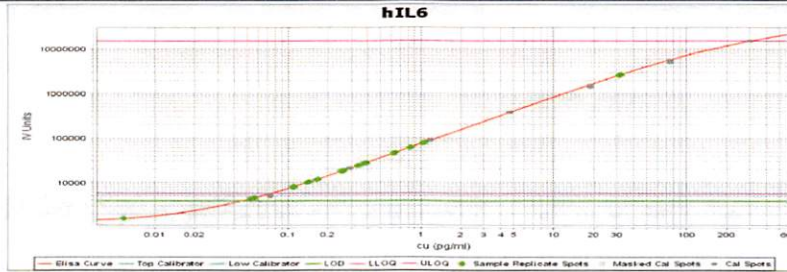
**IFNγ Calibration Curve**



Calibrator Levels (pg/mL)	
1	50
2	13
3	3.13
4	0.781
5	0.195
6	0.049
7	0.012
8	0.000



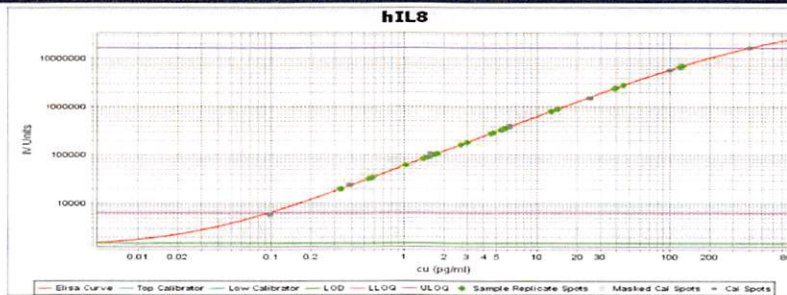
**IL-6 Calibration Curve**



**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

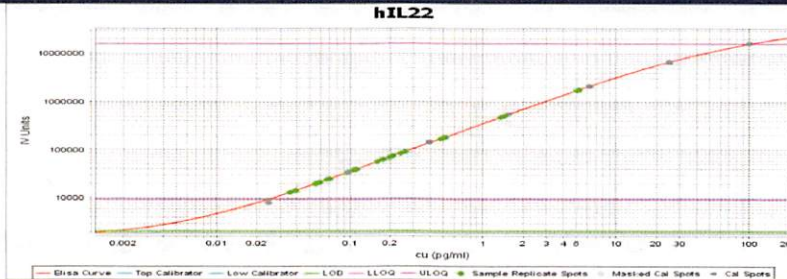
**IL-8 Calibration Curve**



**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

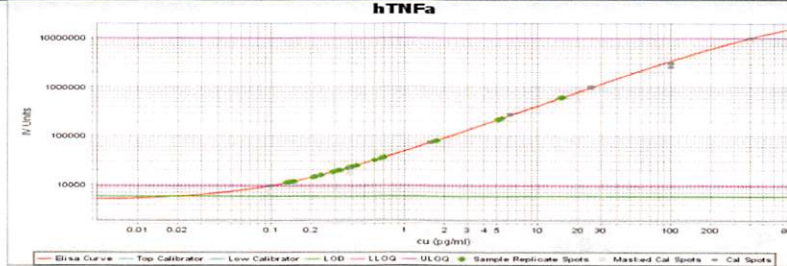
**IL-22 Calibration Curve**



**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

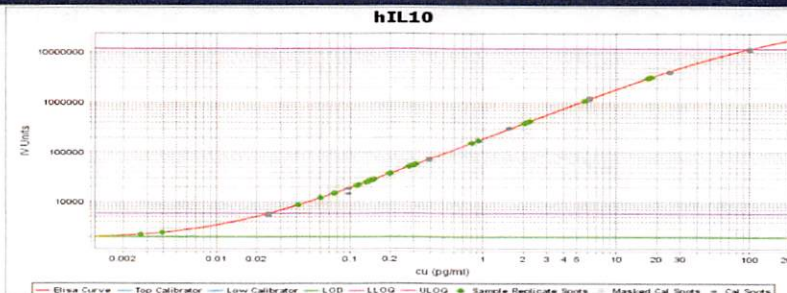
**TNFα Calibration Curve**



**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