

Product Number:	85-0329
Lot Number:	319279
Expiration:	12-Nov-2022
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	319279P
Biotinylated Antibody Rgt	98-0329	319279B
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	193.1	140.6 - 239.0
Panel 2	318935	50.13	44.36 - 73.70
Panel 3	318935	18.13	15.74 - 23.61
Panel 4	318935	5.078	4.400 - 6.601
Panel 5	318935	1.663	1.243 - 2.397

**IL-1 $\beta$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	23.41	16.60 - 33.00
Panel 2	318935	6.828	5.060 - 10.28
Panel 3	318935	2.412	1.780 - 3.430
Panel 4	318935	0.709	0.496 - 1.030
Panel 5	318935	0.252	0.151 - 0.367

**IL-4**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	70.63	53.27 - 82.46
Panel 2	318935	19.49	16.02 - 24.04
Panel 3	318935	7.311	6.011 - 9.016
Panel 4	318935	2.491	1.729 - 2.776
Panel 5	318935	1.192	0.434 - 1.312

**IL-5**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	72.81	54.07 - 121.1
Panel 2	318935	20.56	18.39 - 34.26
Panel 3	318935	7.600	6.909 - 11.42
Panel 4	318935	2.088	2.007 - 3.306
Panel 5	318935	0.784	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	11.86	9.152 - 15.15
Panel 2	318935	3.160	2.766 - 5.014
Panel 3	318935	1.144	1.021 - 1.670
Panel 4	318935	0.347	0.320 - 0.515
Panel 5	318935	0.152	0.142 - 0.223

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	182.1	140.8 - 249.6
Panel 2	318935	76.22	68.32 - 107.0
Panel 3	318935	15.18	14.11 - 21.96
Panel 4	318935	6.902	6.379 - 9.569
Panel 5	318935	2.333	1.951 - 3.380

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	477.5	287.6 - 812.3
Panel 2	318935	125.3	99.01 - 213.4
Panel 3	318935	47.69	33.87 - 79.01
Panel 4	318935	13.67	10.72 - 20.44
Panel 5	318935	4.337	3.706 - 7.504

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	135.5	73.64 - 264.5
Panel 2	318935	20.81	13.35 - 41.63
Panel 3	318935	6.872	6.048 - 11.67
Panel 4	318935	1.803	1.025 - 3.238
Panel 5	318935	0.669	0.478 - 1.052

TNF $\alpha$

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	408.0	213.6 - 599.8
Panel 2	318935	115.5	81.39 - 179.7
Panel 3	318935	42.01	30.07 - 60.48
Panel 4	318935	11.77	9.413 - 16.87
Panel 5	318935	3.888	2.187 - 6.149

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	105.0	86.85 - 130.3
Panel 2	318935	30.60	25.35 - 40.54
Panel 3	318935	10.53	9.246 - 13.87
Panel 4	318935	3.196	2.649 - 4.032
Panel 5	318935	1.186	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 17 Jun 2021

Name

Title

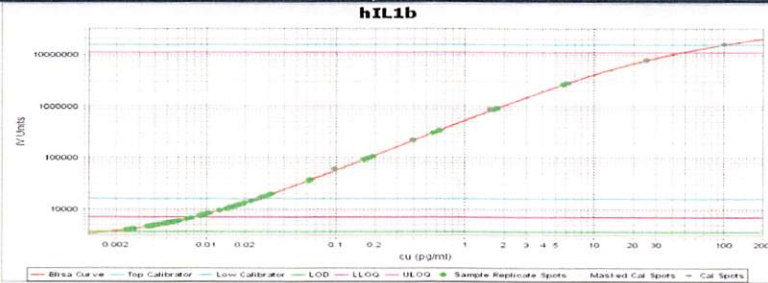
Signature/Date

**IL-12p70 Calibration Curve**



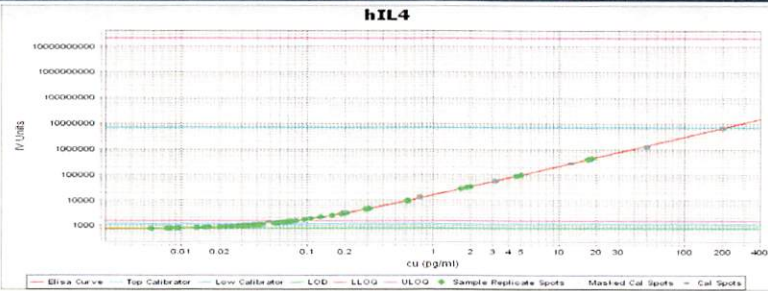
Calibrator Levels (pg/ml)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

**IL-1 $\beta$  Calibration Curve**



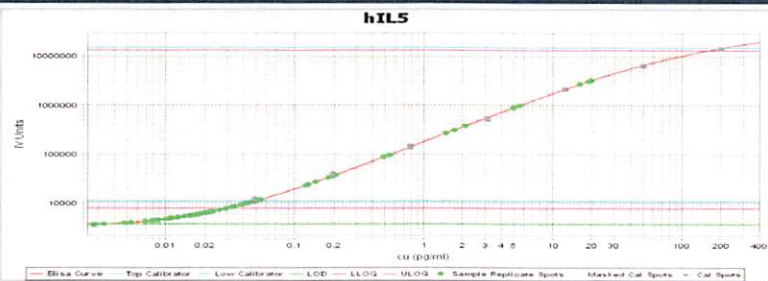
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

**IL-4 Calibration Curve**



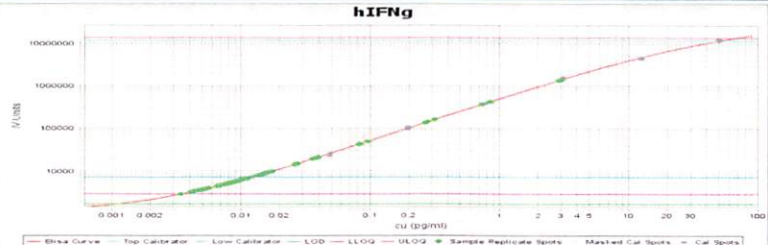
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

**IL-5 Calibration Curve**



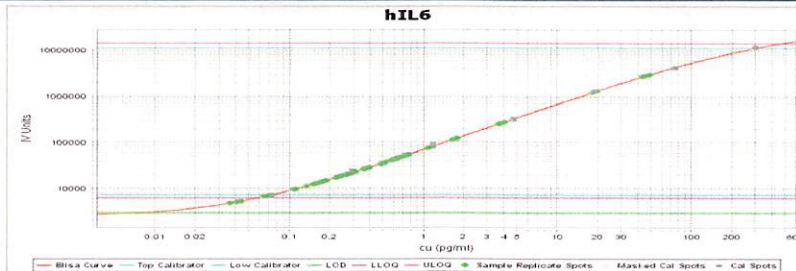
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

**IFN $\gamma$  Calibration Curve**



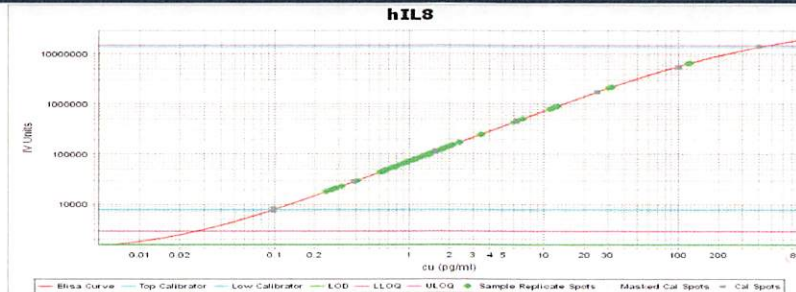
Calibrator Levels (pg/mL)	
1	50
2	12.5
3	3.13
4	0.781
5	0.195
6	0.049
7	0.012
8	0.000

**IL-6 Calibration Curve**



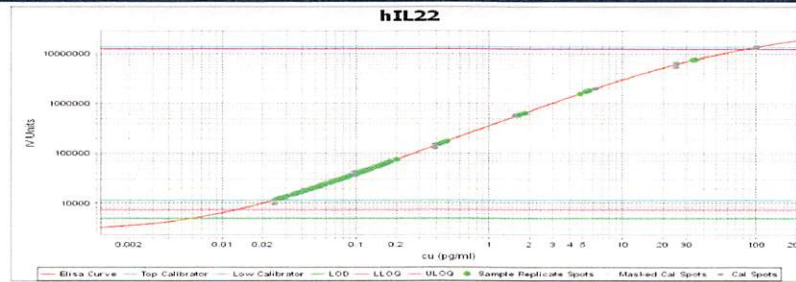
Calibrator Levels (pg/mL)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

**IL-8 Calibration Curve**



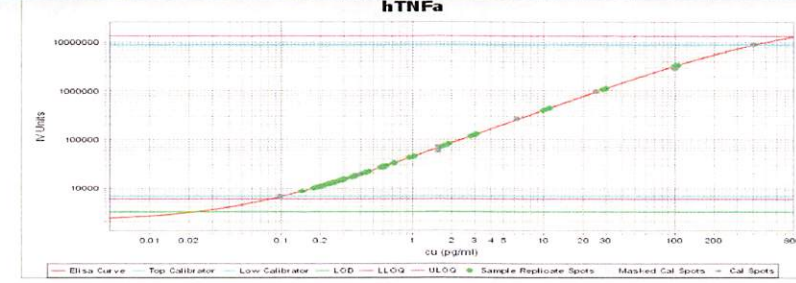
Calibrator Levels (pg/mL)	
1	400
2	100
3	25.00
4	6.250
5	1.563
6	0.391
7	0.098
8	0.000

**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