

Product Number:	85-0329
Lot Number:	319827
Expiration:	23-Dec-2023
Platform(s) :	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	319826P
Biotinylated Antibody Rgt	98-0329	319740B
Sample Diluent 21	1863329	319709
IL-12p70 Calibrator	18910034	319156
IL-1β Calibrator	18910040	319564
IL-4 Calibrator	1858452	319458
IL-5 Calibrator	1858462	318837
IFNγ Calibrator	1858260	319711
IL-6 Calibrator	18910044	319484
IL-8 Calibrator	1858454	319575
IL-22 Calibrator	18910036	319459
TNFα Calibrator	1880012	319247
IL-10 Calibrator	18910043	319215
Wash Buffer (25X) ¹	1860055	N/A
Streptavidin-HRP Reagent ¹	1858263	N/A
Super Signal® Luminol* ¹	1858089	N/A
Super Signal® Peroxide* ¹	1858090	N/A



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

¹ Reagents are not Kit Lot Specific

*SuperSignal® 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 Cytokine1 Kit Instructions (85A-0329).

Release Materials

IL-12p70

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	149.3	131.7 - 203.5
Panel 2	318935	47.90	36.98 - 66.47
Panel 3	318935	16.50	13.40 - 22.07
Panel 4	318935	4.303	3.751 - 6.122
Panel 5	318935	1.432	1.179 - 2.080

IL-1β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	23.65	12.91 - 32.59
Panel 2	318935	7.301	4.217 - 9.740
Panel 3	318935	2.523	1.377 - 3.499
Panel 4	318935	0.684	0.424 - 0.975
Panel 5	318935	0.235	0.137 - 0.343

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	56.36	50.84 - 77.00
Panel 2	318935	16.51	14.18 - 23.93
Panel 3	318935	5.914	5.174 - 8.891
Panel 4	318935	1.674	1.422 - 2.789
Panel 5	318935	0.588	0.410 - 1.241

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	96.67	43.55 - 104.5
Panel 2	318935	29.55	13.38 - 33.50
Panel 3	318935	10.10	4.766 - 11.68
Panel 4	318935	2.901	1.234 - 3.623
Panel 5	318935	1.036	0.470 - 1.263

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

Release Materials

IFN γ

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	10.77	8.245 - 14.41
Panel 2	318935	3.297	2.401 - 4.768
Panel 3	318935	1.134	0.853 - 1.722
Panel 4	318935	0.346	0.271 - 0.505
Panel 5	318935	0.146	0.114 - 0.221

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	166.3	143.3 - 215.0
Panel 2	318935	75.97	62.18 - 101.2
Panel 3	318935	16.03	12.91 - 20.83
Panel 4	318935	6.963	6.002 - 9.003
Panel 5	318935	2.452	1.966 - 3.145

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	455.4	363.4 - 600.3
Panel 2	318935	127.0	91.56 - 189.5
Panel 3	318935	44.77	32.31 - 67.04
Panel 4	318935	12.27	9.930 - 18.44
Panel 5	318935	4.183	3.310 - 6.556

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	72.43	47.37 - 159.2
Panel 2	318935	11.86	6.880 - 28.17
Panel 3	318935	4.062	2.175 - 9.086
Panel 4	318935	1.047	0.686 - 2.605
Panel 5	318935	0.390	0.216 - 0.878

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	305.8	251.0 - 450.4
Panel 2	318935	98.95	69.11 - 155.1
Panel 3	318935	37.00	25.96 - 51.54
Panel 4	318935	9.914	7.461 - 14.87
Panel 5	318935	3.219	2.349 - 4.860

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	85.76	75.65 - 125.4
Panel 2	318935	26.45	21.66 - 39.42
Panel 3	318935	9.389	7.643 - 13.76
Panel 4	318935	2.503	2.344 - 3.891
Panel 5	318935	0.934	0.826 - 1.448

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

Review/Approval

Tony Iem

Quality Control
Supervisor

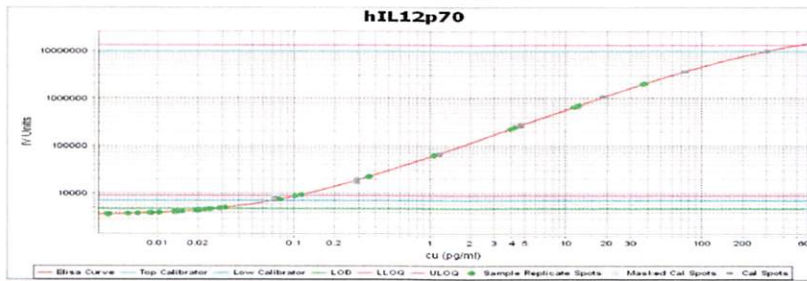
 02/08/2022

Name

Title

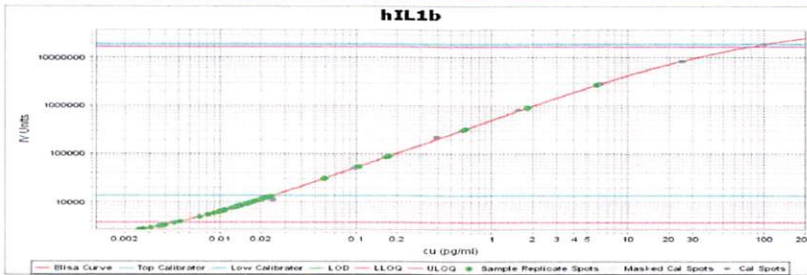
Signature/Date

IL-12p70 Calibration Curve



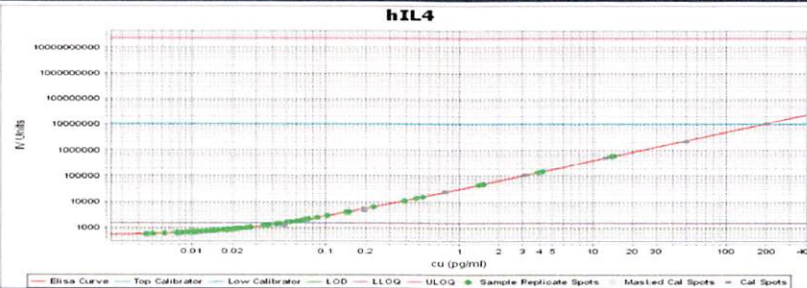
Calibrator Levels (pg/ml)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

IL-1 β Calibration Curve



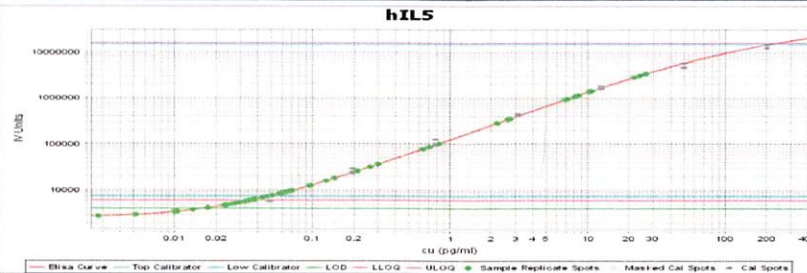
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

IL-4 Calibration Curve



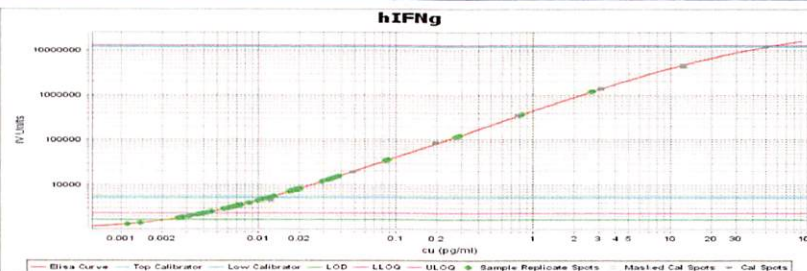
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

IL-5 Calibration Curve



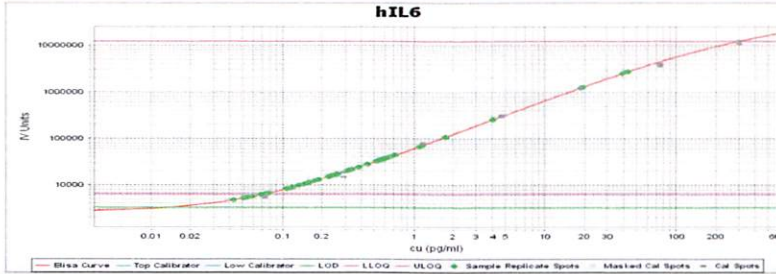
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

IFN γ Calibration Curve



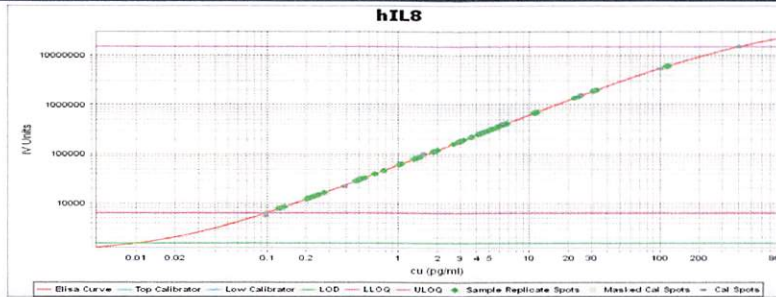
Calibrator Levels (pg/mL)	
1	50
2	12.5
3	3.13
4	0.781
5	0.195
6	0.049
7	0.012
8	0.000

IL-6 Calibration Curve



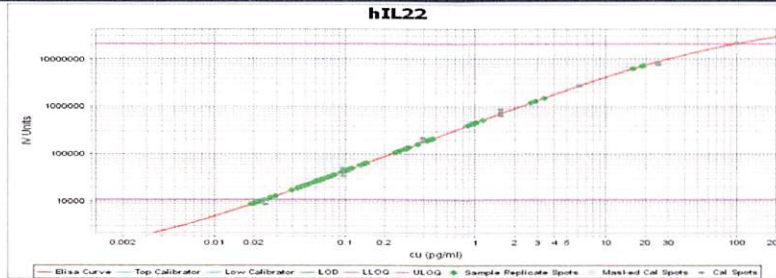
Calibrator Levels (pg/mL)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

IL-8 Calibration Curve



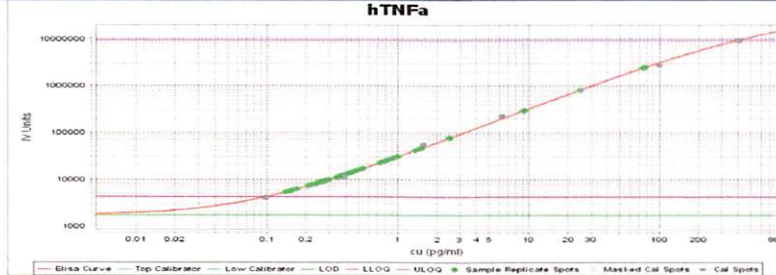
Calibrator Levels (pg/mL)	
1	400
2	100
3	25.00
4	6.250
5	1.563
6	0.391
7	0.098
8	0.000

IL-22 Calibration Curve



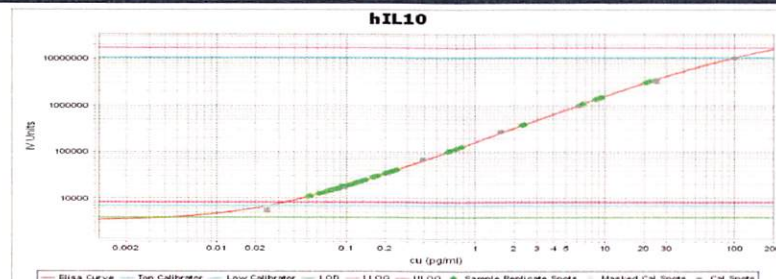
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

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