

Product Number:	85-0329
Lot Number:	318939
Expiration:	26-Dec-2021
Platform(s) :	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	318938P
Biotinylated Antibody Rgt	98-0329	318893B
Sample Diluent 21	1863329	318865
IL-12p70 Calibrator	18910034	318756
IL-1 β Calibrator	18910040	318728
IL-4 Calibrator	1858452	318580
IL-5 Calibrator	1858462	318837
IFN γ Calibrator	1858260	318886
IL-6 Calibrator	18910044	318724
IL-8 Calibrator	1858454	318500
IL-22 Calibrator	18910036	318813
TNF α Calibrator	1880012	318883
IL-10 Calibrator	18910043	318630
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 Cytokine Kit Instructions (85A-0329).

Release Materials

IL-12p70

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	161.3	150.4 - 229.1
Panel 2	318935	53.98	47.22 - 70.84
Panel 3	318935	17.44	15.74 - 23.61
Panel 4	318935	5.349	4.400 - 6.601
Panel 5	318935	1.513	1.358 - 2.282

IL-1 β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	21.76	18.21 - 31.34
Panel 2	318935	7.393	5.580 - 9.76
Panel 3	318935	2.415	1.947 - 3.263
Panel 4	318935	0.755	0.649 - 0.978
Panel 5	318935	0.254	0.172 - 0.345

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	67.30	54.29 - 81.44
Panel 2	318935	21.10	16.02 - 24.04
Panel 3	318935	7.472	6.011 - 9.016
Panel 4	318935	2.165	1.802 - 2.703
Panel 5	318935	0.844	0.522 - 1.224

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	81.79	60.78 - 114.1
Panel 2	318935	24.77	19.98 - 32.68
Panel 3	318935	8.63	7.332 - 11.00
Panel 4	318935	2.636	2.125 - 3.188
Panel 5	318935	0.887	0.681 - 1.205

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	10.35	9.723 - 14.58
Panel 2	318935	3.658	2.991 - 4.789
Panel 3	318935	1.154	1.077 - 1.615
Panel 4	318935	0.364	0.334 - 0.502
Panel 5	318935	0.144	0.146 - 0.219

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	158.1	140.8 - 249.6
Panel 2	318935	78.23	68.32 - 107.0
Panel 3	318935	15.95	14.11 - 21.96
Panel 4	318935	7.236	6.379 - 9.569
Panel 5	318935	2.351	1.951 - 3.380

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	454.9	340.0 - 759.8
Panel 2	318935	143.9	110.40 - 201.9
Panel 3	318935	52.20	38.39 - 74.49
Panel 4	318935	15.33	11.70 - 19.47
Panel 5	318935	5.429	4.086 - 7.124

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	141.4	92.73 - 245.5
Panel 2	318935	25.69	16.18 - 38.80
Panel 3	318935	9.088	6.610 - 11.10
Panel 4	318935	2.092	1.246 - 3.017
Panel 5	318935	0.819	0.535 - 0.994

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	304.2	252.2 - 561.2
Panel 2	318935	104.6	91.10 - 169.9
Panel 3	318935	38.63	33.11 - 57.44
Panel 4	318935	11.18	10.160 - 16.13
Panel 5	318935	3.234	2.584 - 5.753

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	100.4	86.85 - 130.3
Panel 2	318935	31.01	26.35 - 39.53
Panel 3	318935	10.39	9.246 - 13.87
Panel 4	318935	3.166	2.673 - 4.009
Panel 5	318935	1.145	0.955 - 1.480

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

Review/Approval

Shivani Goel

Quality Control
Manager

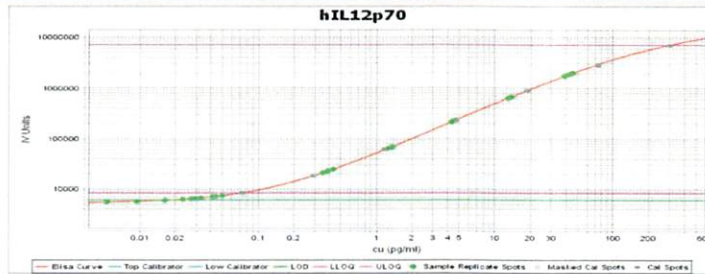
Shivani Goel 02Apr2021

Name

Title

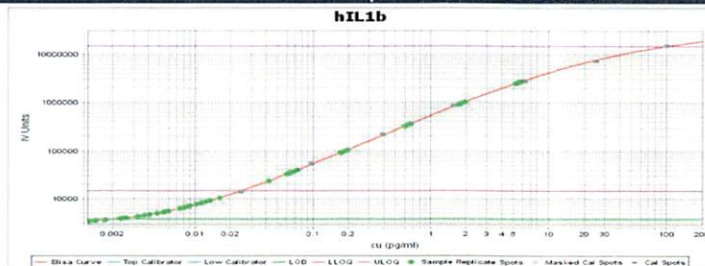
Signature/Date

IL-12p70 Calibration Curve



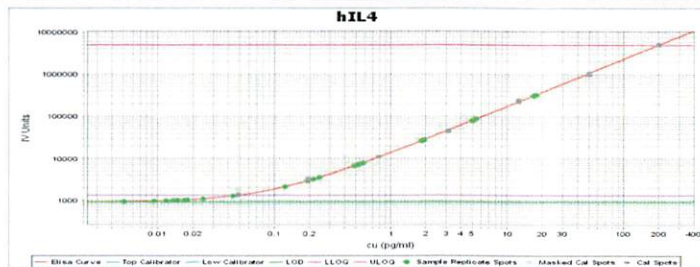
Calibrator Levels (pg/ml)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

IL-1 β Calibration Curve



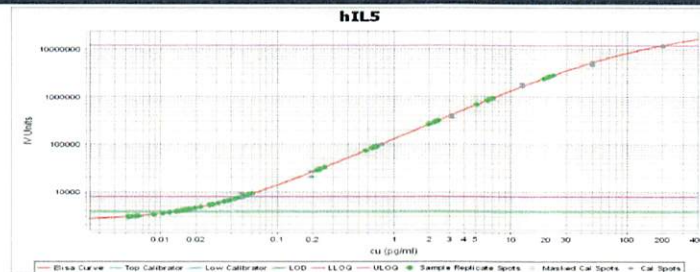
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

IL-4 Calibration Curve



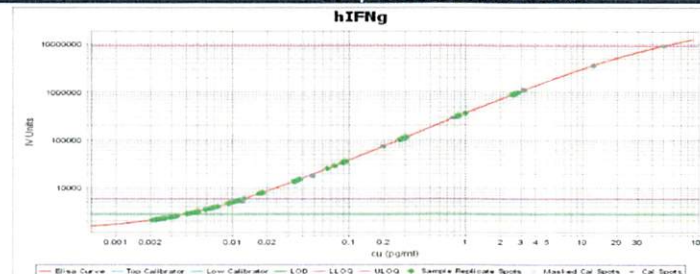
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

IL-5 Calibration Curve



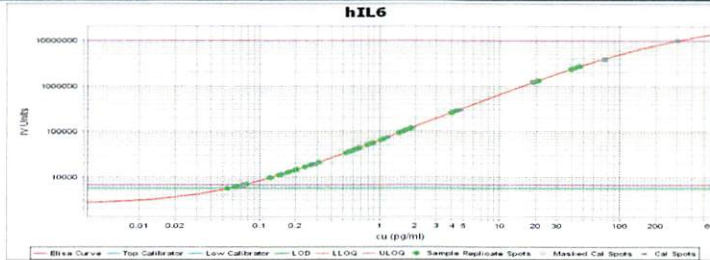
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

IFN γ Calibration Curve



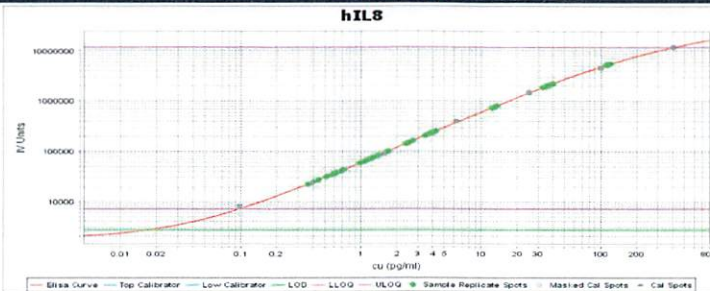
Calibrator Levels (pg/mL)	
1	50
2	13
3	3.13
4	0.781
5	0.195
6	0.049
7	0.012
8	0.000

IL-6 Calibration Curve



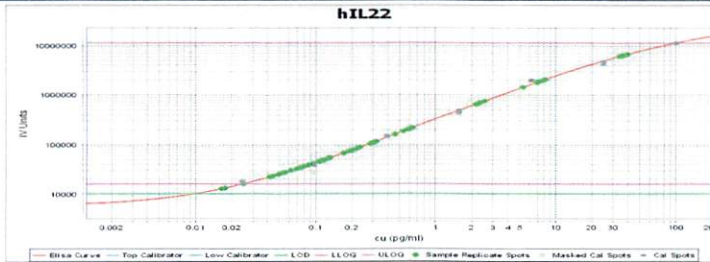
Calibrator Levels (pg/mL)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

IL-8 Calibration Curve



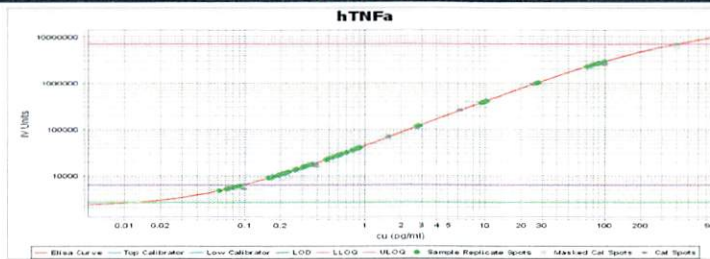
Calibrator Levels (pg/mL)	
1	400
2	100
3	25.00
4	6.250
5	1.563
6	0.391
7	0.098
8	0.000

IL-22 Calibration Curve



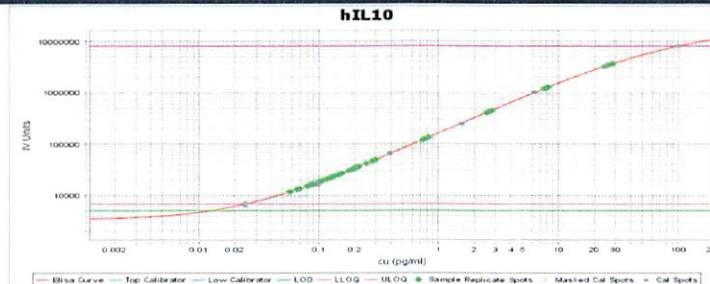
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

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