

Product Number:	85-0329
Lot Number:	319470
Expiration:	20-Oct-2022
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	319470P
Biotinylated Antibody Rgt	98-0329	319470B
Sample Diluent 21	1863329	319333
IL-12p70 Calibrator	18910034	318756
IL-1 β Calibrator	18910040	318728
IL-4 Calibrator	1858452	318480
IL-5 Calibrator	1858462	319223
IFN γ Calibrator	1858260	318898
IL-6 Calibrator	18910044	319173
IL-8 Calibrator	1858454	319011
IL-22 Calibrator	18910036	318813
TNF α Calibrator	1880012	318883
IL-10 Calibrator	18910043	318755
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	192.7	131.8 - 233.4
Panel 2	318935	59.86	39.35 - 72.53
Panel 3	318935	18.61	14.88 - 23.02
Panel 4	318935	5.674	4.159 - 6.457
Panel 5	318935	1.869	1.157 - 2.321

IL-1 β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	23.26	12.01 - 32.00
Panel 2	318935	7.208	3.739 - 9.953
Panel 3	318935	2.283	1.270 - 3.351
Panel 4	318935	0.680	0.387 - 0.986
Panel 5	318935	0.235	0.111 - 0.348

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	72.17	51.07 - 80.47
Panel 2	318935	20.22	14.40 - 24.06
Panel 3	318935	7.078	5.421 - 8.931
Panel 4	318935	2.439	1.585 - 2.879
Panel 5	318935	0.977	0.478 - 1.311

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	60.55	47.13 - 109.0
Panel 2	318935	18.21	15.94 - 32.60
Panel 3	318935	6.370	5.192 - 11.65
Panel 4	318935	1.863	1.649 - 3.323
Panel 5	318935	0.706	0.535 - 1.226

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	11.63	7.193 - 14.98
Panel 2	318935	3.442	1.985 - 4.964
Panel 3	318935	1.183	0.714 - 1.703
Panel 4	318935	0.373	0.249 - 0.513
Panel 5	318935	0.160	0.103 - 0.227

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	182.7	122.5 - 249.4
Panel 2	318935	82.05	57.98 - 109.1
Panel 3	318935	16.63	12.14 - 22.22
Panel 4	318935	7.611	5.891 - 9.589
Panel 5	318935	2.664	1.746 - 3.433

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	504.1	285.9 - 755.9
Panel 2	318935	159.2	89.15 - 204.6
Panel 3	318935	52.77	31.29 - 74.91
Panel 4	318935	15.43	10.10 - 19.60
Panel 5	318935	5.472	3.241 - 7.262

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	132.3	61.61 - 225.7
Panel 2	318935	21.86	11.69 - 35.93
Panel 3	318935	6.576	4.089 - 11.05
Panel 4	318935	1.785	1.243 - 2.815
Panel 5	318935	0.634	0.443 - 0.957

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	379.1	189.4 - 566.2
Panel 2	318935	118.5	66.23 - 174.3
Panel 3	318935	38.73	25.41 - 58.84
Panel 4	318935	11.26	7.756 - 16.68
Panel 5	318935	3.880	2.063 - 5.792

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	106.8	75.66 - 129.7
Panel 2	318935	32.22	21.68 - 40.39
Panel 3	318935	10.36	7.652 - 14.16
Panel 4	318935	3.150	2.387 - 4.063
Panel 5	318935	1.214	0.777 - 1.557

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

Review/Approval

Shivani Goel

Quality Control
Manager

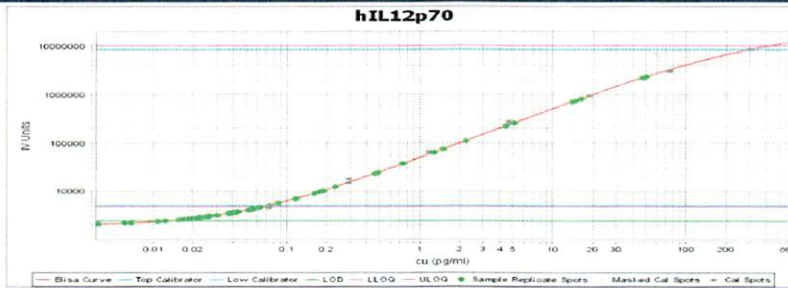
Shivani Goel 29 Sep 2021

Name

Title

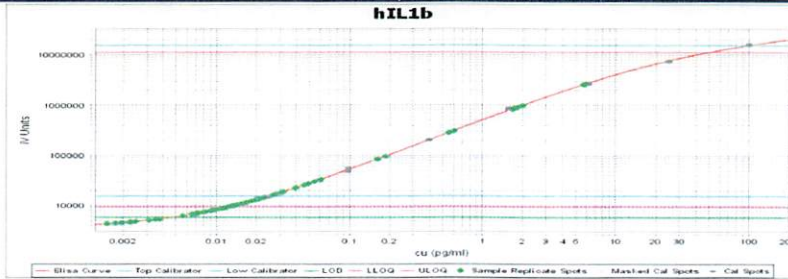
Signature/Date

IL-12p70 Calibration Curve



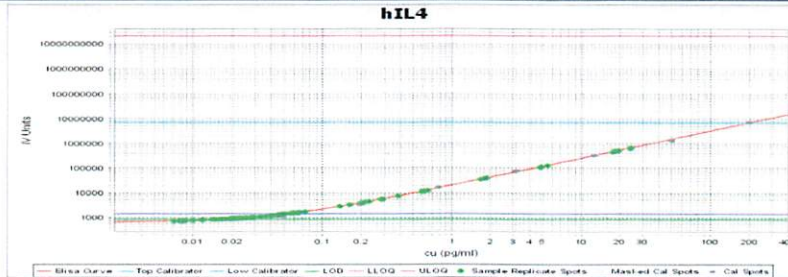
Calibrator Levels (pg/ml)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

IL-1β Calibration Curve



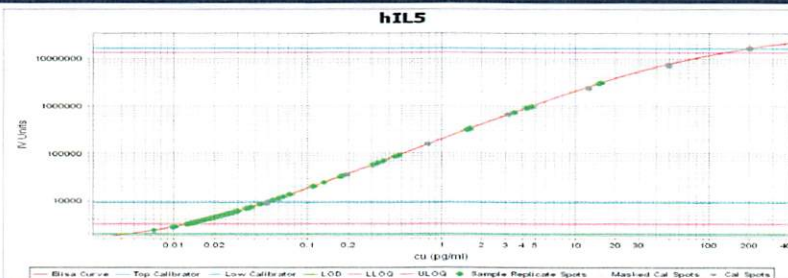
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

IL-4 Calibration Curve



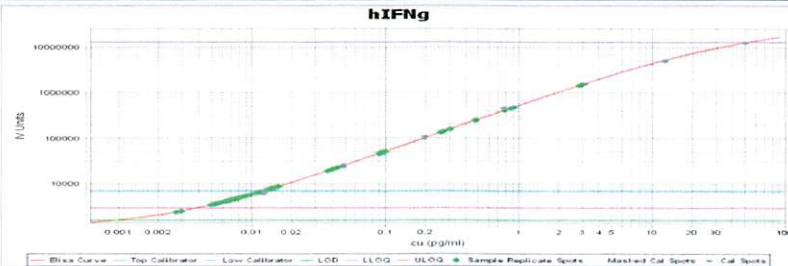
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

IL-5 Calibration Curve



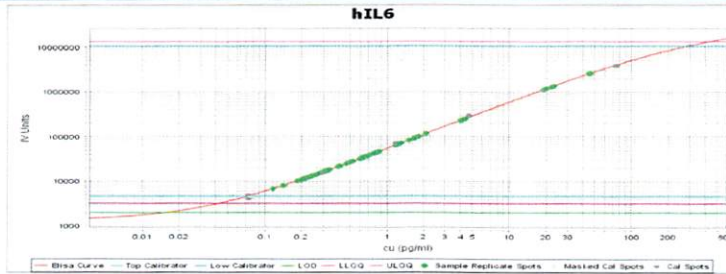
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

IFNγ Calibration Curve



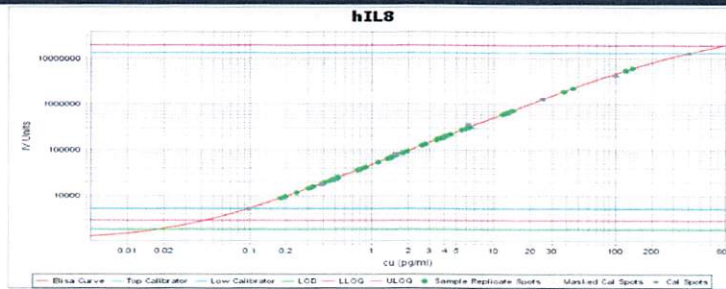
Calibrator Levels (pg/mL)	
1	50
2	12.5
3	3.13
4	0.781
5	0.195
6	0.049
7	0.012
8	0.000

IL-6 Calibration Curve



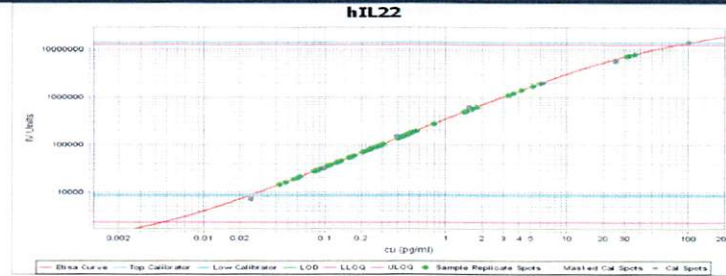
Calibrator Levels (pg/mL)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

IL-8 Calibration Curve



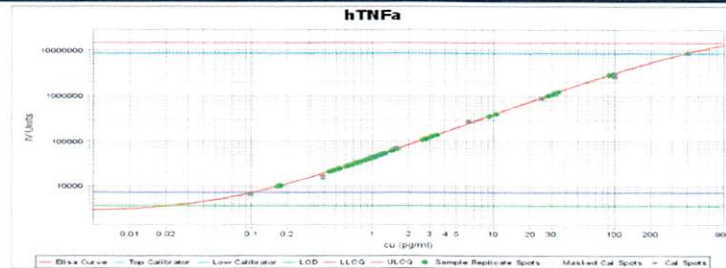
Calibrator Levels (pg/mL)	
1	400
2	100
3	25.00
4	6.250
5	1.563
6	0.391
7	0.098
8	0.000

IL-22 Calibration Curve



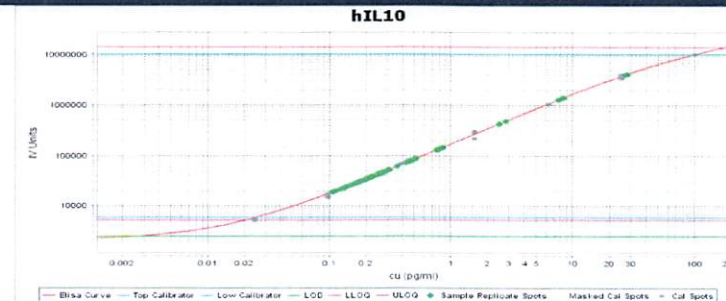
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

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