

Product Number:	85-0329
Lot Number:	319647
Expiration:	12-Nov-2022
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	319647P
Biotinylated Antibody Rgt	98-0329	319647B
Sample Diluent 21	1863329	319579
IL-12p70 Calibrator	18910034	318629
IL-1 β Calibrator	18910040	318728
IL-4 Calibrator	1858452	318480
IL-5 Calibrator	1858462	318837
IFN γ Calibrator	1858260	318631
IL-6 Calibrator	18910044	318724
IL-8 Calibrator	1858454	319011
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

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).



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	318935	185.8	134.8 - 213.6
Panel 2	318935	54.64	36.52 - 68.68
Panel 3	318935	17.37	13.03 - 23.02
Panel 4	318935	5.094	3.581 - 6.569
Panel 5	318935	1.788	0.963 - 2.340

IL-1 β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	23.68	12.01 - 32.00
Panel 2	318935	7.118	3.739 - 9.953
Panel 3	318935	2.588	1.270 - 3.351
Panel 4	318935	0.786	0.387 - 0.986
Panel 5	318935	0.261	0.111 - 0.348

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	66.07	51.07 - 80.47
Panel 2	318935	19.90	14.40 - 24.06
Panel 3	318935	7.420	5.421 - 8.931
Panel 4	318935	2.069	1.585 - 2.879
Panel 5	318935	0.678	0.478 - 1.311

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	89.73	47.13 - 109.0
Panel 2	318935	30.31	15.94 - 32.60
Panel 3	318935	10.66	5.192 - 11.65
Panel 4	318935	3.253	1.649 - 3.323
Panel 5	318935	1.147	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	11.36	7.193 - 14.98
Panel 2	318935	3.813	1.985 - 4.964
Panel 3	318935	1.387	0.714 - 1.703
Panel 4	318935	0.453	0.249 - 0.513
Panel 5	318935	0.186	0.103 - 0.227

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	176.9	122.5 - 249.4
Panel 2	318935	80.91	57.98 - 109.1
Panel 3	318935	17.11	12.14 - 22.22
Panel 4	318935	8.460	5.891 - 9.589
Panel 5	318935	2.730	1.746 - 3.433

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	532.0	285.9 - 755.9
Panel 2	318935	158.5	89.15 - 204.6
Panel 3	318935	56.36	31.29 - 74.91
Panel 4	318935	16.87	10.10 - 19.60
Panel 5	318935	5.557	3.241 - 7.262

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	111.6	61.61 - 225.7
Panel 2	318935	18.84	11.69 - 35.93
Panel 3	318935	6.129	4.089 - 11.05
Panel 4	318935	1.786	1.243 - 2.815
Panel 5	318935	0.582	0.443 - 0.957

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	430.3	189.4 - 566.2
Panel 2	318935	120.6	66.23 - 174.3
Panel 3	318935	41.42	25.41 - 58.84
Panel 4	318935	11.46	7.756 - 16.68
Panel 5	318935	3.917	2.063 - 5.792

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	105.1	75.66 - 129.7
Panel 2	318935	30.41	21.68 - 40.39
Panel 3	318935	10.70	7.652 - 14.16
Panel 4	318935	3.445	2.387 - 4.063
Panel 5	318935	1.254	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 18 Feb 2022

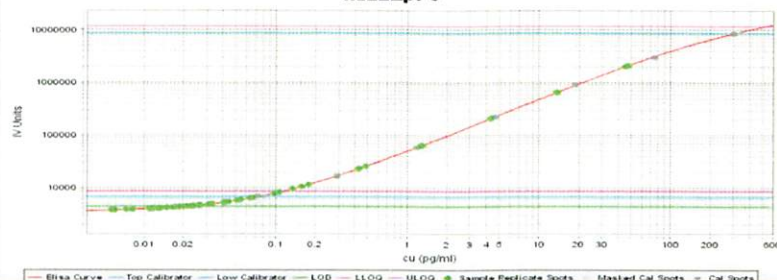
Name

Title

Signature/Date

IL-12p70 Calibration Curve

hIL12p70

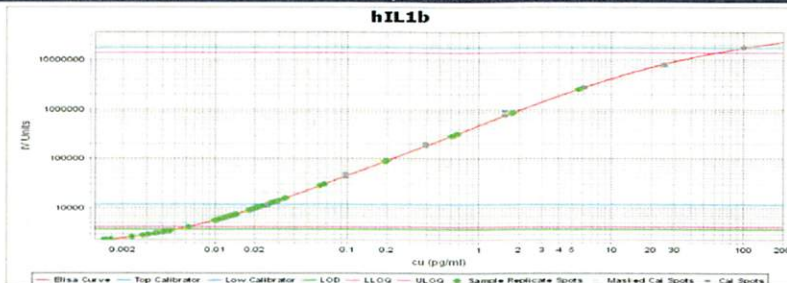


Calibrator Levels (pg/ml)

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

hIL1 β

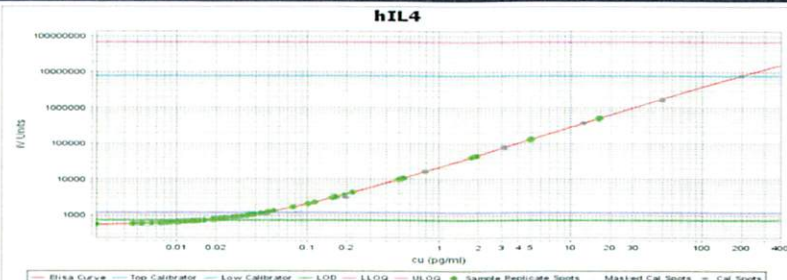


Calibrator Levels (pg/mL)

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

hIL4

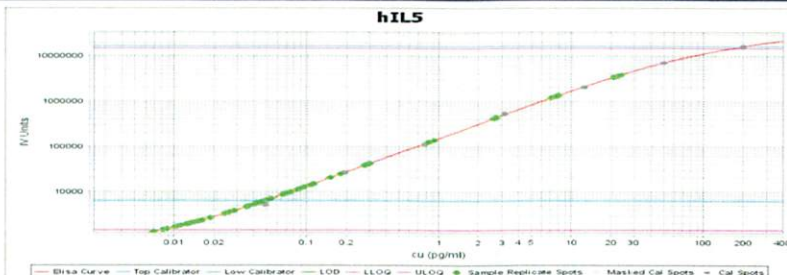


Calibrator Levels (pg/mL)

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

hIL5

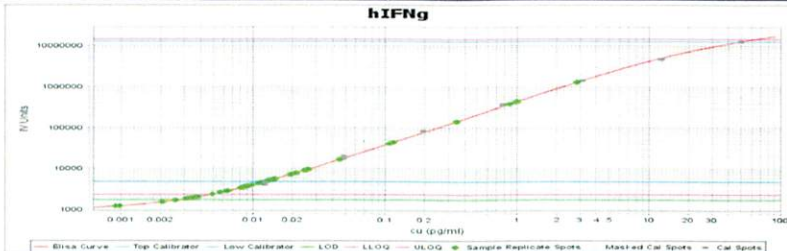


Calibrator Levels (pg/mL)

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

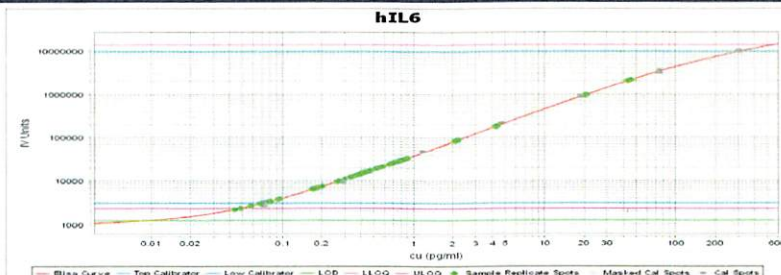
hIFN γ



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

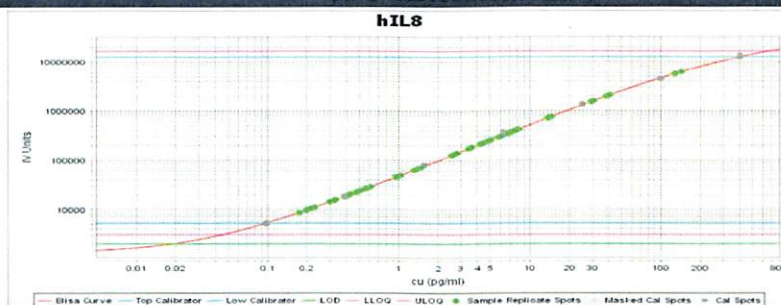
IL-6 Calibration Curve



**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

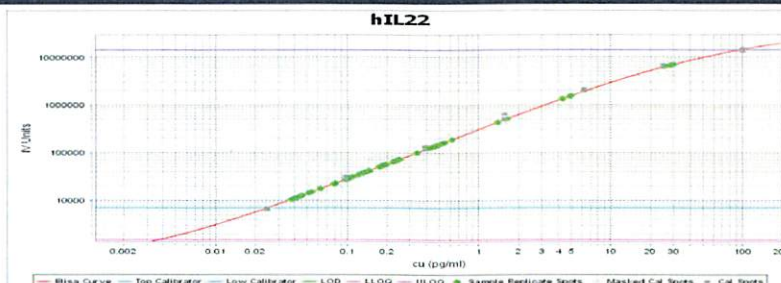
IL-8 Calibration Curve



**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

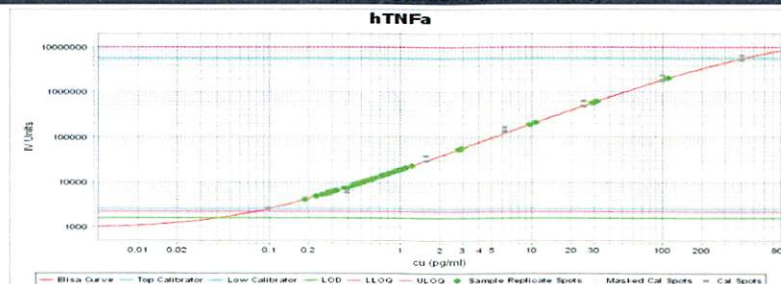
IL-22 Calibration Curve



**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

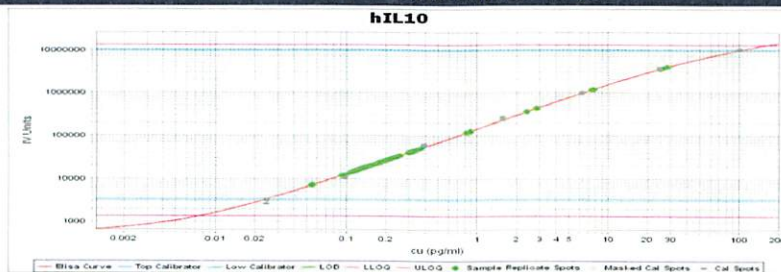
TNFα Calibration Curve



**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