

Product Number:	85-0329
Lot Number:	319620
Expiration:	12-Nov-2022
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	319616P
Biotinylated Antibody Rgt	98-0329	319494B
Sample Diluent 21	1863329	319579
IL-12p70 Calibrator	18910034	319156
IL-1 β Calibrator	18910040	319165
IL-4 Calibrator	1858452	319458
IL-5 Calibrator	1858462	319223
IFN γ Calibrator	1858260	318898
IL-6 Calibrator	18910044	318724
IL-8 Calibrator	1858454	319011
IL-22 Calibrator	18910036	319176
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	142.3	134.8 - 213.6
Panel 2	318935	39.78	36.52 - 68.68
Panel 3	318935	14.94	13.03 - 23.02
Panel 4	318935	4.193	3.581 - 6.569
Panel 5	318935	1.360	0.963 - 2.340

IL-1 β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	16.82	12.01 - 32.00
Panel 2	318935	4.966	3.739 - 9.953
Panel 3	318935	1.864	1.270 - 3.351
Panel 4	318935	0.593	0.387 - 0.986
Panel 5	318935	0.188	0.111 - 0.348

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	54.92	51.07 - 80.47
Panel 2	318935	15.74	14.40 - 24.06
Panel 3	318935	6.181	5.421 - 8.931
Panel 4	318935	1.969	1.585 - 2.879
Panel 5	318935	0.502	0.478 - 1.311

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	67.07	47.13 - 109.0
Panel 2	318935	18.80	15.94 - 32.60
Panel 3	318935	7.373	5.192 - 11.65
Panel 4	318935	2.180	1.649 - 3.323
Panel 5	318935	0.744	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	10.63	7.193 - 14.98
Panel 2	318935	3.103	1.985 - 4.964
Panel 3	318935	1.190	0.714 - 1.703
Panel 4	318935	0.368	0.249 - 0.513
Panel 5	318935	0.135	0.103 - 0.227

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	164.3	122.5 - 249.4
Panel 2	318935	73.01	57.98 - 109.1
Panel 3	318935	15.83	12.14 - 22.22
Panel 4	318935	7.948	5.891 - 9.589
Panel 5	318935	2.347	1.746 - 3.433

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	532.9	285.9 - 755.9
Panel 2	318935	139.0	89.15 - 204.6
Panel 3	318935	50.69	31.29 - 74.91
Panel 4	318935	14.58	10.10 - 19.60
Panel 5	318935	4.441	3.241 - 7.262

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	93.00	61.61 - 225.7
Panel 2	318935	13.87	11.69 - 35.93
Panel 3	318935	4.687	4.089 - 11.05
Panel 4	318935	1.376	1.243 - 2.815
Panel 5	318935	0.428	0.443 - 0.957

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	332.0	189.4 - 566.2
Panel 2	318935	99.32	66.23 - 174.3
Panel 3	318935	38.64	25.41 - 58.84
Panel 4	318935	11.89	7.756 - 16.68
Panel 5	318935	3.317	2.063 - 5.792

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	100.0	75.66 - 129.7
Panel 2	318935	30.13	21.68 - 40.39
Panel 3	318935	11.28	7.652 - 14.16
Panel 4	318935	3.649	2.387 - 4.063
Panel 5	318935	1.187	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

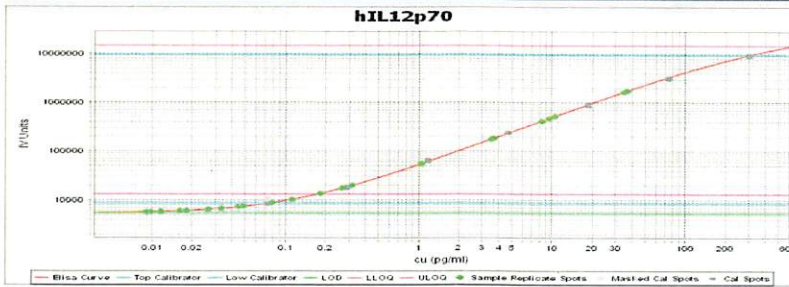
Shivani Goel 28Jan2022

Name

Title

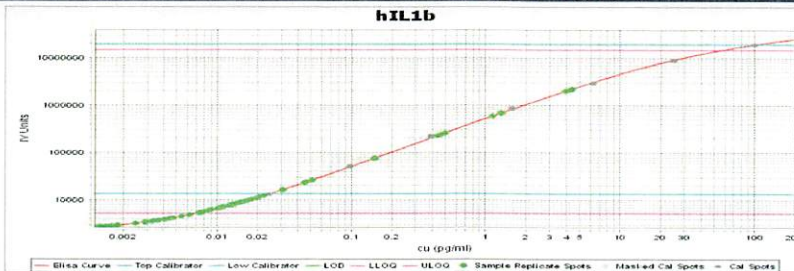
Signature/Date

IL-12p70 Calibration Curve



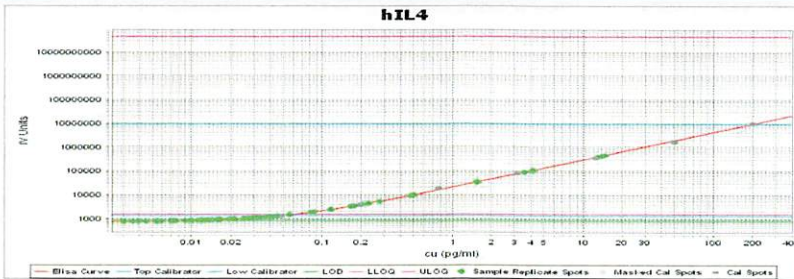
Calibrator Levels (pg/ml)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

IL-1β Calibration Curve



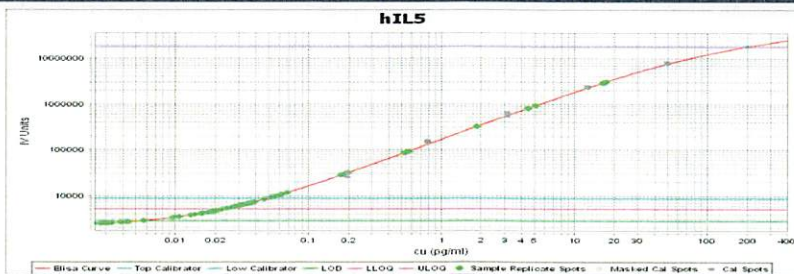
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

IL-4 Calibration Curve



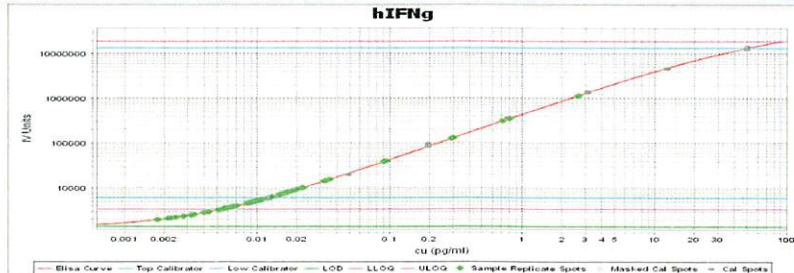
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

IL-5 Calibration Curve

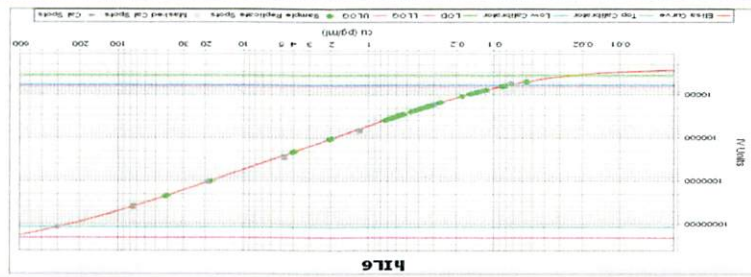


Calibrator Levels (pg/mL)	
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

IFNγ Calibration Curve

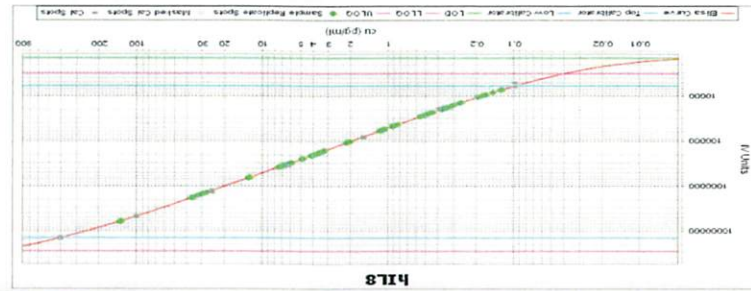


Calibrator Levels (pg/mL)	
1	50
2	12.5
3	3.13
4	0.781
5	0.195
6	0.049
7	0.012
8	0.000



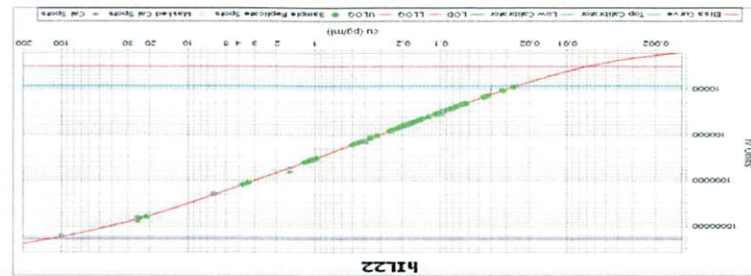
hIL6 Calibration Curve

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



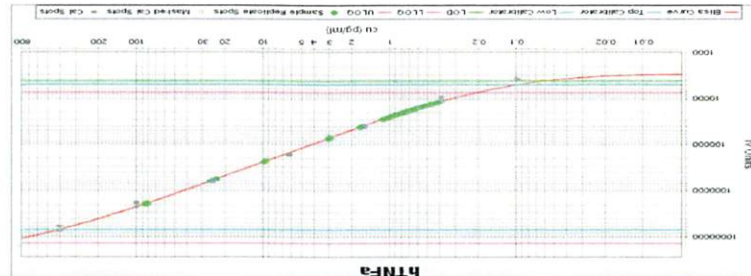
hIL8 Calibration Curve

Callibrator Levels (pg/mL)	
400	1
100	2
25.00	3
6.250	4
1.563	5
0.391	6
0.098	7
0.024	8



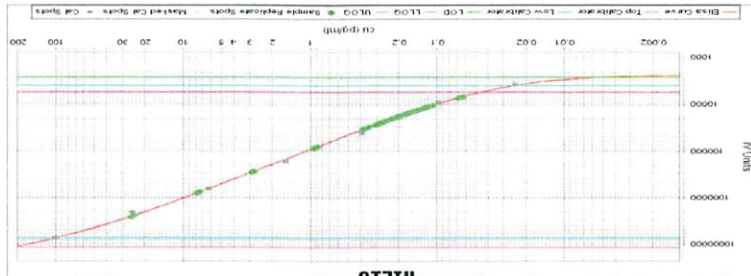
hIL22 Calibration Curve

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



hTNFa Calibration Curve

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



hIL10 Calibration Curve

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