

Product Number:	85-0329
Lot Number:	319836
Expiration:	23-Dec-2023
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	319836P
Biotinylated Antibody Rgt	98-0329	319836B
Sample Diluent 21	1863329	319785
IL-12p70 Calibrator	18910034	319156
IL-1 β Calibrator	18910040	319564
IL-4 Calibrator	1858452	319458
IL-5 Calibrator	1858462	319223
IFN γ Calibrator	1858260	319711
IL-6 Calibrator	18910044	319484
IL-8 Calibrator	1858454	319575
IL-22 Calibrator	18910036	319459
TNF α 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 Cytokine Kit Instructions (85A-0329).

Release Materials

IL-12p70

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	156.6	131.7 - 203.5
Panel 2	318935	45.70	36.98 - 66.47
Panel 3	318935	15.18	13.40 - 22.07
Panel 4	318935	4.148	3.751 - 6.122
Panel 5	318935	1.213	1.179 - 2.080

IL-1 β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	24.98	12.91 - 32.59
Panel 2	318935	7.105	4.217 - 9.740
Panel 3	318935	2.494	1.377 - 3.499
Panel 4	318935	0.697	0.424 - 0.975
Panel 5	318935	0.230	0.137 - 0.343

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	54.48	50.84 - 77.00
Panel 2	318935	15.20	14.18 - 23.93
Panel 3	318935	5.273	5.174 - 8.891
Panel 4	318935	1.564	1.422 - 2.789
Panel 5	318935	0.562	0.410 - 1.241

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	67.35	43.55 - 104.5
Panel 2	318935	17.34	13.38 - 33.50
Panel 3	318935	6.830	4.766 - 11.68
Panel 4	318935	1.419	1.234 - 3.623
Panel 5	318935	0.582	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	10.44	8.245 - 14.41
Panel 2	318935	3.093	2.401 - 4.768
Panel 3	318935	1.100	0.853 - 1.722
Panel 4	318935	0.334	0.271 - 0.505
Panel 5	318935	0.144	0.114 - 0.221

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	180.7	143.3 - 215.0
Panel 2	318935	78.97	62.18 - 101.2
Panel 3	318935	16.08	12.91 - 20.83
Panel 4	318935	7.104	6.002 - 9.003
Panel 5	318935	2.445	1.966 - 3.145

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	413.6	363.4 - 600.3
Panel 2	318935	111.4	91.56 - 189.5
Panel 3	318935	40.98	32.31 - 67.04
Panel 4	318935	11.10	9.930 - 18.44
Panel 5	318935	3.605	3.310 - 6.556

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	81.8	47.37 - 159.2
Panel 2	318935	11.95	6.880 - 28.17
Panel 3	318935	3.859	2.175 - 9.086
Panel 4	318935	1.006	0.686 - 2.605
Panel 5	318935	0.323	0.216 - 0.878

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	336.8	251.0 - 450.4
Panel 2	318935	95.4	69.11 - 155.1
Panel 3	318935	34.76	25.96 - 51.54
Panel 4	318935	9.67	7.461 - 14.87
Panel 5	318935	3.115	2.349 - 4.860

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	100.6	75.65 - 125.4
Panel 2	318935	26.95	21.66 - 39.42
Panel 3	318935	9.21	7.643 - 13.76
Panel 4	318935	2.723	2.344 - 3.891
Panel 5	318935	0.982	0.826 - 1.448

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

Review/Approval

Matthew Koulouris

Quality Control
Supervisor

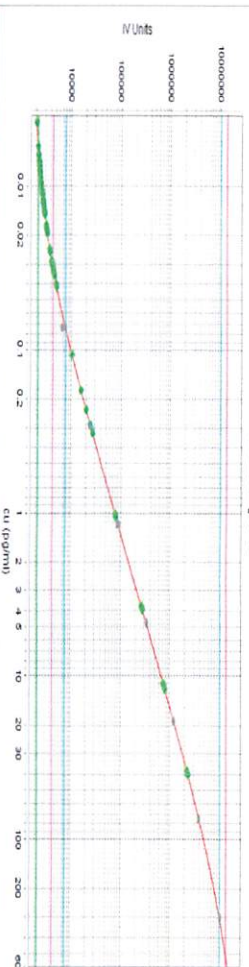
 28 NOV 2022

Name

Title

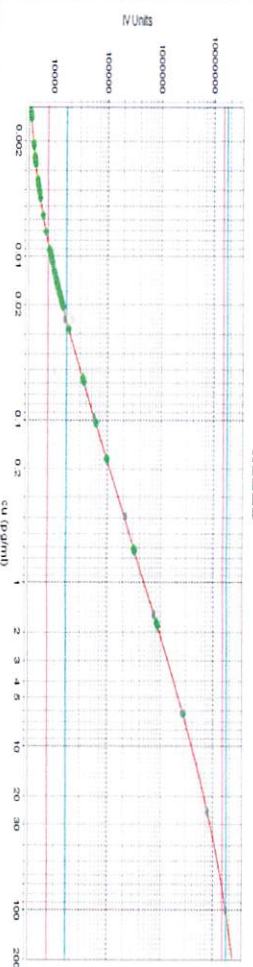
Signature/Date

hIL12p70
hIL12p70 Calibration Curve



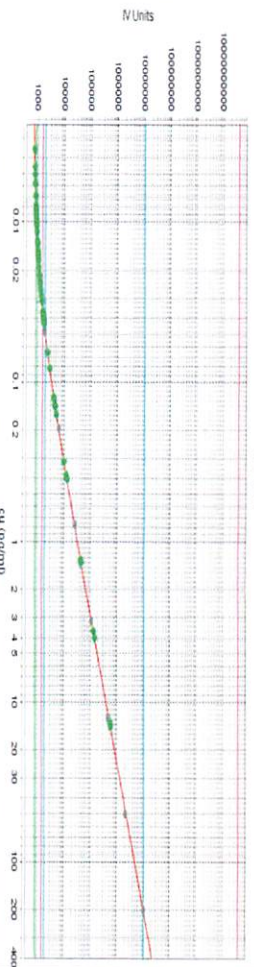
Calibrator Levels (pg/mL)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

hIL1b
hIL1b Calibration Curve



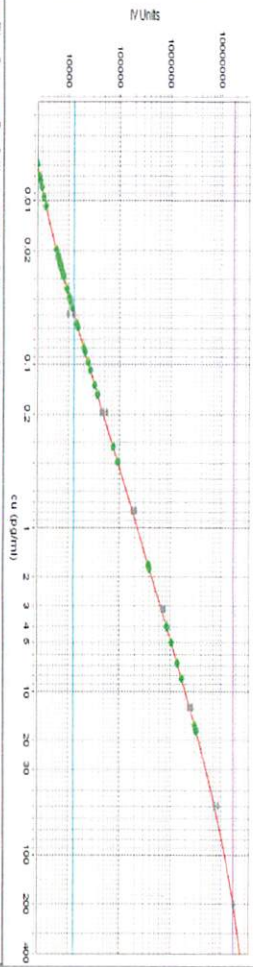
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

hIL4
hIL4 Calibration Curve



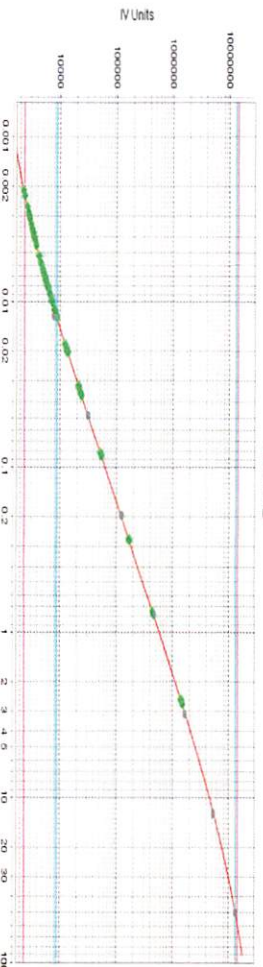
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

hIL5
hIL5 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

hIFNγ
hIFNγ 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

