

Product Number:	85-0329
Lot Number:	318359
Expiration:	11-May-2021
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	318359P
Biotinylated Antibody Rgt	98-0329	318359B
Sample Diluent 21 ¹	1863329	N/A
IL-12p70 Calibrator	18910034	317901
IL-1 β Calibrator	18910040	318067
IL-4 Calibrator	1858452	317999
IL-5 Calibrator	1858462	317984
IFN γ Calibrator	1858260	318097
IL-6 Calibrator	18910044	318089
IL-8 Calibrator	1858454	318342
IL-22 Calibrator	18910036	317907
TNF α Calibrator	1880012	317986
IL-10 Calibrator	18910043	318068
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 Cytokine 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)
Control 1	1-Feb-19	51.2	39.6 - 59.4
Control 2	1-Feb-19	N/A	N/A
Panel 1	1-Feb-19	117	95.3 - 142.9
Panel 2	1-Feb-19	13.4	10.86 - 16.29
Panel 3	1-Feb-19	1.35	0.643 - 2.814

IL-1 β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	15.1	11.09 - 19.74
Control 2	1-Feb-19	0.746	0.418 - 1.074
Panel 1	1-Feb-19	0.560	0.218 - 1.123
Panel 2	1-Feb-19	3.86	2.795 - 5.111
Panel 3	1-Feb-19	16.7	12.02 - 20.84

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	35.4	26.81 - 40.21
Control 2	1-Feb-19	2.15	1.239 - 2.764
Panel 1	1-Feb-19	48.9	35.60 - 53.40
Panel 2	1-Feb-19	5.46	3.819 - 7.482
Panel 3	1-Feb-19	1.249	0.533 - 2.174

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	30.2	23.20 - 35.41
Control 2	1-Feb-19	1.56	1.057 - 2.068
Panel 1	1-Feb-19	71.5	53.92 - 82.17
Panel 2	1-Feb-19	7.67	6.223 - 9.864
Panel 3	1-Feb-19	1.24	0.857 - 1.764

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)
Control 1	1-Feb-19	8.38	5.79 - 9.77
Control 2	1-Feb-19	0.544	0.212 - 0.532
Panel 1	1-Feb-19	14.9	10.3 - 18.4
Panel 2	1-Feb-19	2.09	0.909 - 1.82
Panel 3	1-Feb-19	N/A	N/A

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	42.7	36.3 - 54.6
Control 2	1-Feb-19	2.57	1.67 - 3.26
Panel 1	1-Feb-19	73.6	56.4 - 87.4
Panel 2	1-Feb-19	10.56	7.75 - 13.0
Panel 3	1-Feb-19	1.63	0.975 - 2.52

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	56.4	47.2 - 72.4
Control 2	1-Feb-19	2.84	2.17 - 3.70
Panel 1	1-Feb-19	299	248 - 410
Panel 2	1-Feb-19	11.3	11.1 - 18.7
Panel 3	1-Feb-19	0.99	0.493 - 3.20

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	19.3	13.5 - 20.3
Control 2	1-Feb-19	0.997	0.609 - 1.19
Panel 1	1-Feb-19	36.1	24.4 - 37.0
Panel 2	1-Feb-19	3.59	2.07 - 4.25
Panel 3	1-Feb-19	0.258	0.129 - 0.576

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	49.7	40.5 - 68.9
Control 2	1-Feb-19	2.78	1.54 - 3.52
Panel 1	1-Feb-19	92.7	84.6 - 143
Panel 2	1-Feb-19	25.2	17.9 - 31.4
Panel 3	1-Feb-19	N/A	N/A

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	14.5	10.5 - 19.8
Control 2	1-Feb-19	0.719	0.414 - 1.02
Panel 1	1-Feb-19	0.724	0.478 - 0.971
Panel 2	1-Feb-19	4.31	2.95 - 6.12
Panel 3	1-Feb-19	0.266	0.211 - 0.412

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

Review/Approval

Patrizia Stadler

Associate
Director of QC

P. Stadler

23 Jan 2010

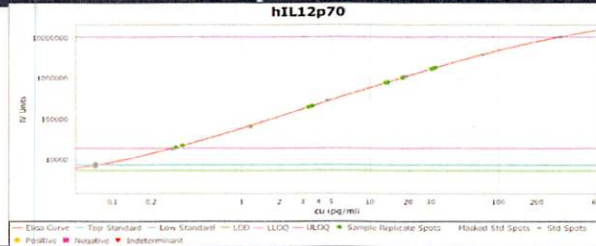
Name

Title

Signature/Date

**CorPlex 10-Plex huAP1 Kit
Certificate of Analysis**

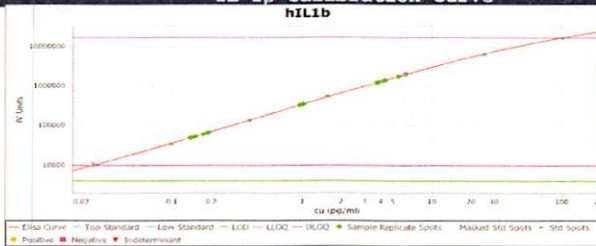
IL-12p70 Calibration Curve



**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

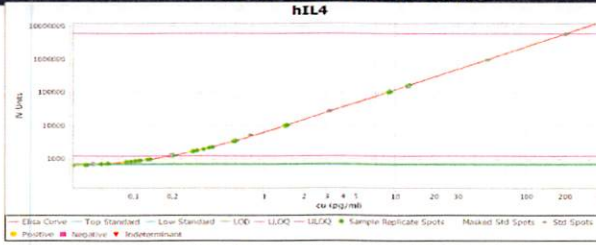
IL-1 β Calibration Curve



**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

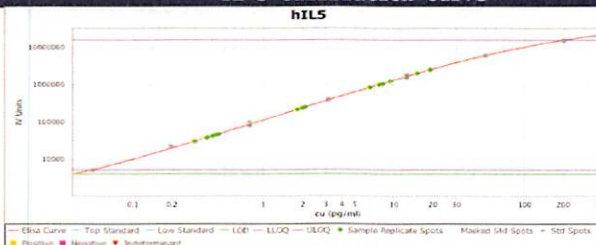
IL-4 Calibration Curve



**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

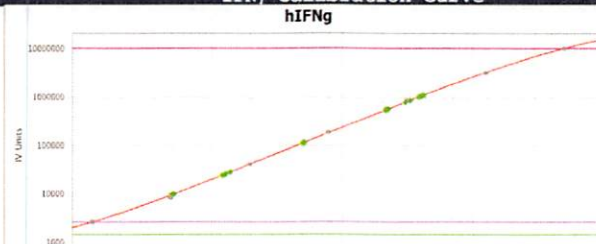
IL-5 Calibration Curve



**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

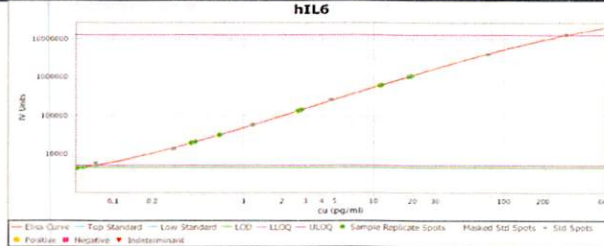


**Calibrator Levels
(pg/mL)**

1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

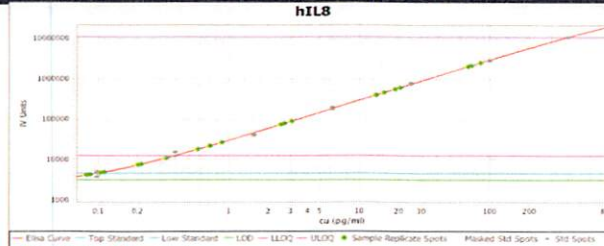
CorPlex 10plex huAP1 Kit
Certificate of Analysis

IL-6 Calibration Curve



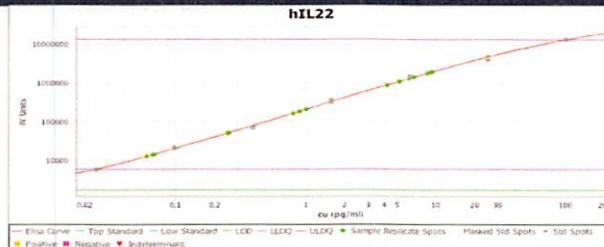
Calibrator Levels (pg/mL)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

IL-8 Calibration Curve



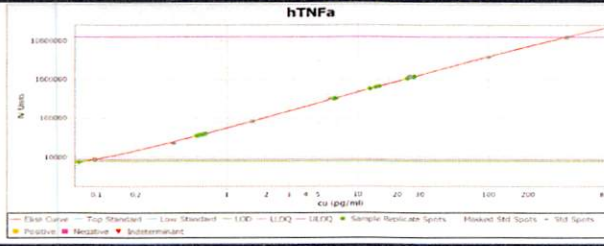
Calibrator Levels (pg/mL)	
1	400
2	100
3	25.00
4	6.250
5	1.563
6	0.391
7	0.098
8	0.000

IL-22 Calibration Curve



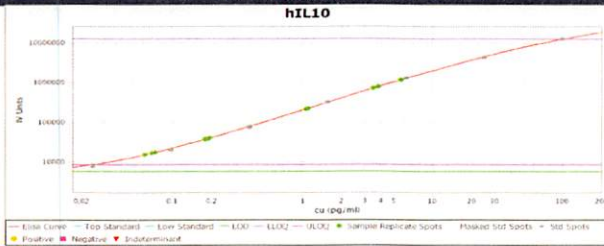
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

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