

Product Number:	85-0329
Lot Number:	318007
Expiration:	13-Aug-2020
Platform(s):	SP-X



Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	318007
Biotinylated Antibody Rgt	98-0329	318007
Sample Diluent 21 ¹	1863329	318027
IL-12p70 Calibrator	18910034	317876
IL-1 β Calibrator	18910040	317793
IL-4 Calibrator	1858452	317726
IL-5 Calibrator	1858462	317984
IFN γ Calibrator	1858260	317863
IL-6 Calibrator	18910044	317985
IL-8 Calibrator	1858454	317862
IL-22 Calibrator	18910036	317907
TNF α Calibrator	1880012	317986
IL-10 Calibrator	18910043	317877
Wash Buffer (25X) ¹	1860055	17512600
Streptavidin-HRP Reagent ¹	1858263	317972
Super Signal [®] Luminol* ¹	1858089	TJ275117
Super Signal [®] Peroxide* ¹	1858090	TJ275118

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 Cytokine1 Kit Instructions (85A-0329).

Release Materials

IL-12p70

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	49.71	37.77 - 56.65
Control 2	1-Feb-19	N/A	N/A
Panel 1	1-Feb-19	121.7	91.26 - 136.9
Panel 2	1-Feb-19	12.49	10.21 - 15.32
Panel 3	1-Feb-19	1.830	0.864 - 2.578

IL-1 β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	16.38	12.13 - 18.29
Control 2	1-Feb-19	0.846	0.535 - 0.984
Panel 1	1-Feb-19	0.727	0.364 - 0.923
Panel 2	1-Feb-19	4.096	3.219 - 4.828
Panel 3	1-Feb-19	17.10	14.03 - 21.04

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	30.32	24.91 - 37.37
Control 2	1-Feb-19	1.775	1.159 - 2.389
Panel 1	1-Feb-19	42.64	33.77 - 50.65
Panel 2	1-Feb-19	5.390	2.891 - 7.442
Panel 3	1-Feb-19	1.630	0.290 - 2.252

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	30.51	24.41 - 36.62
Control 2	1-Feb-19	1.500	1.254 - 1.882
Panel 1	1-Feb-19	72.35	56.42 - 84.63
Panel 2	1-Feb-19	7.553	6.609 - 9.914
Panel 3	1-Feb-19	1.286	1.017 - 1.690

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	7.844	6.094 - 9.142
Control 2	1-Feb-19	0.396	0.226 - 0.491
Panel 1	1-Feb-19	15.78	12.27 - 18.40
Panel 2	1-Feb-19	1.538	1.190 - 1.841
Panel 3	1-Feb-19	1.549	0.652 - 2.383

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	47.68	36.13 - 56.32
Control 2	1-Feb-19	2.282	1.689 - 3.013
Panel 1	1-Feb-19	74.00	57.22 - 85.83
Panel 2	1-Feb-19	10.055	8.131 - 12.71
Panel 3	1-Feb-19	1.833	0.878 - 2.460

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	60.77	38.73 - 83.91
Control 2	1-Feb-19	2.799	1.663 - 4.130
Panel 1	1-Feb-19	350.2	243.0 - 438.9
Panel 2	1-Feb-19	14.75	8.026 - 23.33
Panel 3	1-Feb-19	2.126	0.788 - 3.246

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	17.30	13.23 - 19.84
Control 2	1-Feb-19	0.833	0.550 - 1.109
Panel 1	1-Feb-19	31.56	22.83 - 35.10
Panel 2	1-Feb-19	3.494	2.027 - 4.121
Panel 3	1-Feb-19	0.411	0.078 - 0.580

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	58.31	42.94 - 76.80
Control 2	1-Feb-19	2.207	1.734 - 3.531
Panel 1	1-Feb-19	124.0	92.40 - 146.9
Panel 2	1-Feb-19	25.84	19.15 - 35.96
Panel 3	1-Feb-19	18.34	8.273 - 30.64

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	16.10	11.08 - 20.27
Control 2	1-Feb-19	0.704	0.436 - 0.985
Panel 1	1-Feb-19	0.759	0.448 - 1.079
Panel 2	1-Feb-19	4.827	3.489 - 6.614
Panel 3	1-Feb-19	0.360	0.157 - 0.487

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

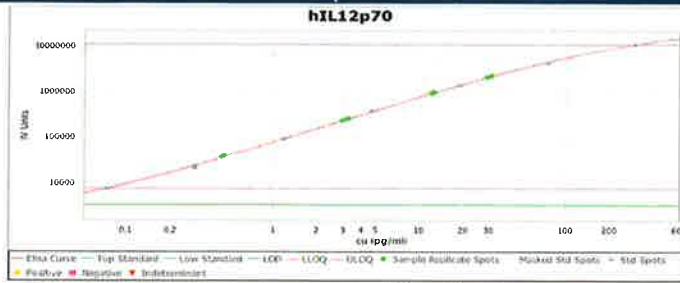
Review/Approval

Kathleen Cornell
Name

QC Manager
Title

Michael L. Dow 28 Mar 2019
Signature/Date

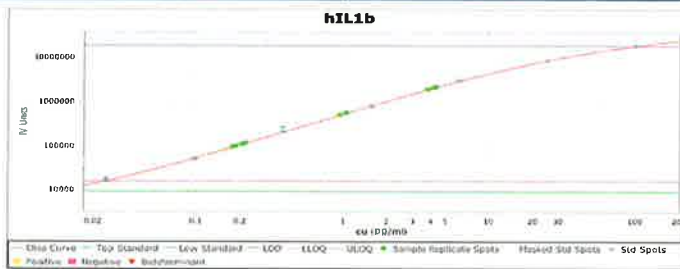
IL-12p70 Calibration Curve



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

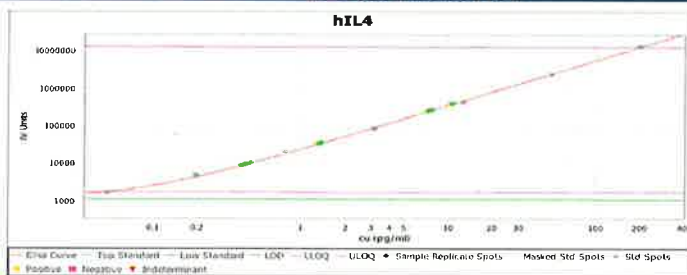
IL-1β Calibration Curve



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

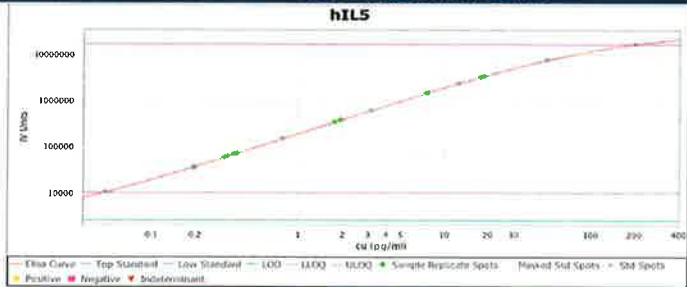
IL-4 Calibration Curve



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

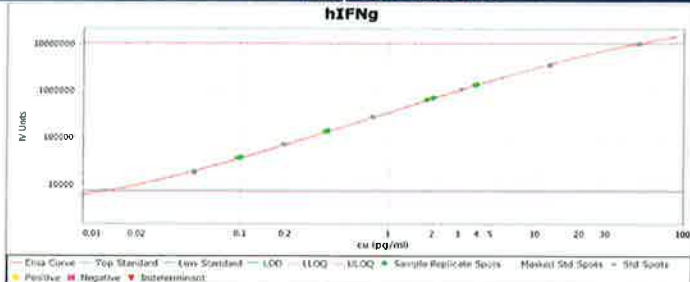
IL-5 Calibration Curve



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

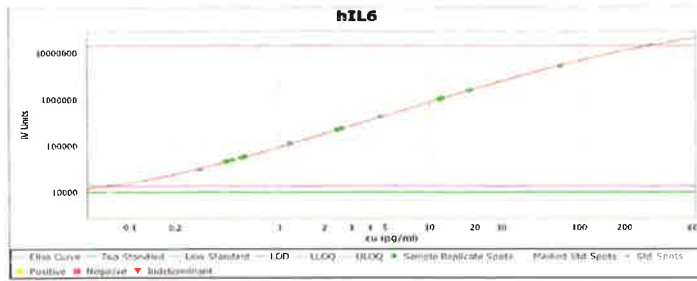
IFNγ Calibration Curve



Calibrator Levels
(pg/mL)

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

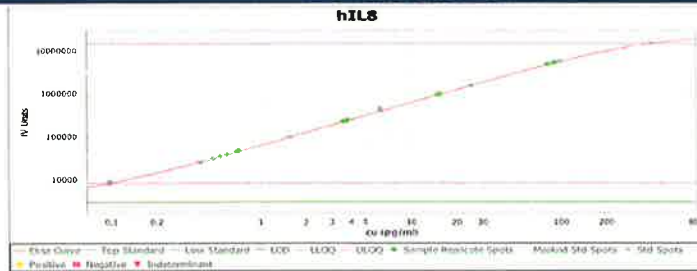
IL-6 Calibration Curve



**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

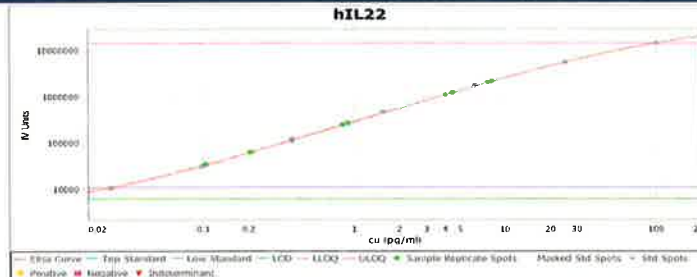
IL-8 Calibration Curve



**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

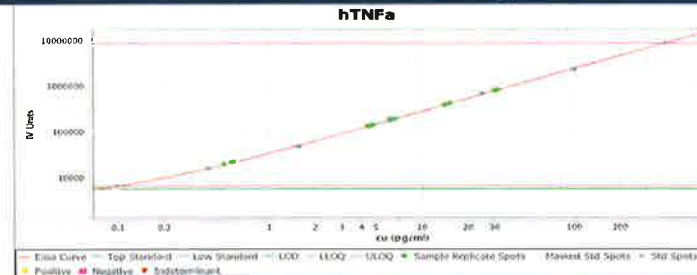
IL-22 Calibration Curve



**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

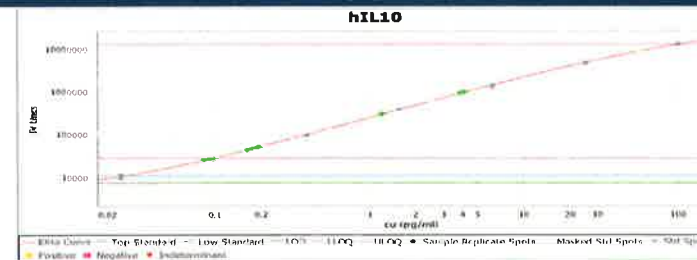
TNF α Calibration Curve



**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