

Product Number:	85-0329
Lot Number:	318607
Expiration:	12-Aug-2021
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	318607P
Biotinylated Antibody Rgt	98-0329	318404B
Sample Diluent 21 ¹	1863329	N/A
IL-12p70 Calibrator	18910034	318571
IL-1 β Calibrator	18910040	318067
IL-4 Calibrator	1858452	318480
IL-5 Calibrator	1858462	318465
IFN γ Calibrator	1858260	318538
IL-6 Calibrator	18910044	318089
IL-8 Calibrator	1858454	318081
IL-22 Calibrator	18910036	318408
TNF α Calibrator	1880012	318436
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)
Panel 1	318641	184	144-216
Panel 2	318641	64.6	48.6-72.9
Panel 3	318641	22.3	16.7-25.0
Panel 4	318641	7.40	5.89-8.84
Panel 5	318641	2.56	1.55-3.13

IL-1 β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	21.2	16.2-24.3
Panel 2	318641	7.37	5.63-8.44
Panel 3	318641	2.48	1.88-2.82
Panel 4	318641	0.827	0.661-0.992
Panel 5	318641	0.301	0.222-0.398

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	79.6	62.3-93.5
Panel 2	318641	26.8	20.2-30.4
Panel 3	318641	9.71	6.89-10.3
Panel 4	318641	3.32	2.34-4.09
Panel 5	318641	1.15	0.712-1.77

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	62.6	53.3-80.9
Panel 2	318641	21.7	18.7-28.0
Panel 3	318641	7.54	5.76-9.49
Panel 4	318641	2.59	2.15-3.58
Panel 5	318641	0.920	0.837-1.29

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	318641	9.38	7.98-12.0
Panel 2	318641	3.44	2.81-4.22
Panel 3	318641	1.21	0.948-1.42
Panel 4	318641	0.484	0.357-0.574
Panel 5	318641	0.188	0.147-0.220

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	125	113-169
Panel 2	318641	4.04	3.55-5.32
Panel 3	318641	1.35	1.10-1.65
Panel 4	318641	0.546	0.426-0.676
Panel 5	318641	0.250	0.089-0.357

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	358	393-589
Panel 2	318641	129.4	121-181
Panel 3	318641	47.3	39.7-59.6
Panel 4	318641	18.0	13.7-22.2
Panel 5	318641	6.14	4.59-6.89

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	20.1	18.3-28.9
Panel 2	318641	6.21	5.11-8.17
Panel 3	318641	1.87	1.65-2.53
Panel 4	318641	0.779	0.624-0.936
Panel 5	318641	0.215	0.190-0.342

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	74.0	53.8-89.2
Panel 2	318641	27.2	19.5-30.7
Panel 3	318641	9.09	5.57-11.0
Panel 4	318641	3.13	1.91-4.36
Panel 5	318641	1.19	0.682-1.86

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	51.4	50.0-75.0
Panel 2	318641	19.0	17.4-26.1
Panel 3	318641	6.62	5.94-8.91
Panel 4	318641	2.31	2.26-3.39
Panel 5	318641	0.815	0.825-1.24

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

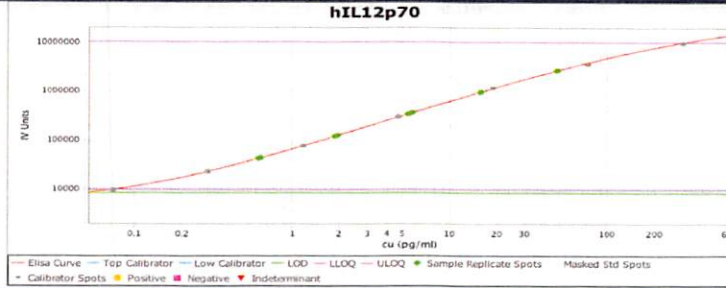
03 Apr 2020

Name

Title

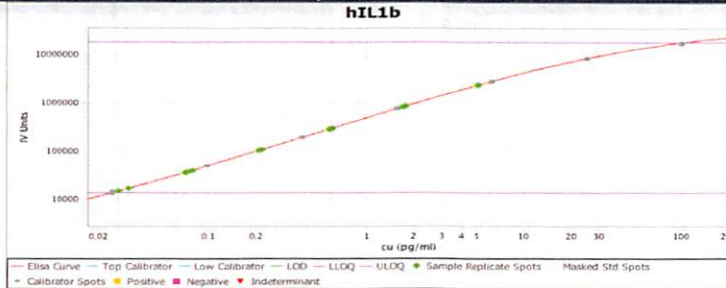
Signature/Date

IL-12p70 Calibration Curve



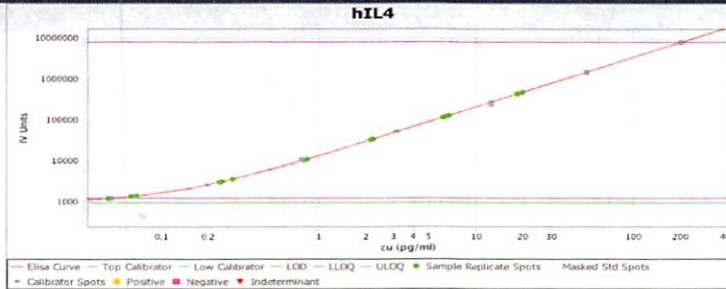
Calibrator Levels (pg/ml)	
1	300
2	75
3	18.8
4	4.69
5	1.17
6	0.293
7	0.073
8	0.000

IL-1 β Calibration Curve



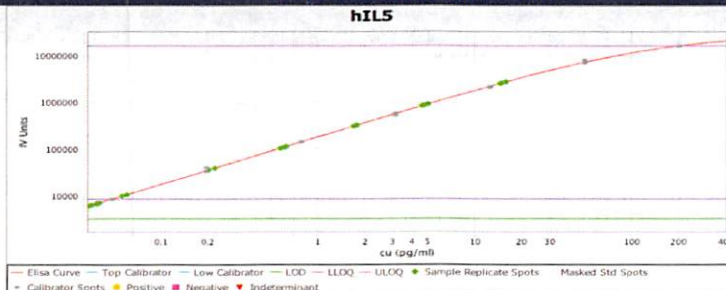
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.56
5	0.391
6	0.098
7	0.024
8	0.000

IL-4 Calibration Curve



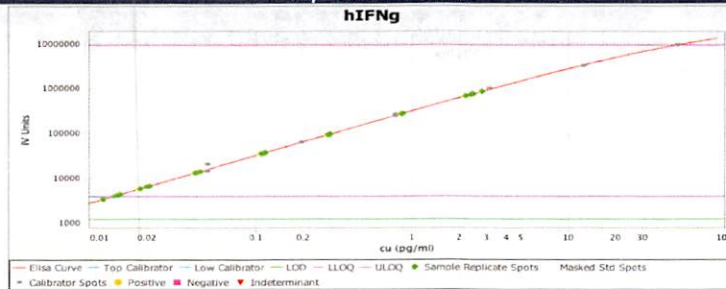
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.5
4	3.13
5	0.781
6	0.195
7	0.049
8	0.000

IL-5 Calibration Curve



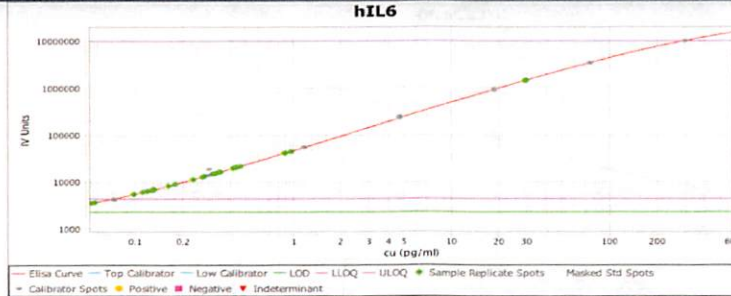
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.5
4	3.13
5	0.781
6	0.195
7	0.049
8	0.000

IFN γ Calibration Curve



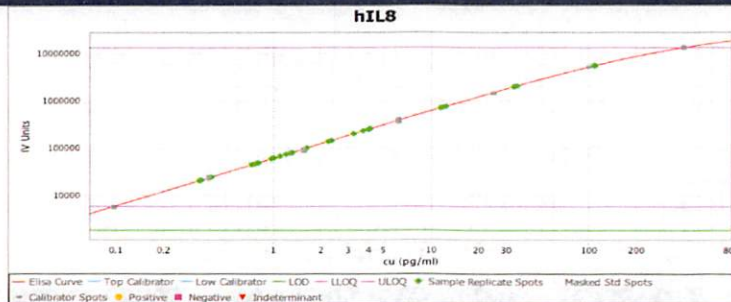
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.5
4	3.13
5	0.781
6	0.195
7	0.049
8	0.000

IL-6 Calibration Curve



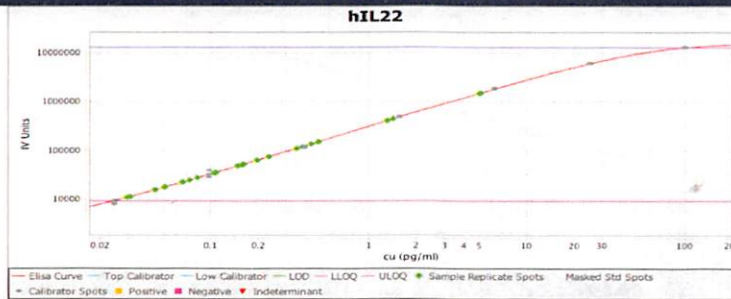
Calibrator Levels (pg/mL)	
1	300
2	75
3	18.8
4	4.69
5	1.17
6	0.293
7	0.073
8	0.000

IL-8 Calibration Curve



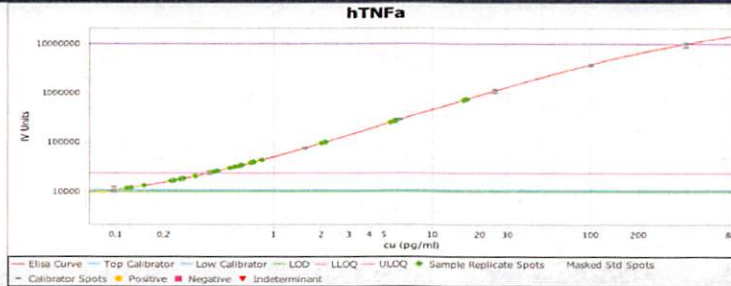
Calibrator Levels (pg/mL)	
1	400
2	100
3	25.0
4	6.25
5	1.56
6	0.391
7	0.098
8	0.000

IL-22 Calibration Curve



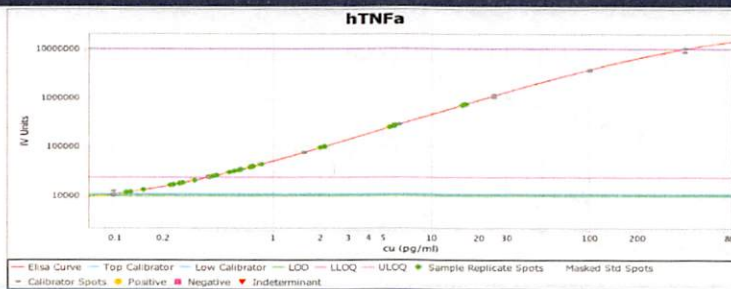
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.56
5	0.391
6	0.098
7	0.024
8	0.000

TNFα Calibration Curve



Calibrator Levels (pg/mL)	
1	400
2	100
3	25.0
4	6.25
5	1.56
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	1.56
5	0.391
6	0.098
7	0.024
8	0.000