

Product Number:	85-0329
Lot Number:	318337
Expiration:	9-Apr-2021
Platform(s):	SP-X



Data below represents results generated on the SP-X Imaging and Analysis System

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	318337P
Biotinylated Antibody Rgt	98-0329	318351B
Sample Diluent 21 ¹	1863329	N/A
IL-12p70 Calibrator	18910034	317901
IL-1β Calibrator	18910040	318067
IL-4 Calibrator	1858452	317999
IL-5 Calibrator	1858462	317984
IFNγ Calibrator	1858260	318097
IL-6 Calibrator	18910044	318089
IL-8 Calibrator	1858454	318081
IL-22 Calibrator	18910036	317907
TNFα Calibrator	1880012	317986
IL-10 Calibrator	18910043	318068
Wash Buffer (25X) ¹	1860055	N/A
Streptavidin-ERP 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).

Release Materials

IL-12p70

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	49.8	39.6 - 59.4
Control 2	1-Feb-19	NA	N/A
Panel 1	1-Feb-19	117	95.3 - 1423
Panel 2	1-Feb-19	12.8	10.9 - 16.3
Panel 3	1-Feb-19	1.40	0.643 - 2.81

IL-1β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	15.5	11.1 - 19.7
Control 2	1-Feb-19	0.770	0.418 - 1.07
Panel 1	1-Feb-19	0.570	0.218 - 1.12
Panel 2	1-Feb-19	4.14	2.80 - 5.11
Panel 3	1-Feb-19	20.1	12.0 - 20.8

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	33.4	26.8 - 40.2
Control 2	1-Feb-19	2.07	1.24 - 2.76
Panel 1	1-Feb-19	43.7	35.6 - 53.4
Panel 2	1-Feb-19	5.10	3.82 - 7.48
Panel 3	1-Feb-19	1.57	0.533 - 2.17

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	31.6	23.2 - 35.4
Control 2	1-Feb-19	1.77	1.06 - 2.07
Panel 1	1-Feb-19	75.1	53.9 - 82.2
Panel 2	1-Feb-19	9.43	6.22 - 9.86
Panel 3	1-Feb-19	1.48	0.857 - 1.76

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	8.00	5.79 - 9.77
Control 2	1-Feb-19	0.400	0.212 - 0.532
Panel 1	1-Feb-19	14.1	10.3 - 18.4
Panel 2	1-Feb-19	1.40	0.909 - 1.82
Panel 3	1-Feb-19	0.250	0.418 - 2.15

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	41.7	36.3 - 54.6
Control 2	1-Feb-19	2.46	1.67 - 3.26
Panel 1	1-Feb-19	78.1	56.4 - 87.4
Panel 2	1-Feb-19	10.3	7.75 - 13.0
Panel 3	1-Feb-19	1.59	0.975 - 2.52

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	50.9	47.2 - 72.4
Control 2	1-Feb-19	2.76	2.17 - 3.70
Panel 1	1-Feb-19	273	248 - 410
Panel 2	1-Feb-19	11.7	11.1 - 18.7
Panel 3	1-Feb-19	1.99	0.493 - 3.20

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	19.0	13.5 - 20.3
Control 2	1-Feb-19	1.01	0.609 - 1.19
Panel 1	1-Feb-19	39.5	24.4 - 37.0
Panel 2	1-Feb-19	3.56	2.07 - 4.25
Panel 3	1-Feb-19	0.280	0.129 - 0.576

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	53.9	40.5 - 68.9
Control 2	1-Feb-19	2.63	1.54 - 3.52
Panel 1	1-Feb-19	98.7	84.6 - 143
Panel 2	1-Feb-19	25.0	17.9 - 31.4
Panel 3	1-Feb-19	2.75	8.42 - 25.0

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	12-Dec-19	14.8	10.5 - 19.8
Control 2	12-Dec-19	0.760	0.414 - 1.02
Panel 1	12-Dec-19	0.740	0.478 - 0.971
Panel 2	12-Dec-19	4.48	2.95 - 6.12
Panel 3	12-Dec-19	0.270	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

78Lodler

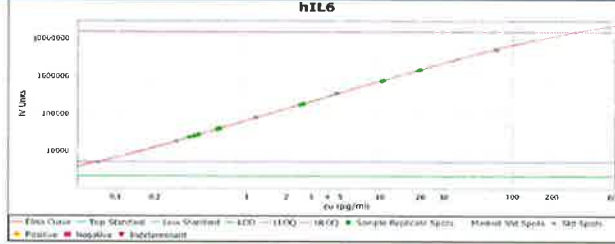
16 MAR 2020

Name

Title

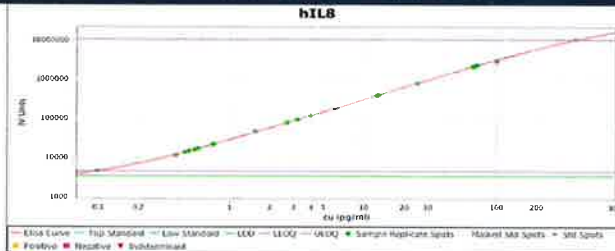
Signature/Date

IL-6 Calibration Curve



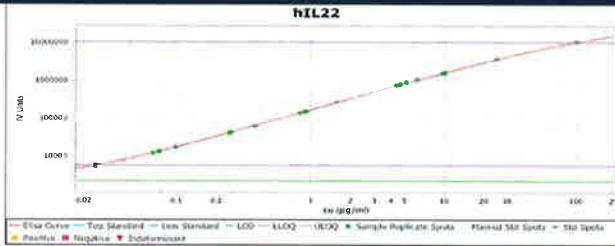
Calibrator Levels (pg/mL)	
1	300
2	75.0
3	18.8
4	4.69
5	1.17
6	0.293
7	0.073
8	0.000

IL-8 Calibration Curve



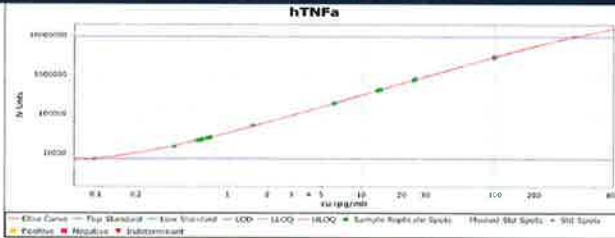
Calibrator Levels (pg/mL)	
1	400
2	100
3	25.0
4	6.25
5	1.56
6	0.391
7	0.098
8	0.000

IL-22 Calibration Curve



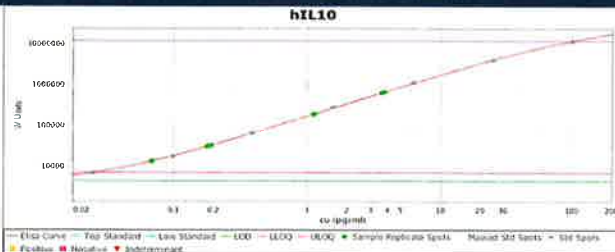
Calibrator Levels (pg/mL)	
1	100
2	25.0
3	6.25
4	1.56
5	0.391
6	0.098
7	0.024
8	0.000

TNFα Calibration Curve



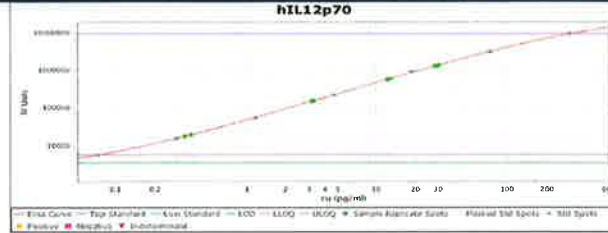
Calibrator Levels (pg/mL)	
1	400
2	100
3	25.0
4	6.25
5	1.56
6	0.391
7	0.098
8	0.000

IL-10 Calibration Curve



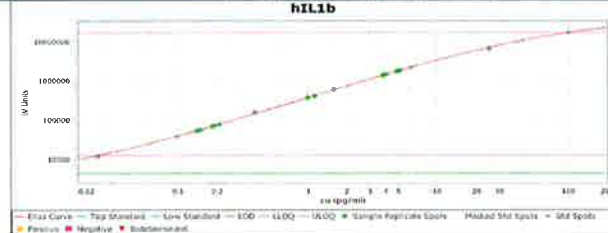
Calibrator Levels (pg/mL)	
1	100
2	25.0
3	6.25
4	1.56
5	0.391
6	0.098
7	0.024
8	0.000

IL-12p70 Calibration Curve



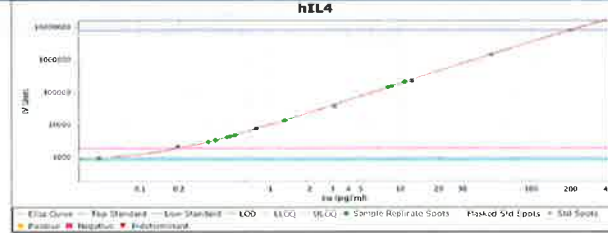
Calibrator Levels (pg/ml)	
1	300
2	75.0
3	18.8
4	4.69
5	1.17
6	0.293
7	0.0732
8	0.000

IL-1 β Calibration Curve



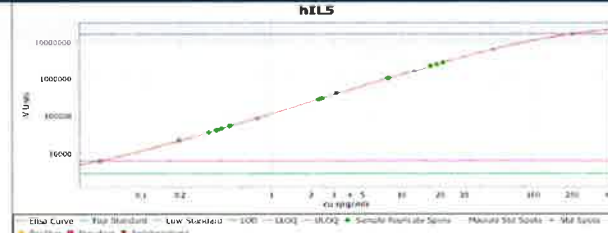
Calibrator Levels (pg/mL)	
1	100
2	25.0
3	6.25
4	1.56
5	0.391
6	0.098
7	0.024
8	0.000

IL-4 Calibration Curve



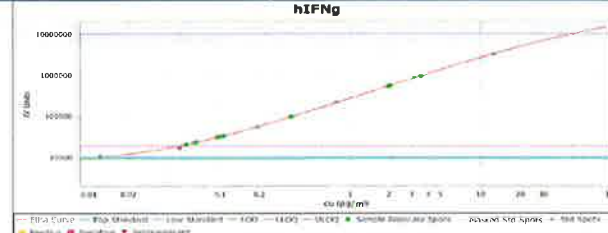
Calibrator Levels (pg/mL)	
1	200
2	50.0
3	12.5
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

IL-5 Calibration Curve



Calibrator Levels (pg/mL)	
1	200
2	50.0
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	50.0
3	12.5
4	3.13
5	0.781
6	0.195
7	0.049
8	0.000