

Product Number:	85-0329
Lot Number:	318086
Expiration:	13-Aug-2020
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	318086
Biotinylated Antibody Rgt	98-0329	318086
Sample Diluent 21 <sup>1</sup>	1863329	318027
IL-12p70 Calibrator	18910034	317901
IL-1β Calibrator	18910040	318067
IL-4 Calibrator	1858452	317999
IL-5 Calibrator	1858462	317984
IFNγ Calibrator	1858260	317863
IL-6 Calibrator	18910044	317985
IL-8 Calibrator	1858454	317862
IL-22 Calibrator	18910036	317907
TNFα Calibrator	1880012	317986
IL-10 Calibrator	18910043	318068
Wash Buffer (25X) <sup>1</sup>	1860055	17512600
Streptavidin-HRP Reagent <sup>1</sup>	1858263	318044
Super Signal <sup>™</sup> Luminol* <sup>1</sup>	1858089	TJ275117
Super Signal <sup>™</sup> Peroxide* <sup>1</sup>	1858090	TJ275118



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

<sup>1</sup> Reagents are not Kit Lot Specific

\*SuperSignal<sup>®</sup> technology is protected by US Patent #6,432,662 and is a registered trademark of ThermoFisher Scie

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	47.31	39.45 - 59.17
Control 2	1-Feb-19	N/A	N/A
Panel 1	1-Feb-19	114.3	94.96 - 142.4
Panel 2	1-Feb-19	13.83	10.87 - 16.30
Panel 3	1-Feb-19	1.471	0.726 - 2.340

**IL-1β**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	13.03	11.16 - 20.25
Control 2	1-Feb-19	0.614	0.492 - 1.060
Panel 1	1-Feb-19	0.535	0.265 - 1.147
Panel 2	1-Feb-19	3.530	3.020 - 5.126
Panel 3	1-Feb-19	14.40	12.42 - 21.26

**IL-4**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	32.51	26.16 - 39.23
Control 2	1-Feb-19	1.718	1.370 - 2.378
Panel 1	1-Feb-19	43.85	34.85 - 52.28
Panel 2	1-Feb-19	4.637	3.660 - 7.771
Panel 3	1-Feb-19	0.873	0.657 - 2.080

**IL-5**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	22.03	20.89 - 38.03
Control 2	1-Feb-19	1.188	1.018 - 2.005
Panel 1	1-Feb-19	54.52	53.78 - 81.53
Panel 2	1-Feb-19	6.577	6.136 - 9.655
Panel 3	1-Feb-19	0.922	0.850 - 1.762

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

Release Materials

IFN $\gamma$

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	6.781	6.302 - 9.193
Control 2	1-Feb-19	0.300	0.209 - 0.517
Panel 1	1-Feb-19	13.31	11.70 - 17.99
Panel 2	1-Feb-19	1.182	1.045 - 1.793
Panel 3	1-Feb-19	0.641	0.237 - 2.576

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	42.46	36.86 - 55.29
Control 2	1-Feb-19	2.158	1.904 - 2.959
Panel 1	1-Feb-19	71.56	58.81 - 88.21
Panel 2	1-Feb-19	9.79	8.376 - 13.08
Panel 3	1-Feb-19	1.381	1.133 - 2.549

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	54.82	48.71 - 73.06
Control 2	1-Feb-19	2.667	2.070 - 3.782
Panel 1	1-Feb-19	316.3	272.4 - 408.5
Panel 2	1-Feb-19	13.11	12.33 - 18.54
Panel 3	1-Feb-19	1.288	0.665 - 3.496

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	15.71	13.45 - 20.17
Control 2	1-Feb-19	0.689	0.621 - 1.127
Panel 1	1-Feb-19	26.93	24.22 - 37.21
Panel 2	1-Feb-19	2.405	2.074 - 4.357
Panel 3	1-Feb-19	0.225	0.101 - 0.649

TNF $\alpha$

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	45.19	39.20 - 68.99
Control 2	1-Feb-19	2.019	1.602 - 3.215
Panel 1	1-Feb-19	93.69	84.72 - 144.1
Panel 2	1-Feb-19	20.27	17.35 - 32.66
Panel 3	1-Feb-19	14.71	9.813 - 24.63

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	10.20	9.305 - 21.29
Control 2	1-Feb-19	0.508	0.395 - 1.014
Panel 1	1-Feb-19	0.565	0.495 - 0.977
Panel 2	1-Feb-19	3.347	3.160 - 6.238
Panel 3	1-Feb-19	0.279	0.214 - 0.422

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

Review/Approval

Patrizia Stadler  
Name

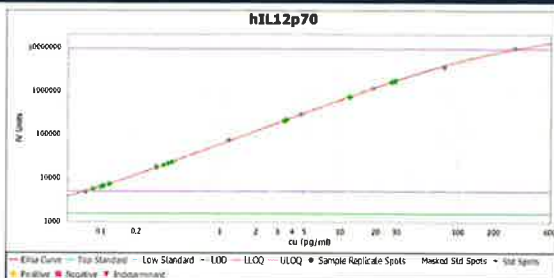
Associate  
Director of QC  
Title

*P Stadler*

16 MAR 2020

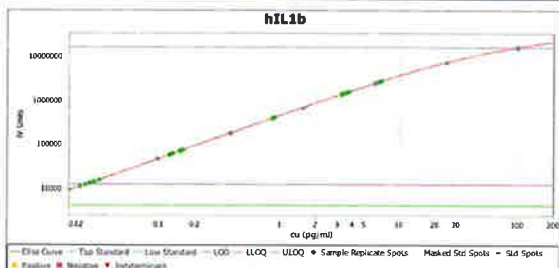
Signature/Date

**IL-12p70 Calibration Curve**



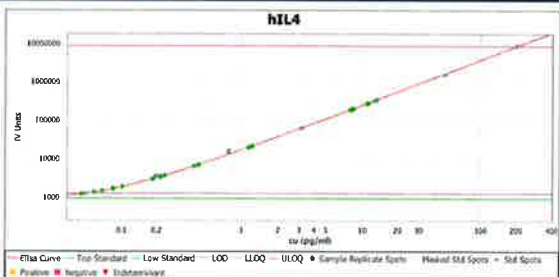
Calibrator Levels (pg/mL)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

**IL-1β Calibration Curve**



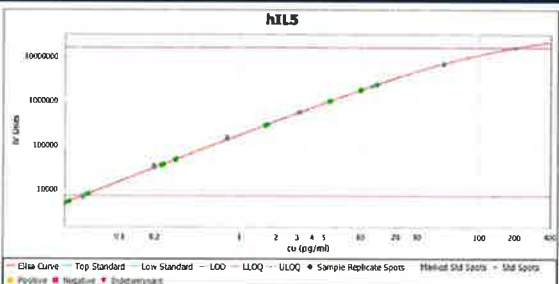
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

**IL-4 Calibration Curve**



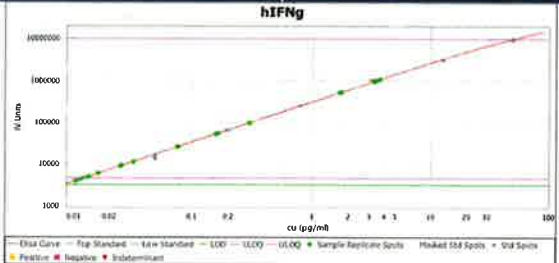
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

**IL-5 Calibration Curve**



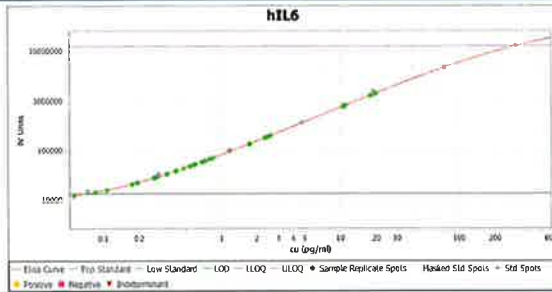
Calibrator Levels (pg/mL)	
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

**IFNγ Calibration Curve**



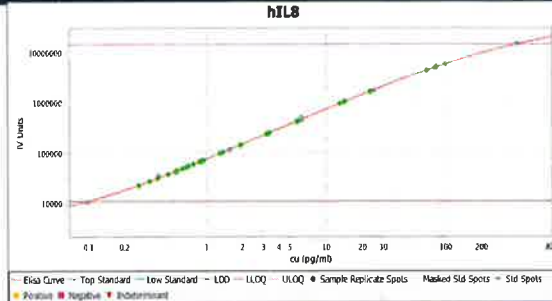
Calibrator Levels (pg/mL)	
1	50
2	12.5
3	3.125
4	0.781
5	0.195
6	0.049
7	0.012
8	0.000

**IL-6 Calibration Curve**



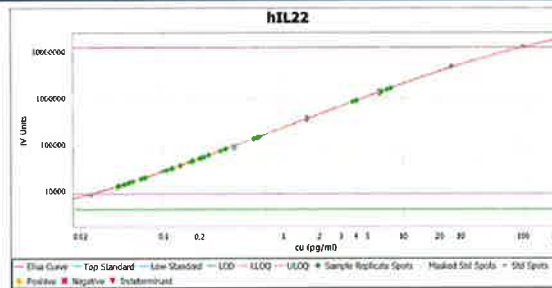
Calibrator Levels (pg/mL)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

**IL-8 Calibration Curve**



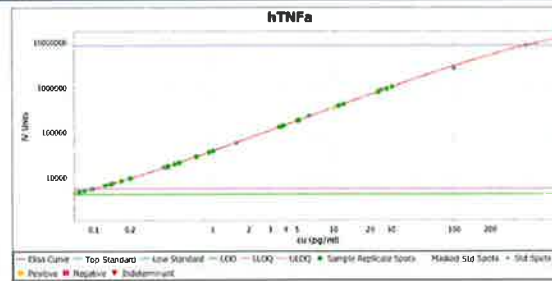
Calibrator Levels (pg/mL)	
1	400
2	100
3	25.00
4	6.250
5	1.563
6	0.391
7	0.098
8	0.000

**IL-22 Calibration Curve**



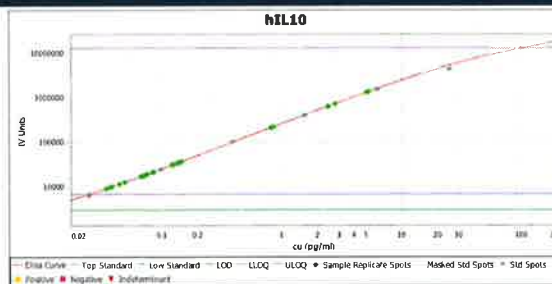
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

**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