

Product Number:	85-0329
Lot Number:	318046
Expiration:	13-Aug-2020
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	318046
Biotinylated Antibody Rgt	98-0329	318046
Sample Diluent 21 <sup>1</sup>	1863329	318027
IL-12p70 Calibrator	18910034	317876
IL-1 $\beta$ Calibrator	18910040	317793
IL-4 Calibrator	1858452	317726
IL-5 Calibrator	1858462	317984
IFN $\gamma$ Calibrator	1858260	317863
IL-6 Calibrator	18910044	317985
IL-8 Calibrator	1858454	317862
IL-22 Calibrator	18910036	317907
TNF $\alpha$ Calibrator	1880012	317986
IL-10 Calibrator	18910043	317877
Wash Buffer (25X) <sup>1</sup>	1860055	17512600
Streptavidin-HRP Reagent <sup>1</sup>	1858263	317962
Super Signal <sup>™</sup> Luminol* <sup>1</sup>	1858089	TJ275117
Super Signal <sup>™</sup> Peroxide* <sup>1</sup>	1858090	TJ275118

<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 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	46.95	37.77 - 56.65
Control 2	1-Feb-19	0.128	N/A
Panel 1	1-Feb-19	107.4	91.26 -136.9
Panel 2	1-Feb-19	14.39	10.21 - 15.32
Panel 3	1-Feb-19	2.099	0.864 - 2.578

IL-1 $\beta$

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	15.88	12.13 - 18.29
Control 2	1-Feb-19	0.823	0.535 - 0.984
Panel 1	1-Feb-19	0.700	0.364 - 0.923
Panel 2	1-Feb-19	4.056	3.219 - 4.828
Panel 3	1-Feb-19	18.56	14.03 - 21.04

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	33.27	24.91 - 37.37
Control 2	1-Feb-19	1.886	1.159 - 2.389
Panel 1	1-Feb-19	44.03	33.77 - 50.65
Panel 2	1-Feb-19	6.864	2.891 - 7.442
Panel 3	1-Feb-19	1.824	0.290 - 2.252

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	31.29	24.41 - 36.62
Control 2	1-Feb-19	1.658	1.254 - 1.882
Panel 1	1-Feb-19	72.28	56.42 - 84.63
Panel 2	1-Feb-19	9.231	6.609 - 9.914
Panel 3	1-Feb-19	1.465	1.017 - 1.690

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	7.441	6.094 - 9.142
Control 2	1-Feb-19	0.387	0.226 - 0.491
Panel 1	1-Feb-19	12.49	12.27 - 18.40
Panel 2	1-Feb-19	1.345	1.190 - 1.841
Panel 3	1-Feb-19	1.573	0.652 - 2.383

**IL-6**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	41.42	36.13 - 56.32
Control 2	1-Feb-19	2.241	1.689 - 3.013
Panel 1	1-Feb-19	67.50	57.22 - 85.83
Panel 2	1-Feb-19	12.010	8.131 - 12.71
Panel 3	1-Feb-19	1.603	0.878 - 2.460

**IL-8**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	58.04	38.73 - 83.91
Control 2	1-Feb-19	2.962	1.663 - 4.130
Panel 1	1-Feb-19	285.7	243.0 - 438.9
Panel 2	1-Feb-19	17.76	8.026 - 23.33
Panel 3	1-Feb-19	2.401	0.788 - 3.246

**IL-22**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	16.53	13.23 - 19.84
Control 2	1-Feb-19	0.867	0.550 - 1.109
Panel 1	1-Feb-19	27.25	22.83 - 35.10
Panel 2	1-Feb-19	3.738	2.027 - 4.121
Panel 3	1-Feb-19	0.469	0.078 - 0.580

**TNF $\alpha$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	50.74	42.94 - 76.80
Control 2	1-Feb-19	2.236	1.734 - 3.531
Panel 1	1-Feb-19	96.2	92.40 - 146.9
Panel 2	1-Feb-19	23.59	19.15 - 35.96
Panel 3	1-Feb-19	20.94	8.273 - 30.64

**IL-10**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	1-Feb-19	15.61	11.08 - 20.27
Control 2	1-Feb-19	0.778	0.436 - 0.985
Panel 1	1-Feb-19	0.725	0.448 - 1.079
Panel 2	1-Feb-19	4.643	3.489 - 6.614
Panel 3	1-Feb-19	0.366	0.157 - 0.487

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

Review/Approval

Patrizia Stadler QC Assoc Director

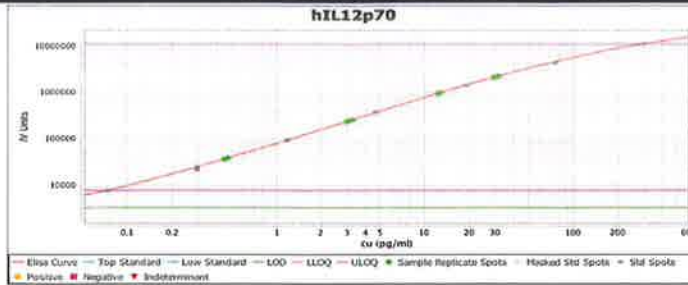
*P. Stadler* 04 April 2019

Name

Title

Signature/Date

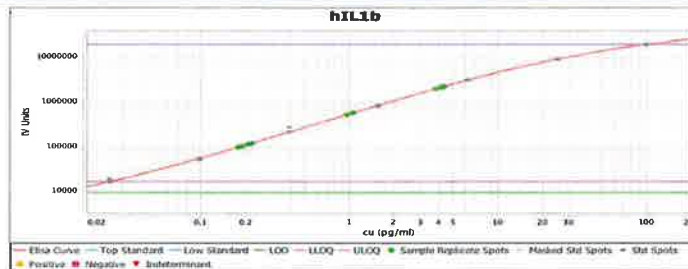
**IL-12p70 Calibration Curve**



**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

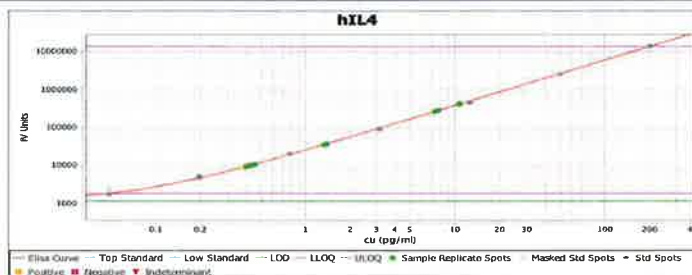
**IL-1β Calibration Curve**



**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

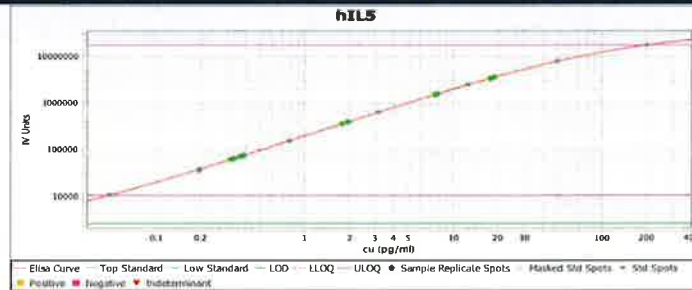
**IL-4 Calibration Curve**



**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

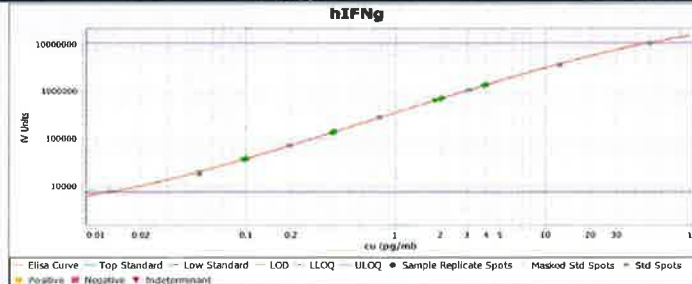
**IL-5 Calibration Curve**



**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

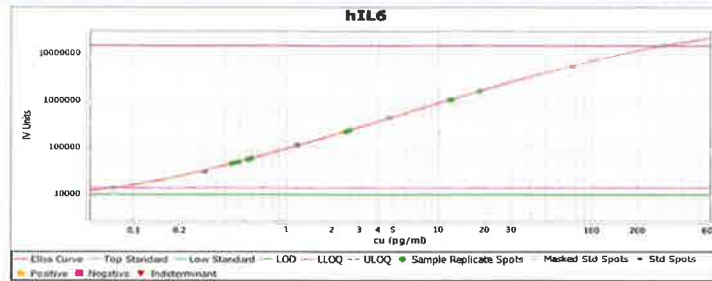
**IFNγ Calibration Curve**



**Calibrator Levels (pg/mL)**

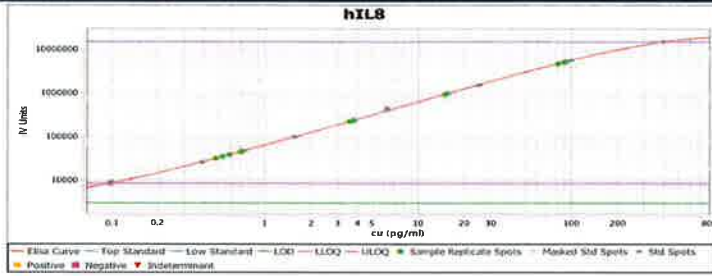
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

**IL-6 Calibration Curve**



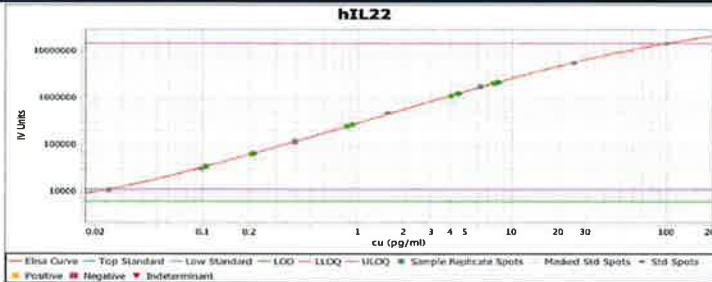
Calibrator Levels (pg/mL)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

**IL-8 Calibration Curve**



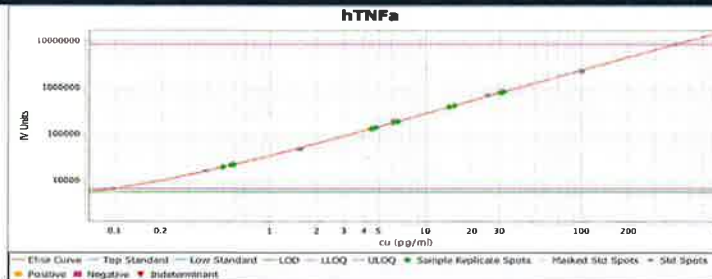
Calibrator Levels (pg/mL)	
1	400
2	100
3	25.00
4	6.250
5	1.563
6	0.391
7	0.098
8	0.000

**IL-22 Calibration Curve**



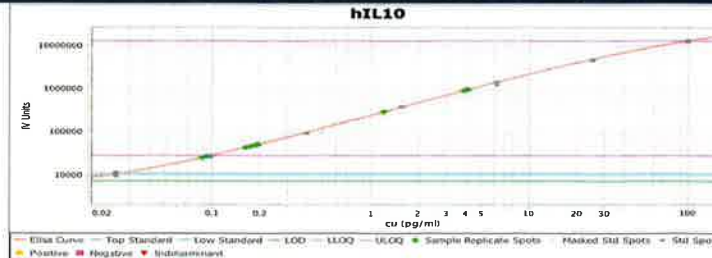
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

**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