

Product Number:	85-0329
Lot Number:	318735
Expiration:	12-Aug-2021
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	318627P
Biotinylated Antibody Rgt	98-0329	318734B
Sample Diluent 21 <sup>1</sup>	1863329	N/A
IL-12p70 Calibrator	18910034	318571
IL-1 $\beta$ Calibrator	18910040	318673
IL-4 Calibrator	1858452	318480
IL-5 Calibrator	1858462	318465
IFN $\gamma$ Calibrator	1858260	318538
IL-6 Calibrator	18910044	318583
IL-8 Calibrator	1858454	318342
IL-22 Calibrator	18910036	318408
TNF $\alpha$ Calibrator	1880012	318436
IL-10 Calibrator	18910043	318630
Wash Buffer (25X) <sup>1</sup>	1860055	N/A
Streptavidin-HRP Reagent <sup>1</sup>	1858263	N/A
Super Signal <sup>®</sup> Luminol* <sup>1</sup>	1858089	N/A
Super Signal <sup>®</sup> Peroxide* <sup>1</sup>	1858090	N/A



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 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)
Panel 1	318641	174.2	74.90	- 312.8
Panel 2	318641	59.37	25.27	- 105.7
Panel 3	318641	21.16	8.118	- 35.35
Panel 4	318641	7.395	2.411	- 12.41
Panel 5	318641	2.559	0.828	- 4.244

**IL-1 $\beta$**

	Lot	Result (pg/mL)	Mean	Range (pg/mL)
Panel 1	318641	23.94	9.353	- 36.93
Panel 2	318641	8.881	3.051	- 12.44
Panel 3	318641	3.073	0.942	- 4.185
Panel 4	318641	1.130	0.288	- 1.506
Panel 5	318641	0.427	0.053	- 0.563

**IL-4**

	Lot	Result (pg/mL)	Mean	Range (pg/mL)
Panel 1	318641	73.36	31.71	- 139.3
Panel 2	318641	26.67	10.51	- 46.13
Panel 3	318641	9.355	3.155	- 16.82
Panel 4	318641	3.835	1.070	- 6.178
Panel 5	318641	1.239	0.416	- 2.046

**IL-5**

	Lot	Result (pg/mL)	Mean	Range (pg/mL)
Panel 1	318641	60.90	23.48	- 97.00
Panel 2	318641	21.73	7.797	- 32.93
Panel 3	318641	7.610	2.673	- 11.72
Panel 4	318641	2.829	0.879	- 4.207
Panel 5	318641	0.987	0.306	- 1.549

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)
Panel 1	318641	10.75	4.077	- 16.34
Panel 2	318641	3.960	1.401	- 5.834
Panel 3	318641	1.388	0.475	- 2.058
Panel 4	318641	0.551	0.163	- 0.798
Panel 5	318641	0.228	0.067	- 0.338

**IL-6**

	Lot	Result (pg/mL)	Mean	Range (pg/mL)
Panel 1	318641	122.2	53.84	- 235.8
Panel 2	318641	4.542	1.795	- 9.227
Panel 3	318641	1.477	0.585	- 3.138
Panel 4	318641	0.599	0.219	- 1.262
Panel 5	318641	0.240	0.214	- 0.437

**IL-8**

	Lot	Result (pg/mL)	Mean	Range (pg/mL)
Panel 1	318641	458.7	175.2	- 729.0
Panel 2	318641	122.4	53.79	- 240.7
Panel 3	318641	43.83	18.28	- 84.37
Panel 4	318641	15.20	5.861	- 29.62
Panel 5	318641	5.398	0.997	- 10.30

**IL-22**

	Lot	Result (pg/mL)	Mean	Range (pg/mL)
Panel 1	318641	10.92	8.132	- 41.07
Panel 2	318641	3.350	2.348	- 11.98
Panel 3	318641	1.100	0.776	- 4.049
Panel 4	318641	0.408	0.274	- 1.445
Panel 5	318641	0.139	0.080	- 0.499

**TNF $\alpha$**

	Lot	Result (pg/mL)	Mean	Range (pg/mL)
Panel 1	318641	75.30	28.83	- 127.9
Panel 2	318641	24.81	10.37	- 45.07
Panel 3	318641	8.39	3.411	- 15.41
Panel 4	318641	3.166	1.115	- 5.442
Panel 5	318641	1.393	0.402	- 2.035

**IL-10**

	Lot	Result (pg/mL)	Mean	Range (pg/mL)
Panel 1	318641	61.79	22.96	- 85.32
Panel 2	318641	22.11	7.491	- 30.02
Panel 3	318641	8.230	2.644	- 10.62
Panel 4	318641	3.254	0.964	- 3.944
Panel 5	318641	1.266	0.339	- 1.486

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*

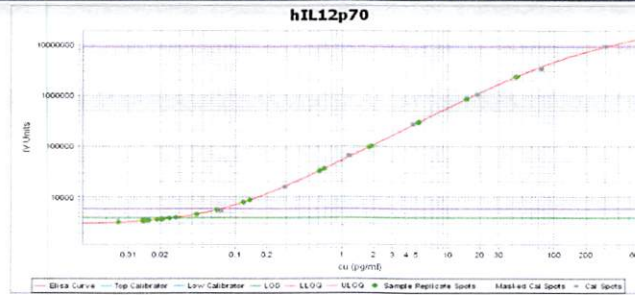
29 May 2020

Name

Title

Signature/Date

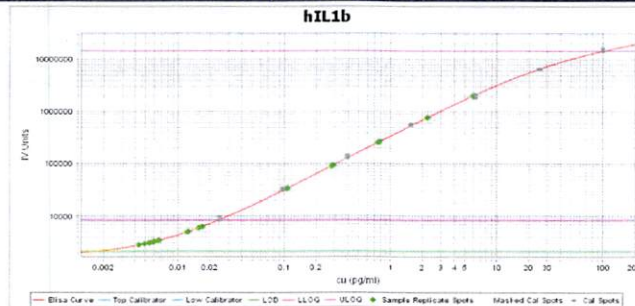
**IL-12p70 Calibration Curve**



**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

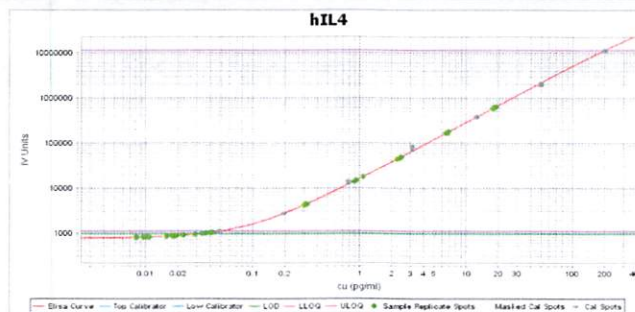
**IL-1β Calibration Curve**



**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

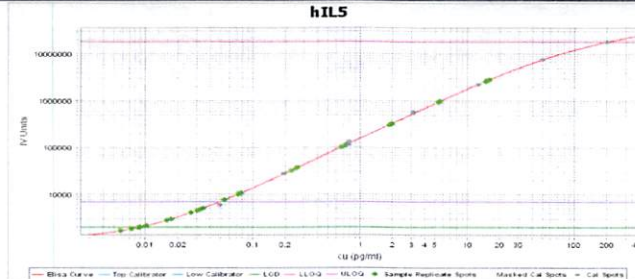
**IL-4 Calibration Curve**



**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

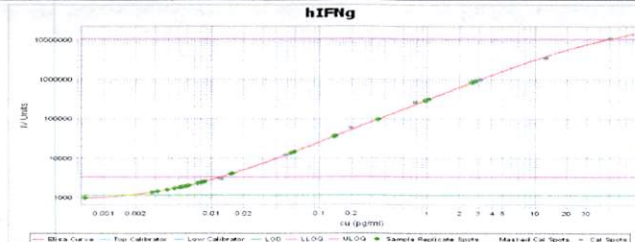
**IL-5 Calibration Curve**



**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

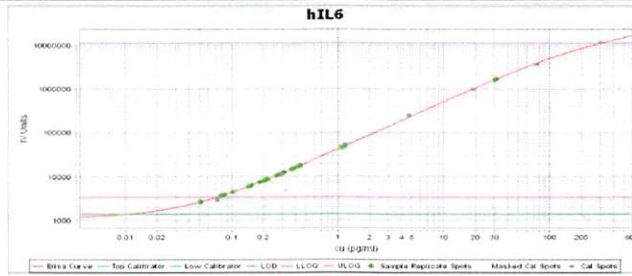
**IFNγ Calibration Curve**



**Calibrator Levels (pg/mL)**

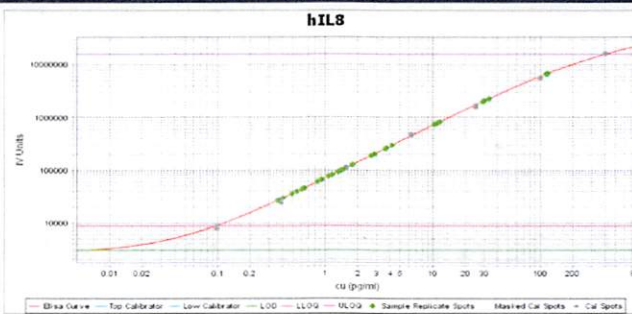
1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

**IL-6 Calibration Curve**



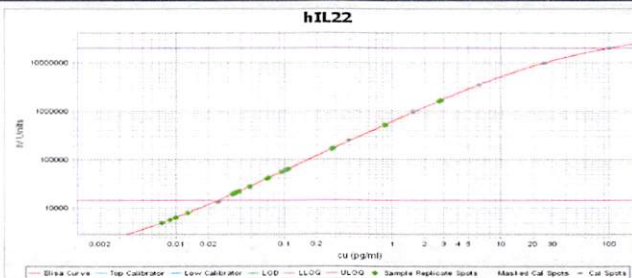
Calibrator Levels (pg/mL)	
1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

**IL-8 Calibration Curve**



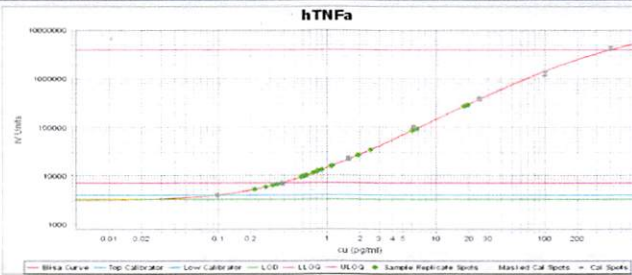
Calibrator Levels (pg/mL)	
1	400
2	100
3	25.00
4	6.250
5	1.563
6	0.391
7	0.098
8	0.000

**IL-22 Calibration Curve**



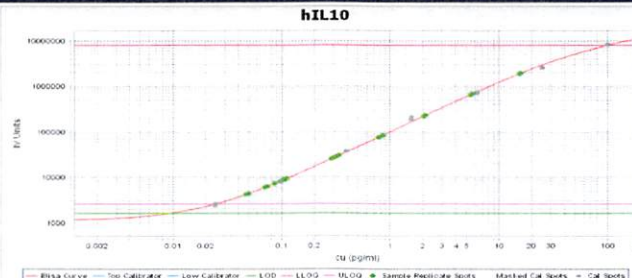
Calibrator Levels (pg/mL)	
1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

**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