

Product Number:	85-0329
Lot Number:	318737
Expiration:	18-Sep-2021
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	318657P
Biotinylated Antibody Rgt	98-0329	318684B
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<sup>1</sup> Kit Instructions (85A-0329).

**Release Materials**

**IL-12p70**

	Lot	Result (pg/mL)	Mean	Range (pg/mL)
Panel 1	318641	160.90	74.90	- 312.8
Panel 2	318641	55.07	25.27	- 105.7
Panel 3	318641	21.88	8.118	- 35.35
Panel 4	318641	7.13	2.411	- 12.41
Panel 5	318641	2.47	0.828	- 4.244

**IL-1 $\beta$**

	Lot	Result (pg/mL)	Mean	Range (pg/mL)
Panel 1	318641	21.2	9.353	- 36.93
Panel 2	318641	7.49	3.051	- 12.44
Panel 3	318641	3.15	0.942	- 4.185
Panel 4	318641	1.040	0.288	- 1.506
Panel 5	318641	0.370	0.053	- 0.563

**IL-4**

	Lot	Result (pg/mL)	Mean	Range (pg/mL)
Panel 1	318641	66.4	31.71	- 139.3
Panel 2	318641	21.4	10.51	- 46.13
Panel 3	318641	8.58	3.155	- 16.82
Panel 4	318641	2.93	1.070	- 6.178
Panel 5	318641	1.03	0.416	- 2.046

**IL-5**

	Lot	Result (pg/mL)	Mean	Range (pg/mL)
Panel 1	318641	58.0	23.48	- 97.00
Panel 2	318641	19.3	7.797	- 32.93
Panel 3	318641	8.09	2.673	- 11.72
Panel 4	318641	2.54	0.879	- 4.207
Panel 5	318641	0.980	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)	an Range (pg/mL)	
Panel 1	318641	9.27	4.077	- 16.34
Panel 2	318641	3.15	1.401	- 5.834
Panel 3	318641	1.29	0.475	- 2.058
Panel 4	318641	0.440	0.163	- 0.798
Panel 5	318641	0.190	0.067	- 0.338

**IL-6**

	Lot	Result (pg/mL)	an Range (pg/mL)	
Panel 1	318641	110	53.84	- 235.8
Panel 2	318641	3.55	1.795	- 9.227
Panel 3	318641	1.25	0.585	- 3.138
Panel 4	318641	0.440	0.219	- 1.262
Panel 5	318641	0.290	0.214	- 0.437

**IL-8**

	Lot	Result (pg/mL)	an Range (pg/mL)	
Panel 1	318641	384	175.2	- 729.0
Panel 2	318641	122.1	53.79	- 240.7
Panel 3	318641	48.6	18.28	- 84.37
Panel 4	318641	15.3	5.861	- 29.62
Panel 5	318641	5.10	0.997	- 10.30

**IL-22**

	Lot	Result (pg/mL)	an Range (pg/mL)	
Panel 1	318641	9.5	8.132	- 41.07
Panel 2	318641	2.68	2.348	- 11.98
Panel 3	318641	0.97	0.776	- 4.049
Panel 4	318641	0.300	0.274	- 1.445
Panel 5	318641	0.110	0.080	- 0.499

**TNF $\alpha$**

	Lot	Result (pg/mL)	an Range (pg/mL)	
Panel 1	318641	55.8	28.83	- 127.9
Panel 2	318641	20.2	10.37	- 45.07
Panel 3	318641	8.16	3.411	- 15.41
Panel 4	318641	2.65	1.115	- 5.442
Panel 5	318641	1.01	0.402	- 2.035

**IL-10**

	Lot	Result (pg/mL)	an Range (pg/mL)	
Panel 1	318641	60.2	22.96	- 85.32
Panel 2	318641	20.8	7.491	- 30.02
Panel 3	318641	8.40	2.644	- 10.62
Panel 4	318641	2.83	0.964	- 3.944
Panel 5	318641	1.090	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*

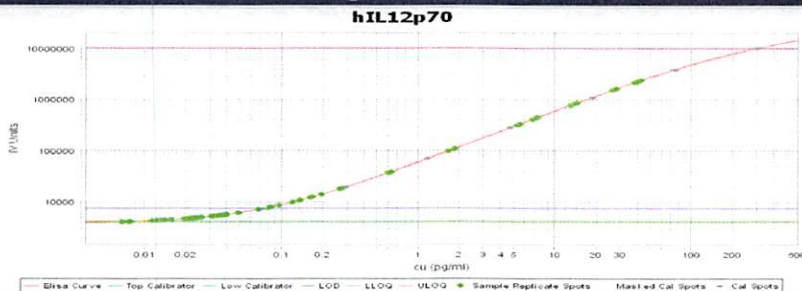
27 May 2020

Name

Title

Signature/Date

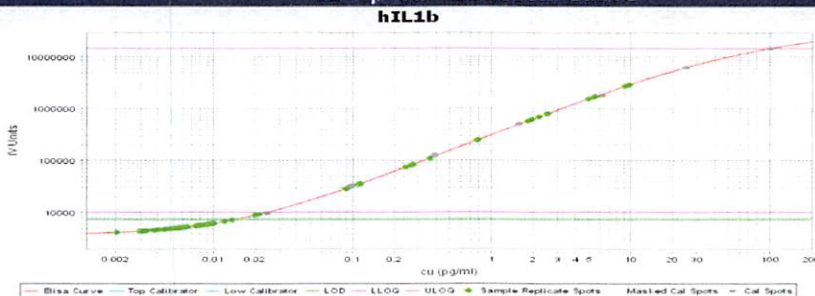
**IL-12p70 Calibration Curve**



**Calibrator Levels (pg/ml)**

1	300
2	75
3	
4	4.69
5	1.17
6	0.293
7	0.073
8	0.000

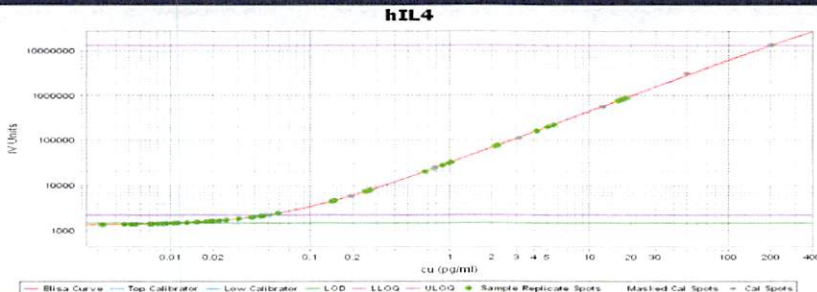
**IL-1β Calibration Curve**



**Calibrator Levels (pg/mL)**

1	100
2	25
3	6.25
4	1.56
5	0.391
6	0.098
7	0.024
8	0.000

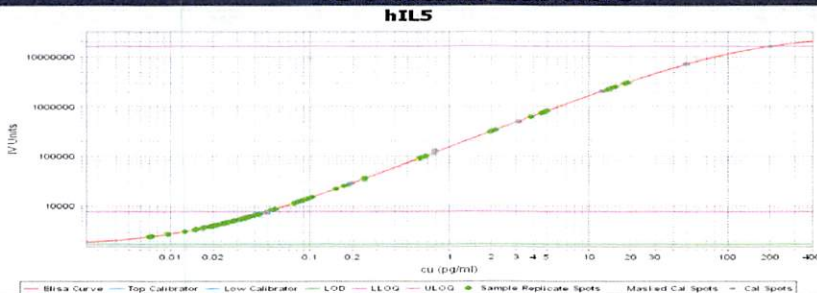
**IL-4 Calibration Curve**



**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

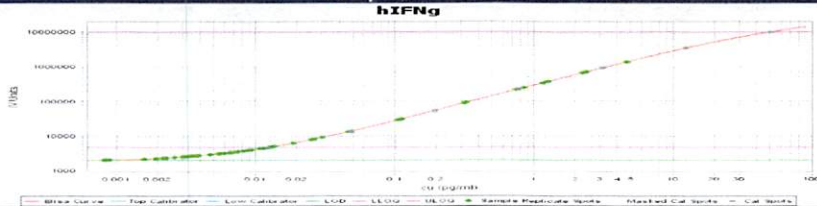
**IL-5 Calibration Curve**



**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

**IFNγ Calibration Curve**

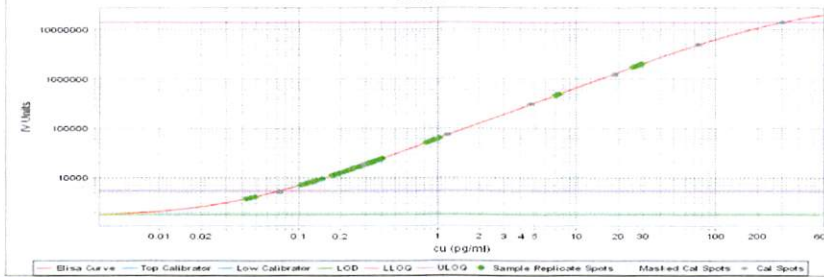


**Calibrator Levels (pg/mL)**

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**

**hIL6**

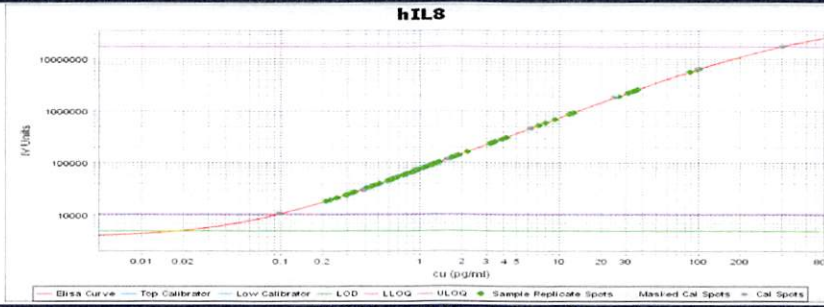


**Calibrator Levels (pg/mL)**

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**

**hIL8**

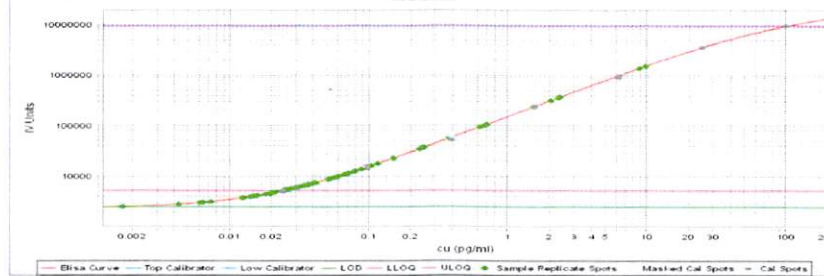


**Calibrator Levels (pg/mL)**

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**

**hIL22**

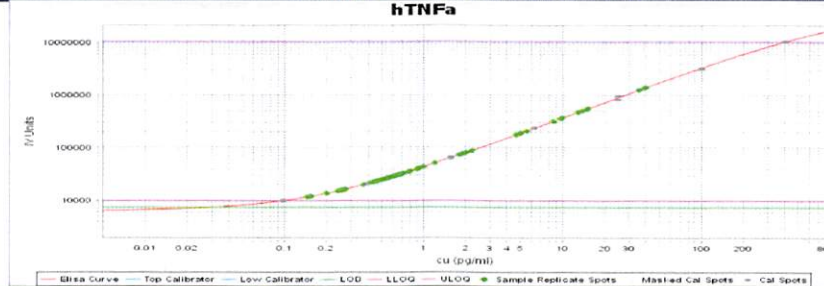


**Calibrator Levels (pg/mL)**

1	100
2	25
3	6.25
4	1.56
5	0.391
6	0.098
7	0.024
8	0.000

**TNFA Calibration Curve**

**hTNFA**

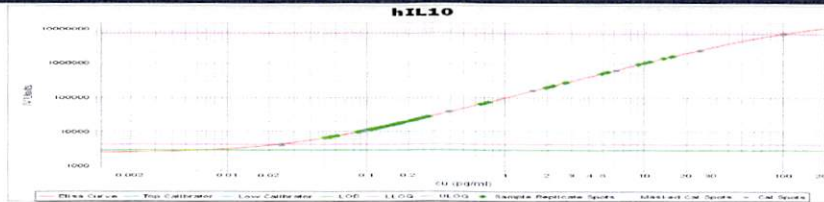


**Calibrator Levels (pg/mL)**

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**

**hIL10**



**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