

<b>Product Number:</b>	85-0329
<b>Lot Number:</b>	504244
<b>Expiration:</b>	5-Mar-2025
<b>Platform(s):</b>	SP-X

CorPlex Cytokine Plate	97-0329	411011P
Biotinylated Antibody Reagent	98-0329	409505B
Sample Diluent 21	1863329	319785
IL-12p70 Calibrator	18910034	319844
IL-1β Calibrator	18910040	319564
IL-4 Calibrator	1858452	318580
IL-5 Calibrator	1858462	318655
IFNγ Calibrator	1858260	319589
IL-6 Calibrator	18910044	319173
IL-8 Calibrator	1858454	319436
IL-22 Calibrator	18910036	319176
TNFα Calibrator	1880012	403812
IL-10 Calibrator	18910043	319842
Wash Buffer (12.5X) <sup>1</sup>	1863332	N/A
Streptavidin-HRP Reagent <sup>1</sup>	1858263	N/A
Super Signal® Luminol <sup>1,2</sup>	1858089	N/A
Super Signal® Peroxide <sup>1,2</sup>	1858090	N/A

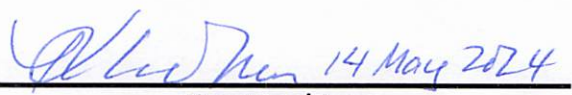


*Data below represents results generated on the SP-X Imaging and Analysis System.*

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

<sup>2</sup> 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 Simoa Complex Cytokine Panel 1 Instruction Book 85-0329 116-7BF-1-AB (KI-0195).

Review/Approval	Kathleen Cornell	Quality Control Supervisor	
	Name	Title	Signature/Date

**Release Materials\***

**IL-12p70**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
Panel 1	319839	147	107 - 186
Panel 2	319839	43.1	33.0 - 51.2
Panel 3	319839	17.7	12.3 - 20.3
Panel 4	319839	4.65	2.59 - 6.31
Panel 5	319839	1.41	0.728 - 2.17

**IL-1β**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
Panel 1	319839	16.4	9.55 - 20.6
Panel 2	319839	4.77	2.82 - 5.78
Panel 3	319839	1.83	1.07 - 2.22
Panel 4	319839	0.562	0.237 - 0.723
Panel 5	319839	0.184	0.081 - 0.271

**IL-4**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
Panel 1	319839	79.1	50.1 - 97.2
Panel 2	319839	24.0	14.5 - 27.5
Panel 3	319839	9.51	5.89 - 10.8
Panel 4	319839	2.72	1.59 - 3.47
Panel 5	319839	1.05	0.553 - 1.34

**IL-5**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
Panel 1	319839	53.4	31.5 - 70.5
Panel 2	319839	18.1	9.01 - 21.6
Panel 3	319839	7.68	3.34 - 8.64
Panel 4	319839	2.34	0.961 - 2.50
Panel 5	319839	0.941	0.370 - 0.968

**IFNγ**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
Panel 1	319839	8.13	5.18 - 14.5
Panel 2	319839	2.68	1.55 - 4.66
Panel 3	319839	1.05	0.710 - 1.83
Panel 4	319839	0.324	0.217 - 0.604
Panel 5	319839	0.150	0.089 - 0.283

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

**Release Materials\***

**IL-6**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	208	149 - 299
<b>Panel 2</b>	319839	55.7	39.7 - 70.6
<b>Panel 3</b>	319839	22.1	14.2 - 29.4
<b>Panel 4</b>	319839	7.76	4.15 - 11.2
<b>Panel 5</b>	319839	2.81	1.32 - 4.52

**IL-8**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	284	183 - 383
<b>Panel 2</b>	319839	75.7	49.8 - 103
<b>Panel 3</b>	319839	27.1	19.9 - 39.3
<b>Panel 4</b>	319839	8.47	5.82 - 12.6
<b>Panel 5</b>	319839	3.16	2.12 - 4.95

**IL-22**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	86.8	36.2 - 101
<b>Panel 2</b>	319839	11.4	6.05 - 15.0
<b>Panel 3</b>	319839	6.41	2.75 - 7.32
<b>Panel 4</b>	319839	1.24	0.565 - 1.49
<b>Panel 5</b>	319839	0.322	0.149 - 0.446

**TNFα**

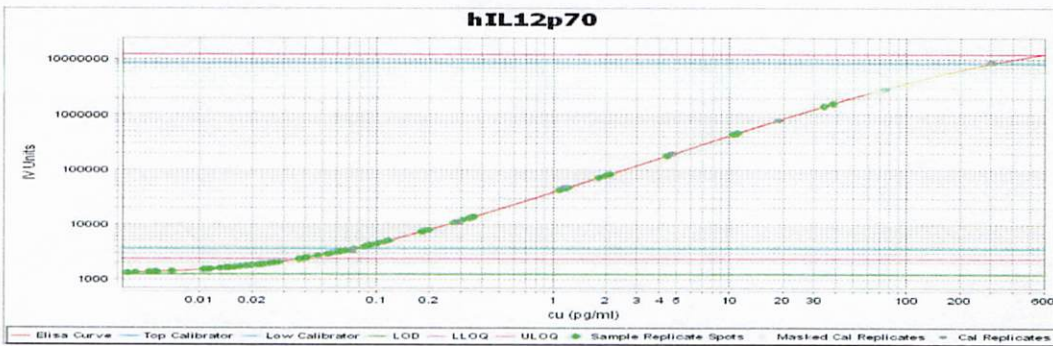
	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	60.2	37.9 - 99.8
<b>Panel 2</b>	319839	17.9	11.8 - 26.5
<b>Panel 3</b>	319839	6.68	4.09 - 10.7
<b>Panel 4</b>	319839	1.35	1.01 - 2.84
<b>Panel 5</b>	319839	0.514	0.369 - 1.09

**IL-10**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	43.8	30.4 - 61.9
<b>Panel 2</b>	319839	13.6	9.02 - 17.8
<b>Panel 3</b>	319839	5.45	3.19 - 7.61
<b>Panel 4</b>	319839	1.57	1.01 - 2.34
<b>Panel 5</b>	319839	0.728	0.409 - 1.02

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

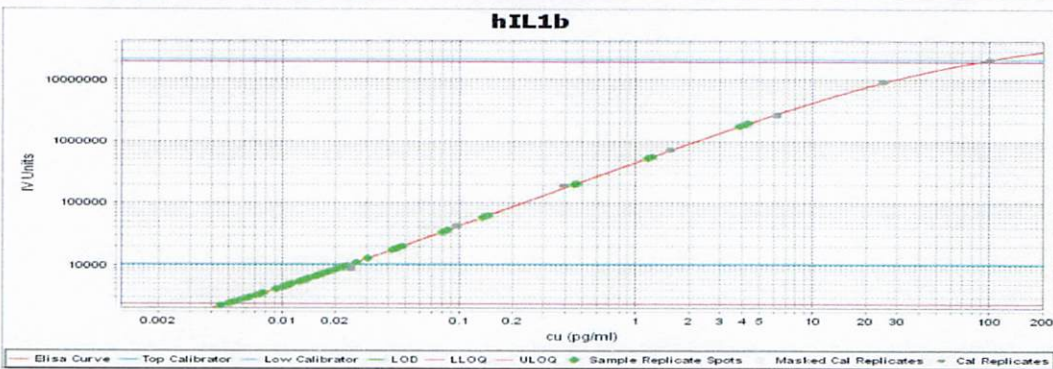
IL-12p70 Calibrator Curve



Calibrator Levels (pg/mL)

A	300
B	75.0
C	18.8
D	4.69
E	1.17
F	0.293
G	0.073
H	0.000

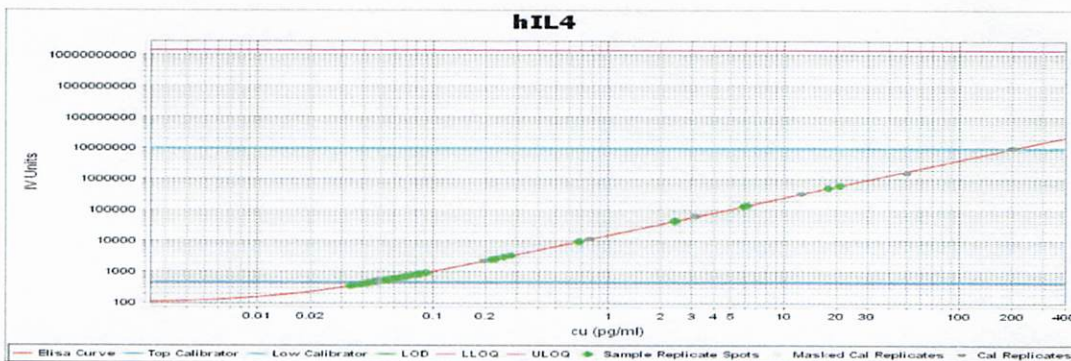
IL-1β Calibrator Curve



Calibrator Levels (pg/mL)

A	100
B	25.0
C	6.25
D	1.56
E	0.391
F	0.098
G	0.024
H	0.000

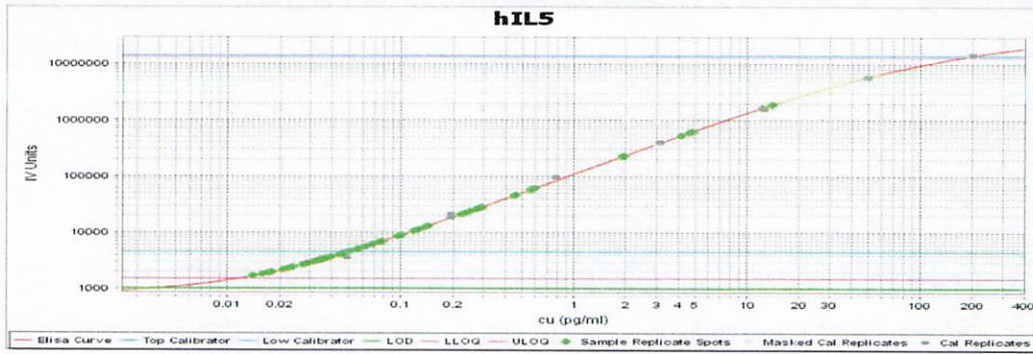
IL-4 Calibrator Curve



Calibrator Levels (pg/mL)

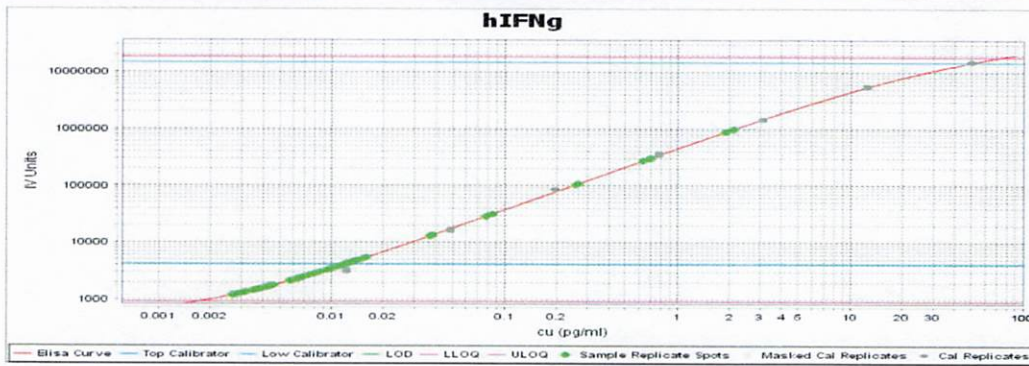
A	200
B	50.0
C	12.5
D	3.13
E	0.781
F	0.195
G	0.049
H	0.000

**IL-5 Calibrator Curve**



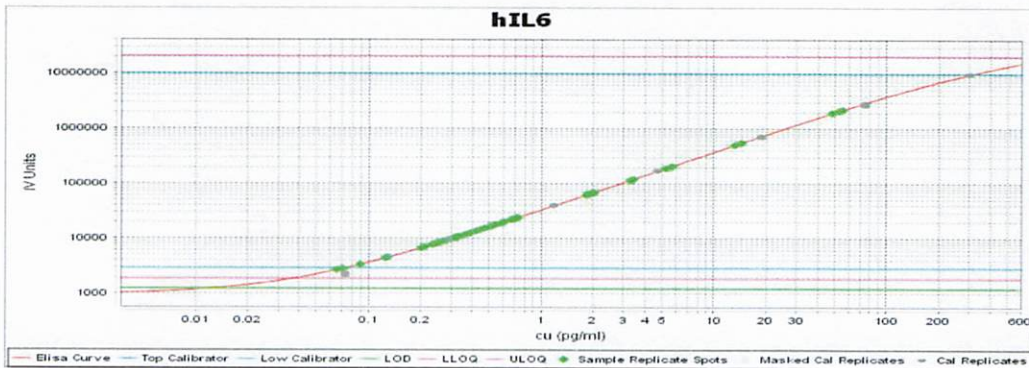
Calibrator Levels (pg/mL)	
A	200
B	50.0
C	12.5
D	3.13
E	0.781
F	0.195
G	0.049
H	0.000

**IFNγ Calibrator Curve**



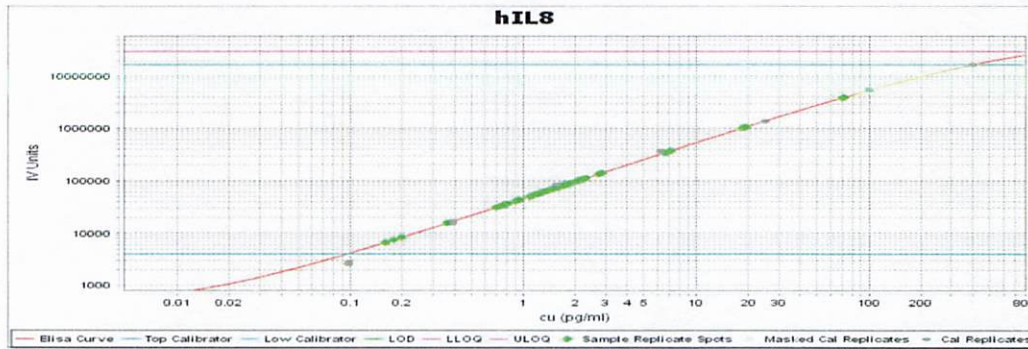
Calibrator Levels (pg/mL)	
A	50.0
B	12.5
C	3.13
D	0.781
E	0.195
F	0.049
G	0.012
H	0.000

**IL-6 Calibrator Curve**



Calibrator Levels (pg/mL)	
A	300
B	75.0
C	18.8
D	4.69
E	1.17
F	0.293
G	0.073
H	0.000

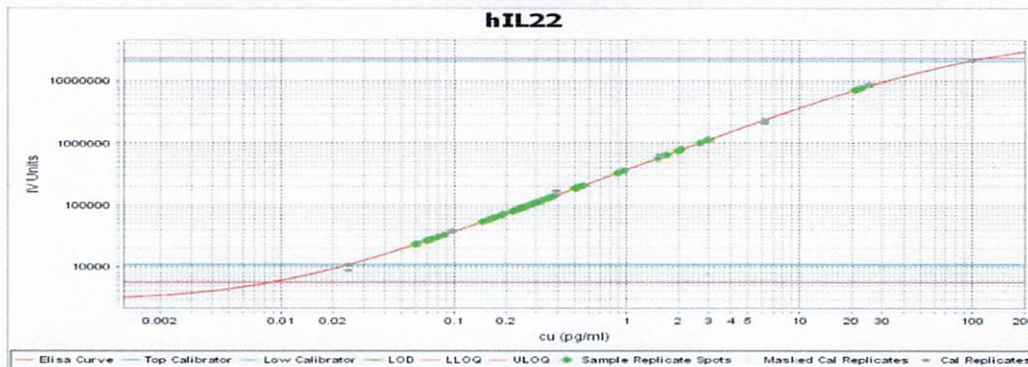
**IL-8 Calibrator Curve**



**Calibrator Levels (pg/mL)**

A	400
B	100
C	25.0
D	6.25
E	1.56
F	0.391
G	0.098
H	0.000

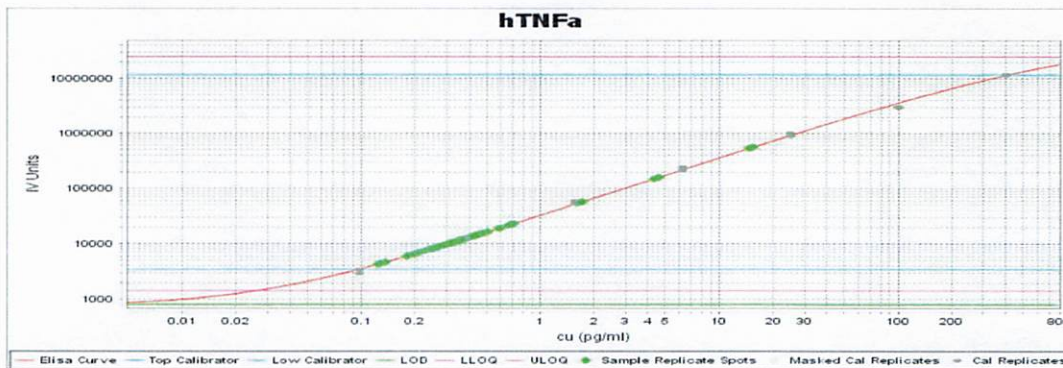
**IL-22 Calibrator Curve**



**Calibrator Levels (pg/mL)**

A	100
B	25.0
C	6.25
D	1.56
E	0.391
F	0.098
G	0.024
H	0.000

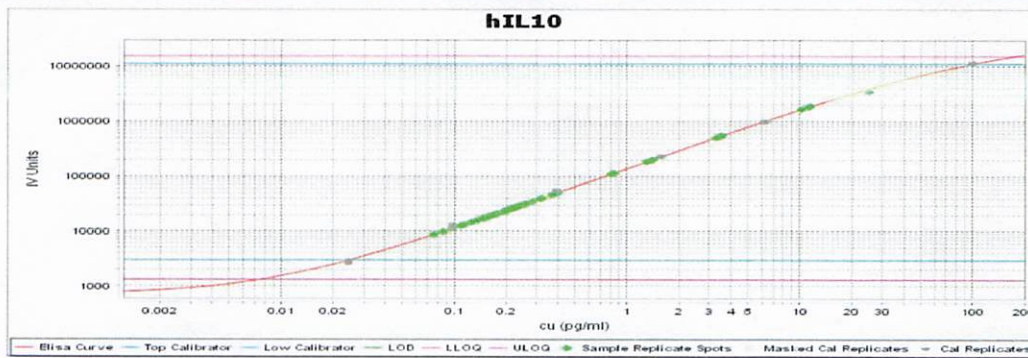
**TNFα Calibrator Curve**



**Calibrator Levels (pg/mL)**

A	400
B	100
C	25.0
D	6.25
E	1.56
F	0.391
G	0.098
H	0.000

**IL-10 Calibrator Curve**



**Calibrator Levels (pg/mL)**

<b>A</b>	100
<b>B</b>	25.0
<b>C</b>	6.25
<b>D</b>	1.56
<b>E</b>	0.391
<b>F</b>	0.098
<b>G</b>	0.024
<b>H</b>	0.000