

<b>Product Number:</b>	85-0329
<b>Lot Number:</b>	504475
<b>Expiration:</b>	26-Mar-2026
<b>Platform(s):</b>	SP-X



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

CorPlex Cytokine Plate	97-0329	433902P
Biotinylated Antibody Reagent	98-0329	427108B
Sample Diluent 21	1863329	405709
IL-12p70 Calibrator	18910034	319844
IL-1β Calibrator	18910040	407211
IL-4 Calibrator	1858452	319438
IL-5 Calibrator	1858462	319838
IFNγ Calibrator	1858260	319843
IL-6 Calibrator	18910044	324303
IL-8 Calibrator	1858454	319436
IL-22 Calibrator	18910036	319459
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

<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 Corplex Cytokine Panel 1 Instruction Book 85-0329 116-7BF-1-AB (KI-0195).

Review/Approval

*Kyle Giroux*

Name

*Senior Document Control Specialist, QA*

Title

*Kyle Giroux 21 Jan 2025*

Signature/Date

**Release Materials\***

**IL-12p70**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	145	107 - 185
<b>Panel 2</b>	319839	41.4	33.4 - 51.2
<b>Panel 3</b>	319839	16.6	12.3 - 20.7
<b>Panel 4</b>	319839	4.65	2.67 - 6.35
<b>Panel 5</b>	319839	1.62	0.760 - 2.18

**IL-1β**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	14.9	9.79 - 20.4
<b>Panel 2</b>	319839	4.32	2.85 - 5.89
<b>Panel 3</b>	319839	1.81	1.06 - 2.31
<b>Panel 4</b>	319839	0.512	0.238 - 0.745
<b>Panel 5</b>	319839	0.192	0.084 - 0.274

**IL-4**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	85.5	50.9 - 96.0
<b>Panel 2</b>	319839	25.9	14.8 - 27.5
<b>Panel 3</b>	319839	10.4	5.94 - 11.0
<b>Panel 4</b>	319839	3.36	1.59 - 3.56
<b>Panel 5</b>	319839	1.61	0.558 - 1.37

**IL-5**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	59.6	31.3 - 73.2
<b>Panel 2</b>	319839	18.0	8.91 - 22.5
<b>Panel 3</b>	319839	8.01	3.45 - 9.05
<b>Panel 4</b>	319839	2.35	0.829 - 2.79
<b>Panel 5</b>	319839	1.04	0.365 - 1.01

**IFNγ**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	11.8	5.23 - 14.5
<b>Panel 2</b>	319839	3.83	1.52 - 4.74
<b>Panel 3</b>	319839	1.69	0.711 - 1.85
<b>Panel 4</b>	319839	0.538	0.216 - 0.614
<b>Panel 5</b>	319839	0.296	0.087 - 0.288

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

**Release Materials\***

**IL-6**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	204	148 - 298
Panel 2	319839	54.4	40.5 - 70.0
Panel 3	319839	21.2	14.6 - 29.2
Panel 4	319839	7.81	4.31 - 11.1
Panel 5	319839	3.18	1.40 - 4.47

**IL-8**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	328	181 - 395
Panel 2	319839	88.4	49.8 - 105
Panel 3	319839	33.2	19.8 - 40.6
Panel 4	319839	10.1	5.79 - 13.0
Panel 5	319839	3.82	2.14 - 5.07

**IL-22**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	49.6	33.7 - 103
Panel 2	319839	7.59	5.99 - 15.1
Panel 3	319839	3.76	2.56 - 7.64
Panel 4	319839	0.802	0.559 - 1.51
Panel 5	319839	0.284	0.148 - 0.449

**TNFα**

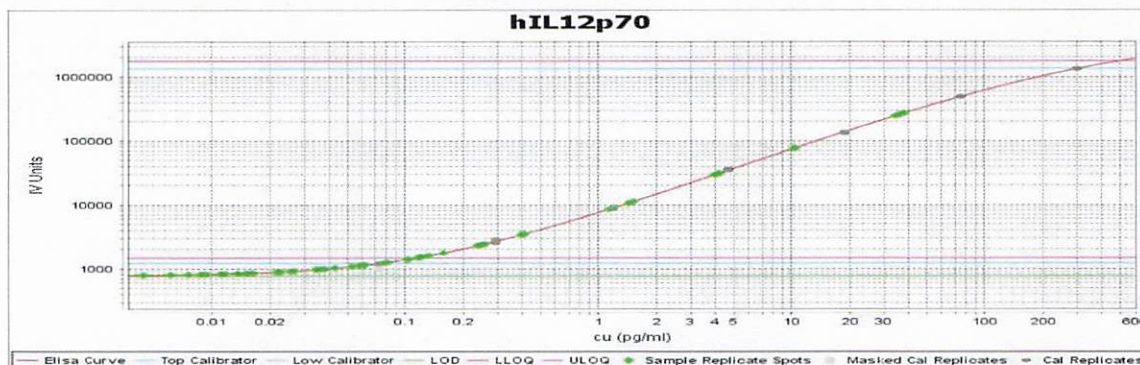
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	80.4	35.7 - 99.4
Panel 2	319839	24.6	12.2 - 26.3
Panel 3	319839	9.75	4.26 - 10.6
Panel 4	319839	2.55	1.05 - 2.81
Panel 5	319839	1.10	0.378 - 1.07

**IL-10**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	48.9	29.2 - 62.2
Panel 2	319839	14.2	9.21 - 17.7
Panel 3	319839	5.67	3.28 - 7.66
Panel 4	319839	1.68	1.02 - 2.39
Panel 5	319839	0.750	0.424 - 1.02

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

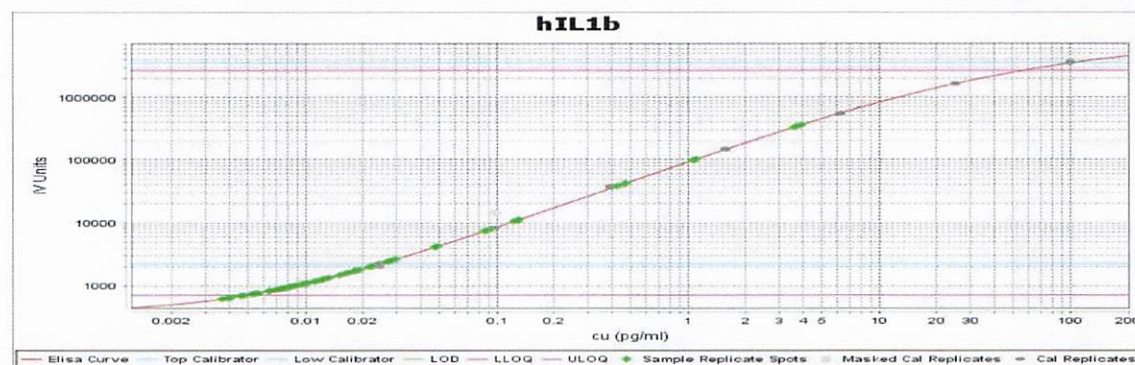
**IL-12p70 Calibrator Curve**



**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

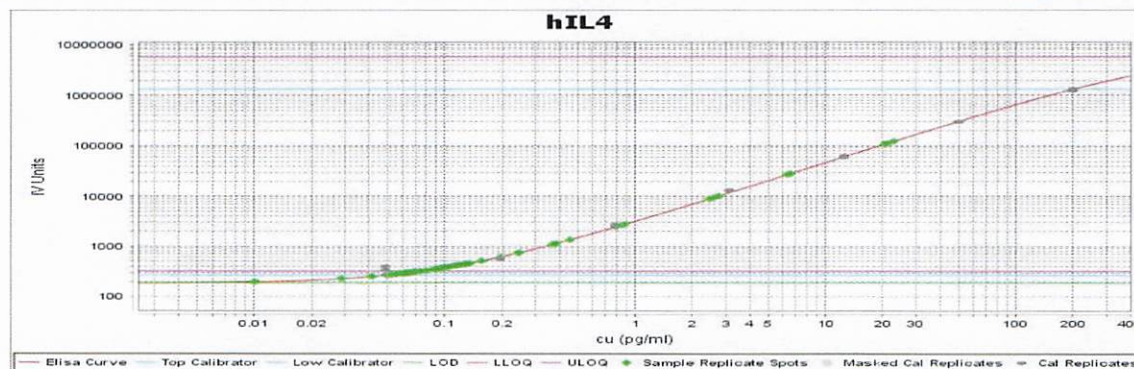
**IL-1β Calibrator Curve**



**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

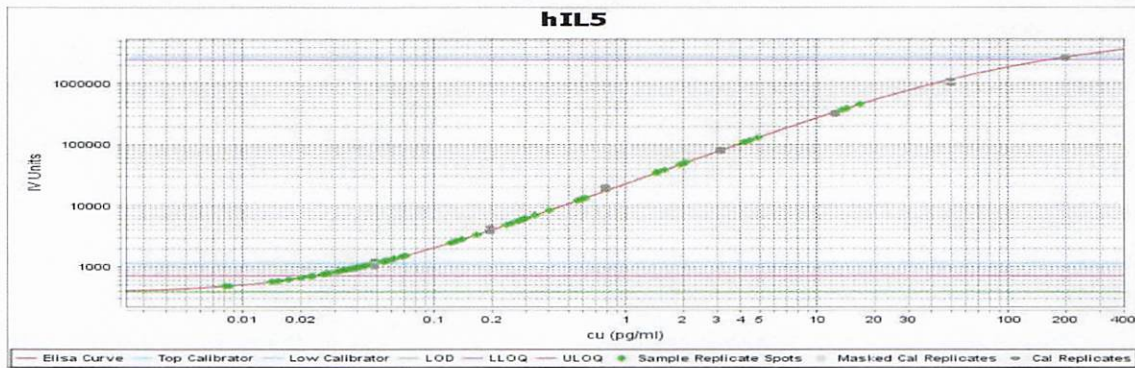
**IL-4 Calibrator Curve**



**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

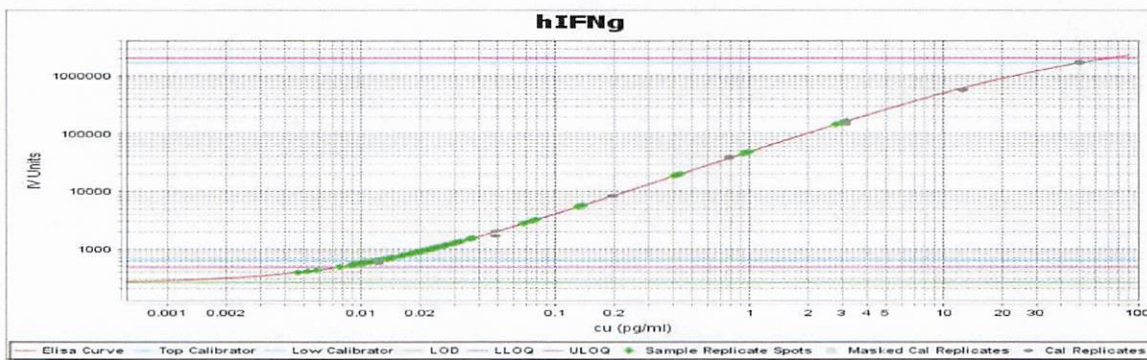
**IL-5 Calibrator Curve**



**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

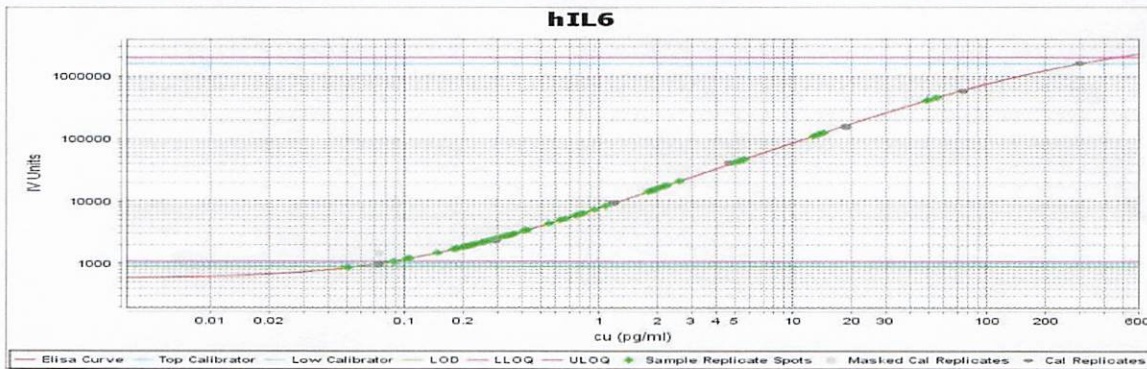
**IFNγ Calibrator Curve**



**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

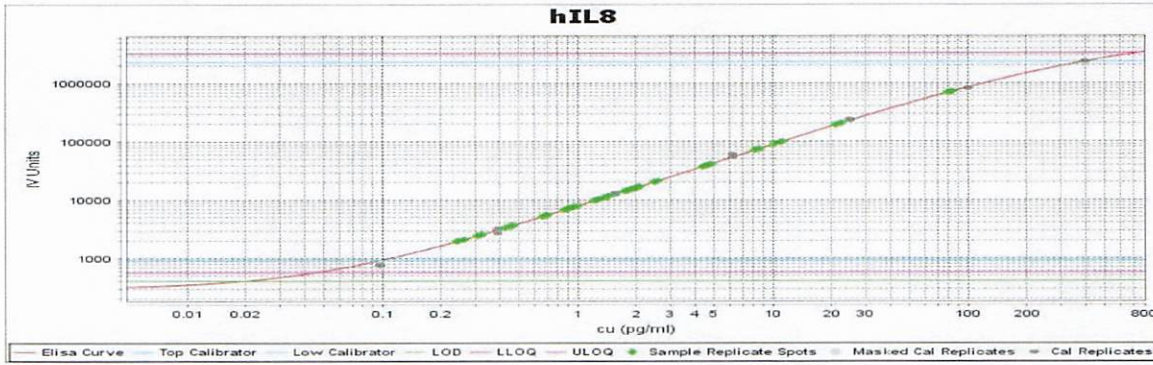
**IL-6 Calibrator Curve**



**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

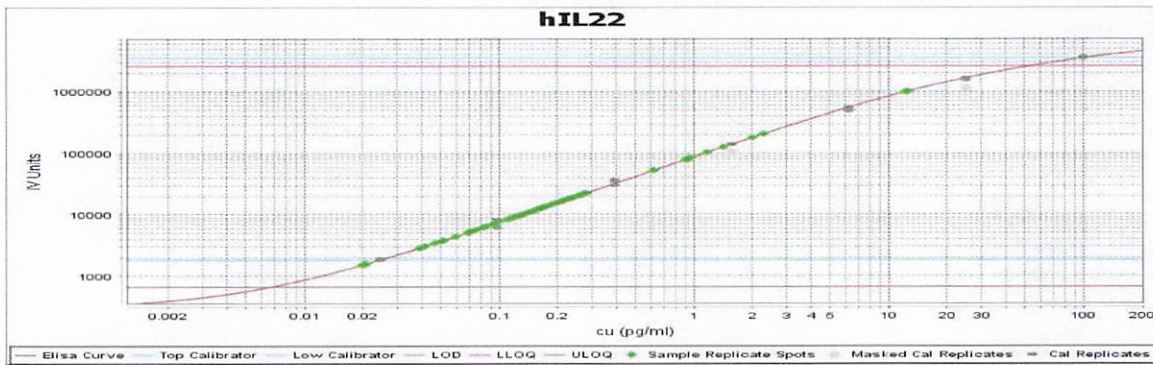
**IL-8 Calibrator Curve**



**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

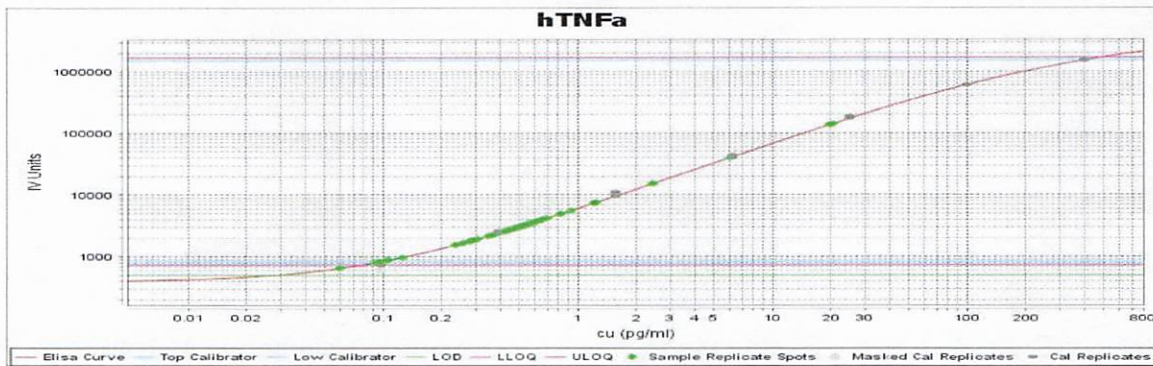
**IL-22 Calibrator Curve**



**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

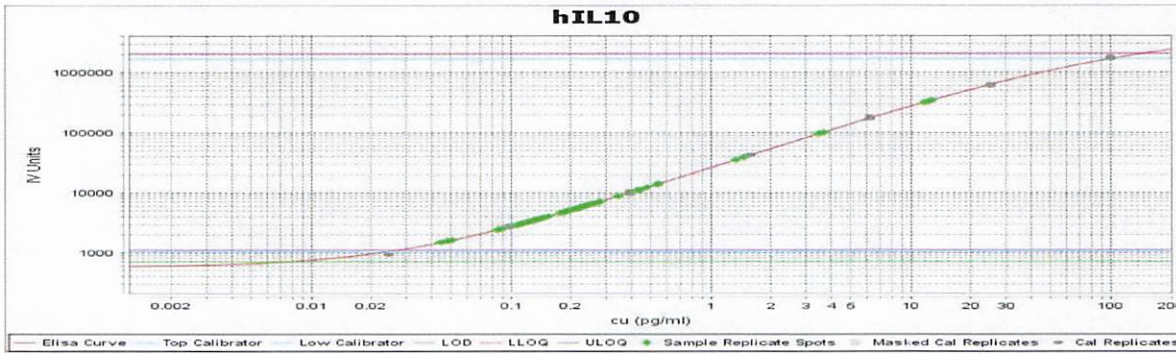
**TNFα Calibrator Curve**



**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)**

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