

<b>Product Number:</b>	85-0329
<b>Lot Number:</b>	504107
<b>Expiration:</b>	5-Mar-2025
<b>Platform(s):</b>	SP-X

Component	Part Number	Lot Number	Control Lot Number
CorPlex Cytokine Plate	97-0329	400507P	N/A
Biotinylated Antibody Reagent	98-0329	334511B	N/A
Sample Diluent 21	1863329	311103	311103A
IL-12p70 Calibrator	18910034	318756	N/A
IL-1β Calibrator	18910040	319165	N/A
IL-4 Calibrator	1858452	318580	N/A
IL-5 Calibrator	1858462	318837	N/A
IFNγ Calibrator	1858260	319843	N/A
IL-6 Calibrator	18910044	319484	N/A
IL-8 Calibrator	1858454	319436	N/A
IL-22 Calibrator	18910036	319176	N/A
TNFα Calibrator	1880012	319460	N/A
IL-10 Calibrator	18910043	318755	N/A
Wash Buffer (12.5X) <sup>1</sup>	1863332	N/A	N/A
Streptavidin-HRP Reagent <sup>1</sup>	1858263	N/A	N/A
Super Signal® Luminol <sup>1,2</sup>	1858089	N/A	N/A
Super Signal® Peroxide <sup>1,2</sup>	1858090	N/A	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

Robert Somogyi  
Name

Quality Control  
Manager  
Title

*[Handwritten Signature]*  
Signature/Date  
31 JAN 2024

**Release Materials\***

**IL-12p70**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	168	107 - 186
Panel 2	319839	47.3	33.0 - 51.2
Panel 3	319839	18.3	12.3 - 20.3
Panel 4	319839	4.57	2.59 - 6.31
Panel 5	319839	1.43	0.728 - 2.17

**IL-1β**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	13.1	9.55 - 20.6
Panel 2	319839	3.76	2.82 - 5.78
Panel 3	319839	1.47	1.07 - 2.22
Panel 4	319839	0.397	0.237 - 0.723
Panel 5	319839	0.146	0.081 - 0.271

**IL-4**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	87.1	50.1 - 97.2
Panel 2	319839	26.3	14.5 - 27.5
Panel 3	319839	10.1	5.89 - 10.8
Panel 4	319839	3.20	1.59 - 3.47
Panel 5	319839	1.09	0.553 - 1.34

**IL-5**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	70.1	31.5 - 70.5
Panel 2	319839	19.3	9.01 - 21.6
Panel 3	319839	7.61	3.34 - 8.64
Panel 4	319839	2.20	0.961 - 2.50
Panel 5	319839	0.771	0.370 - 0.968

**IFNγ**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	12.7	5.18 - 14.5
Panel 2	319839	3.83	1.55 - 4.66
Panel 3	319839	1.57	0.710 - 1.83
Panel 4	319839	0.497	0.217 - 0.604
Panel 5	319839	0.239	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**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	261	149 - 299
Panel 2	319839	63.0	39.7 - 70.6
Panel 3	319839	24.4	14.2 - 29.4
Panel 4	319839	8.49	4.15 - 11.2
Panel 5	319839	3.34	1.32 - 4.52

**IL-8**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	313	183 - 383
Panel 2	319839	89.2	49.8 - 103
Panel 3	319839	32.3	19.9 - 39.3
Panel 4	319839	10.1	5.82 - 12.6
Panel 5	319839	3.57	2.12 - 4.95

**IL-22**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	80.8	36.2 - 101
Panel 2	319839	10.3	6.05 - 15.0
Panel 3	319839	4.77	2.75 - 7.32
Panel 4	319839	0.901	0.565 - 1.49
Panel 5	319839	0.303	0.149 - 0.446

**TNFα**

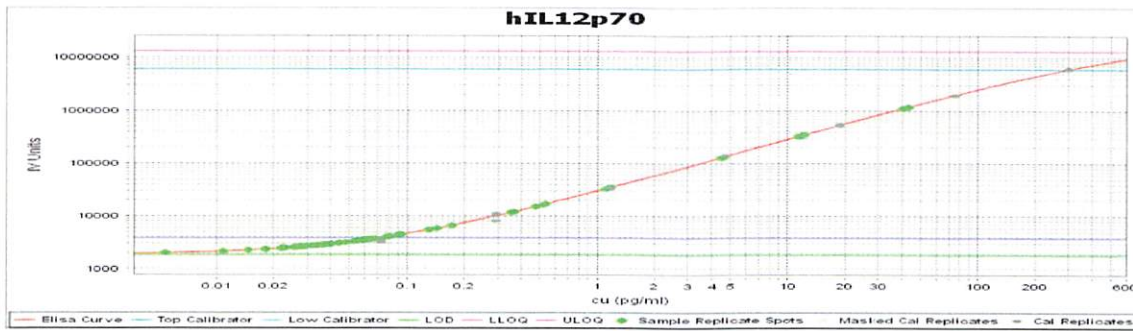
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	66.0	37.9 - 99.8
Panel 2	319839	19.2	11.8 - 26.5
Panel 3	319839	7.22	4.09 - 10.7
Panel 4	319839	1.75	1.01 - 2.84
Panel 5	319839	0.641	0.369 - 1.09

**IL-10**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	62.0	30.4 - 61.9
Panel 2	319839	16.5	9.02 - 17.8
Panel 3	319839	6.18	3.19 - 7.61
Panel 4	319839	1.79	1.01 - 2.34
Panel 5	319839	0.751	0.409 - 1.02

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

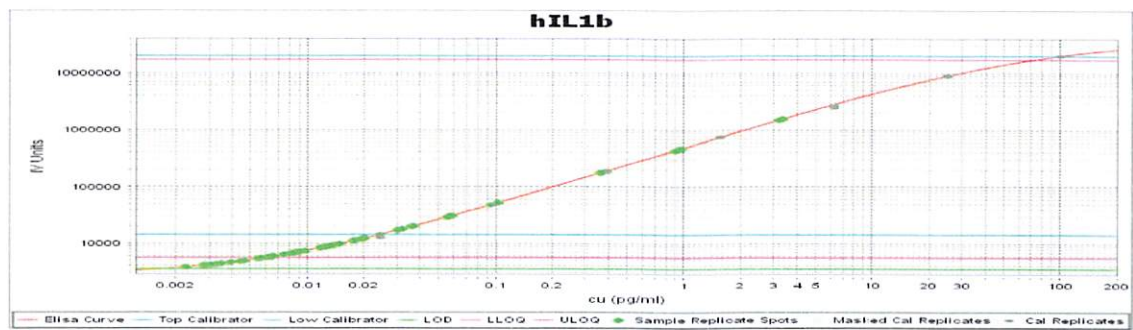
**IL-12p70 Calibrator Curve**



**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

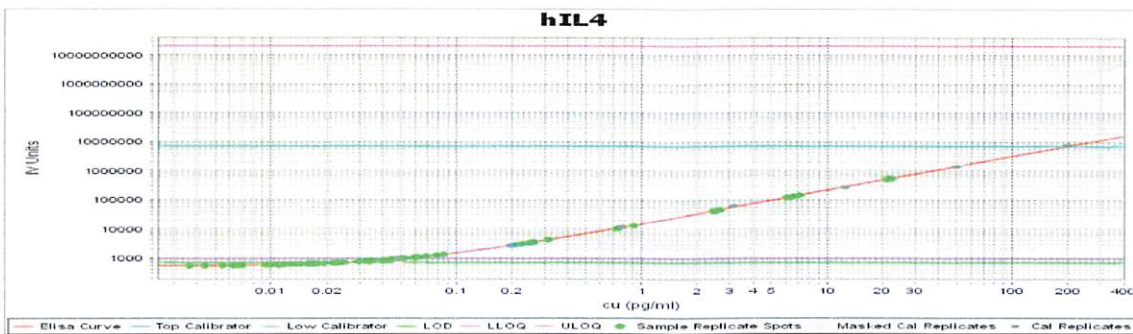
**IL-1β Calibrator Curve**



**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

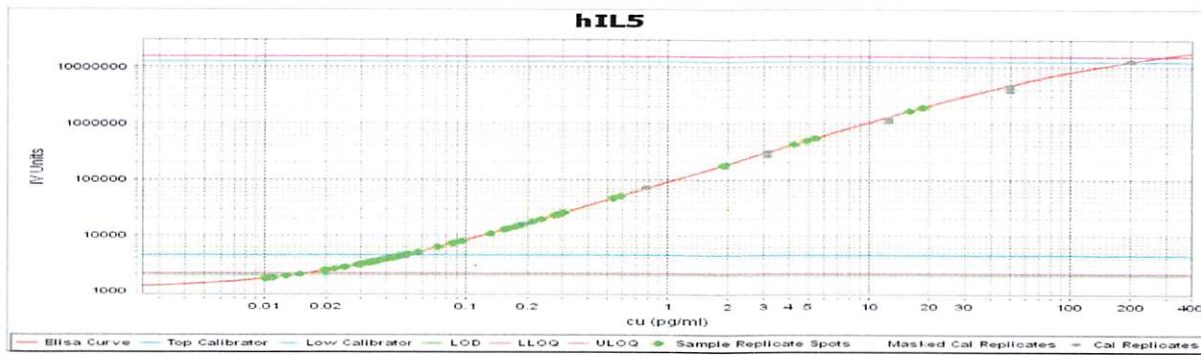
**IL-4 Calibrator Curve**



**Calibrator Levels (pg/mL)**

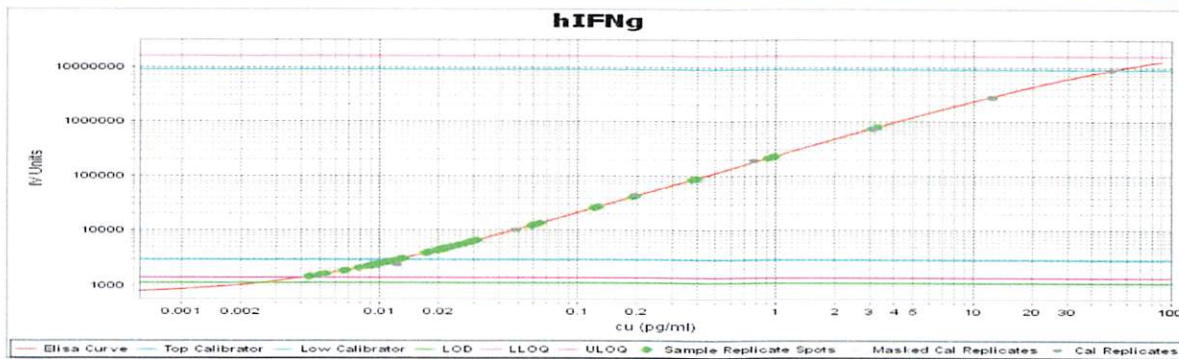
A	200
B	50.0
C	12.5
D	3.13
E	0.781
F	0.195
G	0.049
H	0.000

**IL-5 Calibrator Curve**



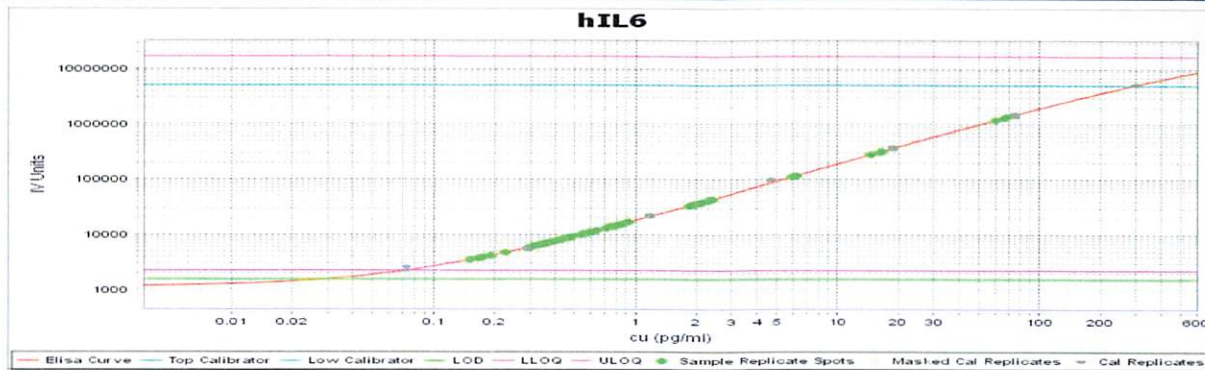
Calibrator Levels (pg/mL)	
A	200
B	50.0
C	12.5
D	3.13
E	0.781
F	0.195
G	0.049
H	0.000

**IFNγ Calibrator Curve**



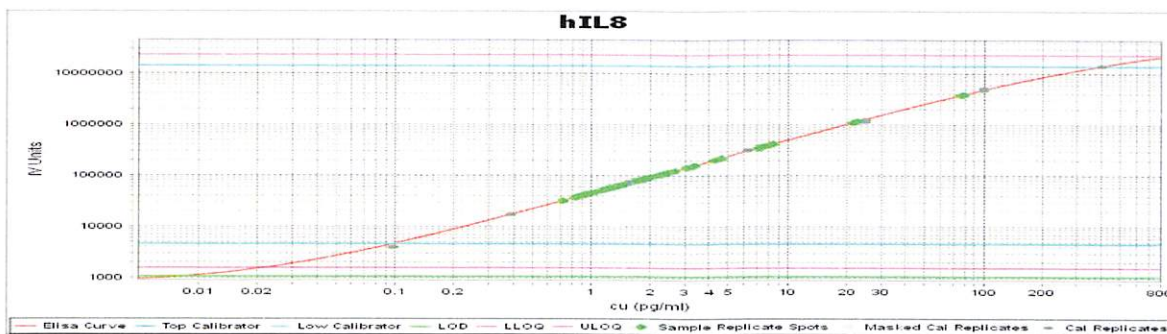
Calibrator Levels (pg/mL)	
A	50.0
B	12.5
C	3.13
D	0.781
E	0.195
F	0.049
G	0.012
H	0.000

**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

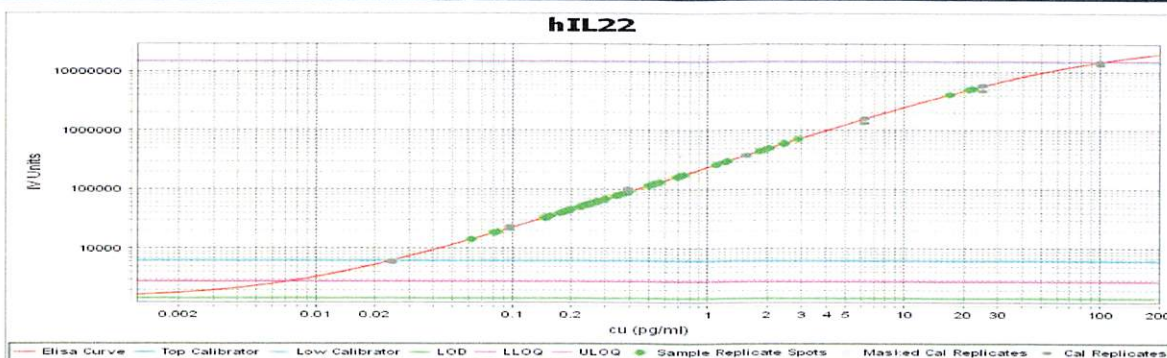
**IL-8 Calibrator Curve**



**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

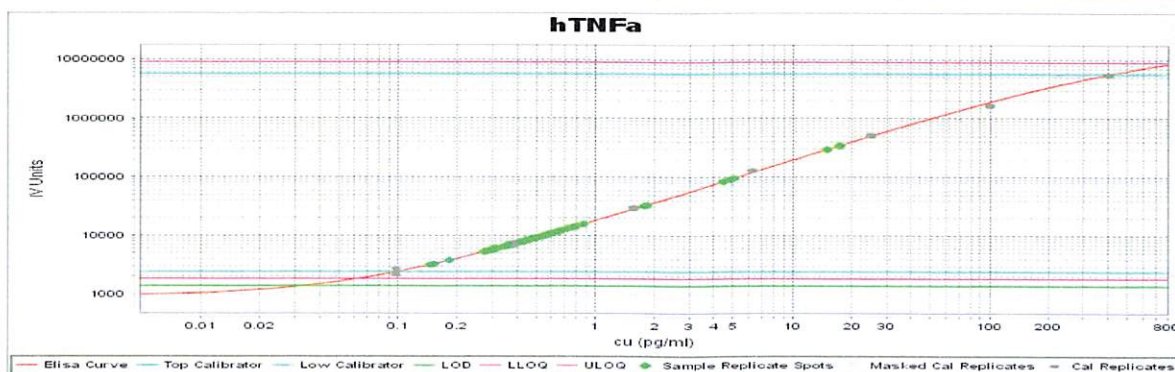
**IL-22 Calibrator Curve**



**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

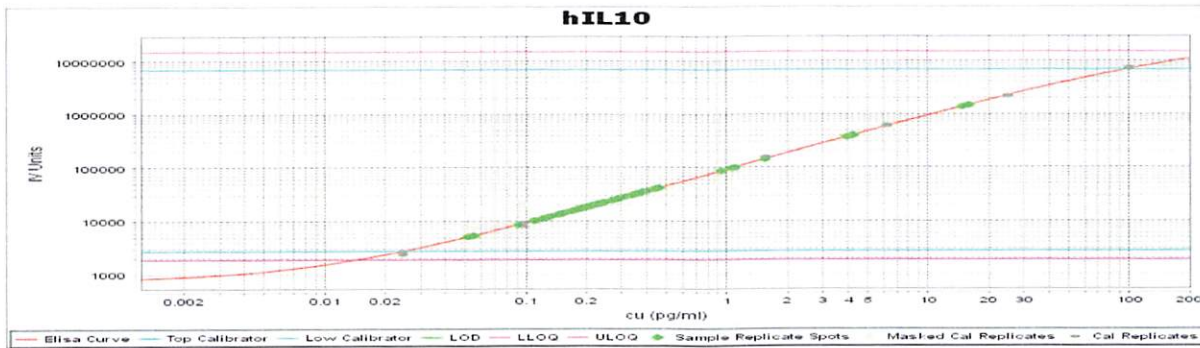
**TNFα Calibrator Curve**



**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)

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