

<b>Product Number:</b>	85-0329
<b>Lot Number:</b>	504015
<b>Expiration:</b>	26-Dec-2024
<b>Platform(s):</b>	SP-X



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

CorPlex Cytokine Plate	97-0329	325510P
Biotinylated Antibody Reagent	98-0329	310302B
Sample Diluent 21	1863329	319785
IL-12p70 Calibrator	18910034	318629
IL-1β Calibrator	18910040	319165
IL-4 Calibrator	1858452	318480
IL-5 Calibrator	1858462	318655
IFNγ Calibrator	1858260	318898
IL-6 Calibrator	18910044	319173
IL-8 Calibrator	1858454	319011
IL-22 Calibrator	18910036	318813
TNFα Calibrator	1880012	319247
IL-10 Calibrator	18910043	318755
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 the Complex Cytokine Panel Kit Instructions (KI-0195).

Review/Approval

Kathleen Cornell  
Name

Quality Control  
Supervisor  
Title

*[Handwritten Signature]* 20 Oct 2023  
Signature/Date



**Release Materials\***

**IL12p70**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	163.3	101.2 - 193.3
Panel 2	319839	46.05	32.24 - 52.86
Panel 3	319839	18.14	12.45 - 20.78
Panel 4	319839	4.669	2.632 - 6.594
Panel 5	319839	1.485	0.764 - 2.272

**IL-1β**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	13.68	10.37 - 21.04
Panel 2	319839	3.714	3.172 - 5.834
Panel 3	319839	1.407	1.201 - 2.331
Panel 4	319839	0.473	0.279 - 0.773
Panel 5	319839	0.134	0.096 - 0.301

**IL-4**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	77.62	45.70 - 99.67
Panel 2	319839	21.81	13.41 - 27.61
Panel 3	319839	9.267	5.612 - 11.10
Panel 4	319839	2.781	1.515 - 3.657
Panel 5	319839	0.738	0.559 - 1.400

**IL-5**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	61.45	27.65 - 80.96
Panel 2	319839	19.36	8.208 - 22.84
Panel 3	319839	8.969	2.848 - 9.877
Panel 4	319839	2.649	1.085 - 2.550
Panel 5	319839	0.919	0.320 - 1.170

**IFNγ**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	9.163	5.622 - 14.47
Panel 2	319839	2.743	1.692 - 4.770
Panel 3	319839	1.142	0.731 - 2.000
Panel 4	319839	0.364	0.215 - 0.692
Panel 5	319839	0.156	0.071 - 0.348

\* 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	223.0	141.9 - 312.1
Panel 2	319839	54.16	39.07 - 73.71
Panel 3	319839	21.60	14.61 - 30.61
Panel 4	319839	7.548	4.493 - 11.67
Panel 5	319839	2.754	1.407 - 4.765

**IL-8**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	416.4	164.0 - 399.7
Panel 2	319839	94.43	45.67 - 107.4
Panel 3	319839	36.35	18.96 - 41.09
Panel 4	319839	10.61	6.105 - 13.08
Panel 5	319839	4.081	2.260 - 5.125

**IL-22**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	88.44	36.22 - 101.9
Panel 2	319839	11.25	6.825 - 14.71
Panel 3	319839	5.525	3.227 - 7.790
Panel 4	319839	1.087	0.646 - 1.509
Panel 5	319839	0.280	0.180 - 0.453

**TNFα**

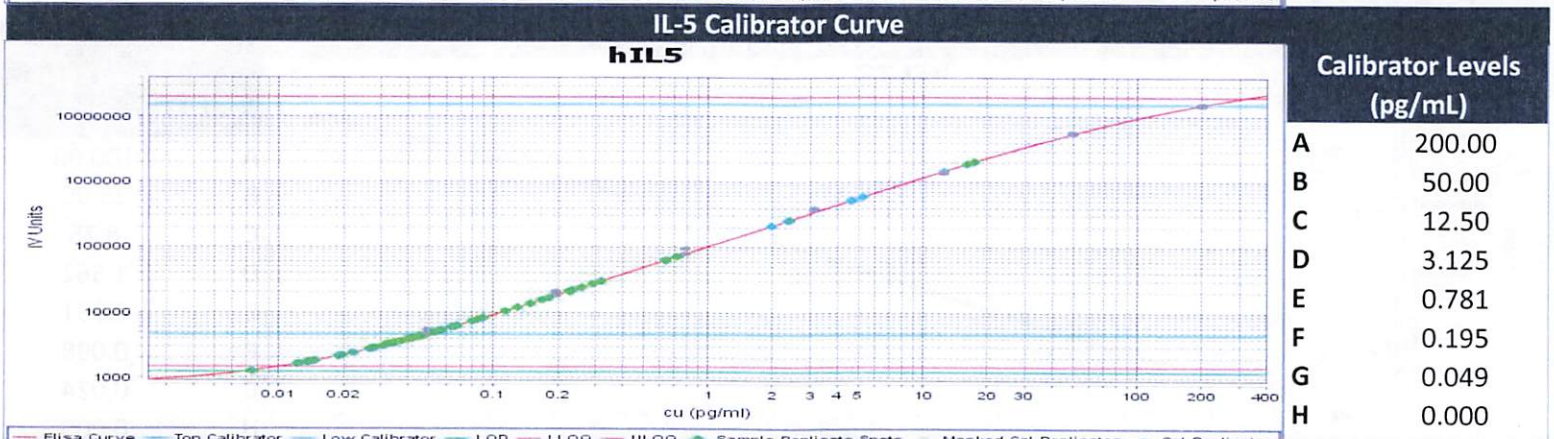
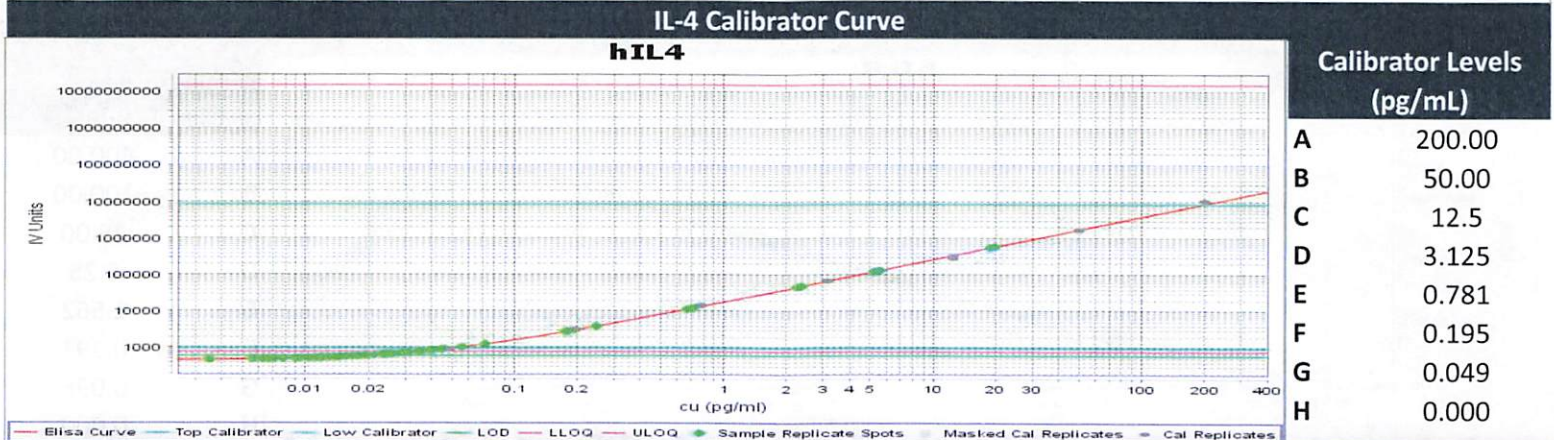
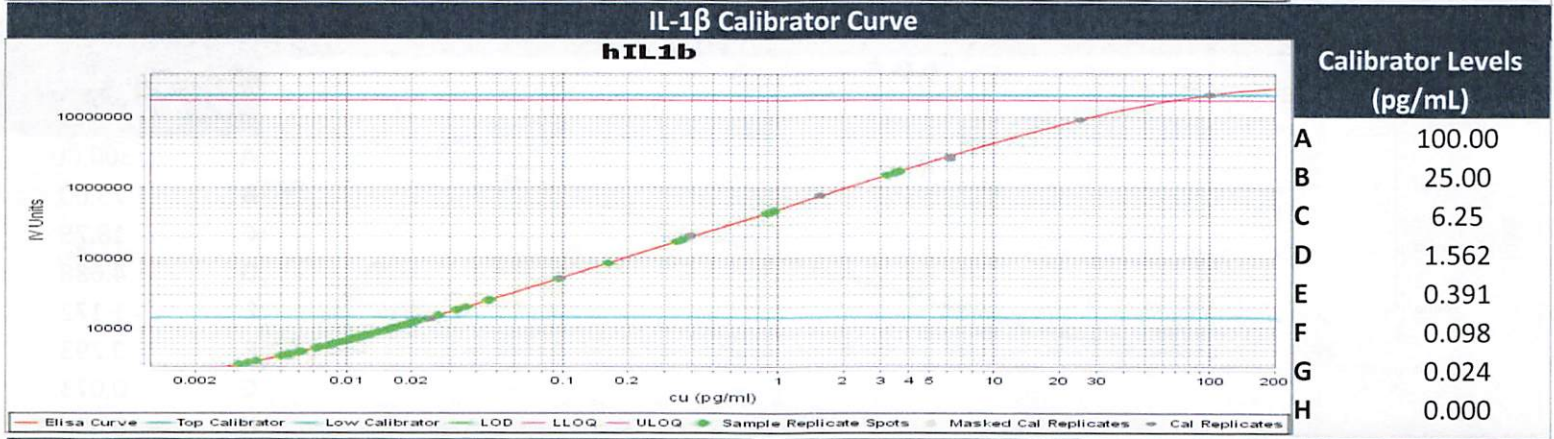
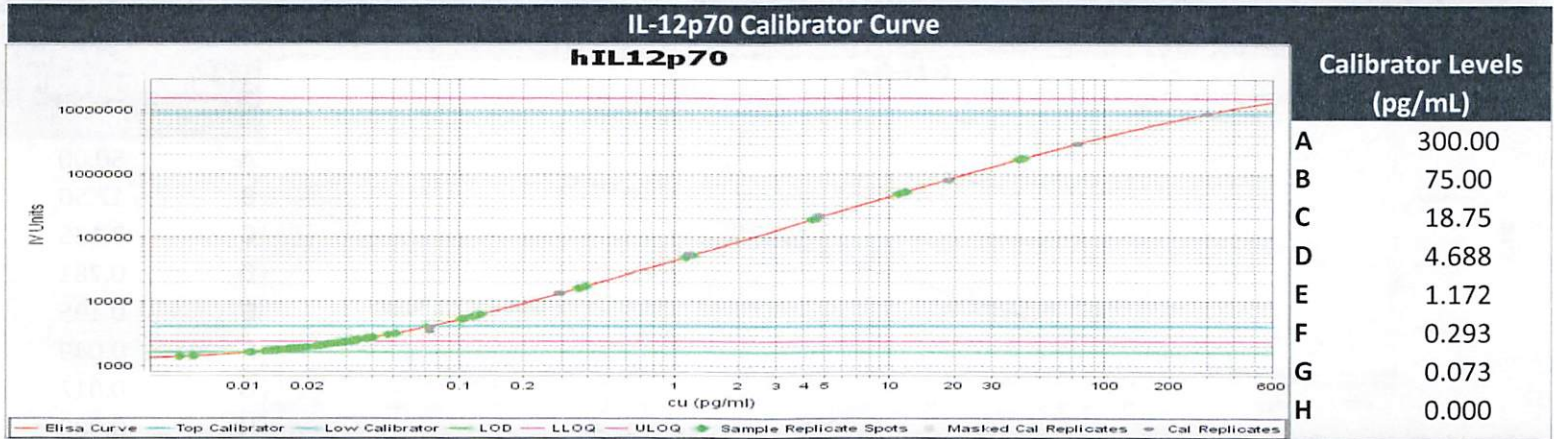
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	64.93	37.50 - 105.1
Panel 2	319839	17.94	12.23 - 27.53
Panel 3	319839	6.640	4.430 - 11.09
Panel 4	319839	1.475	1.213 - 2.913
Panel 5	319839	0.465	0.543 - 1.056

**IL-10**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	53.25	28.83 - 63.34
Panel 2	319839	13.91	9.000 - 18.28
Panel 3	319839	5.496	3.397 - 7.891
Panel 4	319839	1.644	1.141 - 2.386
Panel 5	319839	0.659	0.457 - 1.051

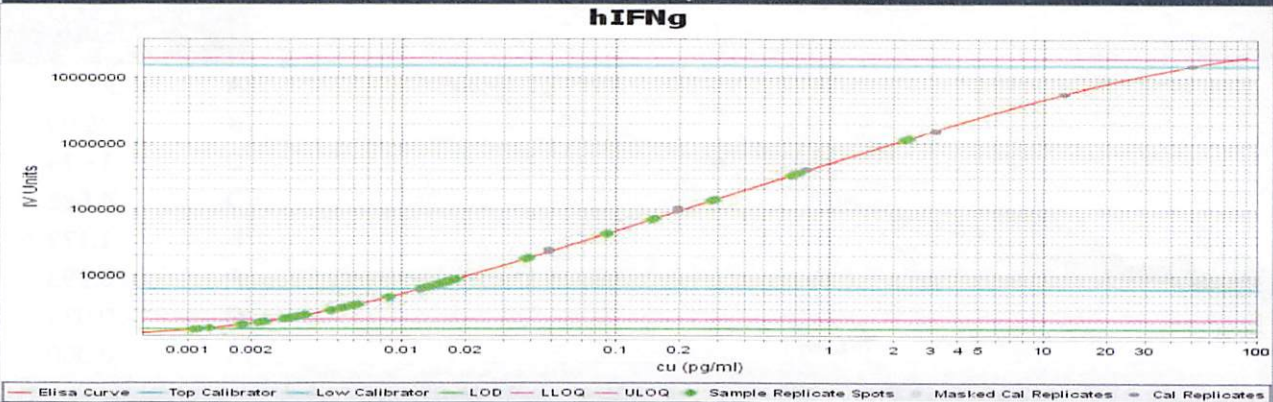
\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.





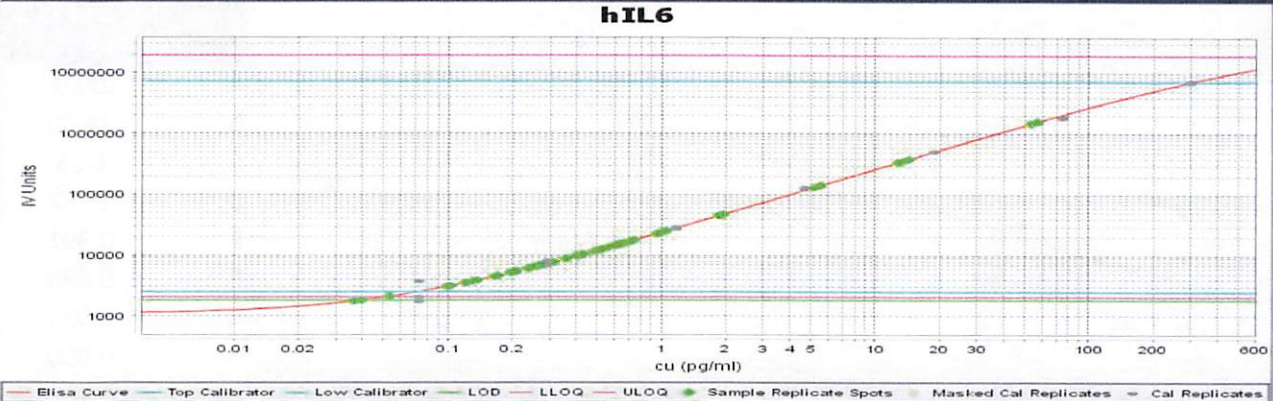


**IFN $\gamma$  Calibrator Curve**



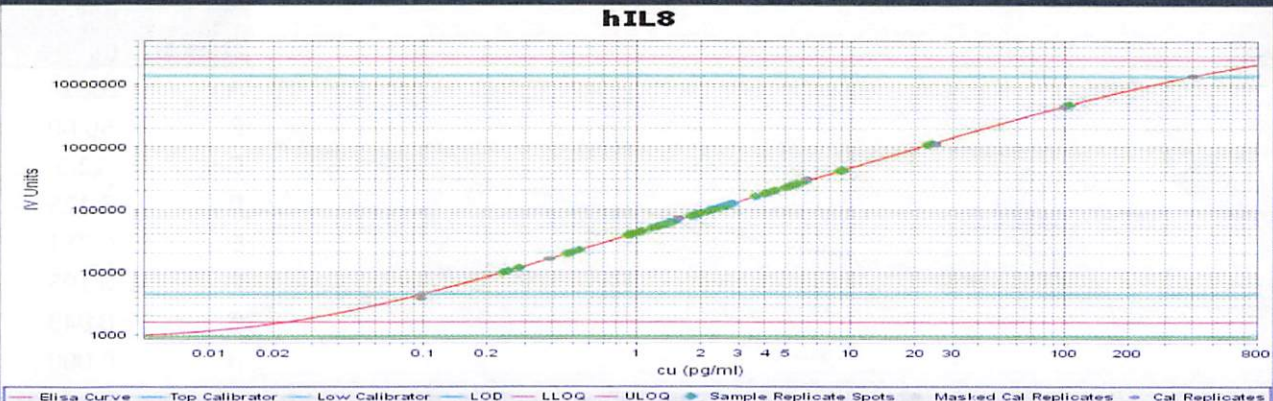
Calibrator Levels (pg/mL)	
A	50.00
B	12.50
C	3.125
D	0.781
E	0.195
F	0.049
G	0.012
H	0.000

**IL-6 Calibrator Curve**



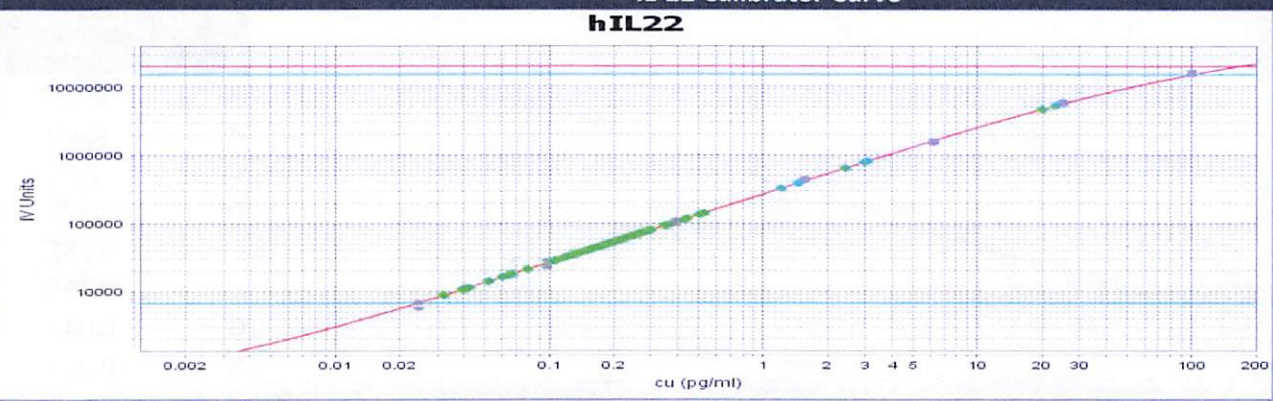
Calibrator Levels (pg/mL)	
A	300.00
B	75.00
C	18.75
D	4.688
E	1.172
F	0.293
G	0.073
H	0.000

**IL-8 Calibrator Curve**



Calibrator Levels (pg/mL)	
A	400.00
B	100.00
C	25.00
D	6.25
E	1.562
F	0.391
G	0.098
H	0.000

**IL-22 Calibrator Curve**

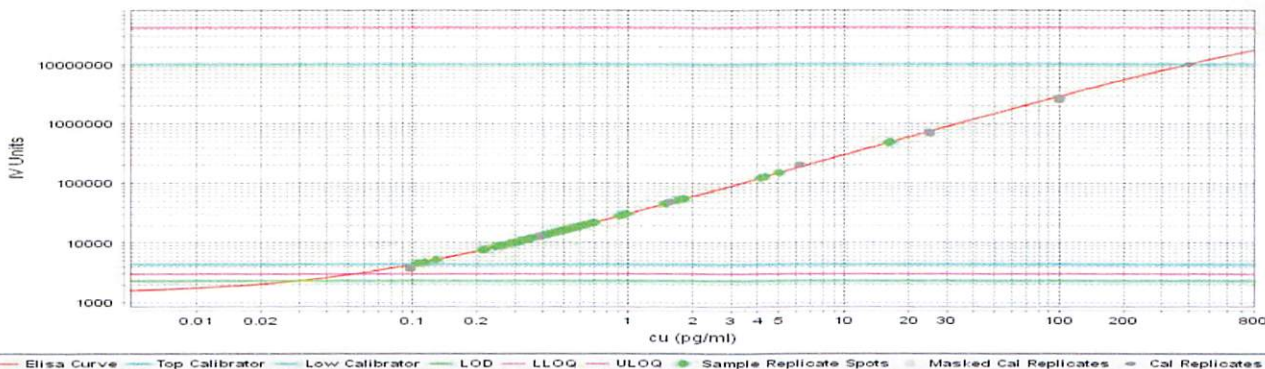


Calibrator Levels (pg/mL)	
A	100.00
B	25.00
C	6.25
D	1.562
E	0.391
F	0.098
G	0.024
H	0.000



**TNFα Calibrator Curve**

**hTNFa**

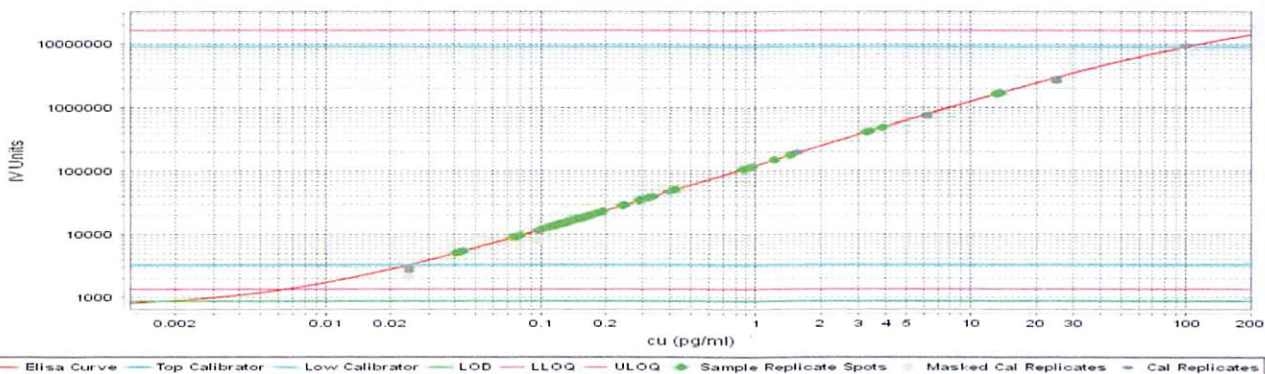


**Calibrator Levels  
(pg/mL)**

A	400.00
B	100.00
C	25.00
D	6.25
E	1.562
F	0.391
G	0.098
H	0.000

**IL-10 Calibrator Curve**

**hIL10**



**Calibrator Levels  
(pg/mL)**

A	100.00
B	25.00
C	6.25
D	1.562
E	0.391
F	0.098
G	0.024
H	0.000