

Product Number:	85-0329
Lot Number:	504068
Expiration:	13-Jan-2025
Platform(s):	SP-X

CorPlex Cytokine Plate	97-0329	325510P
Biotinylated Antibody Reagent	98-0329	332509B
Sample Diluent 21	1863329	319974
IL-12p70 Calibrator	18910034	318756
IL-1β Calibrator	18910040	319165
IL-4 Calibrator	1858452	318480
IL-5 Calibrator	1858462	318655
IFNγ Calibrator	1858260	319437
IL-6 Calibrator	18910044	319173
IL-8 Calibrator	1858454	319436
IL-22 Calibrator	18910036	318813
TNFα Calibrator	1880012	319247
IL-10 Calibrator	18910043	318755
Wash Buffer (12.5X) ¹	1863332	N/A
Streptavidin-HRP Reagent ¹	1858263	N/A
Super Signal® Luminol ^{1,2}	1858089	N/A
Super Signal® Peroxide ^{1,2}	1858090	N/A

¹ Reagents are not Kit Lot Specific

² 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).



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

Review/Approval

Kathleen Cornell

Name

Quality Control
Supervisor

Title

[Signature] 30 Nov 2023

Signature/Date

Release Materials*

IL-12p70

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	152.1	107.0 - 186.2
Panel 2	319839	47.51	32.96 - 51.22
Panel 3	319839	18.30	12.33 - 20.31
Panel 4	319839	4.875	2.590 - 6.308
Panel 5	319839	1.446	0.728 - 2.167

IL-1β

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	11.97	9.553 - 20.59
Panel 2	319839	3.858	2.823 - 5.776
Panel 3	319839	1.490	1.072 - 2.216
Panel 4	319839	0.417	0.237 - 0.723
Panel 5	319839	0.137	0.081 - 0.271

IL-4

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	68.88	50.08 - 97.19
Panel 2	319839	22.54	14.54 - 27.45
Panel 3	319839	8.969	5.892 - 10.84
Panel 4	319839	2.720	1.590 - 3.472
Panel 5	319839	0.818	0.553 - 1.337

IL-5

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	59.47	31.53 - 70.53
Panel 2	319839	18.67	9.014 - 21.60
Panel 3	319839	7.362	3.344 - 8.637
Panel 4	319839	1.738	0.961 - 2.498
Panel 5	319839	0.678	0.370 - 0.968

IFNγ

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	7.320	5.184 - 14.50
Panel 2	319839	2.460	1.553 - 4.656
Panel 3	319839	1.011	0.710 - 1.834
Panel 4	319839	0.314	0.217 - 0.604
Panel 5	319839	0.125	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	201.5	149.3 - 299.3
Panel 2	319839	54.66	39.74 - 70.62
Panel 3	319839	22.57	14.23 - 29.41
Panel 4	319839	7.359	4.151 - 11.20
Panel 5	319839	2.808	1.321 - 4.519

IL-8

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	293.8	182.6 - 382.7
Panel 2	319839	81.98	49.82 - 102.8
Panel 3	319839	30.56	19.87 - 39.34
Panel 4	319839	9.506	5.821 - 12.63
Panel 5	319839	3.412	2.122 - 4.948

IL-22

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	94.39	36.20 - 100.6
Panel 2	319839	13.33	6.046 - 14.95
Panel 3	319839	6.269	2.746 - 7.320
Panel 4	319839	1.140	0.565 - 1.486
Panel 5	319839	0.308	0.149 - 0.446

TNFα

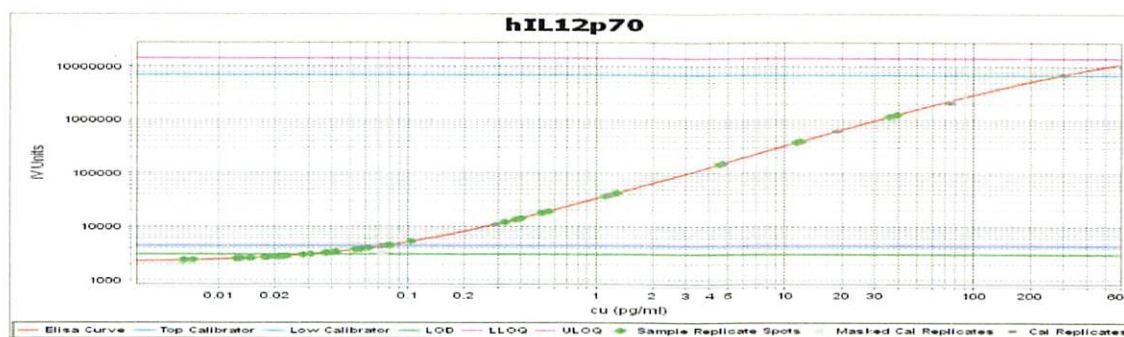
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	69.11	37.89 - 99.76
Panel 2	319839	21.92	11.84 - 26.54
Panel 3	319839	7.767	4.095 - 10.69
Panel 4	319839	1.545	1.007 - 2.838
Panel 5	319839	0.473	0.369 - 1.090

IL-10

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	49.54	30.41 - 61.92
Panel 2	319839	15.08	9.024 - 17.75
Panel 3	319839	6.048	3.193 - 7.609
Panel 4	319839	1.930	1.009 - 2.343
Panel 5	319839	0.765	0.409 - 1.017

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

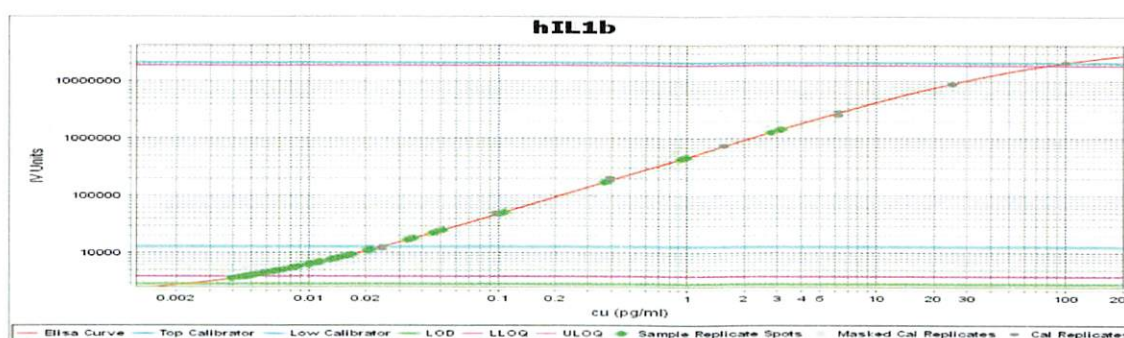
IL-12p70 Calibrator Curve



Calibrator Levels (pg/mL)

A	300.00
B	75.00
C	18.75
D	4.688
E	1.172
F	0.293
G	0.073
H	0.000

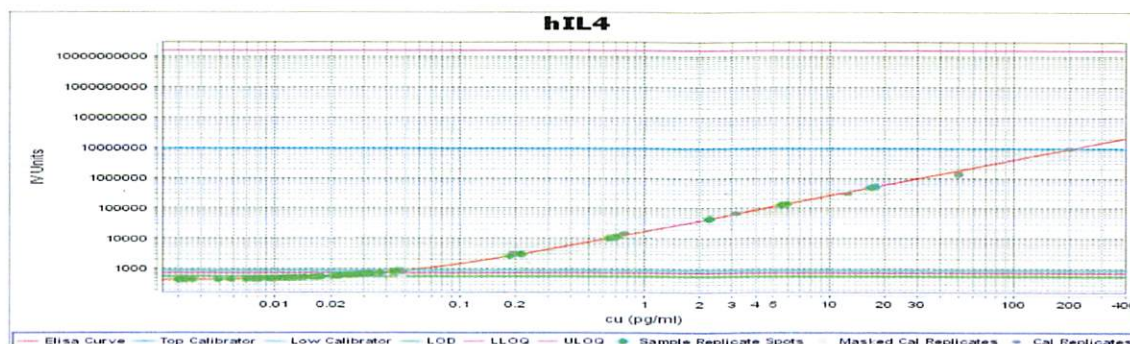
IL-1 β Calibrator Curve



Calibrator Levels (pg/mL)

A	100.00
B	25.00
C	6.250
D	1.562
E	0.391
F	0.098
G	0.024
H	0.000

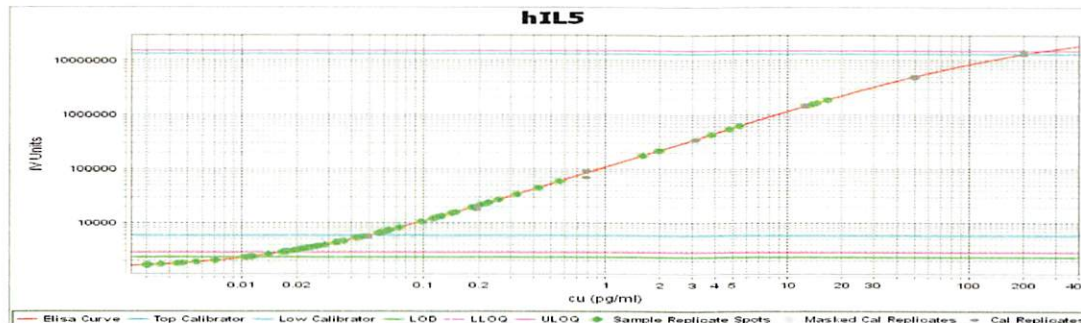
IL-4 Calibrator Curve



Calibrator Levels (pg/mL)

A	200.00
B	50.00
C	12.50
D	3.125
E	0.781
F	0.195
G	0.049
H	0.000

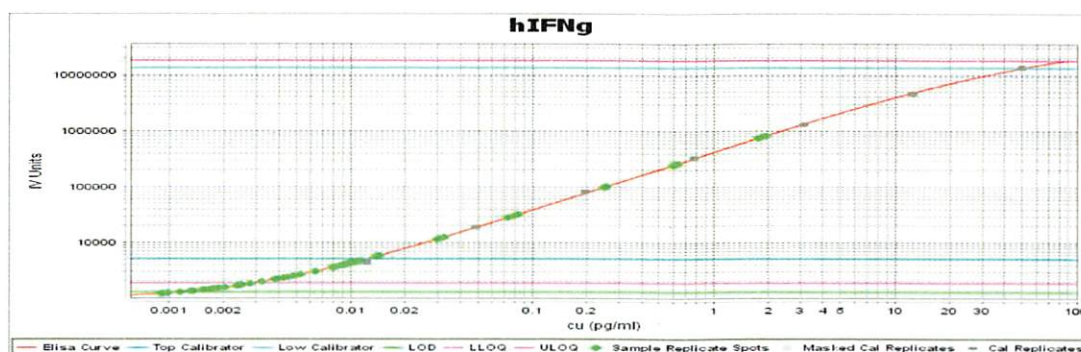
IL-5 Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	200.00
B	50.00
C	12.50
D	3.125
E	0.781
F	0.195
G	0.049
H	0.000

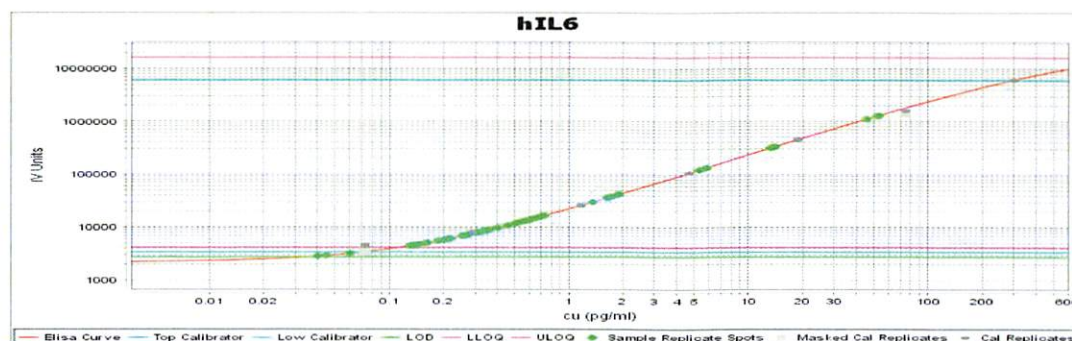
IFNγ Calibrator Curve



**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

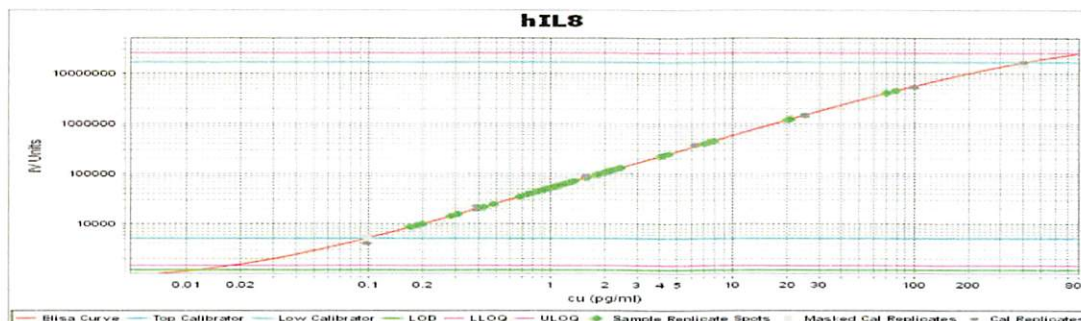
IL-6 Calibrator Curve



**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

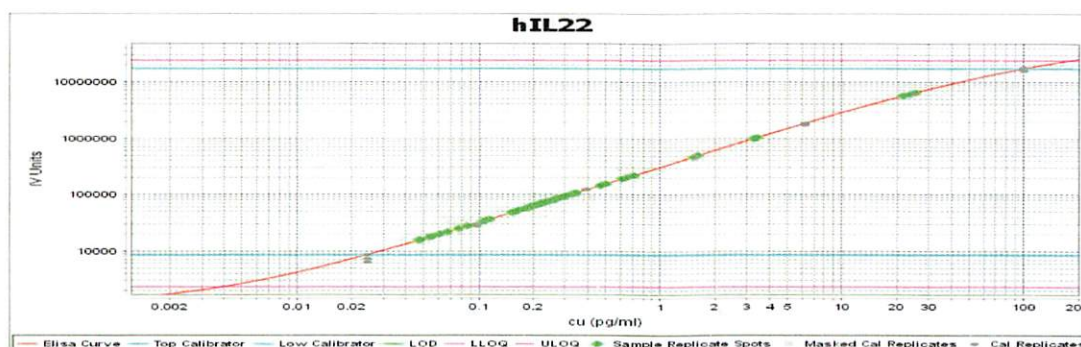
IL-8 Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	400.00
B	100.00
C	25.00
D	6.250
E	1.562
F	0.391
G	0.098
H	0.000

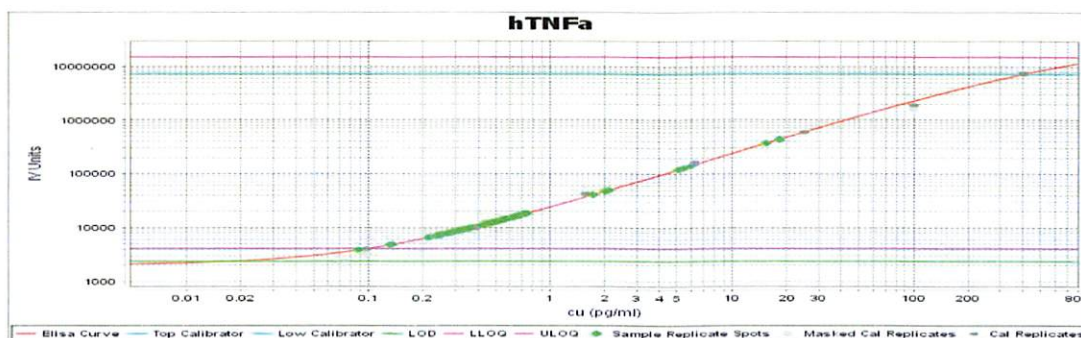
IL-22 Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	100.00
B	25.00
C	6.250
D	1.562
E	0.391
F	0.098
G	0.024
H	0.000

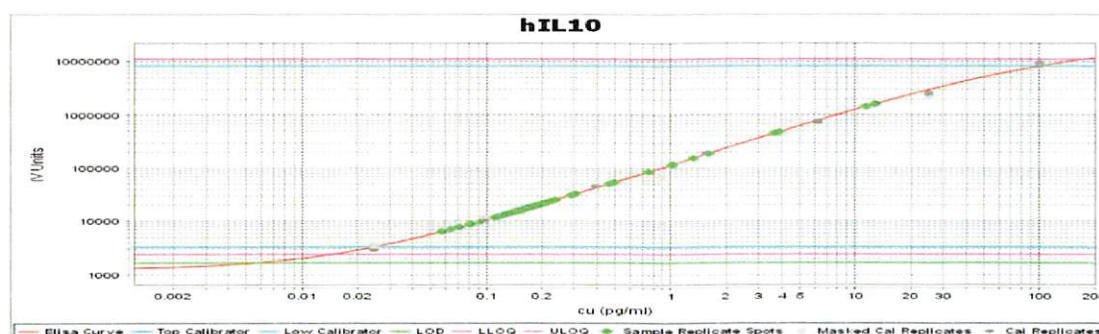
TNFα Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	400.00
B	100.00
C	25.00
D	6.250
E	1.562
F	0.391
G	0.098
H	0.000

IL-10 Calibrator Curve



Calibrator Levels
(pg/mL)

A	100.00
B	25.00
C	6.250
D	1.562
E	0.391
F	0.098
G	0.024
H	0.000