

Product Number:	85-0329
Lot Number:	503984
Expiration:	13-Jan-2025
Platform(s):	SP-X

CorPlex Cytokine Plate	97-0329	326103P
Biotinylated Antibody Reagent	98-0329	326105B
Sample Diluent 21	1863329	319974
IL-12p70 Calibrator	18910034	319844
IL-1β Calibrator	18910040	319165
IL-4 Calibrator	1858452	319016
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	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



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

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

Review/Approval

Kathleen Cornell

Quality Control
Supervisor

[Handwritten Signature] 20 Nov 2023

Name

Title

Signature/Date

Release Materials*

IL-12p70

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	147.3	107.0 - 186.2
Panel 2	319839	41.26	32.96 - 51.22
Panel 3	319839	16.55	12.33 - 20.31
Panel 4	319839	4.596	2.590 - 6.308
Panel 5	319839	1.377	0.728 - 2.167

IL-1β

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	12.53	9.553 - 20.59
Panel 2	319839	3.405	2.823 - 5.776
Panel 3	319839	1.374	1.072 - 2.216
Panel 4	319839	0.358	0.237 - 0.723
Panel 5	319839	0.125	0.081 - 0.271

IL-4

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	80.44	50.08 - 97.19
Panel 2	319839	23.75	14.54 - 27.45
Panel 3	319839	9.764	5.892 - 10.84
Panel 4	319839	2.655	1.590 - 3.472
Panel 5	319839	0.852	0.553 - 1.337

IL-5

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	50.02	31.53 - 70.53
Panel 2	319839	14.49	9.014 - 21.60
Panel 3	319839	6.277	3.344 - 8.637
Panel 4	319839	1.574	0.961 - 2.498
Panel 5	319839	0.593	0.370 - 0.968

IFNγ

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	8.313	5.184 - 14.50
Panel 2	319839	2.487	1.553 - 4.656
Panel 3	319839	1.051	0.710 - 1.834
Panel 4	319839	0.309	0.217 - 0.604
Panel 5	319839	0.132	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	218.6	149.3 - 299.3
Panel 2	319839	54.08	39.74 - 70.62
Panel 3	319839	22.11	14.23 - 29.41
Panel 4	319839	7.213	4.151 - 29.41
Panel 5	319839	2.875	1.321 - 4.519

IL-8

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	277.7	182.6 - 382.7
Panel 2	319839	71.43	49.82 - 102.8
Panel 3	319839	27.62	19.87 - 39.34
Panel 4	319839	8.296	5.821 - 12.63
Panel 5	319839	3.022	2.122 - 4.948

IL-22

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	76.98	36.20 - 100.6
Panel 2	319839	11.11	6.046 - 14.95
Panel 3	319839	5.400	2.746 - 7.320
Panel 4	319839	0.993	0.565 - 1.486
Panel 5	319839	0.297	0.149 - 0.446

TNFα

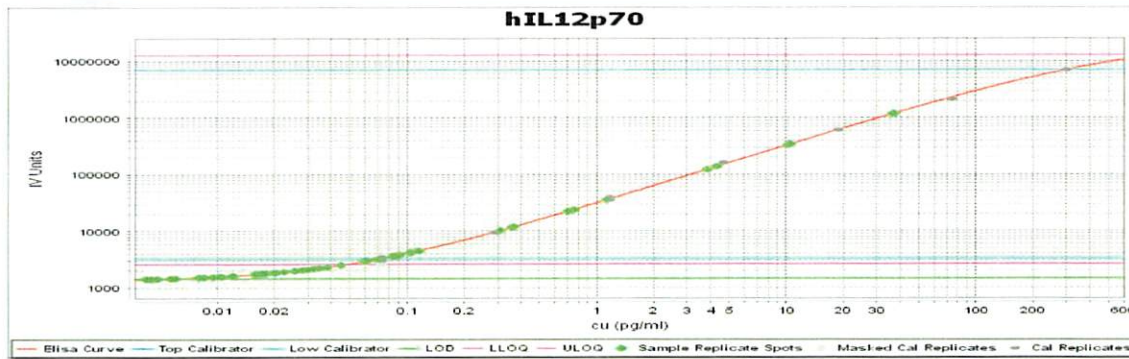
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	69.69	37.89 - 99.76
Panel 2	319839	18.52	11.84 - 26.54
Panel 3	319839	7.663	4.095 - 10.69
Panel 4	319839	1.679	1.007 - 2.838
Panel 5	319839	0.539	0.369 - 1.090

IL-10

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	53.85	30.41 - 61.92
Panel 2	319839	14.27	9.024 - 17.75
Panel 3	319839	5.688	3.193 - 7.609
Panel 4	319839	1.690	1.009 - 2.343
Panel 5	319839	0.655	0.409 - 1.017

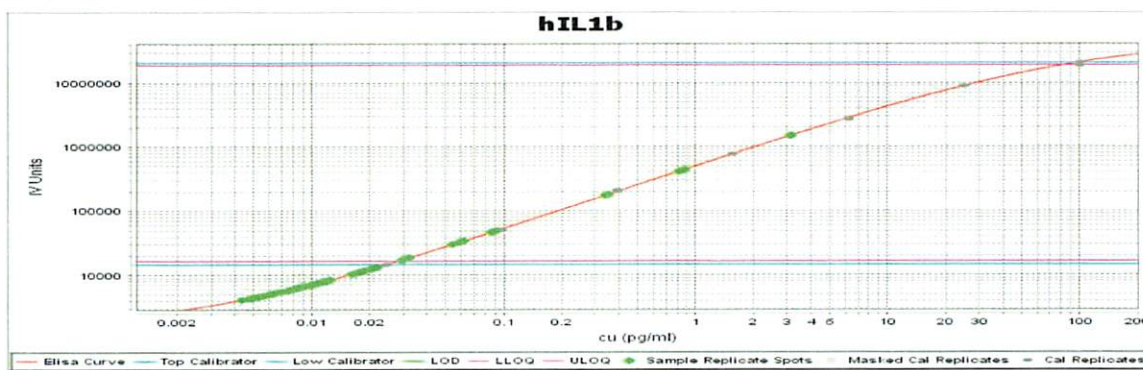
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

IL-12p70 Calibrator Curve



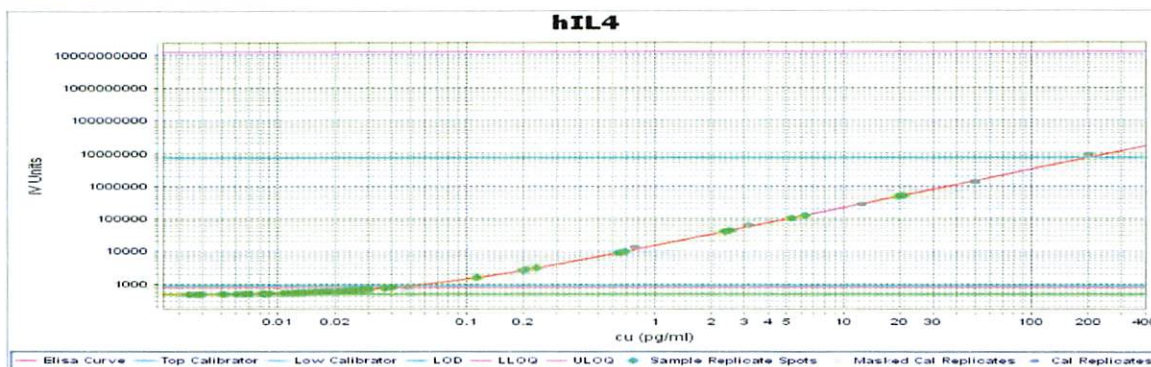
Calibrator Levels (pg/mL)	
A	300.00
B	75.00
C	18.75
D	4.688
E	1.172
F	0.293
G	0.073
H	0.000

IL-1β Calibrator Curve



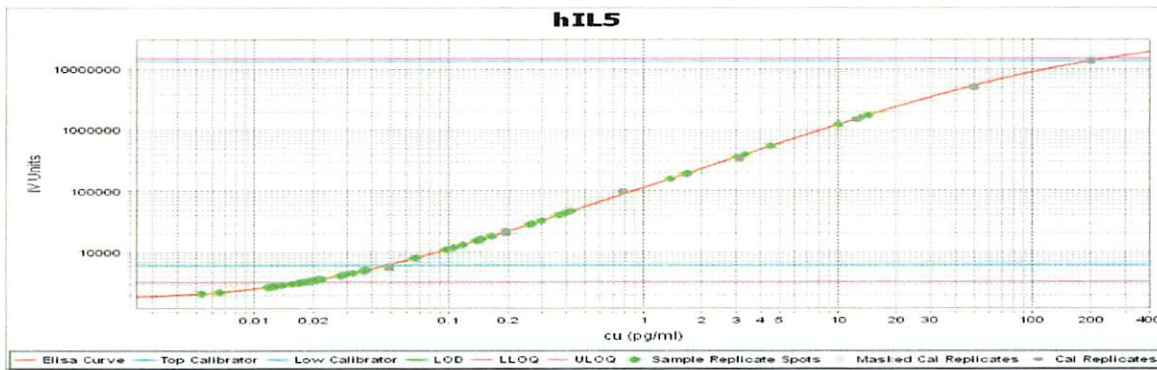
Calibrator Levels (pg/mL)	
A	100.00
B	25.00
C	6.25
D	1.562
E	0.391
F	0.098
G	0.024
H	0.000

IL-4 Calibrator Curve



Calibrator Levels (pg/mL)	
A	200.00
B	50.00
C	12.5
D	3.125
E	0.781
F	0.195
G	0.049
H	0.000

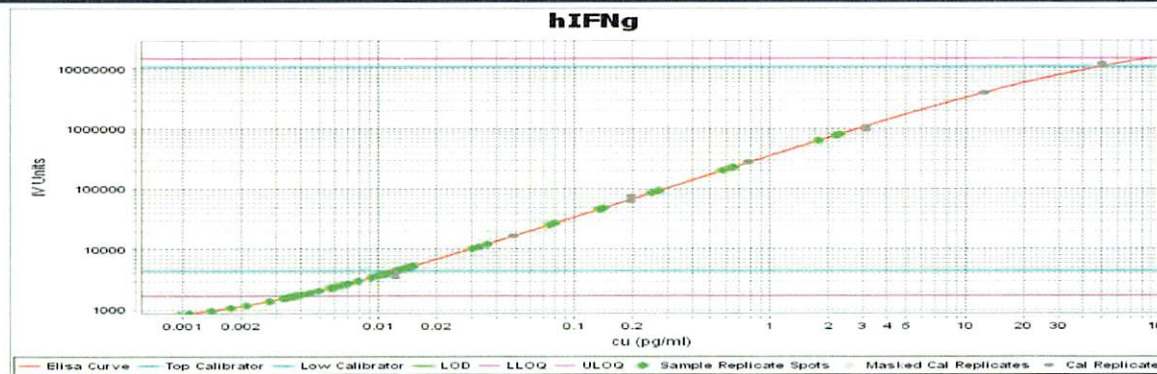
IL-5 Calibrator Curve



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

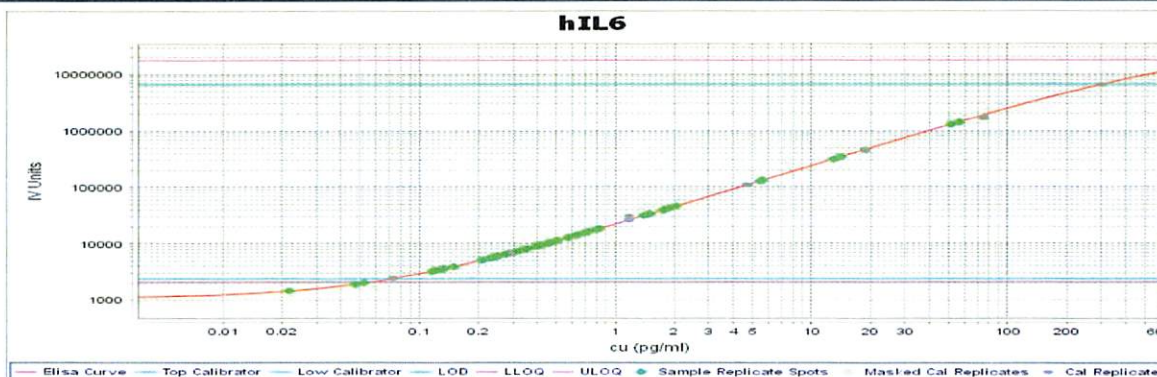
IFN γ Calibrator Curve



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

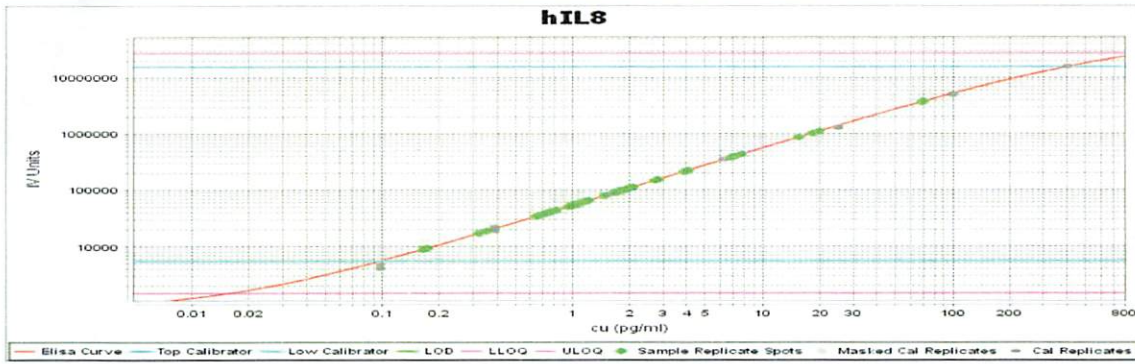
IL-6 Calibrator Curve



Calibrator Levels (pg/mL)

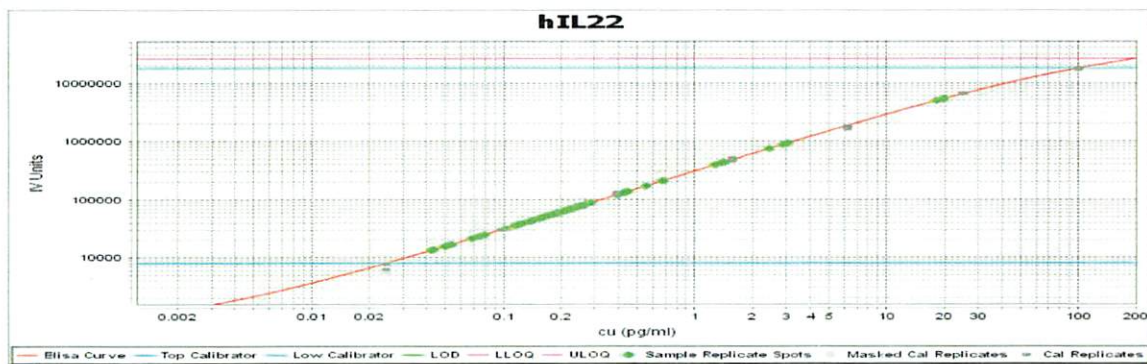
A	300.00
B	75.00
C	18.75
D	4.688
E	1.172
F	0.293
G	0.073
H	0.000

IL-8 Calibrator Curve



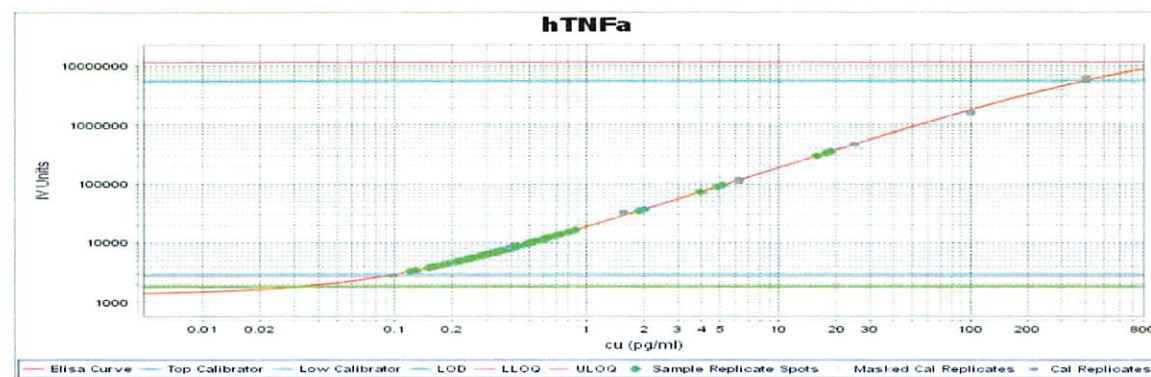
Calibrator Levels (pg/mL)	
A	400.00
B	100.00
C	25.00
D	6.25
E	1.562
F	0.391
G	0.098
H	0.000

IL-22 Calibrator Curve



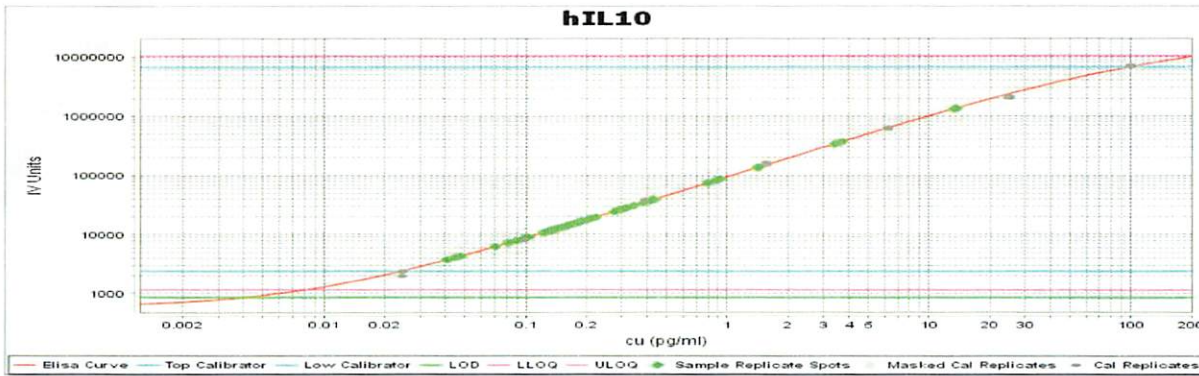
Calibrator Levels (pg/mL)	
A	100.00
B	25.00
C	6.25
D	1.562
E	0.391
F	0.098
G	0.024
H	0.000

TNFα Calibrator Curve



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



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