

Product Number:	85-0329
Lot Number:	504568
Expiration:	26-Mar-2026
Platform(s):	SP-X

Component	Part Number	Lot Number	Control Lot Number
CorPlex Cytokine Plate	97-0329	508406P	N/A
Biotinylated Antibody Reagent	98-0329	502303B	502303AB
Sample Diluent 21	1863329	405709	N/A
IL-12p70 Calibrator	18910034	319844	N/A
IL-1 β Calibrator	18910040	407211	N/A
IL-4 Calibrator	1858452	319438	N/A
IL-5 Calibrator	1858462	319838	N/A
IFN γ Calibrator	1858260	319843	N/A
IL-6 Calibrator	18910044	324303	N/A
IL-8 Calibrator	1858454	319436	N/A
IL-22 Calibrator	18910036	319459	N/A
TNF α Calibrator	1880012	403812	N/A
IL-10 Calibrator	18910043	319842	N/A
Wash Buffer (12.5X) ¹	1863332	N/A	N/A
Streptavidin-HRP Reagent ¹	1858263	N/A	N/A
Super Signal® Luminol ^{1,2}	1858089	N/A	N/A
Super Signal® Peroxide ^{1,2}	1858090	N/A	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 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 16 Apr 2025
Signature/Date

Release Materials*

IL-12p70

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	138	108 - 185
Panel 2	319839	39.7	33.5 - 51.7
Panel 3	319839	15.7	12.4 - 21.1
Panel 4	319839	3.86	2.73 - 6.45
Panel 5	319839	1.25	0.781 - 2.23

IL-1β

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	13.9	9.83 - 20.2
Panel 2	319839	3.69	2.87 - 5.85
Panel 3	319839	1.37	1.07 - 2.32
Panel 4	319839	0.351	0.250 - 0.738
Panel 5	319839	0.131	0.088 - 0.272

IL-4

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	79.1	51.1 - 98.8
Panel 2	319839	23.4	15.2 - 29.5
Panel 3	319839	8.55	6.18 - 11.4
Panel 4	319839	2.19	1.58 - 3.73
Panel 5	319839	1.11	0.531 - 1.50

IL-5

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	48.7	31.1 - 80.5
Panel 2	319839	16.7	8.64 - 23.8
Panel 3	319839	6.61	3.24 - 9.83
Panel 4	319839	2.23	0.78 - 3.01
Panel 5	319839	0.846	0.273 - 1.23

IFNγ

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	11.3	5.46 - 15.5
Panel 2	319839	3.80	1.56 - 4.89
Panel 3	319839	1.73	0.631 - 2.10
Panel 4	319839	0.516	0.200 - 0.670
Panel 5	319839	0.252	0.095 - 0.334

* 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	175	148 - 294
Panel 2	319839	48.0	41.0 - 69.5
Panel 3	319839	19.0	15.0 - 28.9
Panel 4	319839	7.13	4.50 - 11.1
Panel 5	319839	2.65	1.50 - 4.46

IL-8

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	291	189 - 393
Panel 2	319839	78.0	51.3 - 106
Panel 3	319839	28.9	20.3 - 40.9
Panel 4	319839	9.18	6.02 - 13.0
Panel 5	319839	3.38	2.18 - 5.17

IL-22

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	54.2	32.0 - 101
Panel 2	319839	6.94	5.67 - 14.9
Panel 3	319839	3.65	2.42 - 7.55
Panel 4	319839	0.650	0.54 - 1.50
Panel 5	319839	0.226	0.165 - 0.439

TNFα

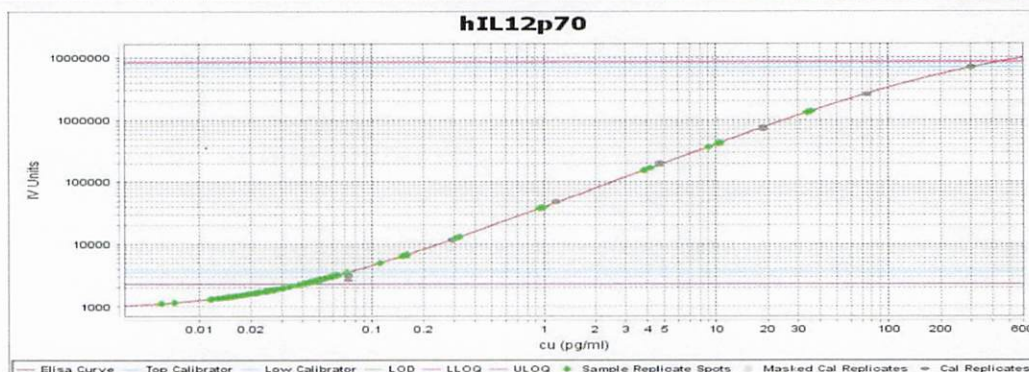
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	74.3	35.8 - 101
Panel 2	319839	24.7	11.6 - 28.3
Panel 3	319839	10.1	4.28 - 11.3
Panel 4	319839	2.89	1.16 - 3.19
Panel 5	319839	1.18	0.413 - 1.34

IL-10

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	43.2	30.3 - 61.7
Panel 2	319839	13.3	9.48 - 17.7
Panel 3	319839	5.45	3.41 - 7.65
Panel 4	319839	1.64	0.944 - 2.59
Panel 5	319839	0.684	0.437 - 1.03

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

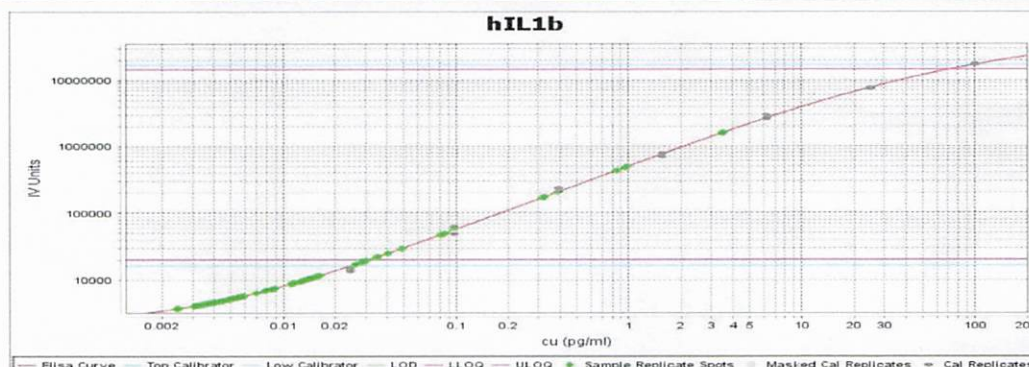
IL-12p70 Calibrator Curve



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

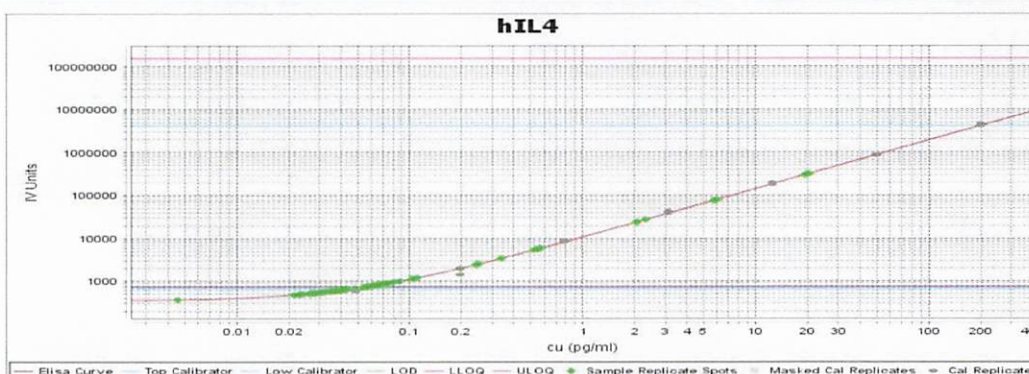
IL-1β Calibrator Curve



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

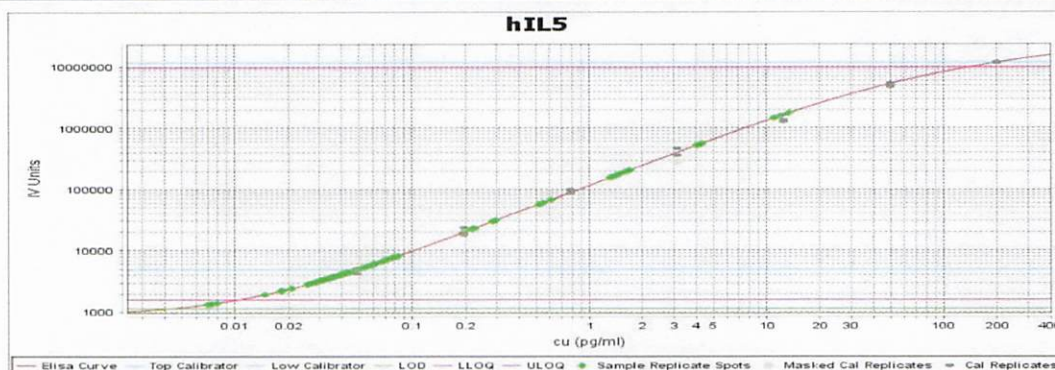
IL-4 Calibrator Curve



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

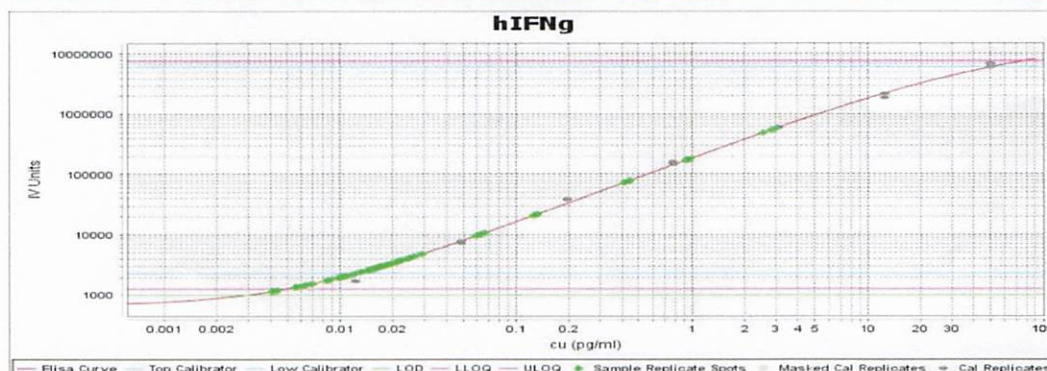
IL-5 Calibrator Curve



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

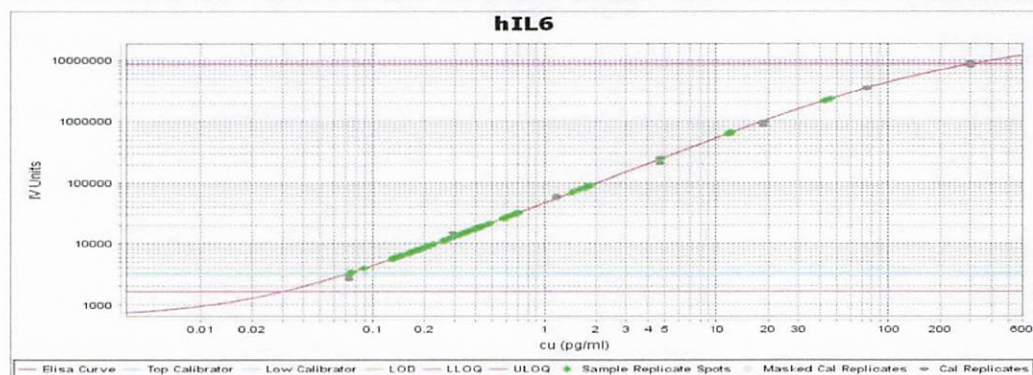
IFNγ Calibrator Curve



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

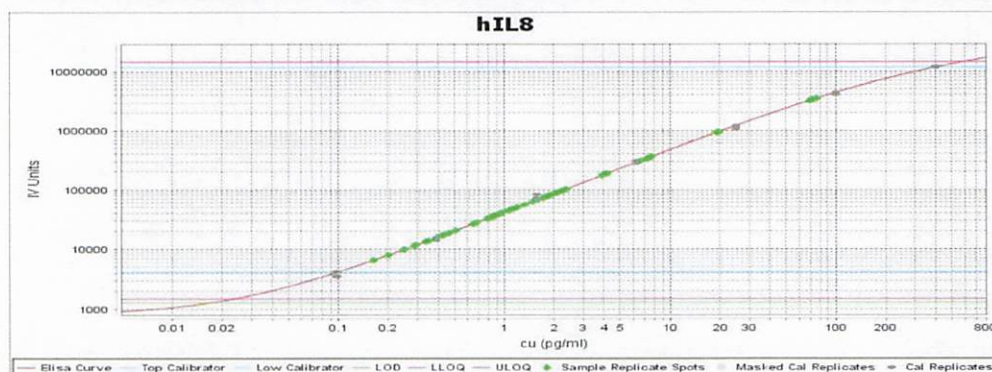
IL-6 Calibrator Curve



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

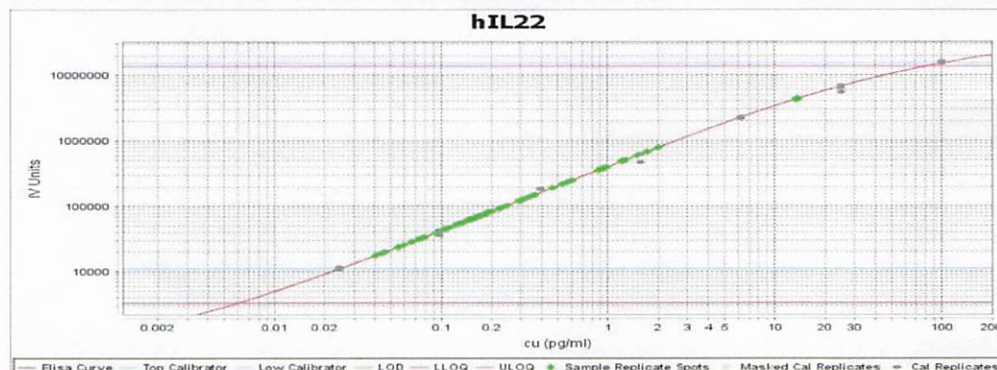
IL-8 Calibrator Curve



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

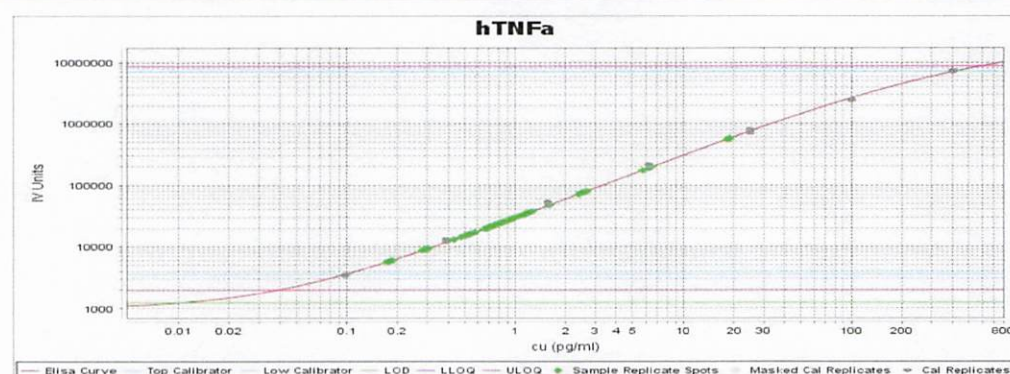
IL-22 Calibrator Curve



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

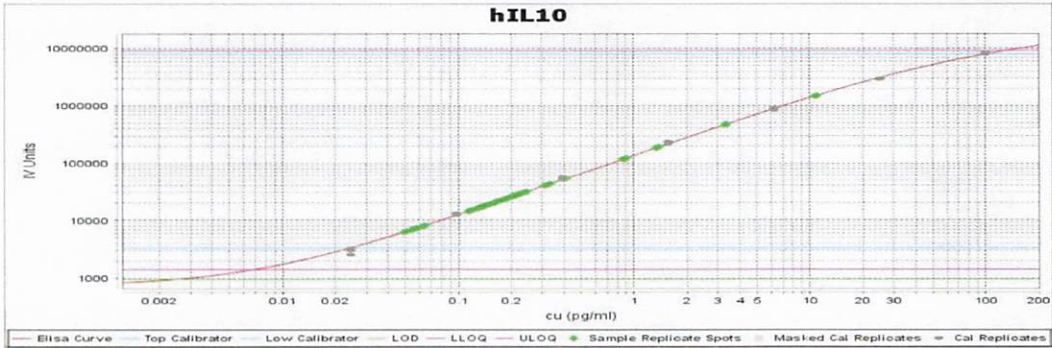
TNFα Calibrator Curve



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