

Product Number:	85-0329
Lot Number:	504715
Expiration:	1-Sep-2026
Platform(s):	SP-X

CorPlex Cytokine Plate	97-0329	523804P
Biotinylated Antibody Reagent	98-0329	524007B
Sample Diluent 21	1863329	505110
IL-12p70 Calibrator	18910034	422511
IL-1β Calibrator	18910040	407211
IL-4 Calibrator	1858452	319458
IL-5 Calibrator	1858462	319838
IFNγ Calibrator	1858260	319843
IL-6 Calibrator	18910044	324303
IL-8 Calibrator	1858454	319575
IL-22 Calibrator	18910036	418312
TNFα Calibrator	1880012	403812
IL-10 Calibrator	18910043	319842
Wash Buffer (12.5X)¹	1863332	N/A
Streptavidin-HRP Reagent¹	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

²SuperSignal[®] 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

Shruti Patel QA Specialist II, QA EP | 29 Sep 2025

Name Title Signature/Date

Release Materials*
IL-12p70

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	142	110 - 185
Panel 2	319839	41.4	33.8 - 51.3
Panel 3	319839	16.5	12.7 - 20.9
Panel 4	319839	4.59	2.87 - 6.32
Panel 5	319839	1.71	0.827 - 2.22

IL-1β

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	12.5	9.78 - 19.9
Panel 2	319839	3.77	2.87 - 5.74
Panel 3	319839	1.49	1.09 - 2.27
Panel 4	319839	0.410	0.258 - 0.720
Panel 5	319839	0.167	0.091 - 0.267

IL-4

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	76.5	52.8 - 98.3
Panel 2	319839	21.9	15.8 - 29.1
Panel 3	319839	8.60	6.35 - 11.4
Panel 4	319839	2.85	1.61 - 3.79
Panel 5	319839	1.49	0.451 - 1.66

IL-5

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	51.6	32.4 - 81.2
Panel 2	319839	18.7	9.15 - 24.0
Panel 3	319839	7.24	3.71 - 9.82
Panel 4	319839	2.50	0.884 - 3.14
Panel 5	319839	0.958	0.321 - 1.20

IFNγ

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	14.6	5.88 - 16.4
Panel 2	319839	4.45	1.70 - 5.12
Panel 3	319839	1.86	0.700 - 2.19
Panel 4	319839	0.601	0.212 - 0.731
Panel 5	319839	0.306	0.078 - 0.385

* 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	207	152 - 288
Panel 2	319839	52.7	41.6 - 68.5
Panel 3	319839	22.0	15.4 - 28.4
Panel 4	319839	8.04	4.73 - 10.9
Panel 5	319839	3.31	1.60 - 4.40

IL-8

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	317	195 - 390
Panel 2	319839	86.9	53.4 - 106
Panel 3	319839	31.1	21.1 - 40.5
Panel 4	319839	10.6	6.27 - 13.0
Panel 5	319839	4.31	2.26 - 5.16

IL-22

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	99.4	32.5 - 115
Panel 2	319839	14.9	6.01 - 15.8
Panel 3	319839	6.60	2.77 - 8.25
Panel 4	319839	1.44	0.527 - 1.66
Panel 5	319839	0.479	0.148 - 0.513

TNFα

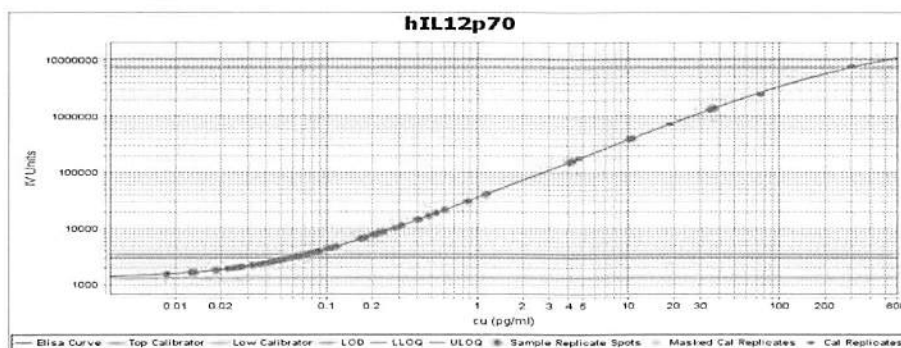
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	91.3	33.8 - 109
Panel 2	319839	28.9	10.6 - 31.5
Panel 3	319839	10.9	3.99 - 12.5
Panel 4	319839	2.98	1.15 - 3.51
Panel 5	319839	1.17	0.414 - 1.51

IL-10

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	57.7	30.7 - 63.4
Panel 2	319839	16.8	9.60 - 18.0
Panel 3	319839	6.91	3.49 - 7.75
Panel 4	319839	2.16	0.987 - 2.58
Panel 5	319839	0.974	0.434 - 1.06

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

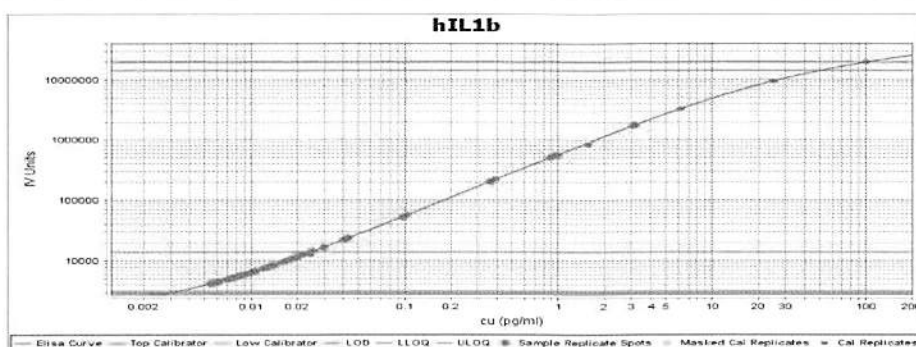
IL-12p70 Calibrator Curve



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

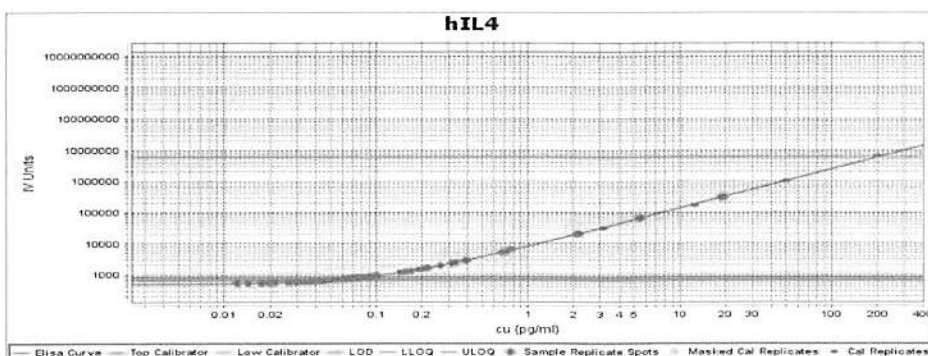
IL-1β Calibrator Curve



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

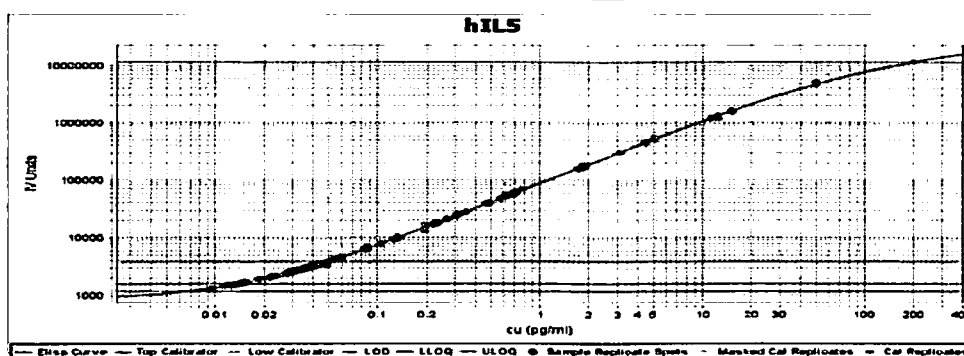
IL-4 Calibrator Curve



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

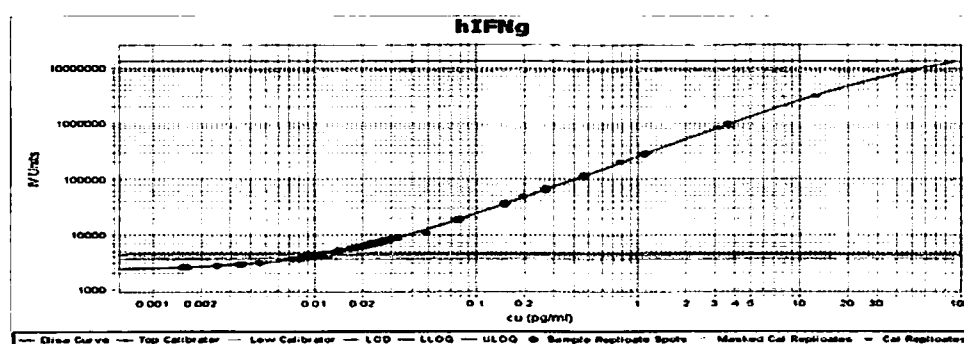
IL-5 Calibrator Curve



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

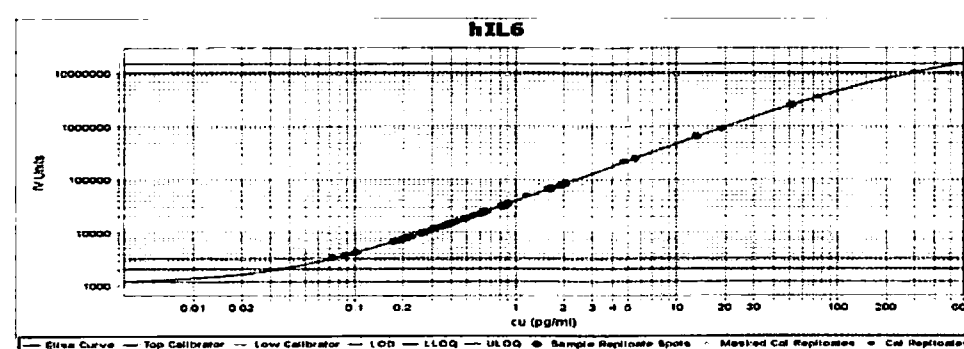
IFNγ Calibrator Curve



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

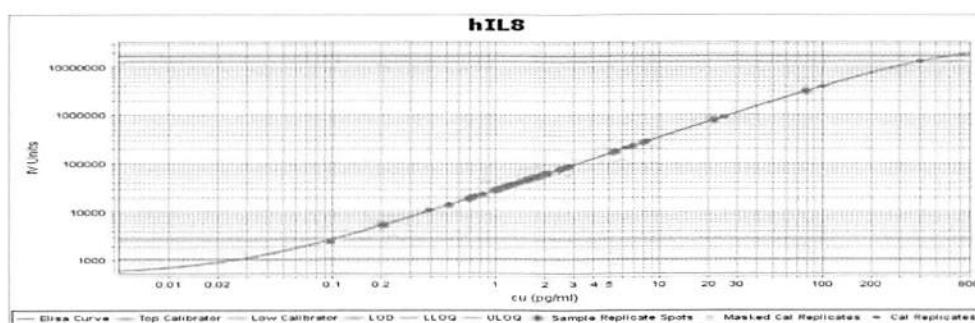
IL-6 Calibrator Curve



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

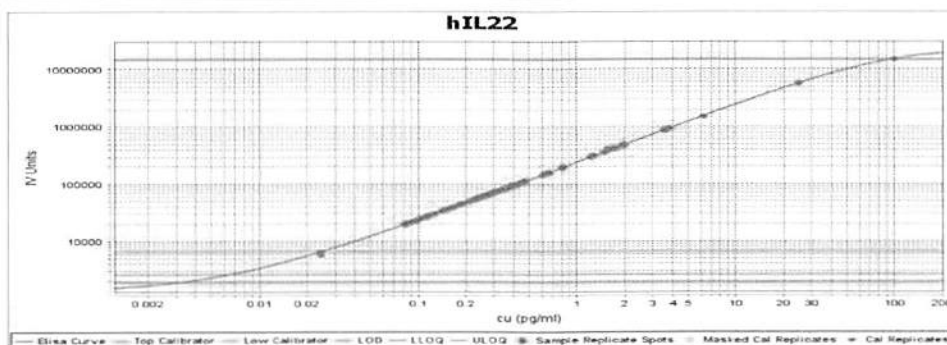
IL-8 Calibrator Curve



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

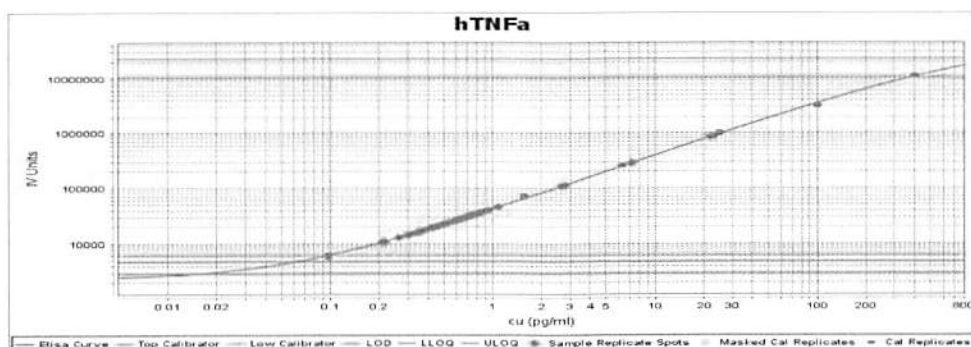
IL-22 Calibrator Curve



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

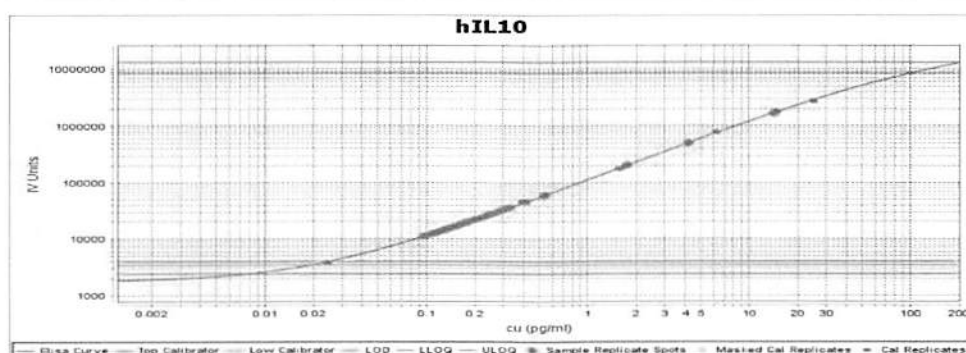
TNFα Calibrator Curve



Calibrator Levels (pg/mL)

A	400
B	100
C	25.0
D	6.25
E	1.562
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