

Product Number:	85-0358
Lot Number:	504338
Expiration:	8-Jun-2025
Platform(s):	SP-X

Component	Part Number	Lot Number	Control Lot Number
CorPlex Cytokine Plate	97-0329	409510P	N/A
Biotinylated Antibody Reagent	98-0329	421304B	N/A
Sample Diluent 21	1863329	311103	311103A
IL-12p70 Calibrator	18910034	318756	N/A
IL-1β Calibrator	18910040	319564	N/A
IL-4 Calibrator	1858452	319016	N/A
IFNγ Calibrator	1858260	319843	N/A
IL-6 Calibrator	18910044	324303	N/A
IL-8 Calibrator	1858454	319436	N/A
TNFα 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 KI-0216 Simoa Corplex Cytokine Panel 1 8-Plex 85-0358.

Review/Approval

Kathleen Cornell

Quality Control Supervisor

Name

Title

Signature/Date

Kathleen Cornell
12 Aug 2024

Release Materials*

IL-12p70

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	130	107 - 186
Panel 2	319839	44.3	33.0 - 51.2
Panel 3	319839	18.6	12.3 - 20.3
Panel 4	319839	5.24	2.59 - 6.31
Panel 5	319839	1.74	0.728 - 2.17

IL-1β

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	14.5	9.55 - 20.6
Panel 2	319839	5.24	2.82 - 5.78
Panel 3	319839	2.11	1.07 - 2.22
Panel 4	319839	0.662	0.237 - 0.723
Panel 5	319839	0.222	0.081 - 0.271

IL-4

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	69.9	50.1 - 97.2
Panel 2	319839	24.1	14.5 - 27.5
Panel 3	319839	9.83	5.89 - 10.8
Panel 4	319839	3.21	1.59 - 3.47
Panel 5	319839	1.12	0.553 - 1.34

IFNγ

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	11.8	5.18 - 14.5
Panel 2	319839	4.20	1.55 - 4.66
Panel 3	319839	1.59	0.710 - 1.83
Panel 4	319839	0.545	0.217 - 0.604
Panel 5	319839	0.252	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	182	149 - 299
Panel 2	319839	55.4	39.7 - 70.6
Panel 3	319839	23.3	14.2 - 29.4
Panel 4	319839	8.89	4.15 - 11.2
Panel 5	319839	3.27	1.32 - 4.52

IL-8

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	319	183 - 383
Panel 2	319839	84.8	49.8 - 103
Panel 3	319839	34.4	19.9 - 39.3
Panel 4	319839	10.9	5.82 - 12.6
Panel 5	319839	4.24	2.12 - 4.95

TNFα

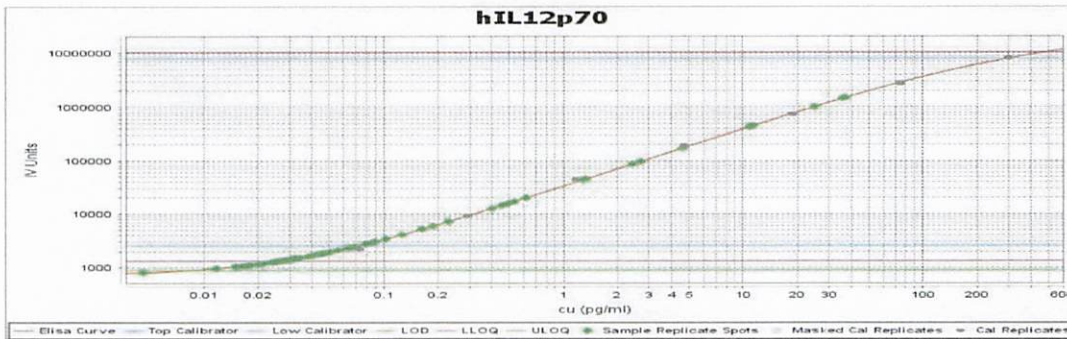
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	46.8	37.9 - 99.8
Panel 2	319839	20.8	11.8 - 26.5
Panel 3	319839	8.25	4.09 - 10.7
Panel 4	319839	2.11	1.01 - 2.84
Panel 5	319839	0.743	0.369 - 1.09

IL-10

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	33.6	30.4 - 61.9
Panel 2	319839	13.6	9.02 - 17.8
Panel 3	319839	5.87	3.19 - 7.61
Panel 4	319839	1.93	1.01 - 2.34
Panel 5	319839	0.794	0.409 - 1.02

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

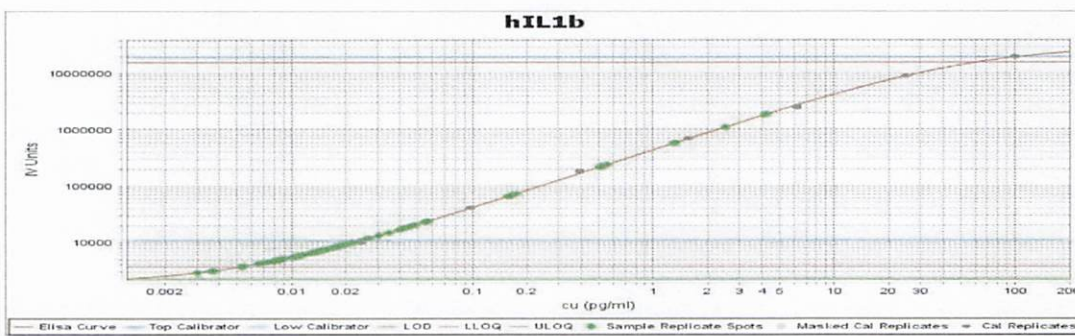
IL-12p70 Calibrator Curve



**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

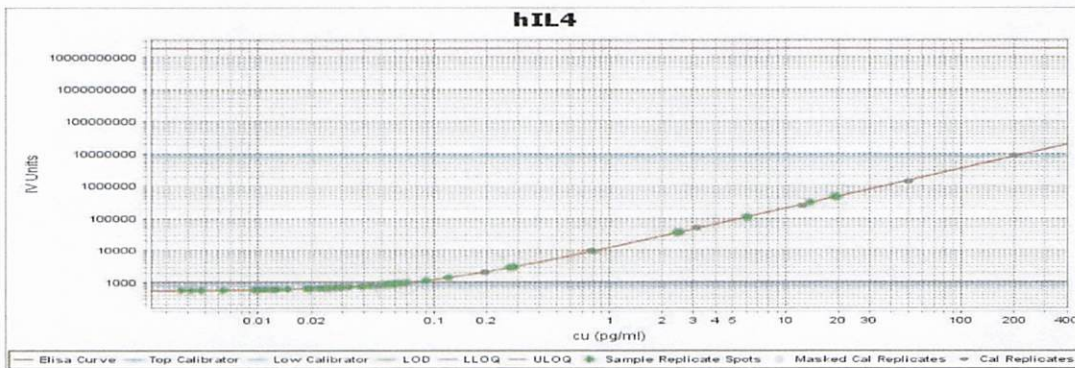
IL-1β Calibrator Curve



**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

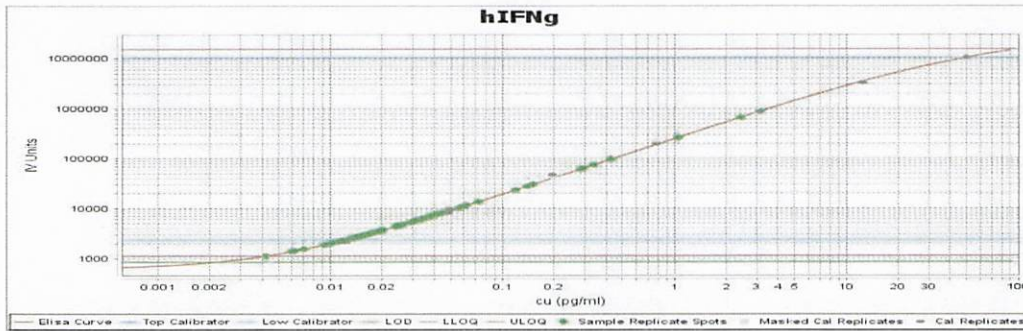
IL-4 Calibrator Curve



**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

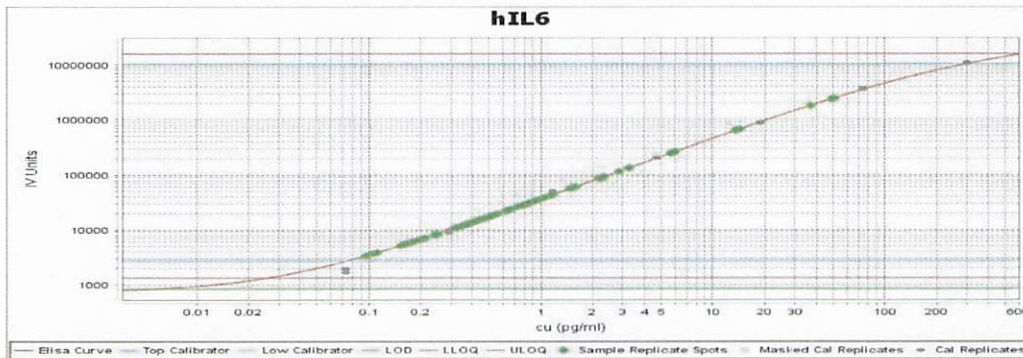
IFN γ Calibrator Curve



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

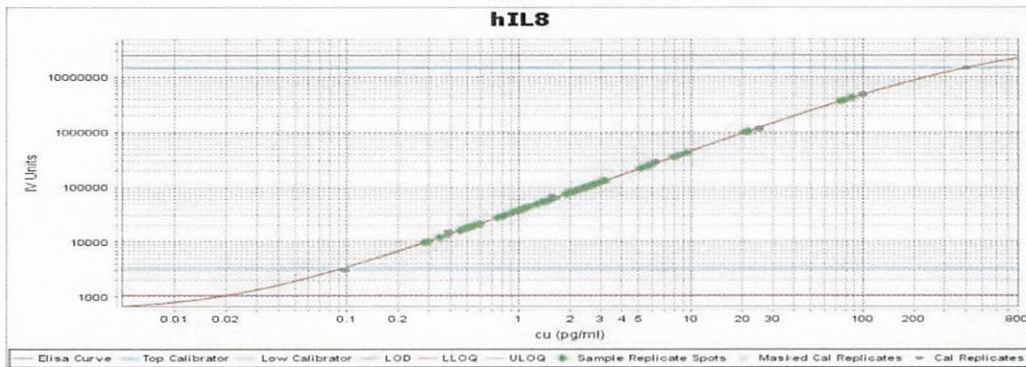
IL-6 Calibrator Curve



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

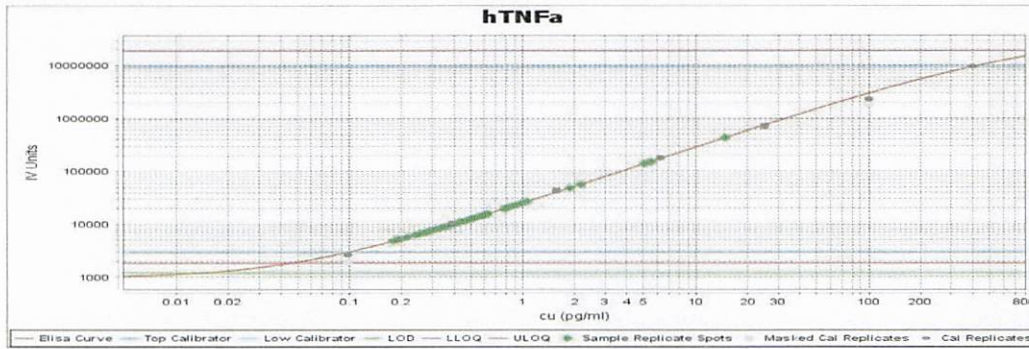
IL-8 Calibrator Curve



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

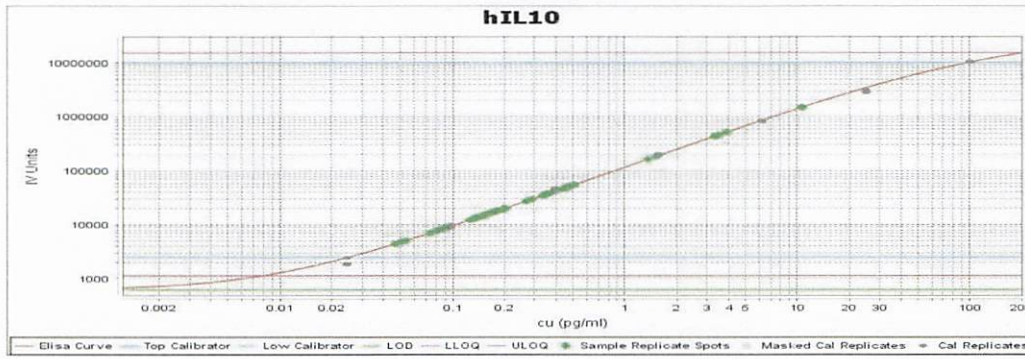
TNFα Calibrator Curve



**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