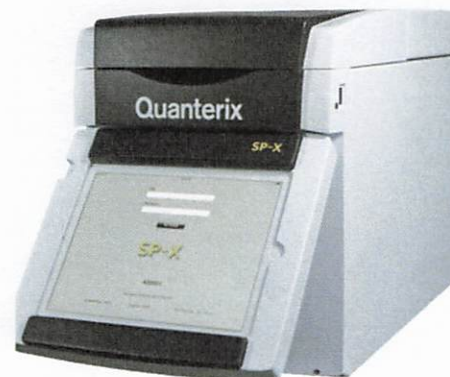


Product Number:	85-0329
Lot Number:	504299
Expiration:	5-Jul-2025
Platform(s):	SP-X

CorPlex Cytokine Plate	97-0329	417205P
Biotinylated Antibody Reagent	98-0329	417908B
Sample Diluent 21	1863329	319974
IL-12p70 Calibrator	18910034	319156
IL-1β Calibrator	18910040	407211
IL-4 Calibrator	1858452	319016
IL-5 Calibrator	1858462	318837
IFNγ Calibrator	1858260	319589
IL-6 Calibrator	18910044	324303
IL-8 Calibrator	1858454	319436
IL-22 Calibrator	18910036	319176
TNFα Calibrator	1880012	403812
IL-10 Calibrator	18910043	319842
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 Corplex Cytokine Panel 1 Instruction Book 85-0329 116-7BF-1-AB (KI-0195).

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title

Kathleen Cornell 23 Jul 2024
Signature/Date

Release Materials*

IL-12p70

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	169	107 - 186
Panel 2	319839	50.1	33.0 - 51.2
Panel 3	319839	19.4	12.3 - 20.3
Panel 4	319839	5.35	2.59 - 6.31
Panel 5	319839	1.63	0.728 - 2.17

IL-1β

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	16.8	9.55 - 20.6
Panel 2	319839	5.02	2.82 - 5.78
Panel 3	319839	1.91	1.07 - 2.22
Panel 4	319839	0.512	0.237 - 0.723
Panel 5	319839	0.172	0.081 - 0.271

IL-4

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	79.7	50.1 - 97.2
Panel 2	319839	27.3	14.5 - 27.5
Panel 3	319839	9.32	5.89 - 10.8
Panel 4	319839	2.86	1.59 - 3.47
Panel 5	319839	0.992	0.553 - 1.34

IL-5

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	64.2	31.5 - 70.5
Panel 2	319839	21.5	9.01 - 21.6
Panel 3	319839	6.97	3.34 - 8.64
Panel 4	319839	2.30	0.961 - 2.50
Panel 5	319839	0.832	0.370 - 0.968

IFNγ

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	9.31	5.18 - 14.5
Panel 2	319839	2.94	1.55 - 4.66
Panel 3	319839	1.22	0.710 - 1.83
Panel 4	319839	0.378	0.217 - 0.604
Panel 5	319839	0.169	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	251	149 - 299
Panel 2	319839	67.4	39.7 - 70.6
Panel 3	319839	25.4	14.2 - 29.4
Panel 4	319839	8.83	4.15 - 11.2
Panel 5	319839	3.20	1.32 - 4.52

IL-8

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	344	183 - 383
Panel 2	319839	97.3	49.8 - 103
Panel 3	319839	37.0	19.9 - 39.3
Panel 4	319839	11.2	5.82 - 12.6
Panel 5	319839	4.50	2.12 - 4.95

IL-22

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	118	36.2 - 101
Panel 2	319839	14.0	6.05 - 15.0
Panel 3	319839	6.60	2.75 - 7.32
Panel 4	319839	1.12	0.565 - 1.49
Panel 5	319839	0.359	0.149 - 0.446

TNFα

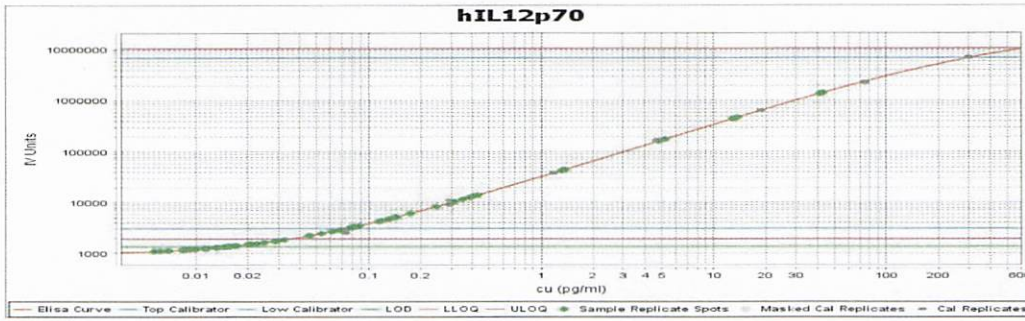
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	93.2	37.9 - 99.8
Panel 2	319839	28.9	11.8 - 26.5
Panel 3	319839	9.92	4.09 - 10.7
Panel 4	319839	2.61	1.01 - 2.84
Panel 5	319839	1.01	0.369 - 1.09

IL-10

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	58.8	30.4 - 61.9
Panel 2	319839	16.0	9.02 - 17.8
Panel 3	319839	7.08	3.19 - 7.61
Panel 4	319839	2.16	1.01 - 2.34
Panel 5	319839	0.928	0.409 - 1.02

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

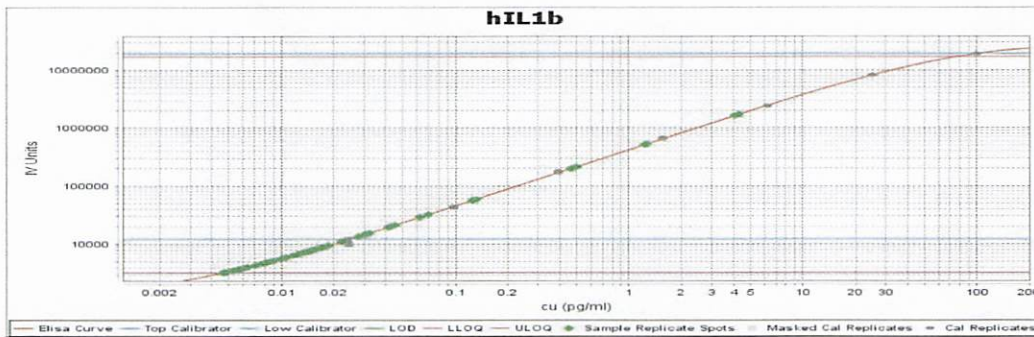
IL-12p70 Calibrator Curve



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

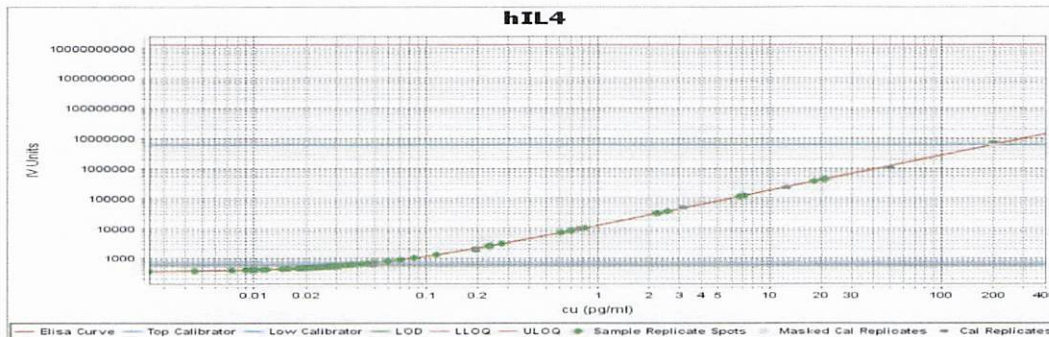
IL-1β Calibrator Curve



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

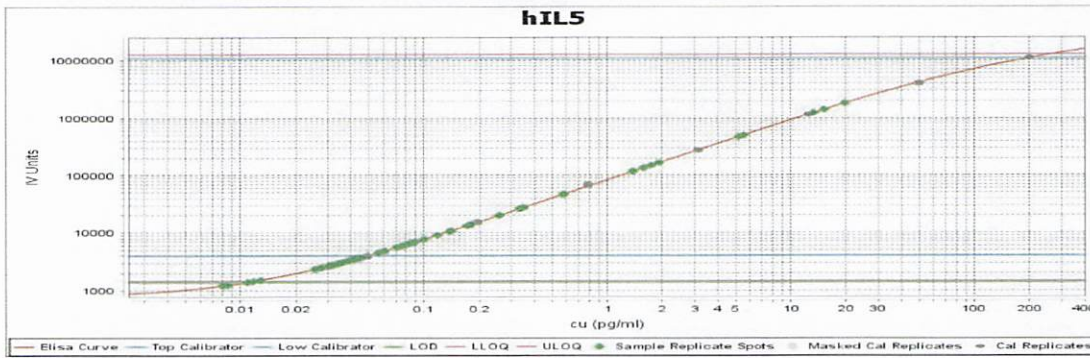
IL-4 Calibrator Curve



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

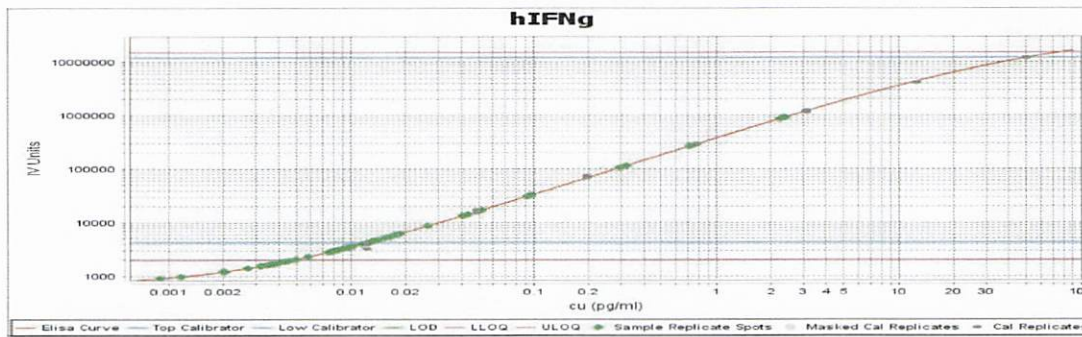
IL-5 Calibrator Curve



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

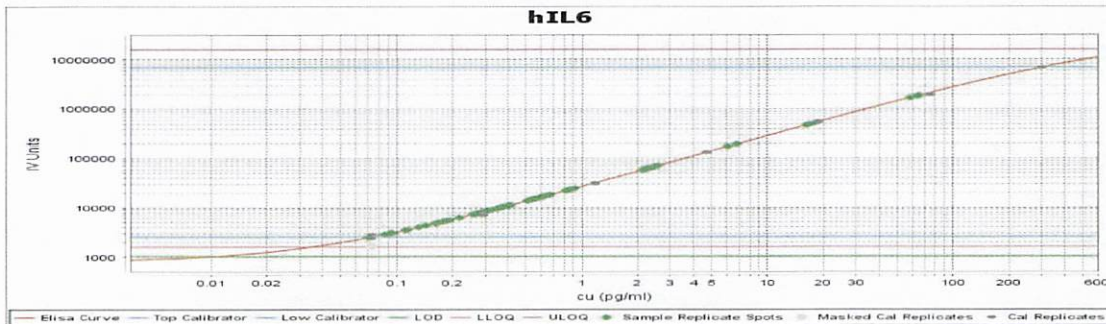
IFNγ Calibrator Curve



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

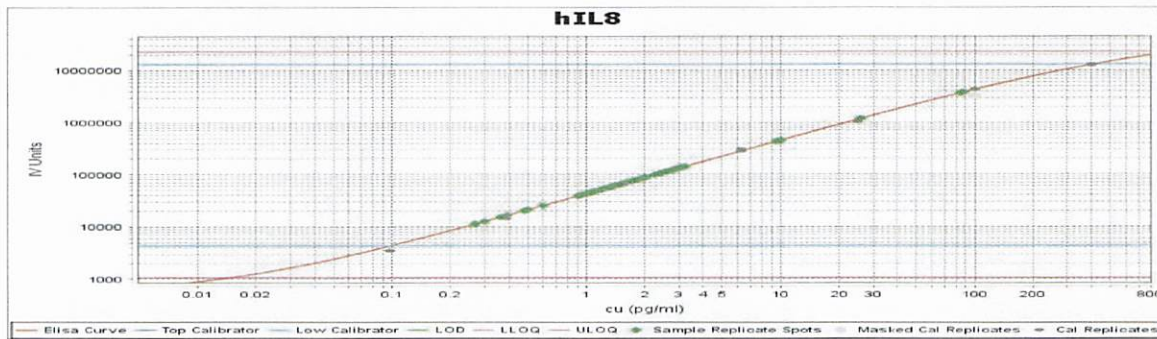
IL-6 Calibrator Curve



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

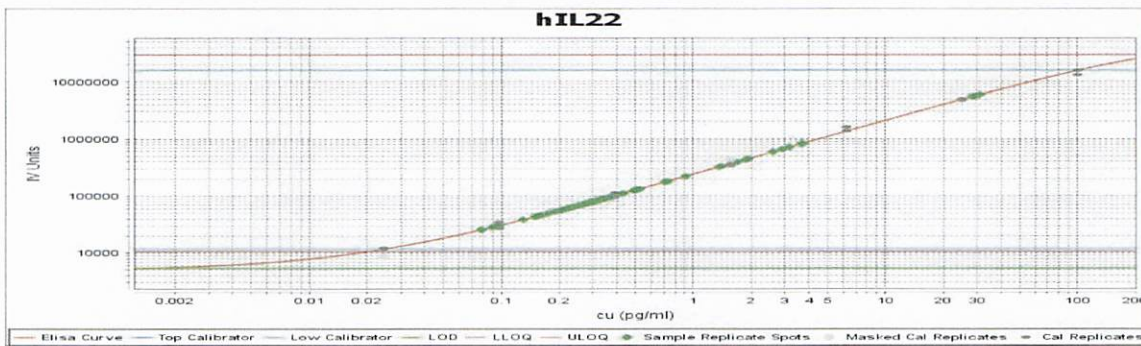
IL-8 Calibrator Curve



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

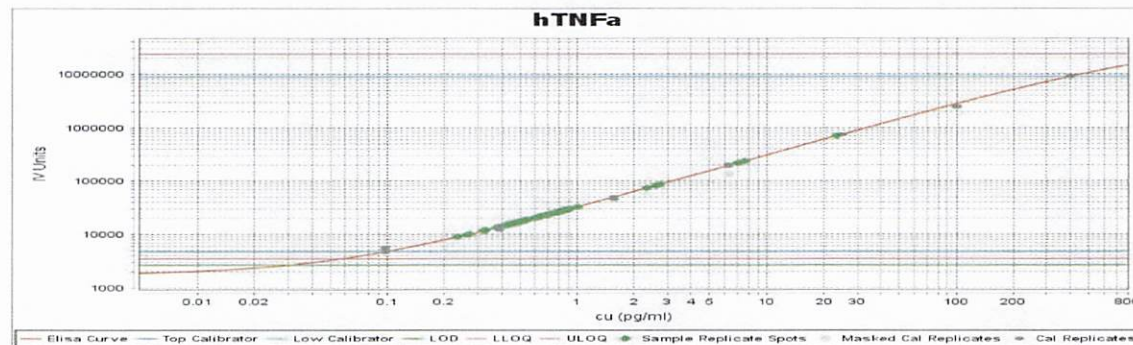
IL-22 Calibrator Curve



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

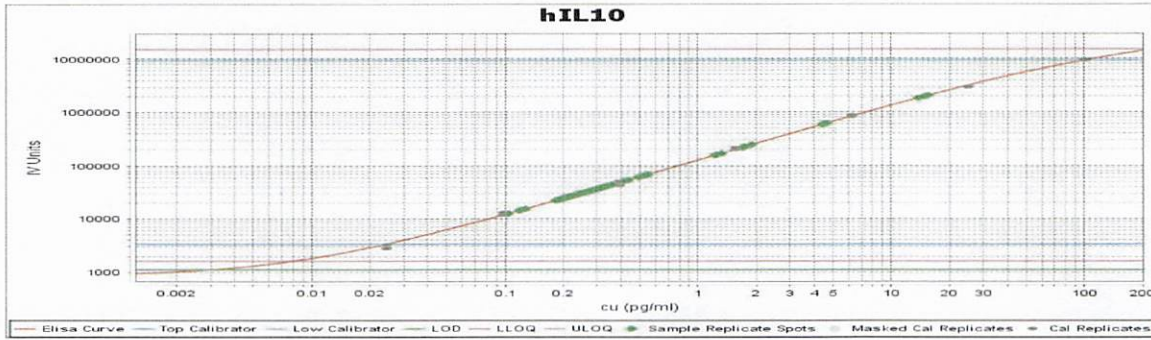
TNFα Calibrator Curve



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