

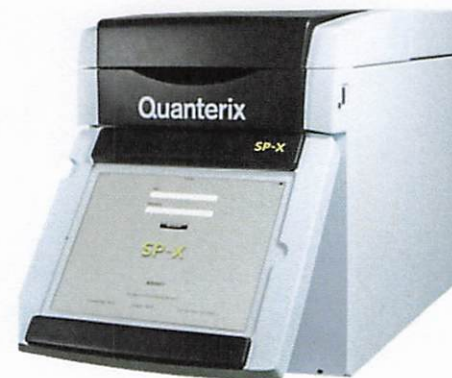
Product Number:	85-0329
Lot Number:	504588
Expiration:	30-Apr-2025
Platform(s):	SP-X

CorPlex Cytokine Plate	97-0329	407512P
Biotinylated Antibody Reagent	98-0329	408108B
Sample Diluent 21	1863329	403020
IL-12p70 Calibrator	18910034	319844
IL-1β Calibrator	18910040	319165
IL-4 Calibrator	1858452	319016
IL-5 Calibrator	1858462	319838
IFNγ Calibrator	1858260	319843
IL-6 Calibrator	18910044	324303
IL-8 Calibrator	1858454	319436
IL-22 Calibrator	18910036	319176
TNFα Calibrator	1880012	319460
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

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).



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

Review/Approval

Kyle Giroux Senior Document Control Specialist, QA [Signature] 16Apr2025
Name Title Signature/Date

Release Materials*
IL-12p70

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	165	108 - 185
Panel 2	319839	44.6	33.5 - 51.7
Panel 3	319839	17.6	12.4 - 21.1
Panel 4	319839	4.55	2.73 - 6.45
Panel 5	319839	1.42	0.781 - 2.23

IL-1β

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	14.5	9.83 - 20.2
Panel 2	319839	3.93	2.87 - 5.85
Panel 3	319839	1.53	1.07 - 2.32
Panel 4	319839	0.430	0.250 - 0.738
Panel 5	319839	0.138	0.088 - 0.272

IL-4

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	83.5	51.1 - 98.8
Panel 2	319839	23.1	15.2 - 29.5
Panel 3	319839	8.78	6.18 - 11.4
Panel 4	319839	2.44	1.58 - 3.73
Panel 5	319839	0.733	0.531 - 1.50

IL-5

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	70.3	31.1 - 80.5
Panel 2	319839	19.9	8.64 - 23.8
Panel 3	319839	8.05	3.24 - 9.83
Panel 4	319839	2.36	0.78 - 3.01
Panel 5	319839	0.831	0.273 - 1.23

IFNγ

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	11.7	5.46 - 15.5
Panel 2	319839	3.44	1.56 - 4.89
Panel 3	319839	1.36	0.631 - 2.10
Panel 4	319839	0.417	0.200 - 0.670
Panel 5	319839	0.184	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	228	148 - 294
Panel 2	319839	57.6	41.0 - 69.5
Panel 3	319839	21.3	15.0 - 28.9
Panel 4	319839	7.40	4.50 - 11.1
Panel 5	319839	2.75	1.50 - 4.46

IL-8

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	277	189 - 393
Panel 2	319839	84.9	51.3 - 106
Panel 3	319839	30.2	20.3 - 40.9
Panel 4	319839	9.57	6.02 - 13.0
Panel 5	319839	3.26	2.18 - 5.17

IL-22

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	109	32.0 - 101
Panel 2	319839	10.9	5.67 - 14.9
Panel 3	319839	6.53	2.42 - 7.55
Panel 4	319839	0.848	0.540 - 1.50
Panel 5	319839	0.329	0.165 - 0.439

TNFα

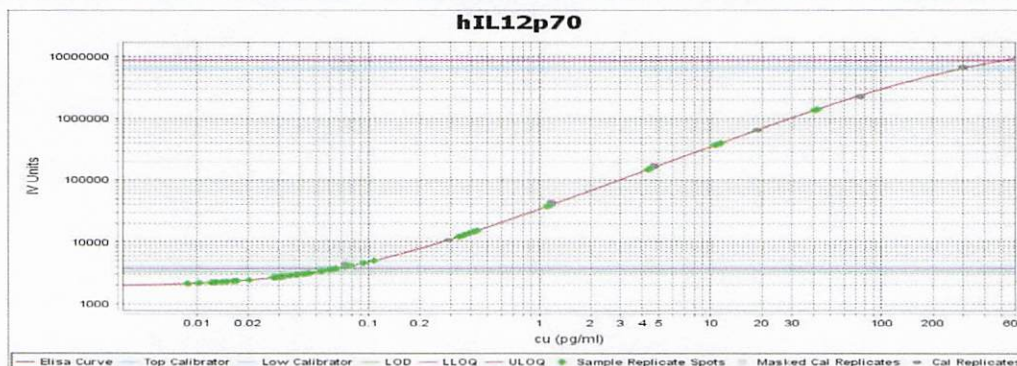
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	64.0	35.8 - 101
Panel 2	319839	17.7	11.6 - 28.3
Panel 3	319839	6.20	4.28 - 11.3
Panel 4	319839	1.50	1.16 - 3.19
Panel 5	319839	0.603	0.413 - 1.34

IL-10

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	51.3	30.3 - 61.7
Panel 2	319839	14.1	9.48 - 17.7
Panel 3	319839	5.33	3.41 - 7.65
Panel 4	319839	1.56	0.944 - 2.59
Panel 5	319839	0.634	0.437 - 1.03

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

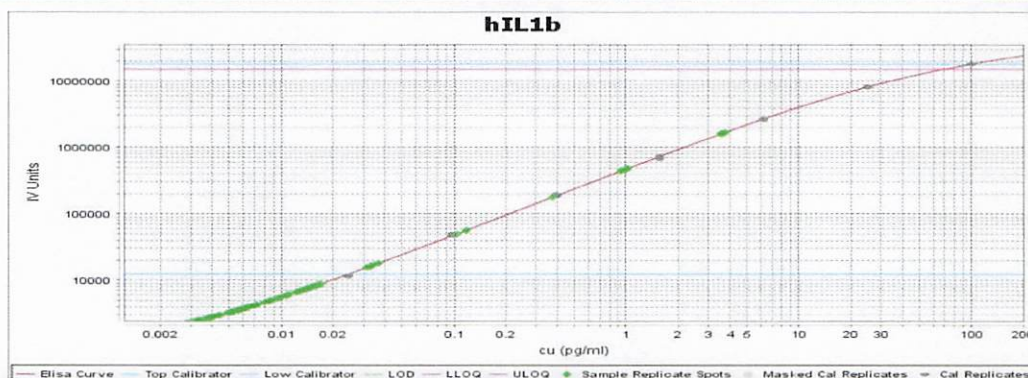
IL-12p70 Calibrator Curve



**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

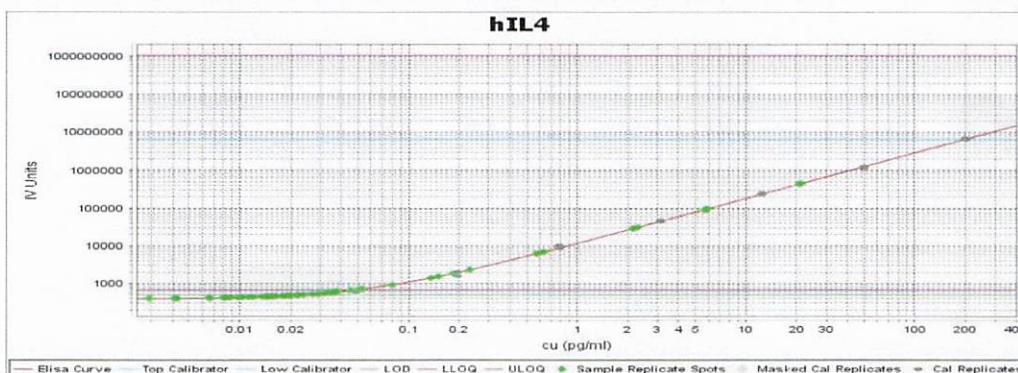
IL-1 β Calibrator Curve



**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

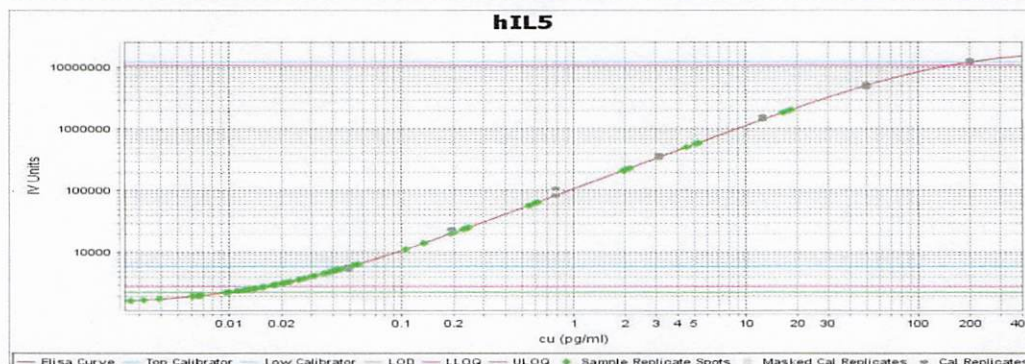
IL-4 Calibrator Curve



**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

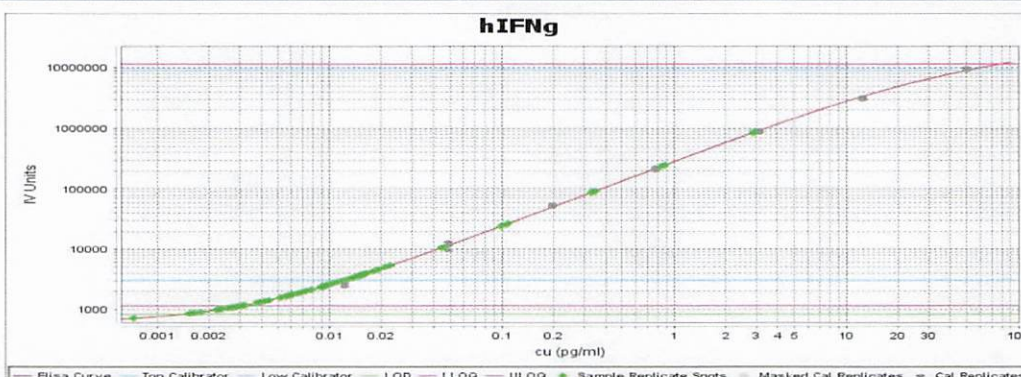
IL-5 Calibrator Curve



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

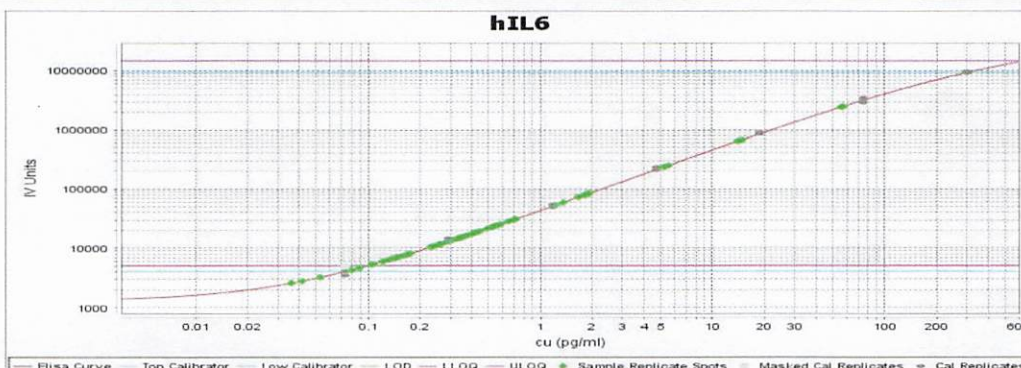
IFNγ Calibrator Curve



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

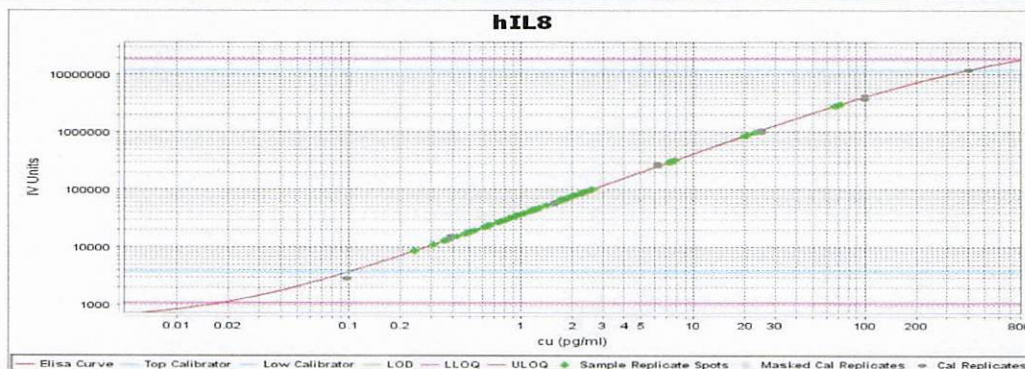
IL-6 Calibrator Curve



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

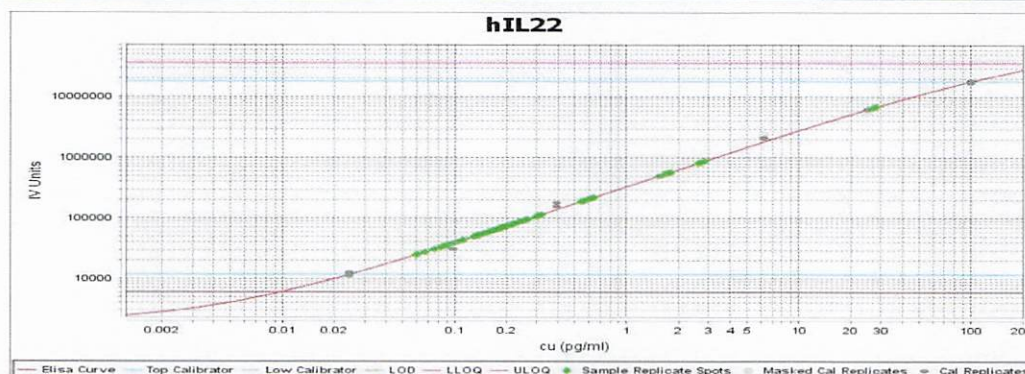
IL-8 Calibrator Curve



**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

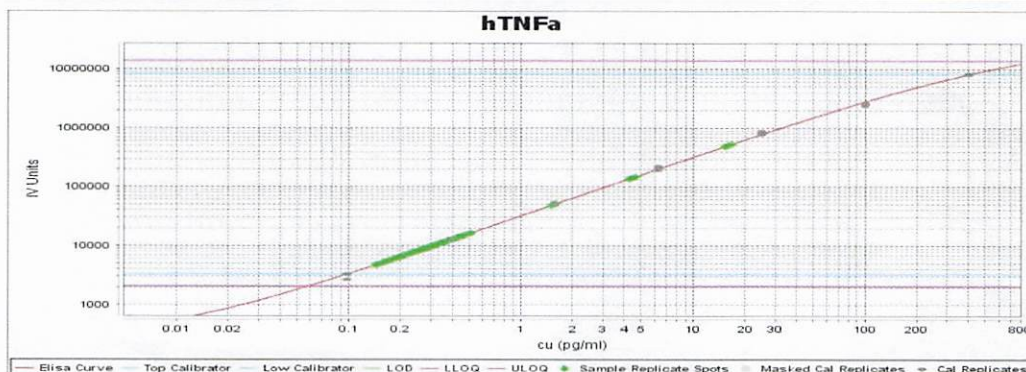
IL-22 Calibrator Curve



**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

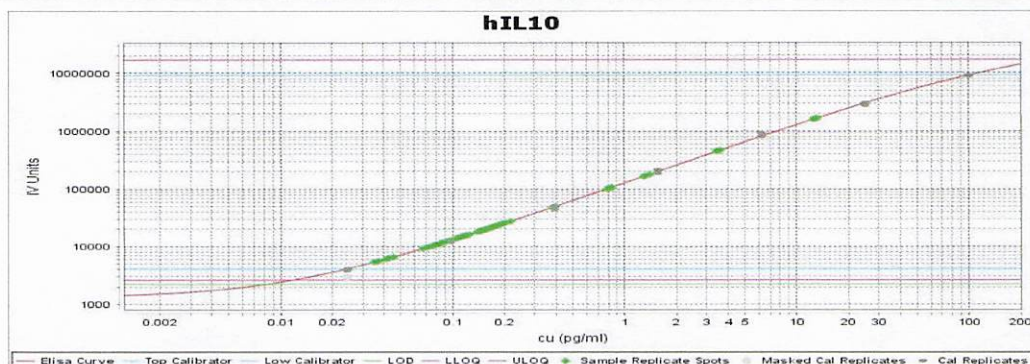
TNFα Calibrator Curve



**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