

Product Number:	85-0329
Lot Number:	504418
Expiration:	8-Apr-2025
Platform(s):	SP-X



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

CorPlex Cytokine Plate	97-0329	427107P
Biotinylated Antibody Reagent	98-0329	427108B
Sample Diluent 21	1863329	405709
IL-12p70 Calibrator	18910034	319844
IL-1β Calibrator	18910040	407211
IL-4 Calibrator	1858452	319016
IL-5 Calibrator	1858462	318655
IFNγ Calibrator	1858260	319843
IL-6 Calibrator	18910044	324303
IL-8 Calibrator	1858454	319436
IL-22 Calibrator	18910036	319459
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

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

[Signature] 27 Nov 2024
Signature/Date

Release Materials*

IL-12p70

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	153	107 - 185
Panel 2	319839	40.8	33.4 - 51.2
Panel 3	319839	16.5	12.3 - 20.7
Panel 4	319839	4.45	2.67 - 6.35
Panel 5	319839	1.52	0.760 - 2.18

IL-1β

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	14.8	9.79 - 20.4
Panel 2	319839	3.79	2.85 - 5.89
Panel 3	319839	1.66	1.06 - 2.31
Panel 4	319839	0.459	0.238 - 0.745
Panel 5	319839	0.197	0.084 - 0.274

IL-4

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	75.2	50.9 - 96.0
Panel 2	319839	20.4	14.8 - 27.5
Panel 3	319839	8.44	5.94 - 11.0
Panel 4	319839	3.04	1.59 - 3.56
Panel 5	319839	1.60	0.558 - 1.37

IL-5

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	54.4	31.3 - 73.2
Panel 2	319839	19.4	8.91 - 22.5
Panel 3	319839	7.09	3.45 - 9.05
Panel 4	319839	2.48	0.829 - 2.79
Panel 5	319839	0.826	0.365 - 1.01

IFNγ

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	10.8	5.23 - 14.5
Panel 2	319839	3.03	1.52 - 4.74
Panel 3	319839	1.40	0.711 - 1.85
Panel 4	319839	0.453	0.216 - 0.614
Panel 5	319839	0.249	0.087 - 0.288

* 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	200	148 - 298
Panel 2	319839	49.9	40.5 - 70.0
Panel 3	319839	19.7	14.6 - 29.2
Panel 4	319839	7.46	4.31 - 11.1
Panel 5	319839	2.69	1.40 - 4.47

IL-8

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	290	181 - 395
Panel 2	319839	78.9	49.8 - 105
Panel 3	319839	30.3	19.8 - 40.6
Panel 4	319839	8.83	5.79 - 13.0
Panel 5	319839	3.44	2.14 - 5.07

IL-22

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	72.2	33.7 - 103
Panel 2	319839	8.65	5.99 - 15.1
Panel 3	319839	4.89	2.56 - 7.64
Panel 4	319839	0.856	0.559 - 1.51
Panel 5	319839	0.292	0.148 - 0.449

TNFα

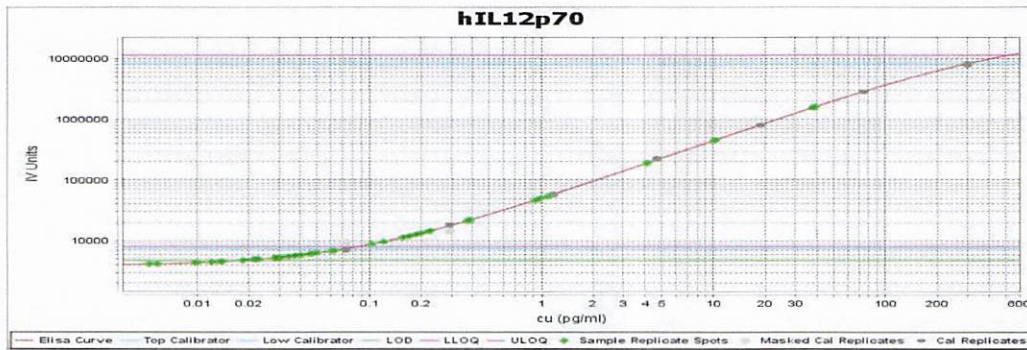
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	66.8	35.7 - 99.4
Panel 2	319839	18.1	12.2 - 26.3
Panel 3	319839	7.24	4.26 - 10.6
Panel 4	319839	1.96	1.05 - 2.81
Panel 5	319839	0.905	0.378 - 1.07

IL-10

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	57.0	29.2 - 62.2
Panel 2	319839	15.3	9.21 - 17.7
Panel 3	319839	5.80	3.28 - 7.66
Panel 4	319839	1.73	1.02 - 2.39
Panel 5	319839	0.723	0.424 - 1.02

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

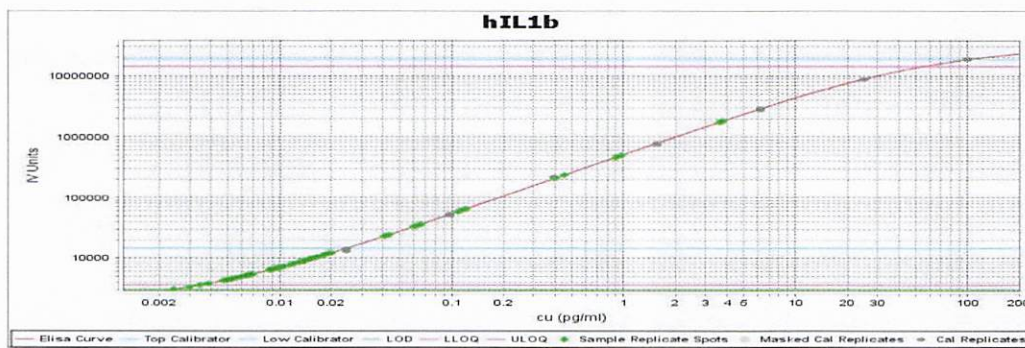
IL-12p70 Calibrator Curve



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

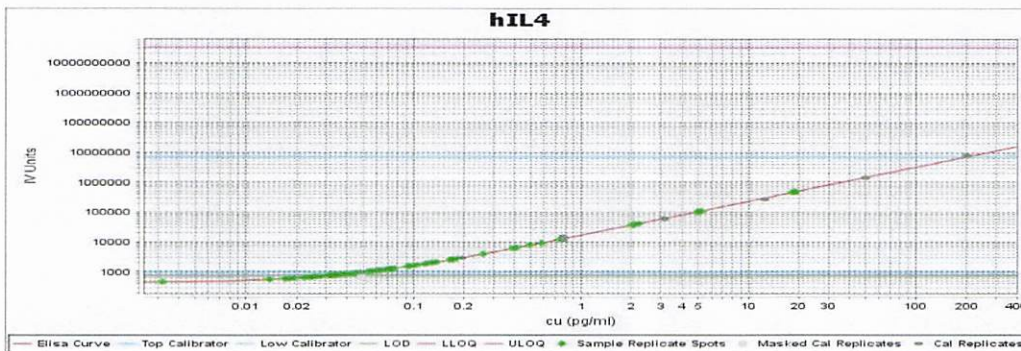
IL-1β Calibrator Curve



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

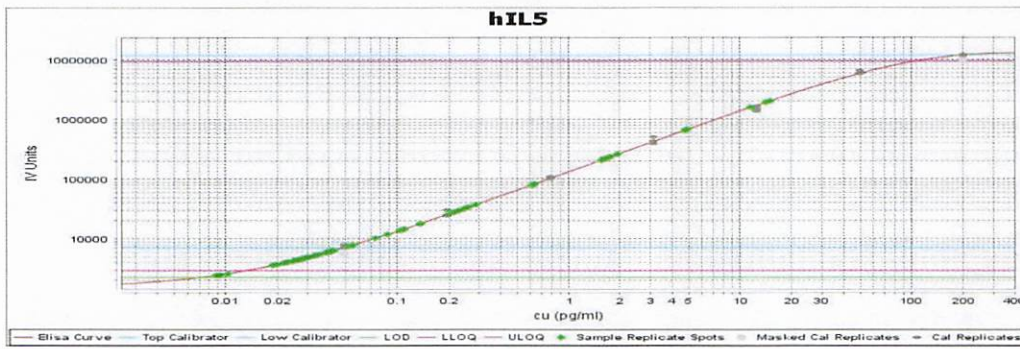
IL-4 Calibrator Curve



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

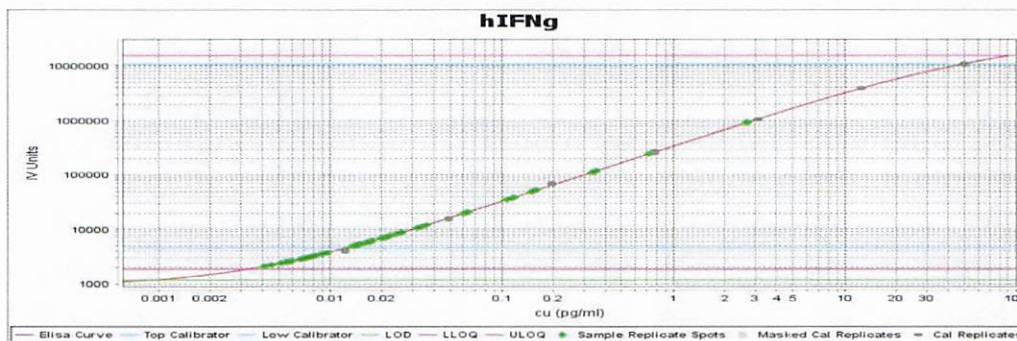
IL-5 Calibrator Curve



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

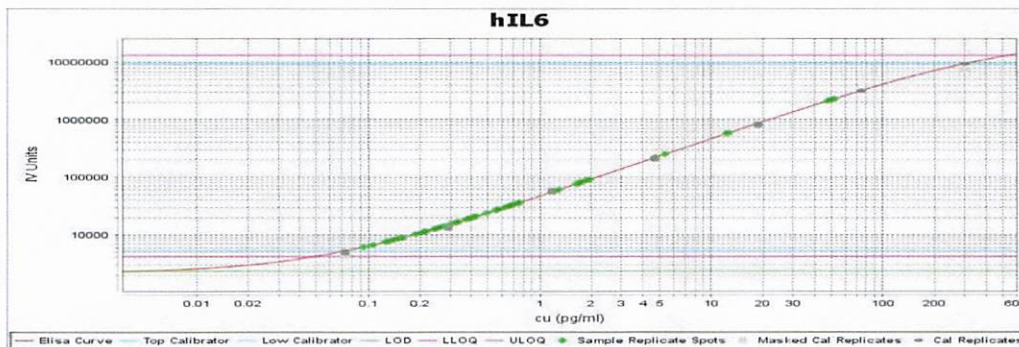
IFNγ Calibrator Curve



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

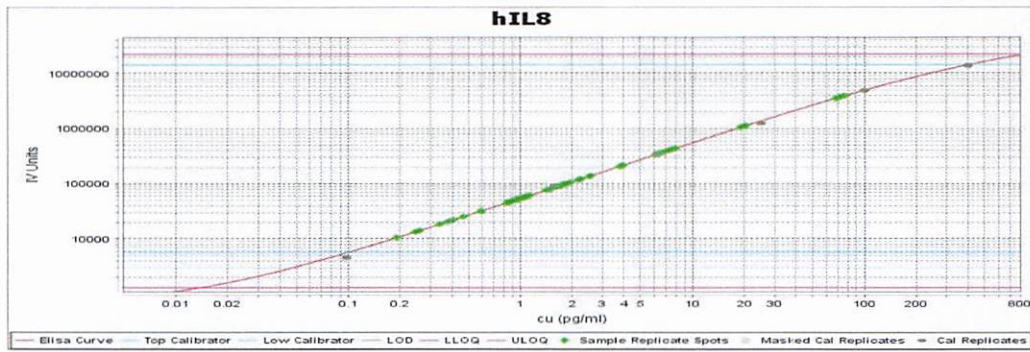
IL-6 Calibrator Curve



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

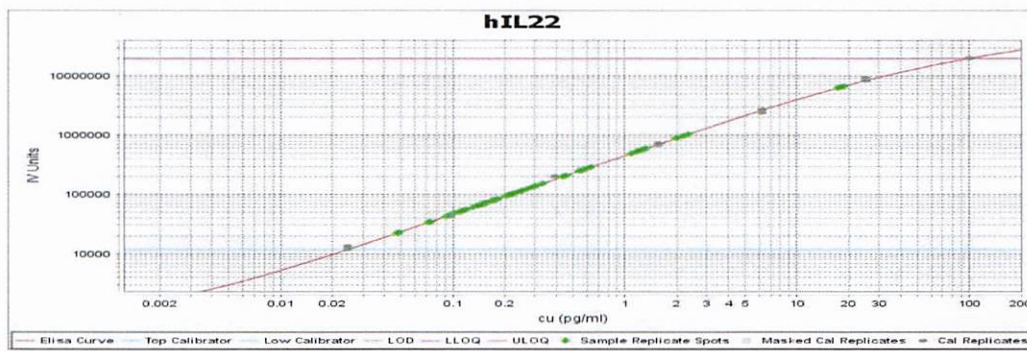
IL-8 Calibrator Curve



**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

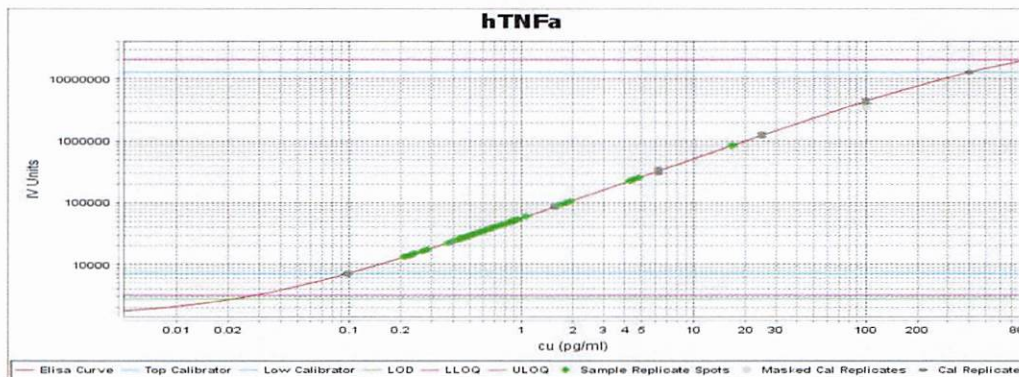
IL-22 Calibrator Curve



**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

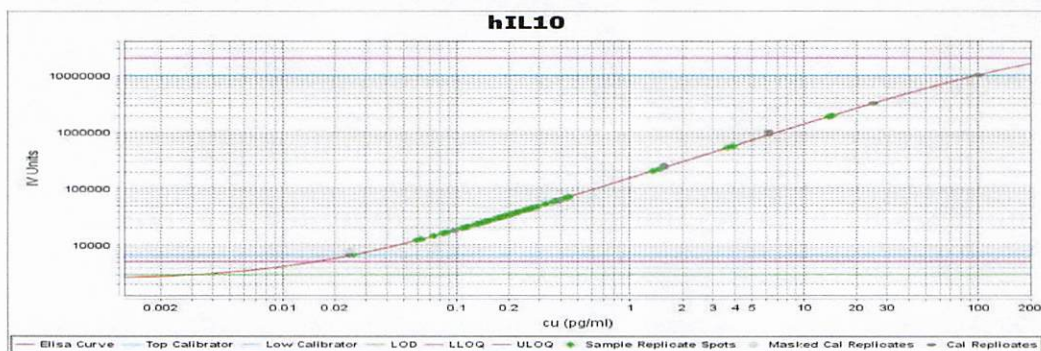
TNFα Calibrator Curve



**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