

Product Number:	85-0363
Lot Number:	504061
Expiration:	13-Jan-2025
Platform(s):	SP-X

CorPlex Cytokine Plate	97-0329	326103P
Biotinylated Antibody Reagent	98-0329	326105B
Sample Diluent 21	1863329	319974
IFNγ Calibrator	1858260	319589
IL-6 Calibrator	18910044	319173
IL-8 Calibrator	1858454	319436
TNFα Calibrator	1880012	319247
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



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 4-Plex Kit Instructions (KI-0220).

Review/Approval

Kathleen Cornell

Quality Control
Supervisor

[Handwritten Signature] 20 Nov 2023

Name

Title

Signature/Date

Release Materials*

IFN γ

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	8.313	5.184 - 14.50
Panel 2	319839	2.487	1.553 - 4.656
Panel 3	319839	1.051	0.710 - 1.834
Panel 4	319839	0.309	0.217 - 0.604
Panel 5	319839	0.132	0.089 - 0.283

IL-6

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	218.6	149.3 - 299.3
Panel 2	319839	54.08	39.74 - 70.62
Panel 3	319839	22.11	14.23 - 29.41
Panel 4	319839	7.213	4.151 - 11.20
Panel 5	319839	2.875	1.321 - 4.519

IL-8

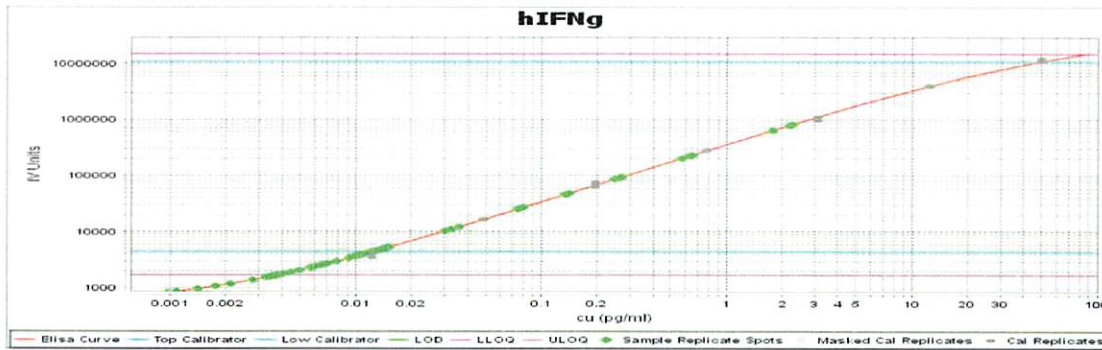
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	277.7	182.6 - 382.7
Panel 2	319839	71.43	49.82 - 102.8
Panel 3	319839	27.62	19.87 - 39.34
Panel 4	319839	8.296	5.821 - 12.63
Panel 5	319839	3.022	2.122 - 4.948

TNF α

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	69.69	37.89 - 99.76
Panel 2	319839	18.52	11.84 - 26.54
Panel 3	319839	7.663	4.095 - 10.69
Panel 4	319839	1.679	1.007 - 2.838
Panel 5	319839	0.539	0.369 - 1.090

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

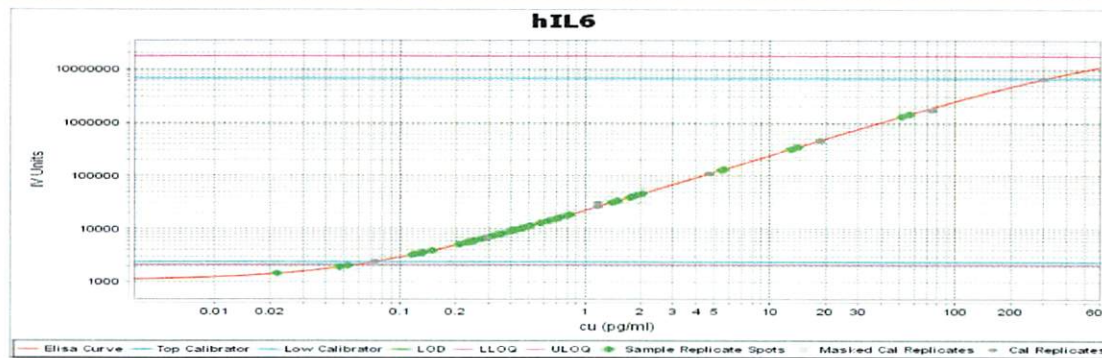
IFN γ Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	50.00
B	12.50
C	3.125
D	0.781
E	0.195
F	0.049
G	0.012
H	0.000

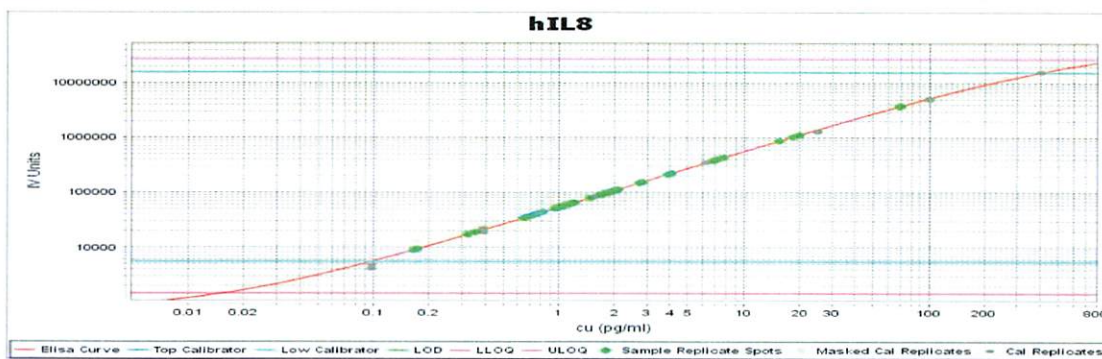
IL-6 Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	300.00
B	75.00
C	18.75
D	4.688
E	1.172
F	0.293
G	0.073
H	0.000

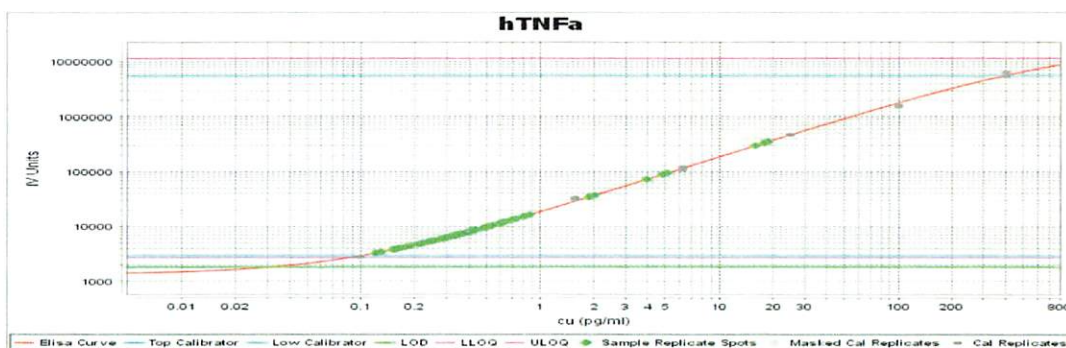
IL-8 Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	400.00
B	100.00
C	25.00
D	6.250
E	1.562
F	0.391
G	0.098
H	0.000

TNFα Calibrator Curve



Calibrator Levels (pg/mL)

A	400.00
B	100.00
C	25.00
D	6.250
E	1.562
F	0.391
G	0.098
H	0.000