

Product Number:	85-0473
Lot Number:	504195
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
IFNγ Calibrator	1858260	319843
IL-6 Calibrator	18910044	324303
IL-8 Calibrator	1858454	319436
TNFα Calibrator	1880012	319460
IL-10 Calibrator	18910043	319842
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 5-Plex Kit Instructions (KI-0258).

Review/Approval	Kathleen Cornell	Quality Control Supervisor		28 Mar 2024
	Name	Title	Signature/Date	

Release Materials*

IFN γ

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	10.1	5.18 - 14.5
Panel 2	319839	2.99	1.55 - 4.66
Panel 3	319839	1.26	0.710 - 1.83
Panel 4	319839	0.393	0.217 - 0.604
Panel 5	319839	0.154	0.089 - 0.283

IL-6

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	199	149 - 299
Panel 2	319839	49.7	39.7 - 70.6
Panel 3	319839	19.4	14.2 - 29.4
Panel 4	319839	7.16	4.15 - 11.2
Panel 5	319839	2.55	1.32 - 4.52

IL-8

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	255	183 - 383
Panel 2	319839	75.1	49.8 - 103
Panel 3	319839	30.1	19.9 - 39.3
Panel 4	319839	9.74	5.82 - 12.6
Panel 5	319839	3.60	2.12 - 4.95

TNF α

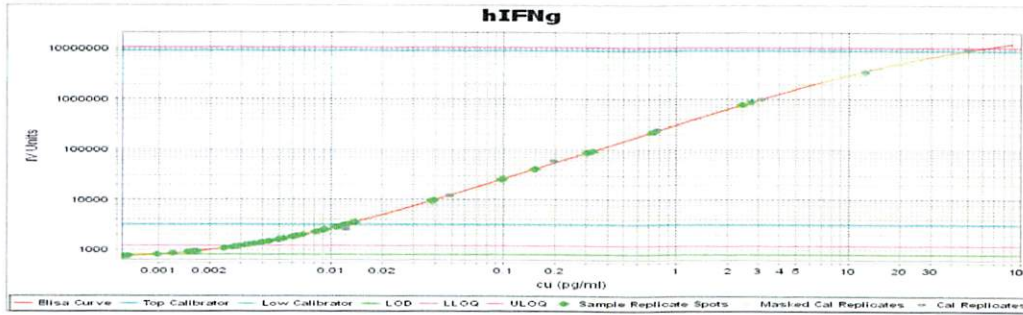
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	62.1	37.9 - 99.8
Panel 2	319839	19.4	11.8 - 26.5
Panel 3	319839	6.78	4.09 - 10.7
Panel 4	319839	1.61	1.01 - 2.84
Panel 5	319839	0.505	0.369 - 1.09

IL-10

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	50.8	30.4 - 61.9
Panel 2	319839	14.6	9.02 - 17.8
Panel 3	319839	5.84	3.19 - 7.61
Panel 4	319839	1.75	1.01 - 2.34
Panel 5	319839	0.718	0.409 - 1.02

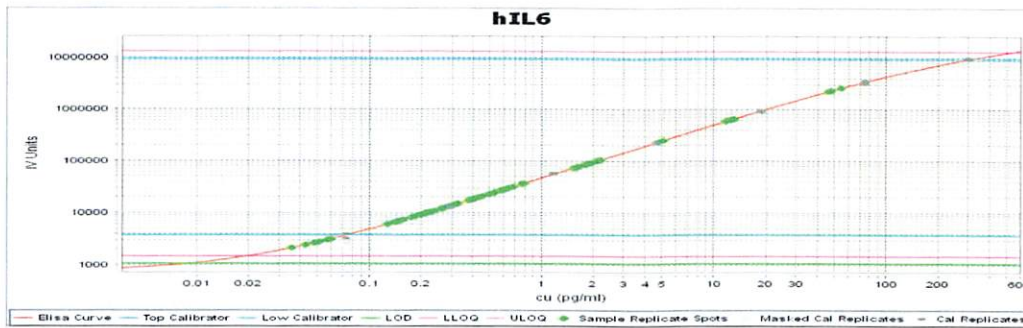
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

IFN γ Calibrator Curve



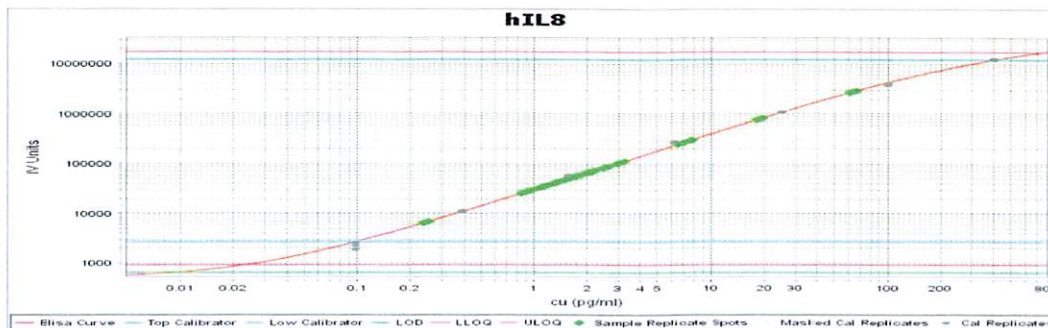
Calibrator Levels (pg/mL)	
A	50.0
B	12.5
C	3.13
D	0.781
E	0.195
F	0.049
G	0.012
H	0.000

IL-6 Calibrator Curve



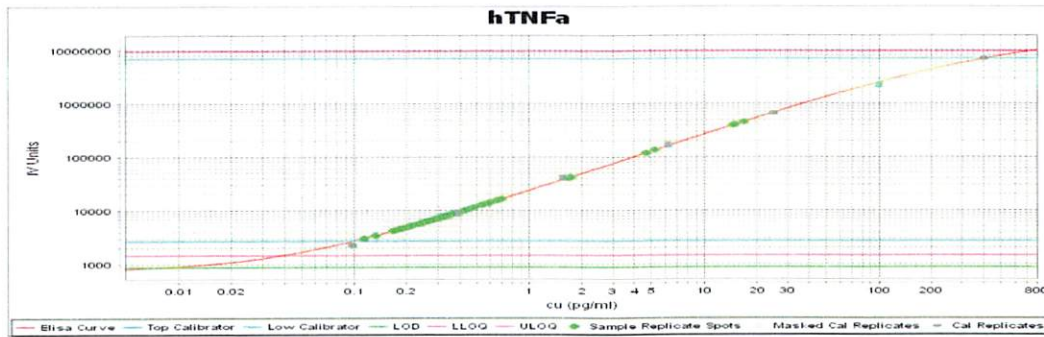
Calibrator Levels (pg/mL)	
A	300
B	75.0
C	18.8
D	4.69
E	1.17
F	0.293
G	0.073
H	0.000

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

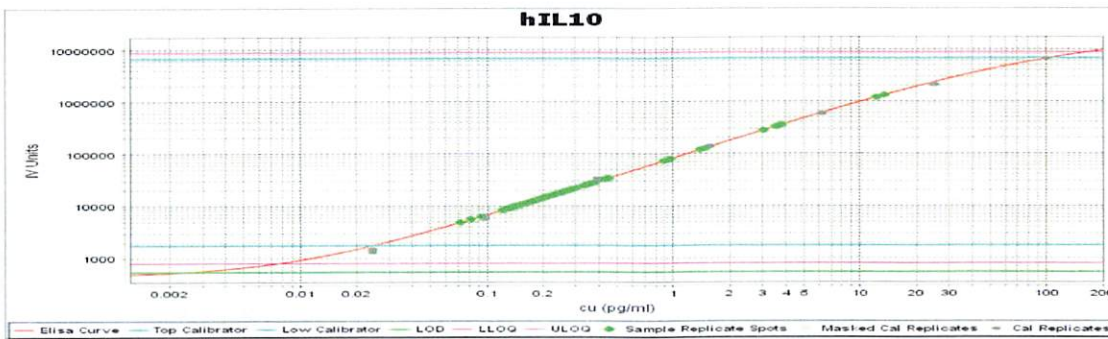
TNF α Calibrator Curve



**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