

Product Number:	85-0353
Lot Number:	504193
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-1 β Calibrator	18910040	319165
IL-6 Calibrator	18910044	324303
IL-8 Calibrator	1858454	319436
TNF α 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



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 85-0353 116-6A2-1-AB (KI-0251).

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title

[Signature] 28 Mar 2024
Signature/Date

Release Materials*

IL-1 β

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	12.3	9.55 - 20.6
Panel 2	319839	3.42	2.82 - 5.78
Panel 3	319839	1.31	1.07 - 2.22
Panel 4	319839	0.372	0.237 - 0.723
Panel 5	319839	0.114	0.081 - 0.271

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

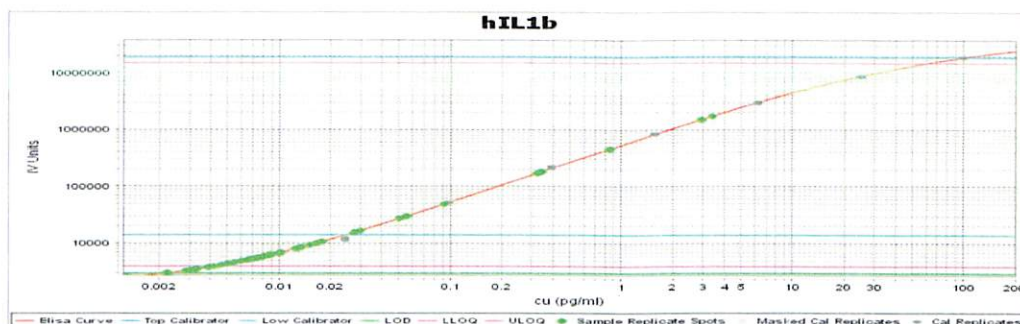
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	62.1	37.9 - 99.8
Panel 2	319839	19.4	11.8 - 26.5
Panel 3	319839	6.78	4.09 - 10.7
Panel 4	319839	1.61	1.01 - 2.84
Panel 5	319839	0.505	0.369 - 1.09

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.

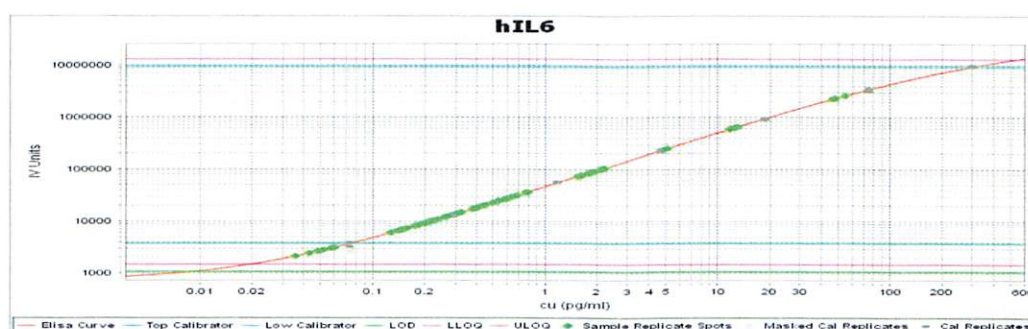
IL-1 β Calibrator Curve



**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

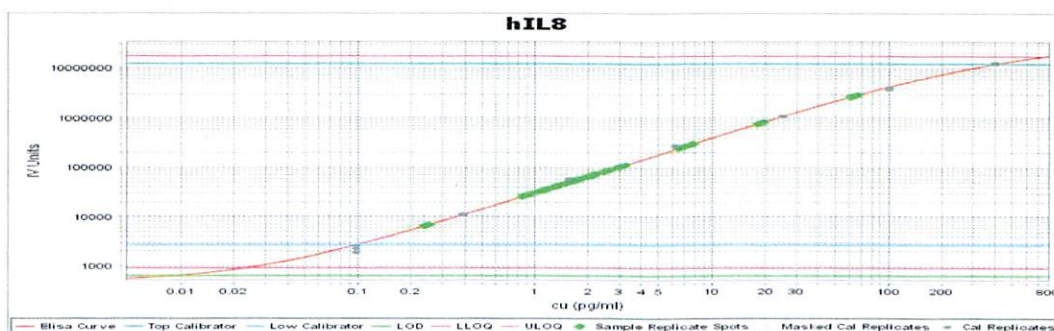
IL-6 Calibrator Curve



**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

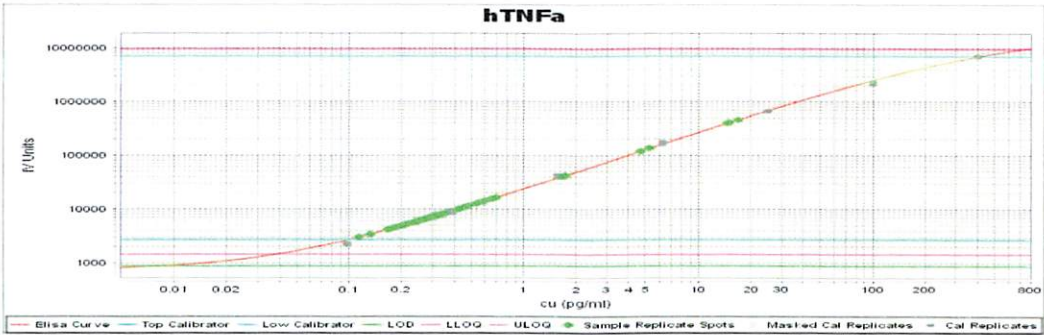
IL-8 Calibrator Curve



**Calibrator Levels
(pg/mL)**

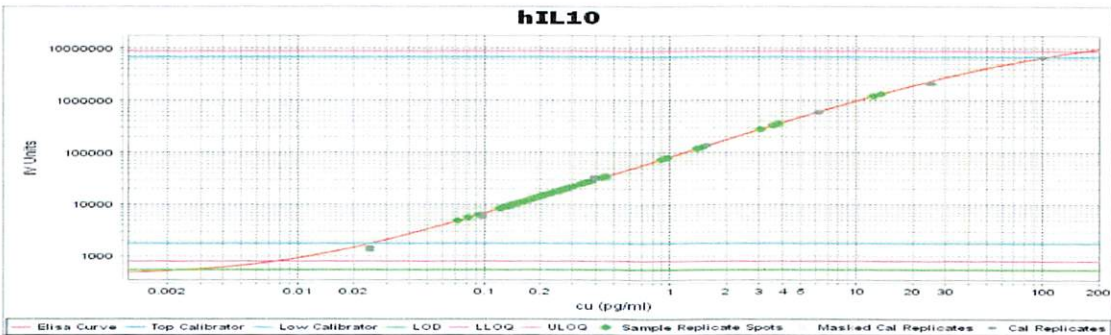
A	400
B	100
C	25.0
D	6.25
E	1.56
F	0.391
G	0.098
H	0.000

TNFα Calibrator Curve



Calibrator Levels (pg/mL)	
A	400
B	100
C	25.0
D	6.25
E	1.562
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