

Product Number:	85-0356
Lot Number:	504194
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
IFN γ Calibrator	1858260	319843
TNF α Calibrator	1880012	319460
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 KI-0260 Simoa Corplex Cytokine Panel 1 3 Plex Instruction Book 85-0356.



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

Review/Approval

Kathleen Cornell

Name

Quality Control
Supervisor

Title

Kathleen Cornell 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

IFN γ

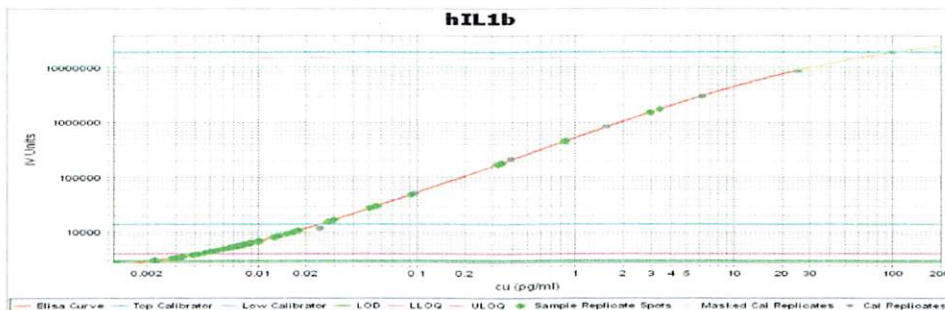
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	10.1	5.18 - 14.5
Panel 2	319839	2.99	1.55 - 4.66
Panel 3	319839	1.26	0.710 - 1.83
Panel 4	319839	0.393	0.217 - 0.604
Panel 5	319839	0.154	0.089 - 0.283

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

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

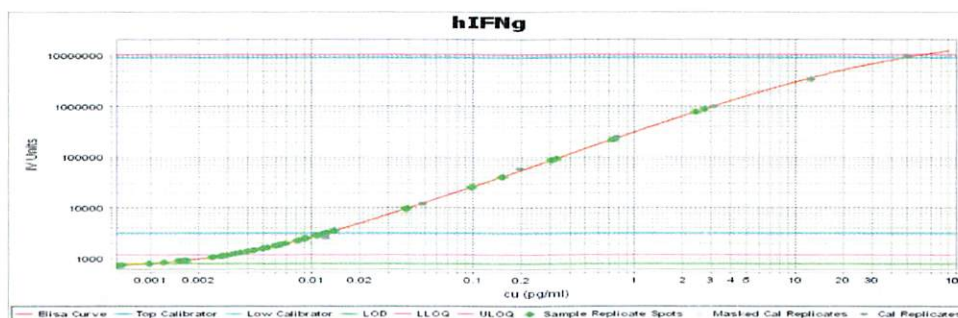
IL-1 β Calibrator Curve



**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

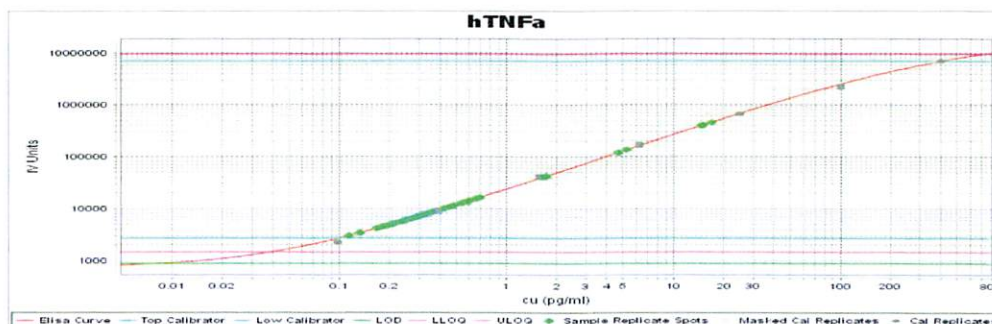
IFN γ Calibrator Curve



**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

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