

Product Number:	85-0339
Lot Number:	503981
Expiration:	26-Dec-2024
Platform(s):	SP-X



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

CorPlex Cytokine Plate	97-0329	325105P
Biotinylated Antibody Reagent	98-0329	310302B
Sample Diluent 21	1863329	319785
IL-6 Calibrator	18910044	319173
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

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 the Corplex Cytokine Panel Kit Instructions (KI-0229).

Release Materials*			
IL-6			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	211.1	141.9 - 312.1
Panel 2	319839	53.67	39.07 - 73.71
Panel 3	319839	20.55	14.61 - 30.61
Panel 4	319839	7.080	4.493 - 11.67
Panel 5	319839	2.550	1.407 - 4.765
TNFα			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	64.79	37.50 - 105.1
Panel 2	319839	18.49	12.23 - 27.53
Panel 3	319839	6.900	4.430 - 11.09
Panel 4	319839	1.650	1.213 - 2.913
Panel 5	319839	0.544	0.543 - 1.056

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

Review/Approval

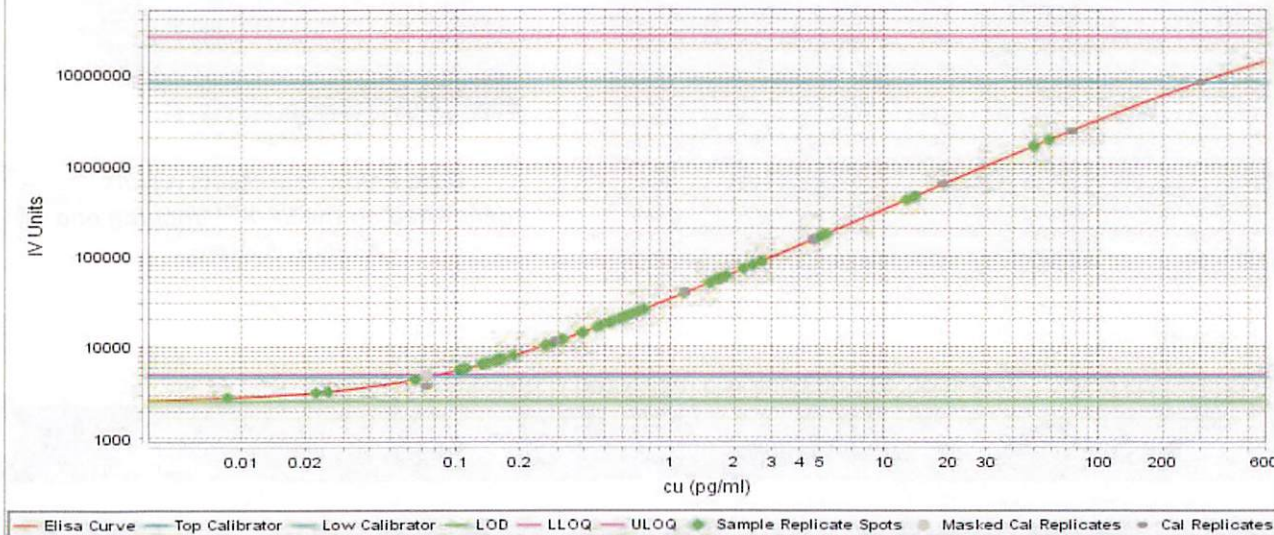
Kathleen Cornell
Name

Quality Control
Supervisor
Title

[Signature] 15 Sep 2023
Signature/Date

IL-6 Calibrator Curve

hIL6

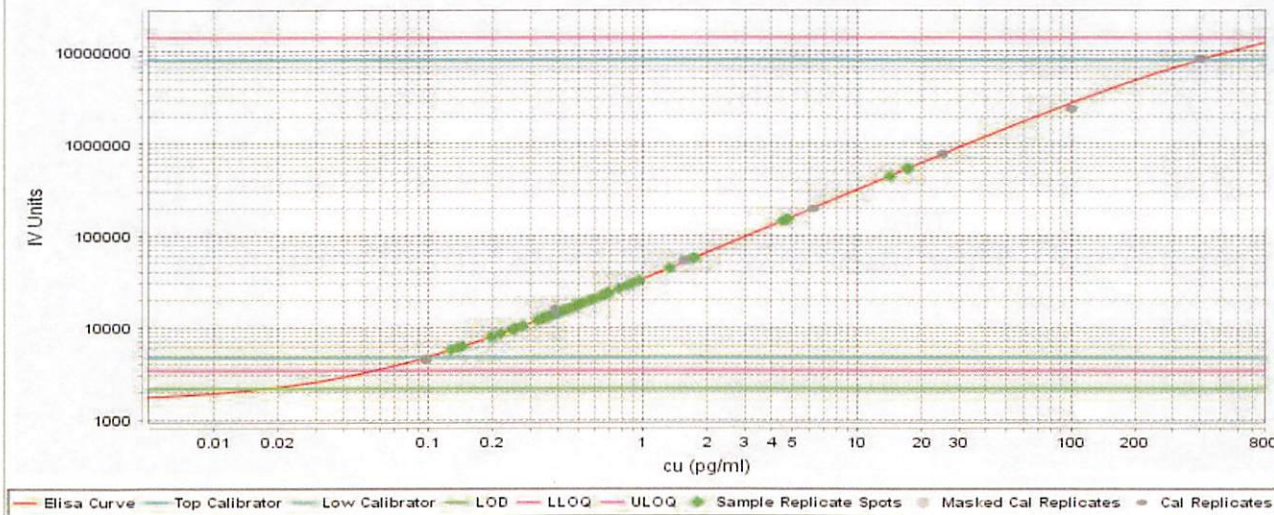


**Calibrator Levels
(pg/mL)**

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

TNFα Calibrator Curve

hTNFa



**Calibrator Levels
(pg/mL)**

A	400.00
B	100.00
C	25.00
D	6.25
E	1.562
F	0.391
G	0.098
H	0.000