

Product Number:	85-0339
Lot Number:	504491
Expiration:	26-Mar-2026
Platform(s):	SP-X



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

CorPlex Cytokine Plate	97-0329	433902P
Biotinylated Antibody Reagent	98-0329	427108B
Sample Diluent 21	1863329	405709
IL-6 Calibrator	18910044	324303
TNFα Calibrator	1880012	403812
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 KI-0229 Complex Cytokine Panel 1 2-Plex Kit 85-0339.

Release Materials*			
IL-6			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	204	148 - 298
Panel 2	319839	54.4	40.5 - 70.0
Panel 3	319839	21.2	14.6 - 29.2
Panel 4	319839	7.81	4.31 - 11.1
Panel 5	319839	3.18	1.40 - 4.47
TNFα			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	80.4	35.7 - 99.4
Panel 2	319839	24.6	12.2 - 26.3
Panel 3	319839	9.75	4.26 - 10.6
Panel 4	319839	2.55	1.05 - 2.81
Panel 5	319839	1.10	0.378 - 1.07

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

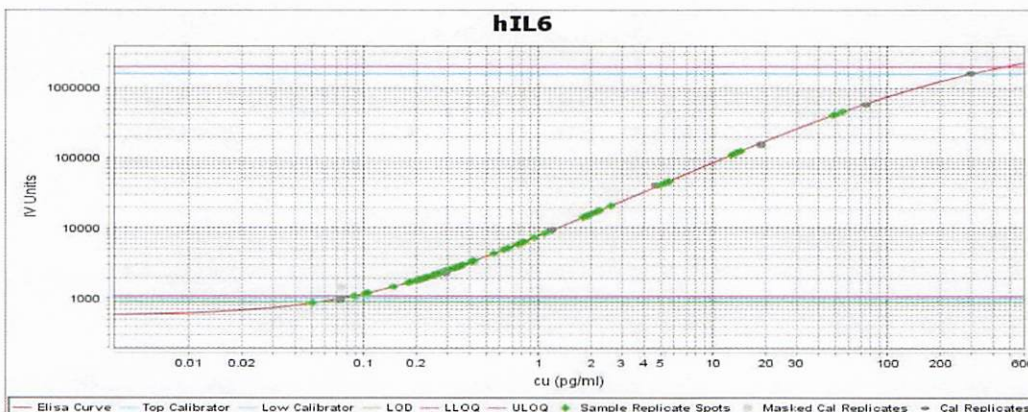
Review/Approval

Kyle Giroux
Name

21 Jan 2025
Senior S Document Control Specialist, QA
Title

[Signature] 21 Jan 2025
Signature/Date

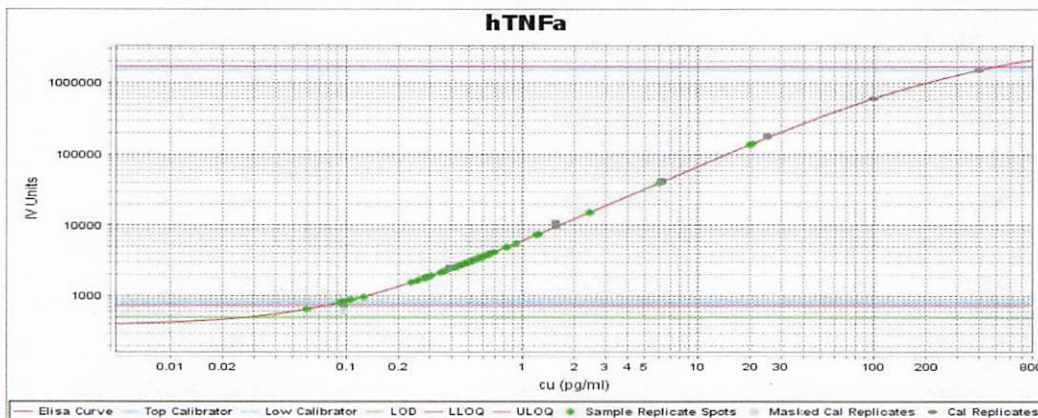
IL-6 Calibrator Curve



**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

TNFα 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