

Product Number:	85-0353
Lot Number:	504085
Expiration:	13-Jan-2025
Platform(s):	SP-X



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

CorPlex Cytokine Plate	97-0329	325510P
Biotinylated Antibody Reagent	98-0329	332509B
Sample Diluent 21	1863329	319974
IL-1β Calibrator	18910040	319165
IL-6 Calibrator	18910044	319173
IL-8 Calibrator	1858454	319436
TNFα Calibrator	1880012	319247
IL-10 Calibrator	18910043	318755
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 Simoa Complex Cytokine Panel 1

Instruction Book 85-0353 (KI-0251)

Review/Approval

Kathleen Cornell

Quality Control
Supervisor

Name

Title

[Handwritten Signature] 11 DEC 2023

Signature/Date

Release Materials*

IL-1 β

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	11.97	9.553 - 20.59
Panel 2	319839	3.858	2.823 - 5.776
Panel 3	319839	1.490	1.072 - 2.216
Panel 4	319839	0.417	0.237 - 0.723
Panel 5	319839	0.137	0.081 - 0.271

IL-6

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	201.5	149.3 - 299.3
Panel 2	319839	54.66	39.74 - 70.62
Panel 3	319839	22.57	14.23 - 29.41
Panel 4	319839	7.359	4.151 - 11.20
Panel 5	319839	2.808	1.321 - 4.519

IL-8

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	293.8	182.6 - 382.7
Panel 2	319839	81.98	49.82 - 102.8
Panel 3	319839	30.56	19.87 - 39.34
Panel 4	319839	9.506	5.821 - 12.63
Panel 5	319839	3.412	2.122 - 4.948

TNF α

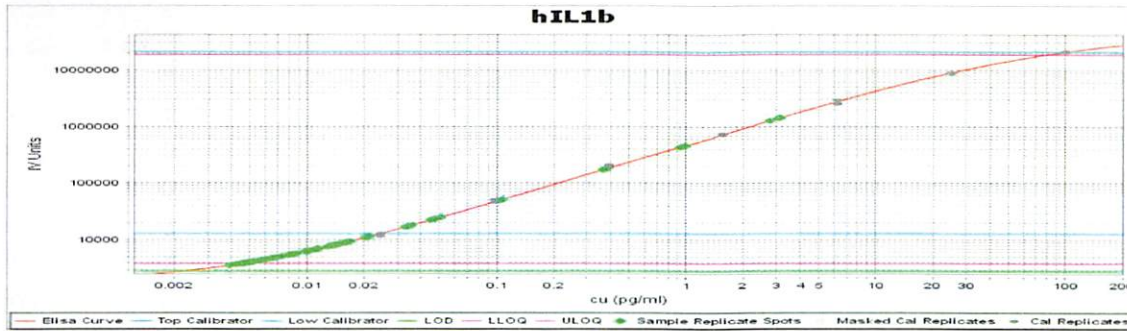
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	69.11	37.89 - 99.76
Panel 2	319839	21.92	11.84 - 26.54
Panel 3	319839	7.767	4.095 - 10.69
Panel 4	319839	1.545	1.007 - 2.838
Panel 5	319839	0.473	0.369 - 1.090

IL-10

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	49.54	30.41 - 61.92
Panel 2	319839	15.08	9.024 - 17.75
Panel 3	319839	6.048	3.193 - 7.609
Panel 4	319839	1.930	1.009 - 2.343
Panel 5	319839	0.765	0.409 - 1.017

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

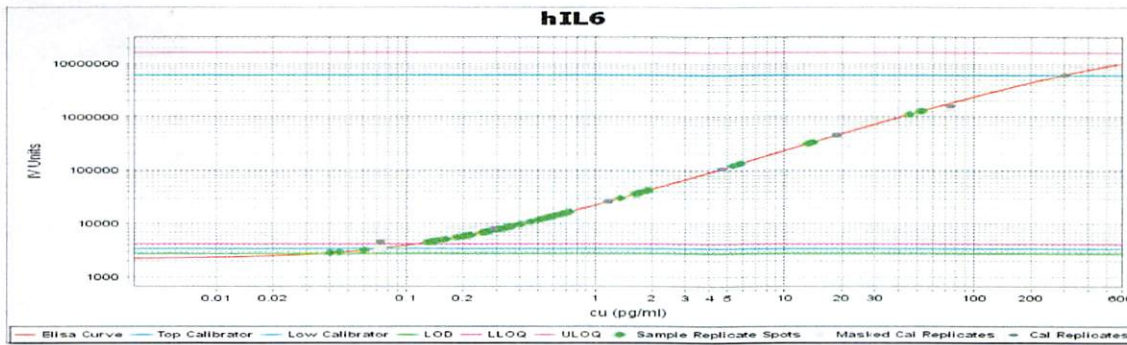
IL-1β Calibrator Curve



Calibrator Levels (pg/mL)

A	100.00
B	25.00
C	6.250
D	1.562
E	0.391
F	0.098
G	0.024
H	0.000

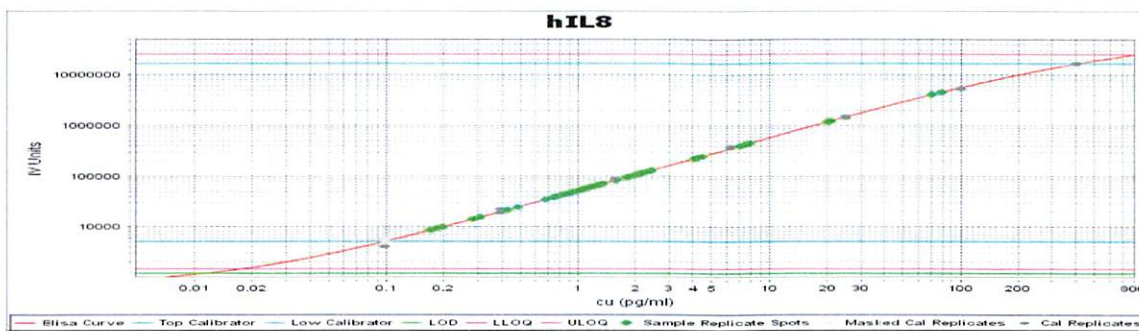
IL-6 Calibrator Curve



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

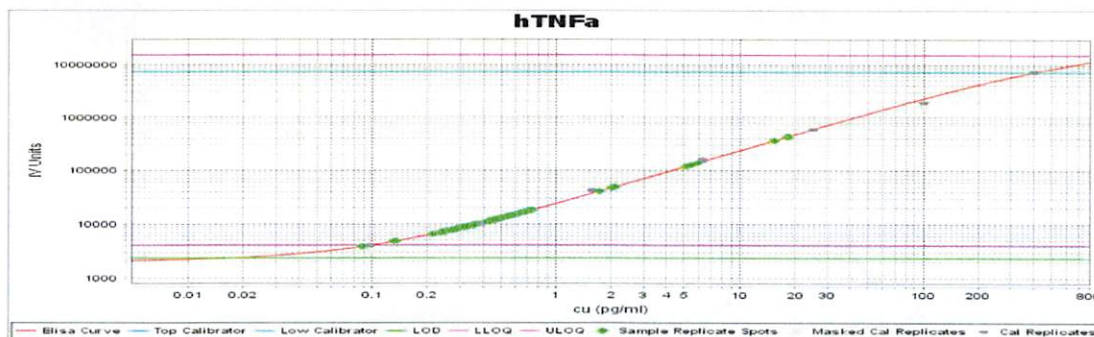
IL-8 Calibrator Curve



Calibrator Levels (pg/mL)

A	400.00
B	100.00
C	25.00
D	6.250
E	1.562
F	0.391
G	0.098
H	0.000

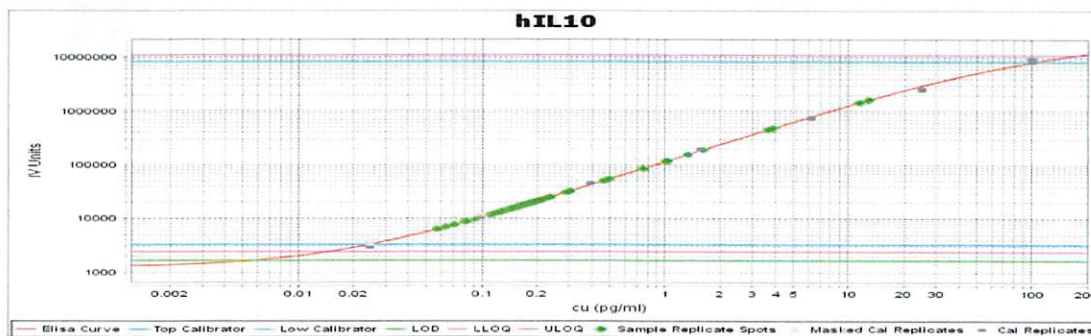
TNFα Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	400.00
B	100.00
C	25.00
D	6.250
E	1.562
F	0.391
G	0.098
H	0.000

IL-10 Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	100.00
B	25.00
C	6.250
D	1.562
E	0.391
F	0.098
G	0.024
H	0.000