

Product Number:	85-0412
Lot Number:	504073
Expiration:	13-Jan-2025
Platform(s):	SP-X

CorPlex Cytokine Plate	97-0329	325510P
Biotinylated Antibody Reagent	98-0329	332509B
Sample Diluent 21	1863329	319974
IL-6 Calibrator	18910044	319173
IL-8 Calibrator	1858454	319436
IL-22 Calibrator	18910036	318813
IL-10 Calibrator	18910043	318755
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



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 4-Plex Kit Instructions (KI-0220).

Review/Approval

Kathleen Cornell

Quality Control
Supervisor

Name

Title

Kathleen Cornell 01 Dec 2023
Signature/Date

Release Materials*

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

IL-22

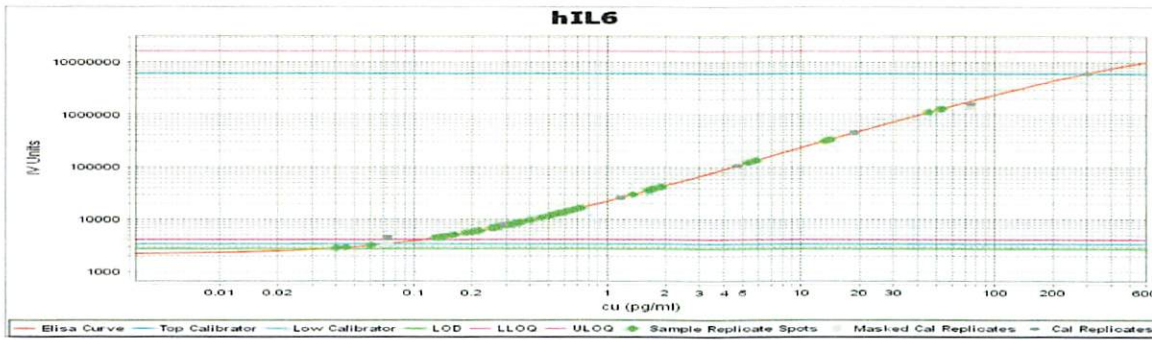
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	94.39	36.20 - 100.6
Panel 2	319839	13.33	6.046 - 14.95
Panel 3	319839	6.269	2.746 - 7.320
Panel 4	319839	1.140	0.565 - 1.486
Panel 5	319839	0.308	0.149 - 0.446

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

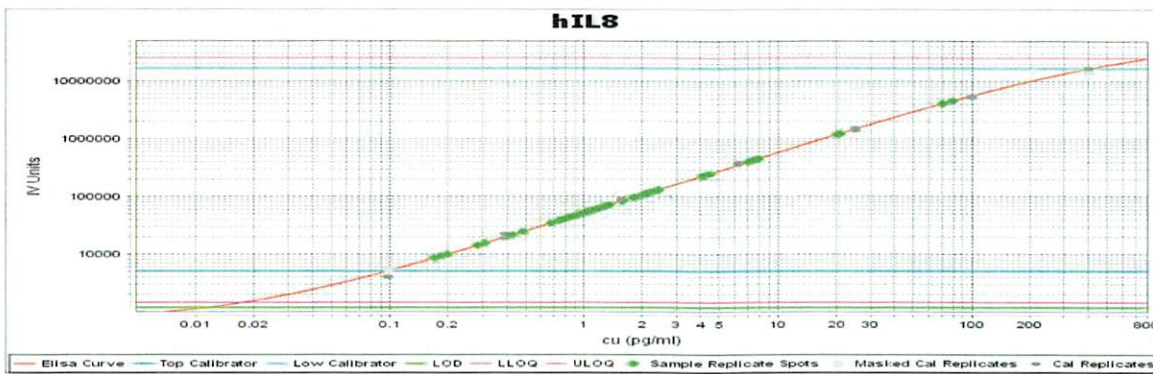
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

IL-6 Calibrator Curve



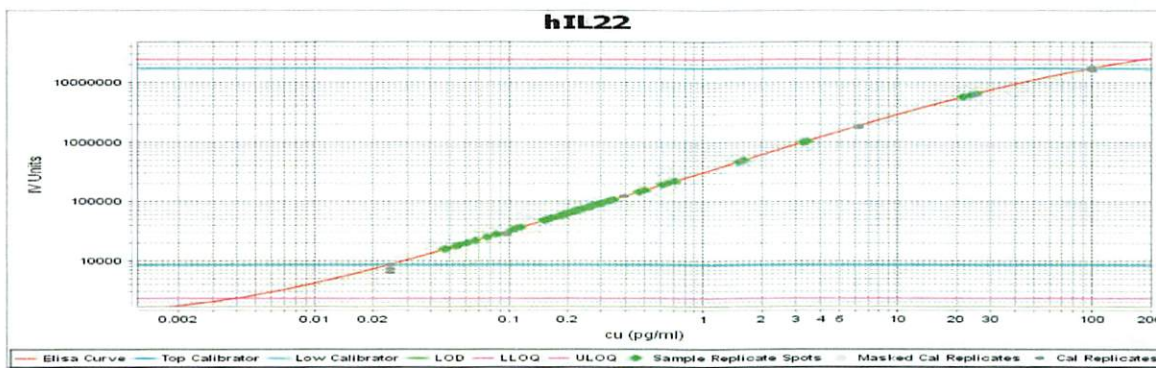
Calibrator Levels (pg/mL)	
A	300.00
B	75.00
C	18.75
D	4.688
E	1.172
F	0.293
G	0.073
H	0.000

IL-8 Calibrator Curve



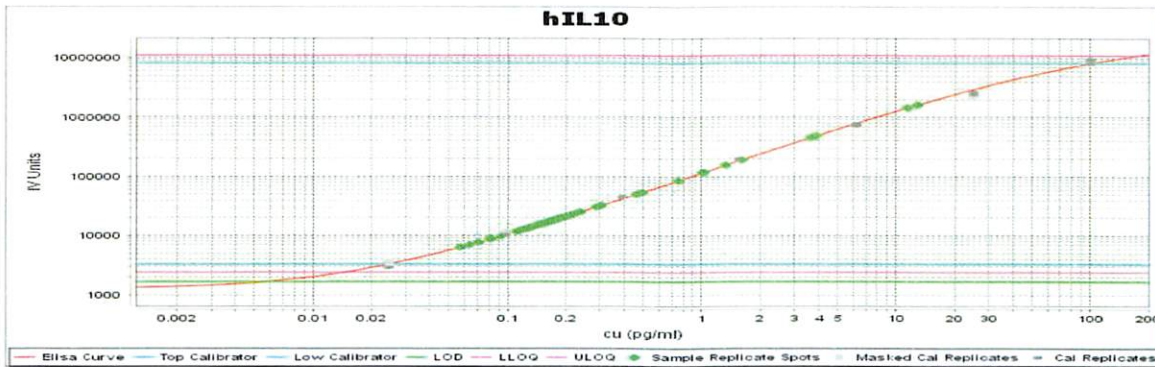
Calibrator Levels (pg/mL)	
A	400.00
B	100.00
C	25.00
D	6.250
E	1.562
F	0.391
G	0.098
H	0.000

IL-22 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-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