

Product Number:	85-0472
Lot Number:	503962
Expiration:	19-Feb-2024
Platform(s):	SP-X

CorPlex Cytokine Plate	97-0329	319986P
Biotinylated Antibody Reagent	98-0329	319352B
Sample Diluent 21	1863329	319785
IL-12p70 Calibrator	18910034	318571
IL-1β Calibrator	18910040	318673
IFNγ Calibrator	1858260	318886
IL-6 Calibrator	18910044	319173
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



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 the Corplex Cytokine Panel Kit Instructions (KI-0195).

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title

[Handwritten Signature]
Signature/Date
22 Aug 2023

Release Materials*

IL12p70

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	167.2	101.2 - 193.3
Panel 2	319839	46.67	32.24 - 52.86
Panel 3	319839	19.02	12.45 - 20.78
Panel 4	319839	5.041	2.632 - 6.594
Panel 5	319839	1.768	0.764 - 2.272

IL-1β

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	19.25	10.37 - 21.04
Panel 2	319839	5.504	3.172 - 5.834
Panel 3	319839	2.217	1.201 - 2.331
Panel 4	319839	0.640	0.279 - 0.773
Panel 5	319839	0.245	0.096 - 0.301

IFNγ

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	8.842	5.622 - 14.47
Panel 2	319839	2.547	1.692 - 4.770
Panel 3	319839	1.096	0.731 - 2.000
Panel 4	319839	0.339	0.215 - 0.692
Panel 5	319839	0.171	0.071 - 0.348

IL-6

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	213.1	141.9 - 312.1
Panel 2	319839	53.36	39.07 - 73.71
Panel 3	319839	21.61	14.61 - 30.61
Panel 4	319839	7.383	4.493 - 11.67
Panel 5	319839	3.124	1.407 - 4.765

TNFα

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	74.26	37.50 - 105.1
Panel 2	319839	20.70	12.23 - 27.53
Panel 3	319839	7.756	4.430 - 11.09
Panel 4	319839	1.930	1.213 - 2.913
Panel 5	319839	0.772	0.543 - 1.056

IL-10

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	54.53	28.83 - 63.34
Panel 2	319839	14.45	9.000 - 18.28
Panel 3	319839	5.760	3.397 - 7.891
Panel 4	319839	1.747	1.141 - 2.386
Panel 5	319839	0.791	0.457 - 1.051

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

IL-12p70 Calibrator Curve

hIL12p70

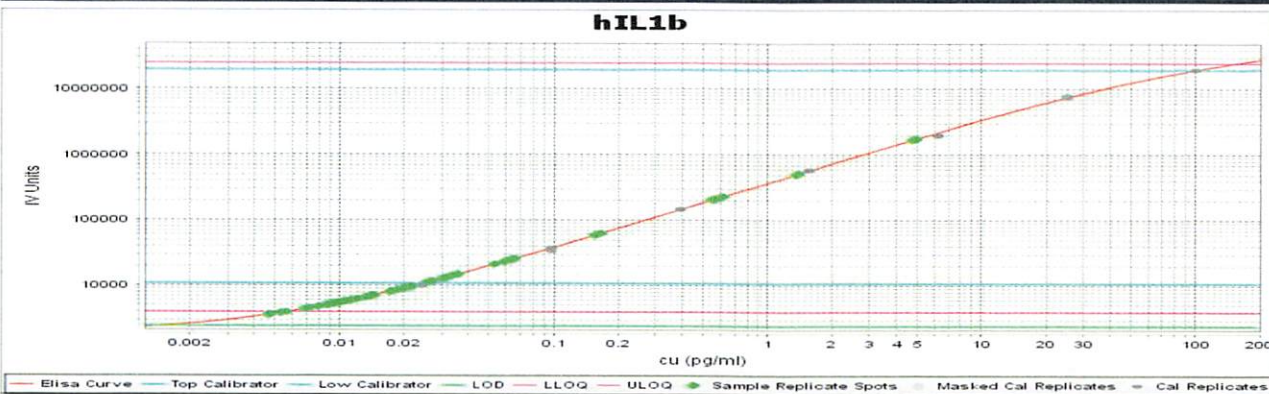


Calibrator Levels (pg/mL)

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-1β Calibrator Curve

hIL1b

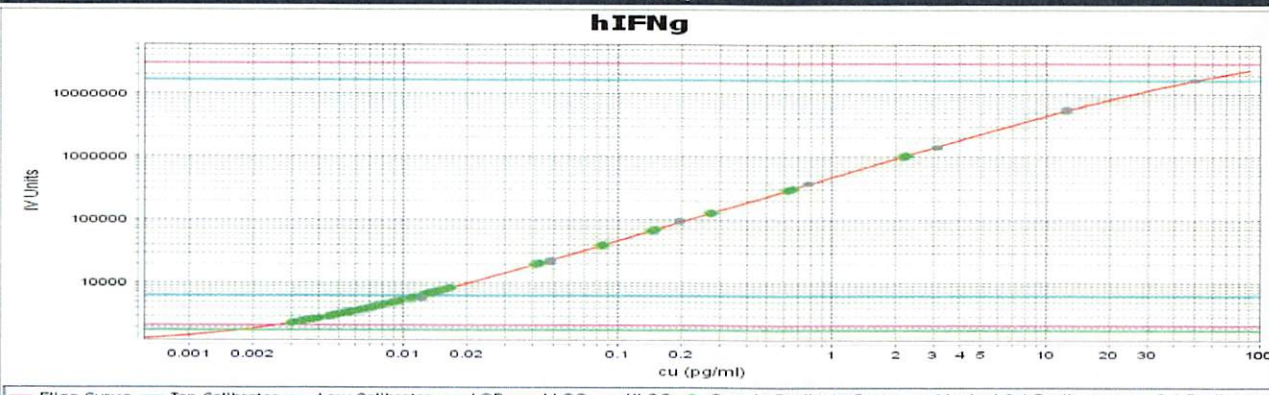


Calibrator Levels (pg/mL)

A	100.00
B	25.00
C	6.25
D	1.562
E	0.391
F	0.098
G	0.024
H	0.000

IFNγ Calibrator Curve

hIFNγ

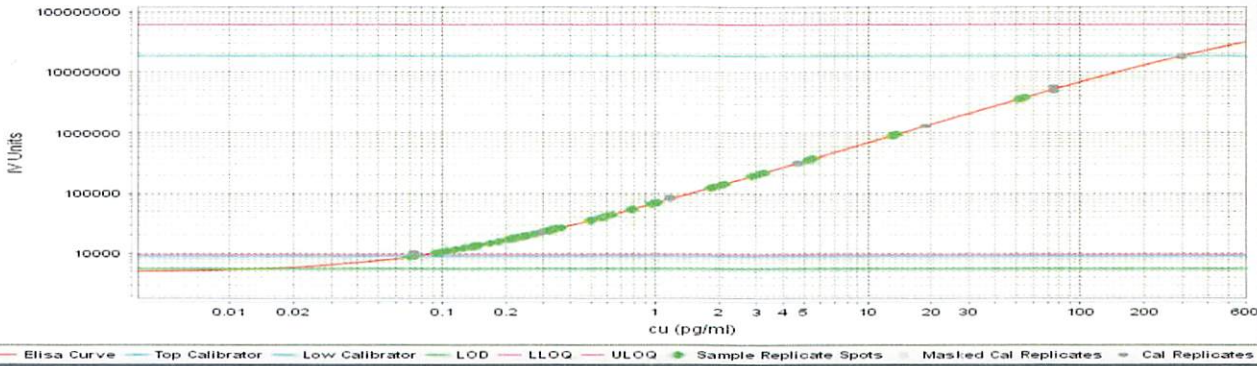


Calibrator Levels (pg/mL)

A	50.00
B	12.50
C	3.125
D	0.781
E	0.195
F	0.049
G	0.012
H	0.000

IL-6 Calibrator Curve

hIL6

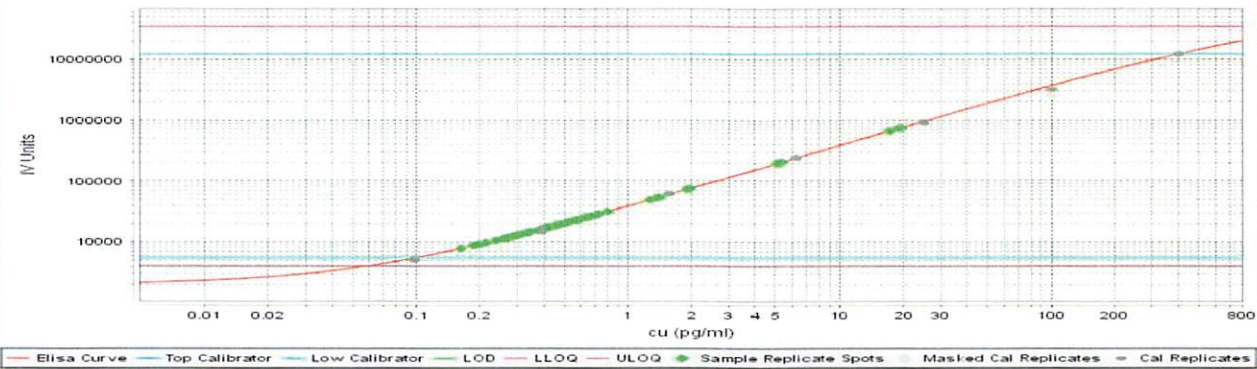


Calibrator Levels
(pg/mL)

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

IL-6 Calibrator Curve

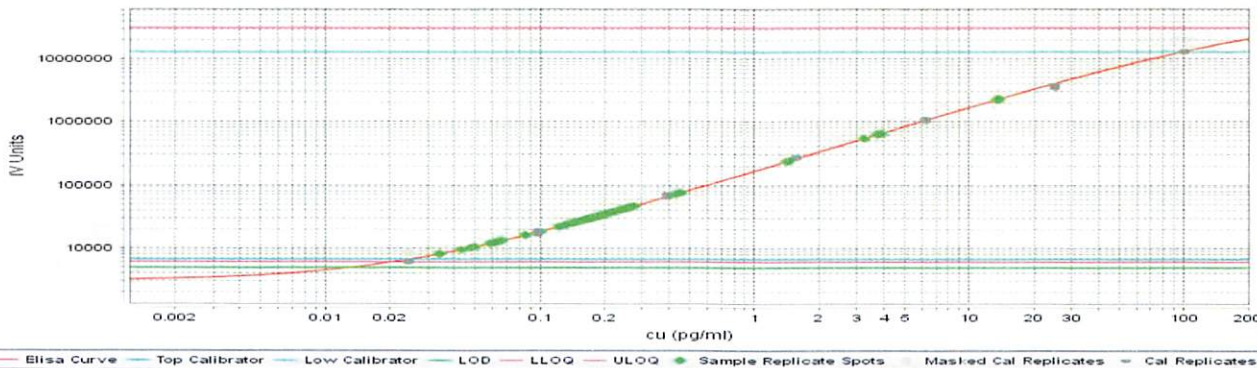
hTNFa



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

hIL10



Calibrator Levels
(pg/mL)

A	100.00
B	25.00
C	6.25
D	1.562
E	0.391
F	0.098
G	0.024
H	0.000