

<b>Product Number:</b>	85-0358
<b>Lot Number:</b>	503982
<b>Expiration:</b>	26-Dec-2024
<b>Platform(s):</b>	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-12p70 Calibrator	18910034	318756
IL-1β Calibrator	18910040	319165
IL-4 Calibrator	1858452	319016
IFNγ Calibrator	1858260	319437
IL-6 Calibrator	18910044	319173
IL-8 Calibrator	1858454	319436
TNFα Calibrator	1880012	319247
IL-10 Calibrator	18910043	319842
Wash Buffer (12.5X) <sup>1</sup>	1863332	N/A
Streptavidin-HRP Reagent <sup>1</sup>	1858263	N/A
Super Signal® Luminol <sup>1,2</sup>	1858089	N/A
Super Signal® Peroxide <sup>1,2</sup>	1858090	N/A

<sup>1</sup> Reagents are not Kit Lot Specific

<sup>2</sup> 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-0216).

Review/Approval

Kathleen Cornell  
Name

Quality Control  
Supervisor  
Title

*[Handwritten Signature]* 15 Sep 2023  
Signature/Date

**Release Materials\***

**IL12p70**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
Panel 1	319839	154.0	101.2 - 193.3
Panel 2	319839	44.19	32.24 - 52.86
Panel 3	319839	17.35	12.45 - 20.78
Panel 4	319839	4.340	2.632 - 6.594
Panel 5	319839	1.410	0.764 - 2.272

**IL-1β**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
Panel 1	319839	12.00	10.37 - 21.04
Panel 2	319839	3.470	3.172 - 5.834
Panel 3	319839	1.350	1.201 - 2.331
Panel 4	319839	0.420	0.279 - 0.773
Panel 5	319839	0.130	0.096 - 0.301

**IL-4**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
Panel 1	319839	68.31	45.70 - 99.67
Panel 2	319839	21.87	13.41 - 27.61
Panel 3	319839	8.110	5.612 - 11.10
Panel 4	319839	2.360	1.515 - 3.657
Panel 5	319839	0.860	0.559 - 1.400

**IFNγ**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
Panel 1	319839	8.920	5.622 - 14.47
Panel 2	319839	2.670	1.692 - 4.770
Panel 3	319839	1.130	0.731 - 2.000
Panel 4	319839	0.370	0.215 - 0.692
Panel 5	319839	0.170	0.071 - 0.348

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

**Release Materials\***

**IL-6**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
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

**IL-8**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
Panel 1	319839	305.6	164.0 - 399.7
Panel 2	319839	81.55	45.67 - 107.4
Panel 3	319839	31.28	18.96 - 41.09
Panel 4	319839	9.020	6.105 - 13.08
Panel 5	319839	3.400	2.260 - 5.125

**TNFα**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
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

**IL-10**

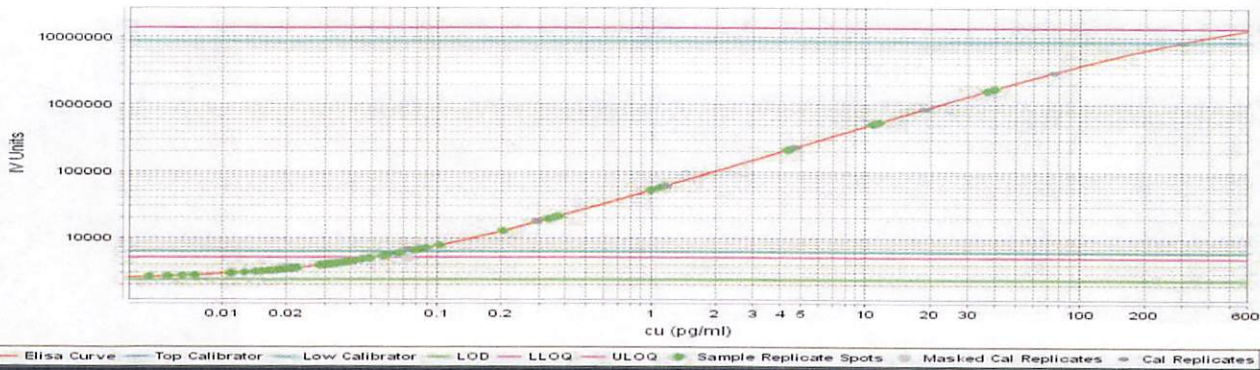
	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
Panel 1	319839	49.18	28.83 - 63.34
Panel 2	319839	12.72	9.000 - 18.28
Panel 3	319839	4.820	3.397 - 7.891
Panel 4	319839	1.460	1.141 - 2.386
Panel 5	319839	0.600	0.457 - 1.051

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



**IL-12p70 Calibrator Curve**

**hIL12p70**

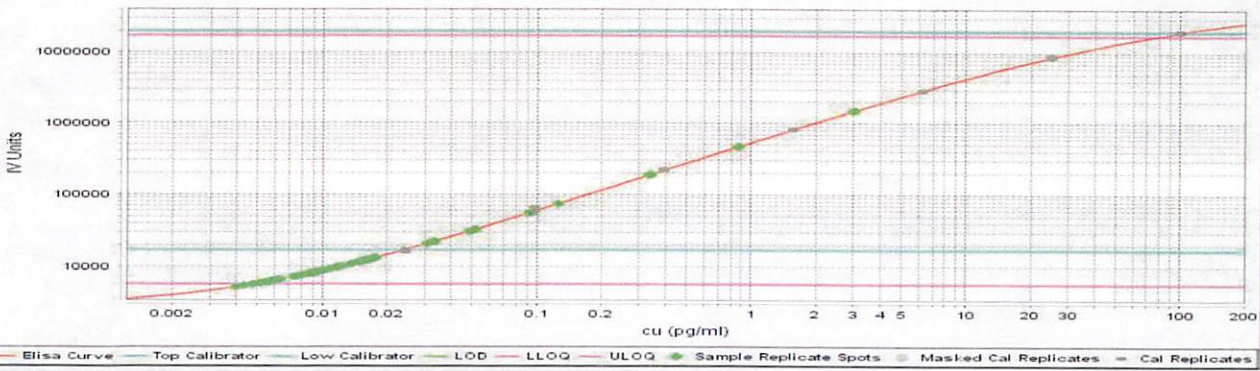


**Calibrator Levels  
(pg/mL)**

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**

**hIL1β**

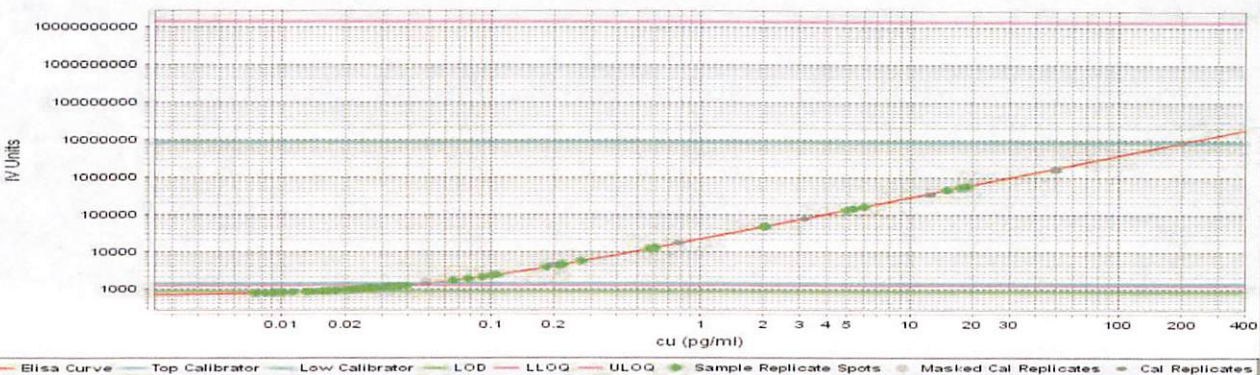


**Calibrator Levels  
(pg/mL)**

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

**IL-4 Calibrator Curve**

**hIL4**

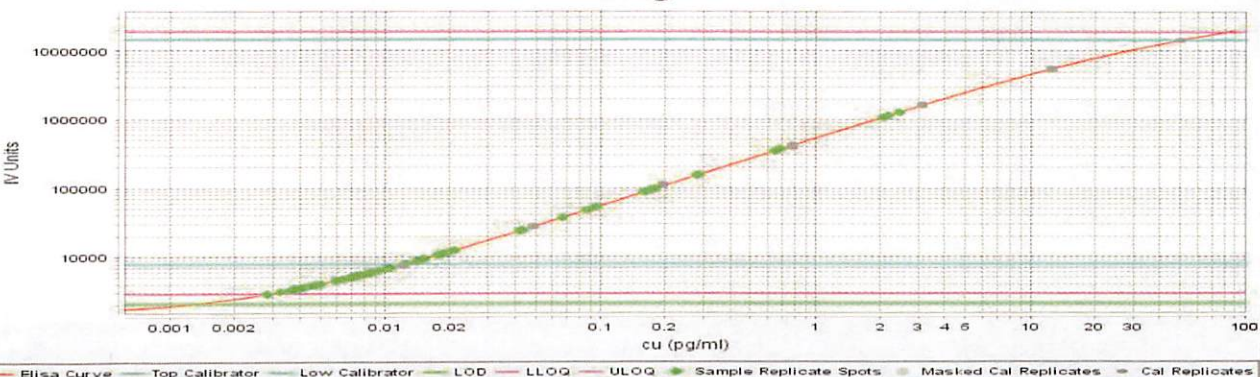


**Calibrator Levels  
(pg/mL)**

A	200.00
B	50.00
C	12.5
D	3.125
E	0.781
F	0.195
G	0.049
H	0.000

**IFNγ Calibrator Curve**

**hIFNγ**



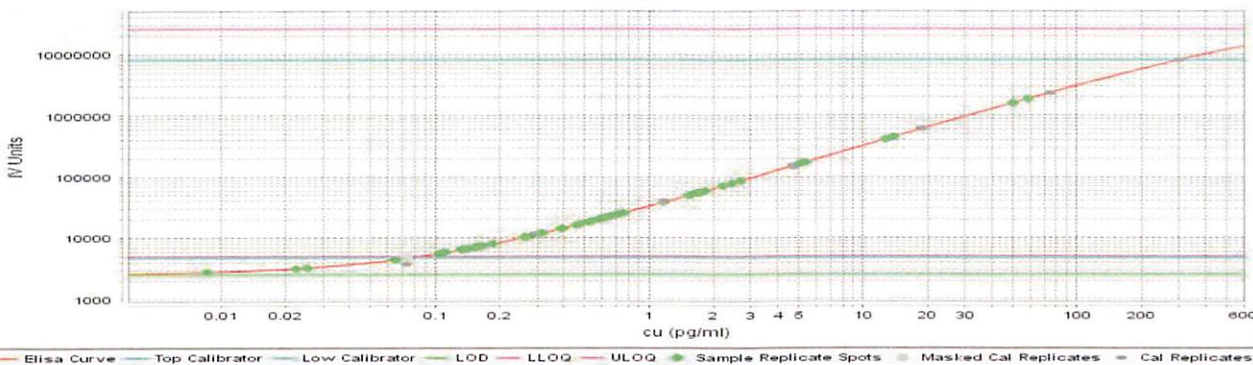
**Calibrator Levels  
(pg/mL)**

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

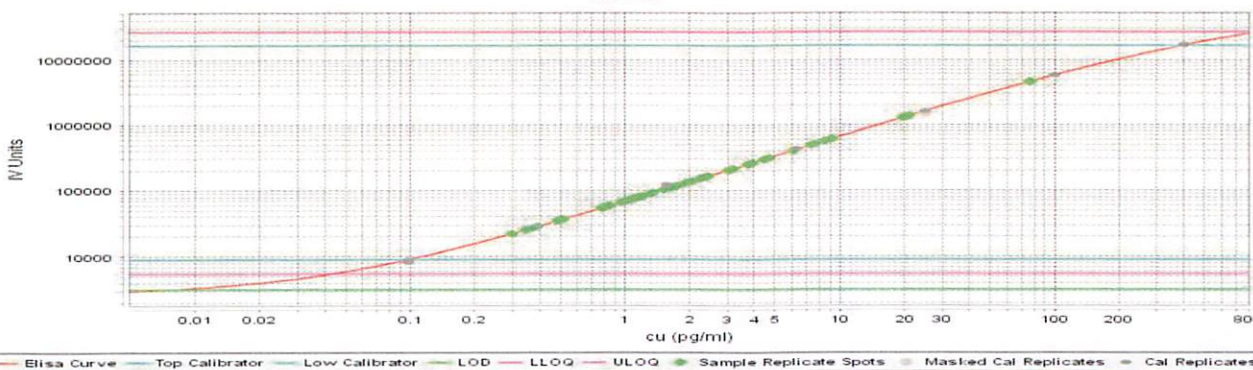


Calibrator Levels (pg/mL)

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-8 Calibrator Curve

hIL8

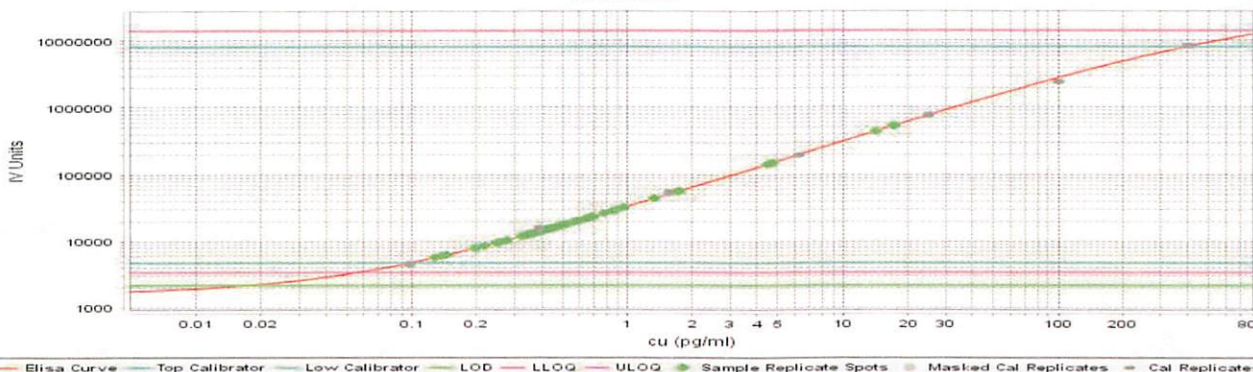


Calibrator Levels (pg/mL)

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

TNFα Calibrator Curve

hTNFα

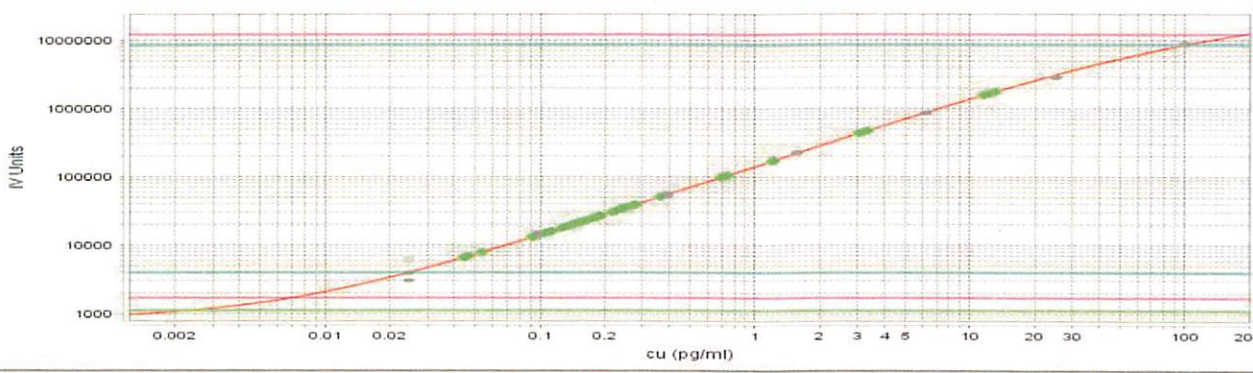


Calibrator Levels (pg/mL)

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

IL-10 Calibrator Curve

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