

<b>Product Number:</b>	85-0422
<b>Lot Number:</b>	504374
<b>Expiration:</b>	8-Jun-2025
<b>Platform(s):</b>	SP-X

Component	Part Number	Lot Number	Control Lot Number
CorPlex Cytokine Plate	97-0329	409510P	N/A
Biotinylated Antibody Reagent	98-0329	421304B	N/A
Sample Diluent 21	1863329	311103	311103A
IL-12p70 Calibrator	18910034	318756	N/A
IL-1β Calibrator	18910040	319564	N/A
IL-6 Calibrator	18910044	324303	N/A
IL-8 Calibrator	1858454	319436	N/A
TNFα Calibrator	1880012	403812	N/A
IL-10 Calibrator	18910043	319842	N/A
Wash Buffer (12.5X) <sup>1</sup>	1863332	N/A	N/A
Streptavidin-HRP Reagent <sup>1</sup>	1858263	N/A	N/A
Super Signal® Luminol <sup>1,2</sup>	1858089	N/A	N/A
Super Signal® Peroxide <sup>1,2</sup>	1858090	N/A	N/A



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

<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 Simoa Corplex Cytokine Panel 1 6-Plex Instruction Book 85-0422 116-6A3-1-AB (KI-0261)

Release Materials*			
IL-12p70			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	130	107 - 186
Panel 2	319839	44.3	33.0 - 51.2
Panel 3	319839	18.6	12.3 - 20.3
Panel 4	319839	5.24	2.59 - 6.31
Panel 5	319839	1.74	0.728 - 2.17

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

Review/Approval

*Kyle Giroux*

*Senior Document Control Specialist, QA*

*[Signature] 03Sep2024*

Name

Title

Signature/Date

**Release Materials\***

**IL-1 $\beta$**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
Panel 1	319839	14.5	9.55 - 20.6
Panel 2	319839	5.24	2.82 - 5.78
Panel 3	319839	2.11	1.07 - 2.22
Panel 4	319839	0.662	0.237 - 0.723
Panel 5	319839	0.222	0.081 - 0.271

**IL-6**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
Panel 1	319839	182	149 - 299
Panel 2	319839	55.4	39.7 - 70.6
Panel 3	319839	23.3	14.2 - 29.4
Panel 4	319839	8.89	4.15 - 11.2
Panel 5	319839	3.27	1.32 - 4.52

**IL-8**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
Panel 1	319839	319	183 - 383
Panel 2	319839	84.8	49.8 - 103
Panel 3	319839	34.4	19.9 - 39.3
Panel 4	319839	10.9	5.82 - 12.6
Panel 5	319839	4.24	2.12 - 4.95

**TNF $\alpha$**

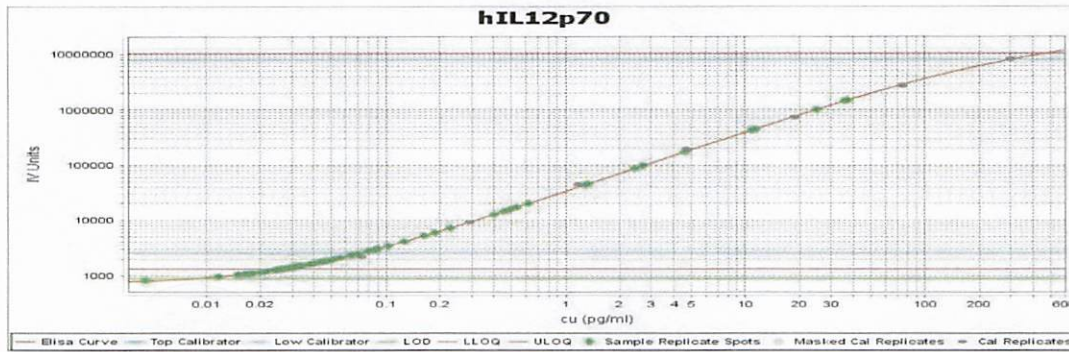
	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
Panel 1	319839	46.8	37.9 - 99.8
Panel 2	319839	20.8	11.8 - 26.5
Panel 3	319839	8.25	4.09 - 10.7
Panel 4	319839	2.11	1.01 - 2.84
Panel 5	319839	0.743	0.369 - 1.09

**IL-10**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
Panel 1	319839	33.6	30.4 - 61.9
Panel 2	319839	13.6	9.02 - 17.8
Panel 3	319839	5.87	3.19 - 7.61
Panel 4	319839	1.93	1.01 - 2.34
Panel 5	319839	0.794	0.409 - 1.02

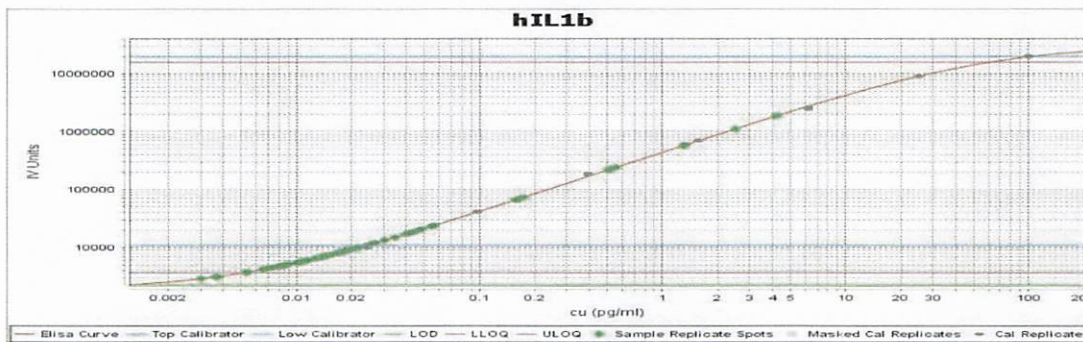
\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

**IL-12p70 Calibrator Curve**



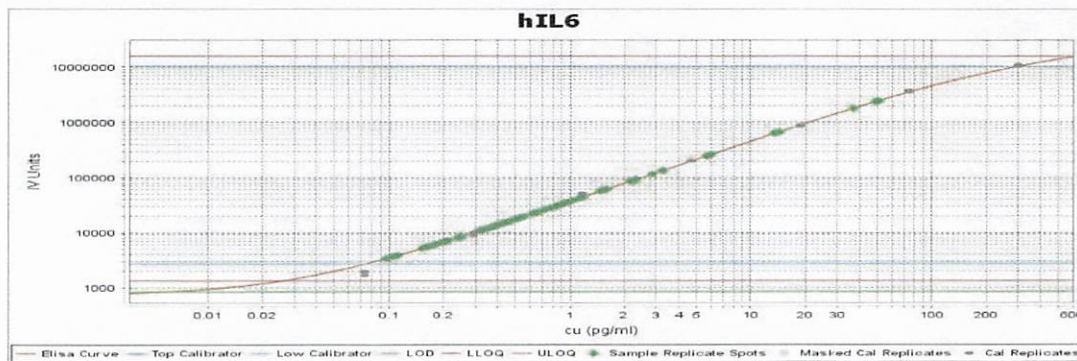
Calibrator Levels (pg/mL)	
A	300
B	75.0
C	18.8
D	4.69
E	1.17
F	0.293
G	0.073
H	0.000

**IL-1β Calibrator Curve**



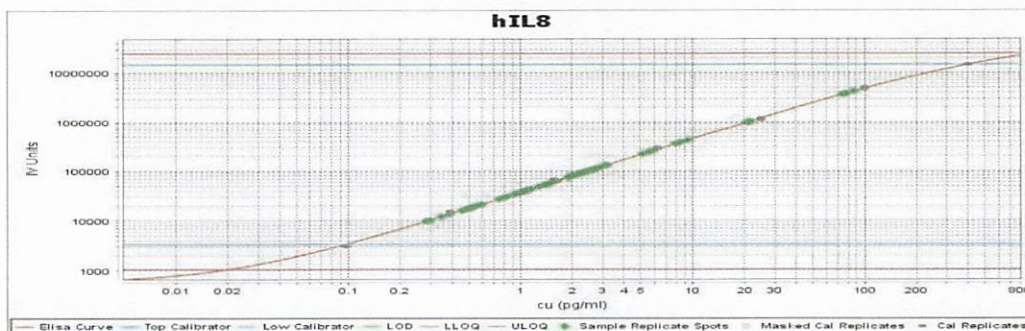
Calibrator Levels (pg/mL)	
A	100
B	25.0
C	6.25
D	1.56
E	0.391
F	0.098
G	0.024
H	0.000

**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

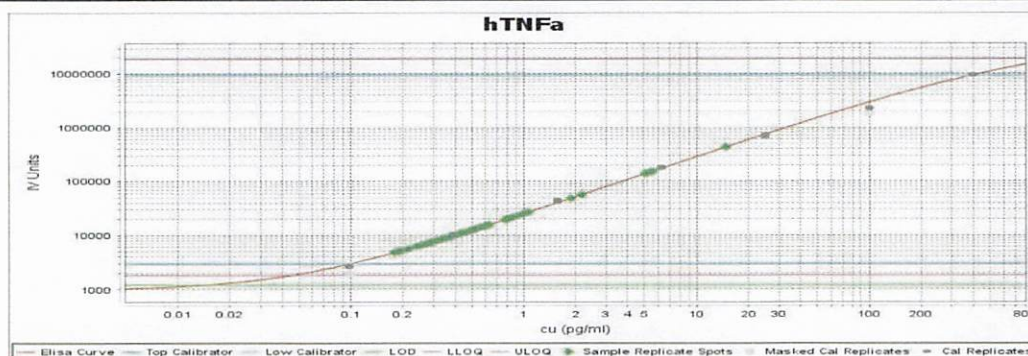
**IL-8 Calibrator Curve**



**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

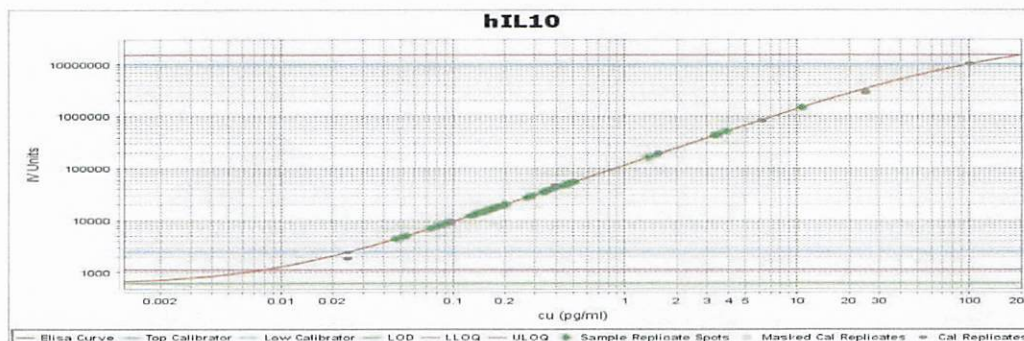
**TNFα Calibrator Curve**



**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

**IL-10 Calibrator Curve**



**Calibrator Levels (pg/mL)**

A	100
B	25.0
C	6.25
D	1.56
E	0.391
F	0.098
G	0.024
H	0.000