

<b>Product Number:</b>	85-0353
<b>Lot Number:</b>	504069
<b>Expiration:</b>	13-Jan-2025
<b>Platform(s):</b>	SP-X



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

<b>CorPlex Cytokine Plate</b>	97-0329	326103P
<b>Biotinylated Antibody Reagent</b>	98-0329	326105B
<b>Sample Diluent 21</b>	1863329	319974
<b>IL-1β Calibrator</b>	18910040	319165
<b>IL-6 Calibrator</b>	18910044	319173
<b>IL-8 Calibrator</b>	1858454	319436
<b>TNFα Calibrator</b>	1880012	319247
<b>IL-10 Calibrator</b>	18910043	318755
<b>Wash Buffer (12.5X)<sup>1</sup></b>	1863332	N/A
<b>Streptavidin-HRP Reagent<sup>1</sup></b>	1858263	N/A
<b>Super Signal® Luminol<sup>1,2</sup></b>	1858089	N/A
<b>Super Signal® Peroxide<sup>1,2</sup></b>	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 Simoa Corplex Cytokine Panel 1

Instruction Book 85-0353 (KI-0251)

Review/Approval	Kathleen Cornell	Quality Control Supervisor	
	Name	Title	Signature/Date

**Release Materials\***

**IL-1 $\beta$**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	12.53	9.553 - 20.59
Panel 2	319839	3.405	2.823 - 5.776
Panel 3	319839	1.374	1.072 - 2.216
Panel 4	319839	0.358	0.237 - 0.723
Panel 5	319839	0.125	0.081 - 0.271

**IL-6**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	218.6	149.3 - 299.3
Panel 2	319839	54.08	39.74 - 70.62
Panel 3	319839	22.11	14.23 - 29.41
Panel 4	319839	7.213	4.151 - 11.20
Panel 5	319839	2.875	1.321 - 4.519

**IL-8**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	277.7	182.6 - 382.7
Panel 2	319839	71.43	49.82 - 102.8
Panel 3	319839	27.62	19.87 - 39.34
Panel 4	319839	8.296	5.821 - 12.63
Panel 5	319839	3.022	2.122 - 4.948

**TNF $\alpha$**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	69.69	37.89 - 99.76
Panel 2	319839	18.52	11.84 - 26.54
Panel 3	319839	7.663	4.095 - 10.69
Panel 4	319839	1.679	1.007 - 2.838
Panel 5	319839	0.539	0.369 - 1.090

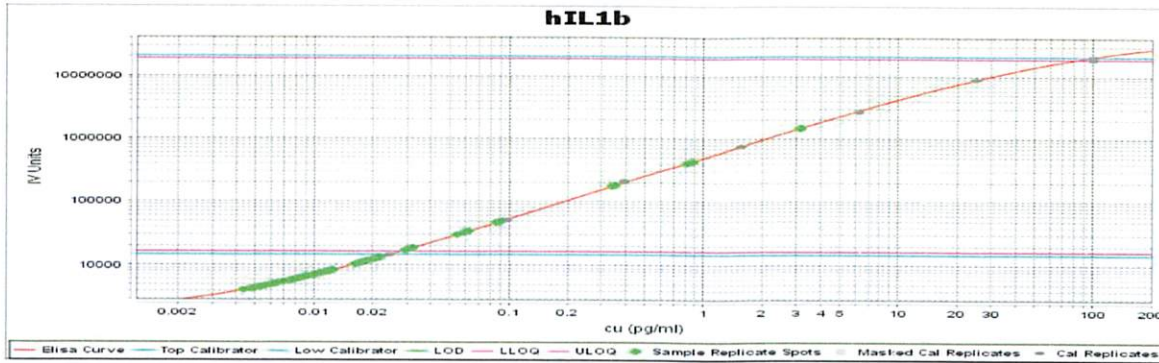
**IL-10**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	53.85	30.41 - 61.92
Panel 2	319839	14.27	9.024 - 17.75
Panel 3	319839	5.688	3.193 - 7.609
Panel 4	319839	1.690	1.009 - 2.343
Panel 5	319839	0.655	0.409 - 1.017

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



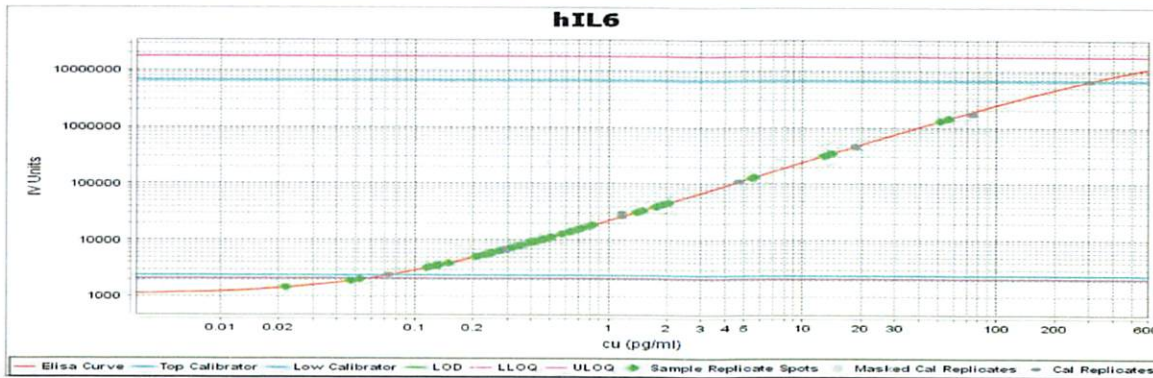
**IL-1 $\beta$  Calibrator Curve**



**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

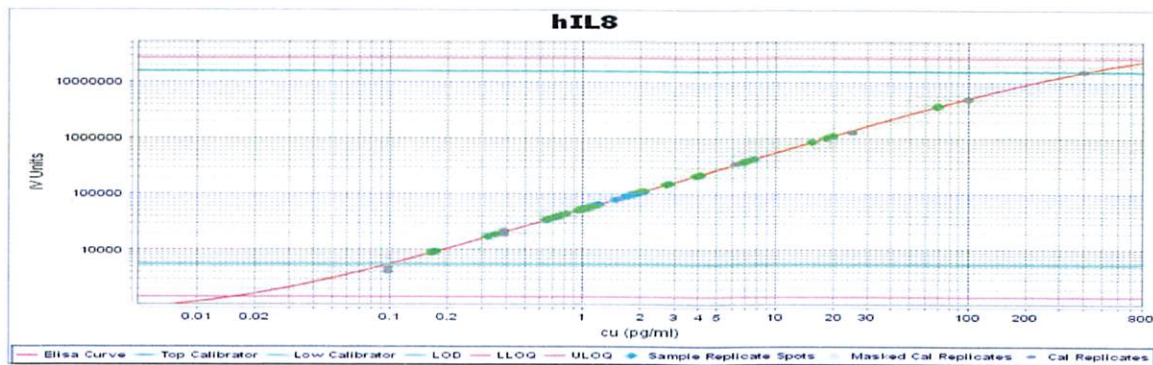
**IL-6 Calibrator Curve**



**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**

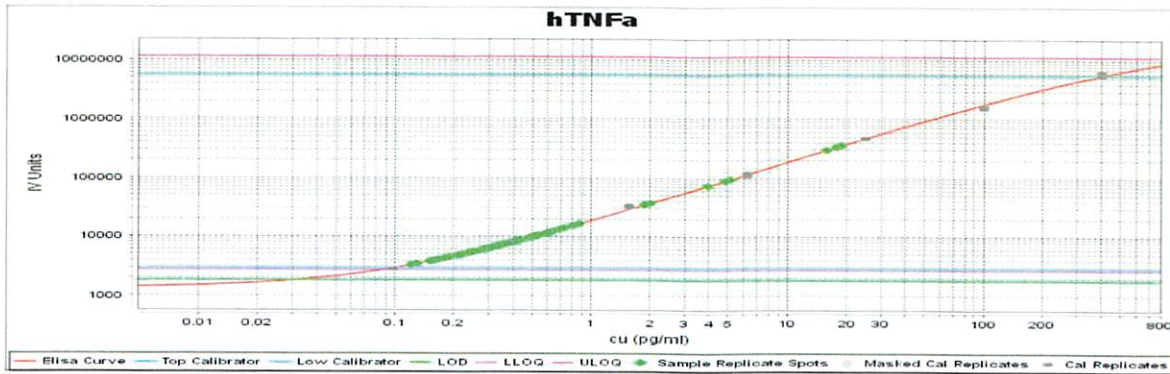


**Calibrator Levels (pg/mL)**

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

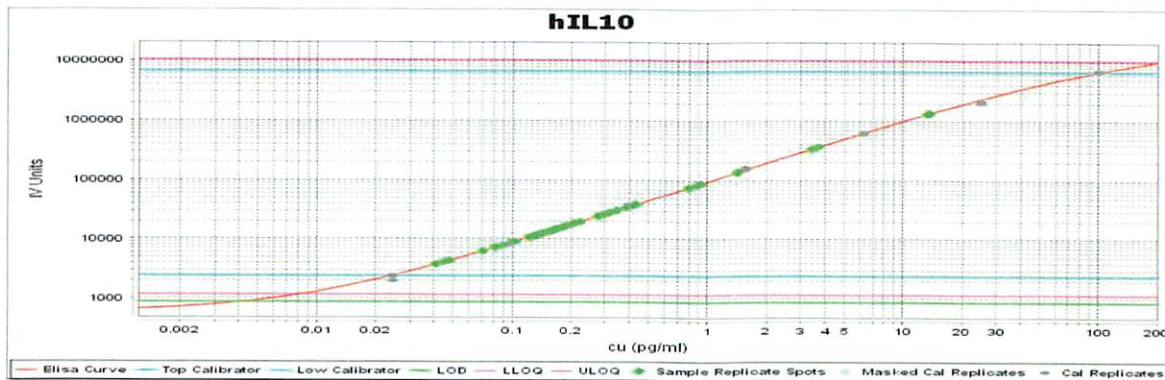


**TNFα Calibrator Curve**



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