

<b>Product Number:</b>	85-0412
<b>Lot Number:</b>	504245
<b>Expiration:</b>	18-Jun-2025
<b>Platform(s):</b>	SP-X

<b>CorPlex Cytokine Plate</b>	97-0329	411011P
<b>Biotinylated Antibody Reagent</b>	98-0329	409505B
<b>Sample Diluent 21</b>	1863329	319785
<b>IL-6 Calibrator</b>	18910044	319173
<b>IL-8 Calibrator</b>	1858454	319436
<b>IL-22 Calibrator</b>	18910036	319176
<b>IL-10 Calibrator</b>	18910043	319842
<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

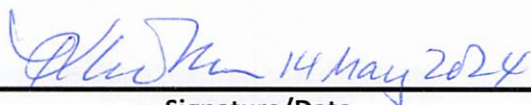


*Data below represents results generated on the SP-X Imaging and Analysis System.*

<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 Instruction Book 85-0353 116-6A2-1-AB (KI-0200).

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

**Release Materials\***

**IL-6**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	208	149 - 299
<b>Panel 2</b>	319839	55.7	39.7 - 70.6
<b>Panel 3</b>	319839	22.1	14.2 - 29.4
<b>Panel 4</b>	319839	7.76	4.15 - 11.2
<b>Panel 5</b>	319839	2.81	1.32 - 4.52

**IL-8**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	284	183 - 383
<b>Panel 2</b>	319839	75.7	49.8 - 103
<b>Panel 3</b>	319839	27.1	19.9 - 39.3
<b>Panel 4</b>	319839	8.47	5.82 - 12.6
<b>Panel 5</b>	319839	3.16	2.12 - 4.95

**IL-22**

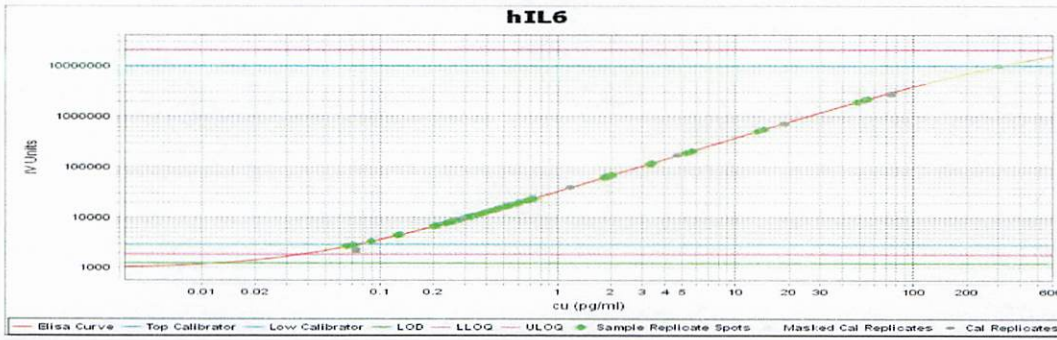
	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	86.8	36.2 - 101
<b>Panel 2</b>	319839	11.4	6.05 - 15.0
<b>Panel 3</b>	319839	6.41	2.75 - 7.32
<b>Panel 4</b>	319839	1.24	0.565 - 1.49
<b>Panel 5</b>	319839	0.322	0.149 - 0.446

**IL-10**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	43.8	30.4 - 61.9
<b>Panel 2</b>	319839	13.6	9.02 - 17.8
<b>Panel 3</b>	319839	5.45	3.19 - 7.61
<b>Panel 4</b>	319839	1.57	1.01 - 2.34
<b>Panel 5</b>	319839	0.728	0.409 - 1.02

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

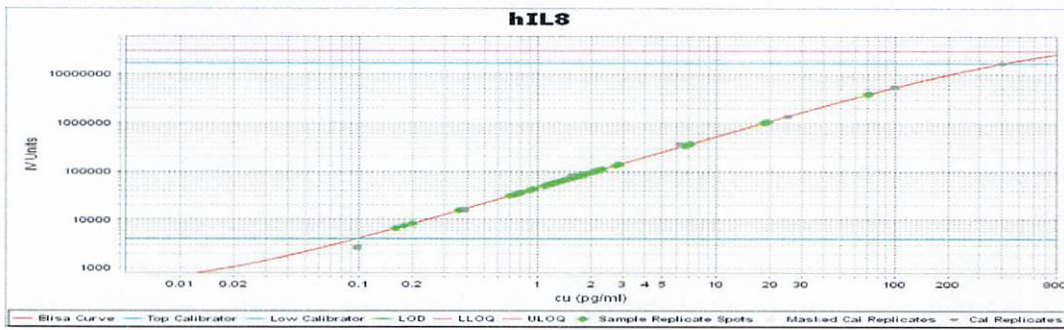
**IL-6 Calibrator Curve**



**Calibrator Levels  
(pg/mL)**

A	300
B	75.0
C	18.8
D	4.69
E	1.17
F	0.293
G	0.073
H	0.000

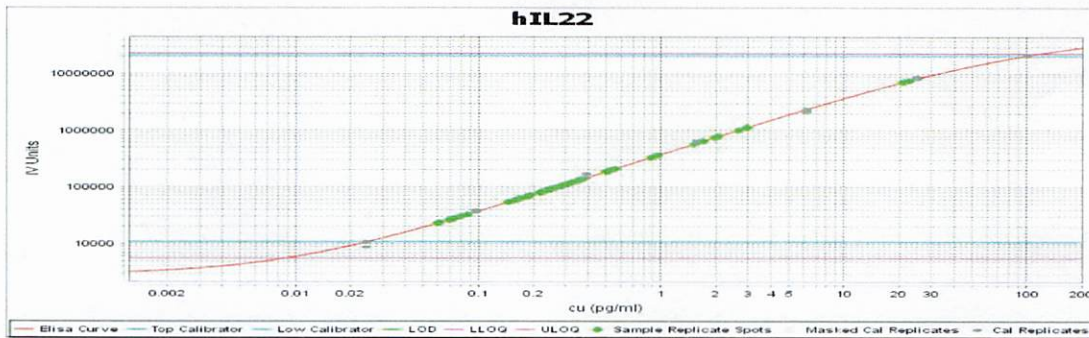
**IL-8 Calibrator Curve**



**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

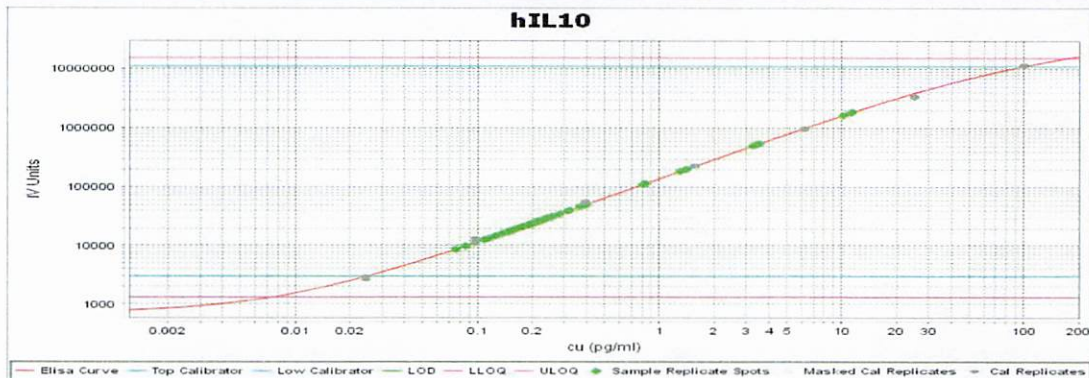
**IL-22 Calibrator Curve**



**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-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