

<b>Product Number:</b>	85-0412
<b>Lot Number:</b>	504428
<b>Expiration:</b>	14-Aug-2025
<b>Platform(s):</b>	SP-X

<b>CorPlex Cytokine Plate</b>	97-0329	426910P
<b>Biotinylated Antibody Reagent</b>	98-0329	422012B
<b>Sample Diluent 21</b>	1863329	403020
<b>IL-6 Calibrator</b>	18910044	324303
<b>IL-8 Calibrator</b>	1858454	319436
<b>IL-22 Calibrator</b>	18910036	319459
<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 Complex Cytokine Panel 1 Instruction Book 85-0412 116-5A0-1-AB (KI-0200).

Review/Approval

*Kyle Grison* Senior Document Control Specialist, QA  
 Name Title

*[Signature]* 30 Oct 2024  
 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	205	148 - 298
<b>Panel 2</b>	319839	53.2	40.5 - 70.0
<b>Panel 3</b>	319839	19.8	14.6 - 29.2
<b>Panel 4</b>	319839	7.01	4.31 - 11.1
<b>Panel 5</b>	319839	2.54	1.40 - 4.47

**IL-8**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	256	181 - 395
<b>Panel 2</b>	319839	75.0	49.8 - 105
<b>Panel 3</b>	319839	26.1	19.8 - 40.6
<b>Panel 4</b>	319839	7.14	5.79 - 13.0
<b>Panel 5</b>	319839	2.33	2.14 - 5.07

**IL-22**

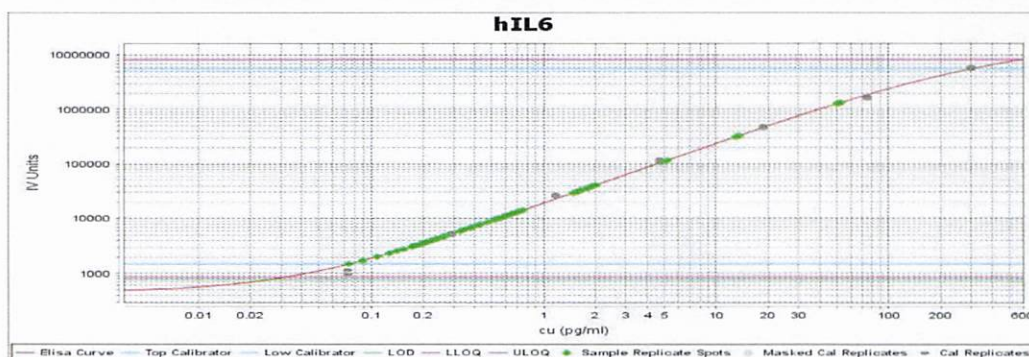
	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	54.6	33.7 - 103
<b>Panel 2</b>	319839	7.98	5.99 - 15.1
<b>Panel 3</b>	319839	4.33	2.56 - 7.64
<b>Panel 4</b>	319839	0.774	0.56 - 1.51
<b>Panel 5</b>	319839	0.222	0.148 - 0.449

**IL-10**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	39.8	29.2 - 62.2
<b>Panel 2</b>	319839	12.3	9.21 - 17.7
<b>Panel 3</b>	319839	4.74	3.28 - 7.66
<b>Panel 4</b>	319839	1.38	1.02 - 2.39
<b>Panel 5</b>	319839	0.554	0.424 - 1.02

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

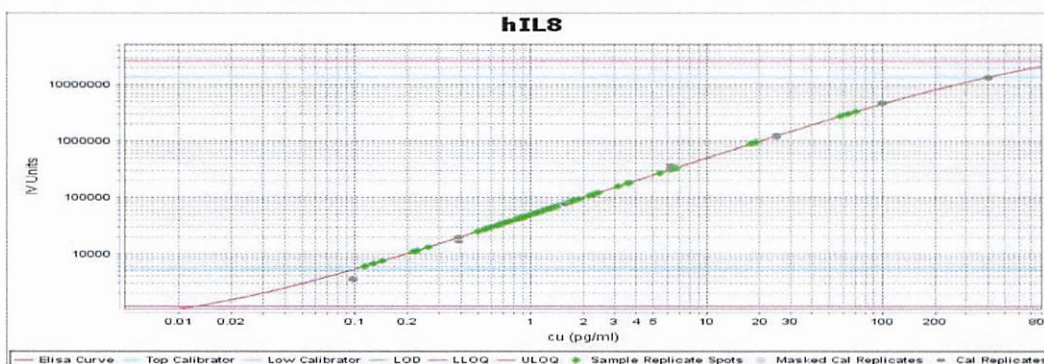
**IL-6 Calibrator Curve**



**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

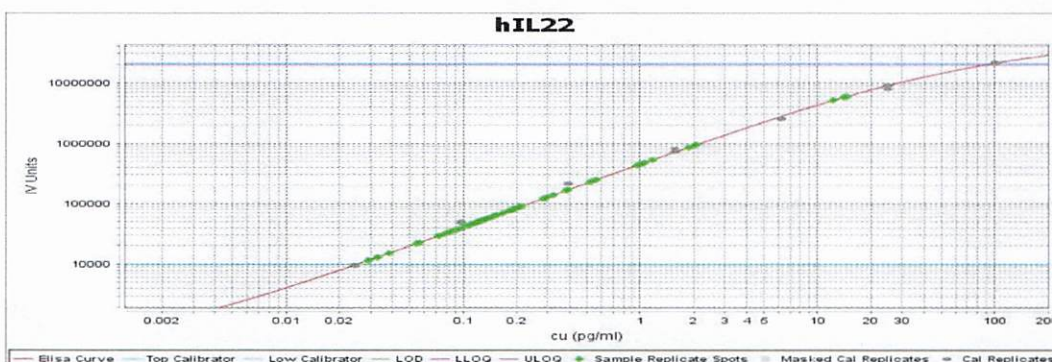
**IL-8 Calibrator Curve**



**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

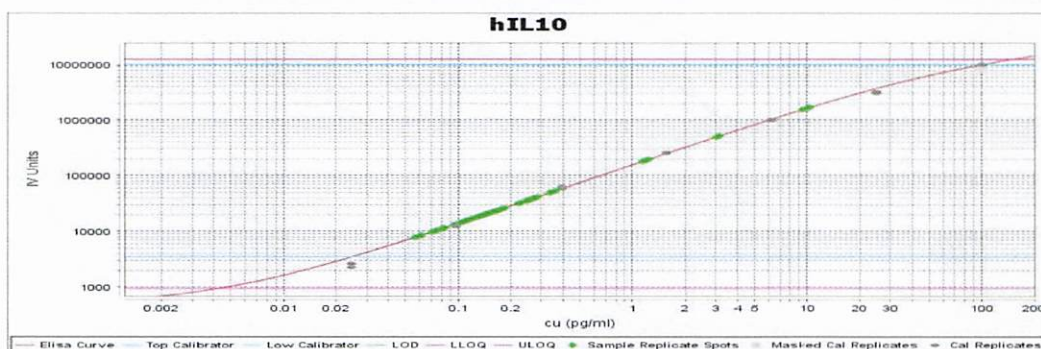
**IL-22 Calibrator Curve**



**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