

<b>Product Number:</b>	85-0412
<b>Lot Number:</b>	504392
<b>Expiration:</b>	14-Aug-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-6 Calibrator	18910044	324303	N/A
IL-8 Calibrator	1858454	319436	N/A
IL-22 Calibrator	18910036	319459	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 Complex Cytokine Panel 1 Instruction Book 85-0412 116-5A0-1-AB (KI-0200).

Review/Approval

*Kyle Giroux*

Name

*Senior Document Control Specialist, QA*

Title

*[Signature] 26 Sep 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	182	148 - 298
<b>Panel 2</b>	319839	55.4	40.5 - 70.0
<b>Panel 3</b>	319839	23.3	14.6 - 29.2
<b>Panel 4</b>	319839	8.89	4.31 - 11.1
<b>Panel 5</b>	319839	3.27	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	319	181 - 395
<b>Panel 2</b>	319839	84.8	49.8 - 105
<b>Panel 3</b>	319839	34.4	19.8 - 40.6
<b>Panel 4</b>	319839	10.9	5.79 - 13.0
<b>Panel 5</b>	319839	4.24	2.14 - 5.07

**IL-22**

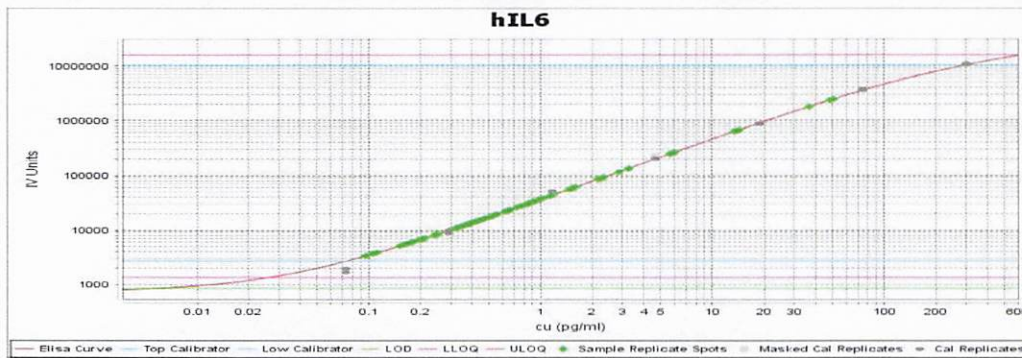
	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	46.8	33.7 - 103
<b>Panel 2</b>	319839	8.81	5.99 - 15.1
<b>Panel 3</b>	319839	4.23	2.56 - 7.64
<b>Panel 4</b>	319839	0.877	0.56 - 1.51
<b>Panel 5</b>	319839	0.237	0.148 - 0.449

**IL-10**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319839	33.6	29.2 - 62.2
<b>Panel 2</b>	319839	13.6	9.21 - 17.7
<b>Panel 3</b>	319839	5.87	3.28 - 7.66
<b>Panel 4</b>	319839	1.93	1.02 - 2.39
<b>Panel 5</b>	319839	0.794	0.424 - 1.02

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

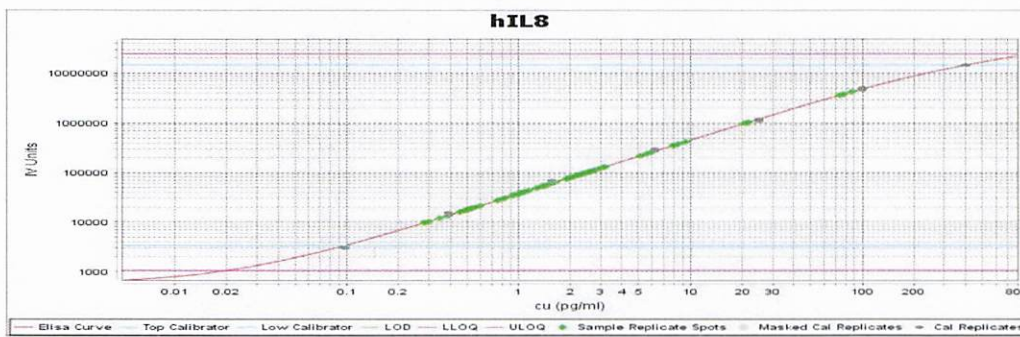
**IL-6 Calibrator Curve**



**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

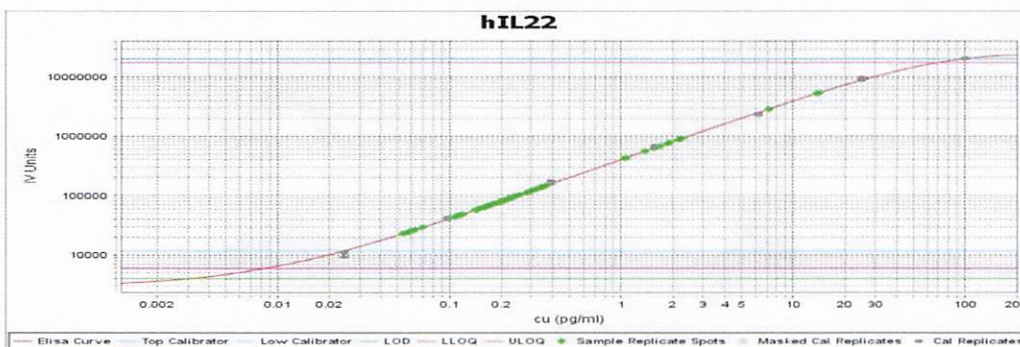
**IL-8 Calibrator Curve**



**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

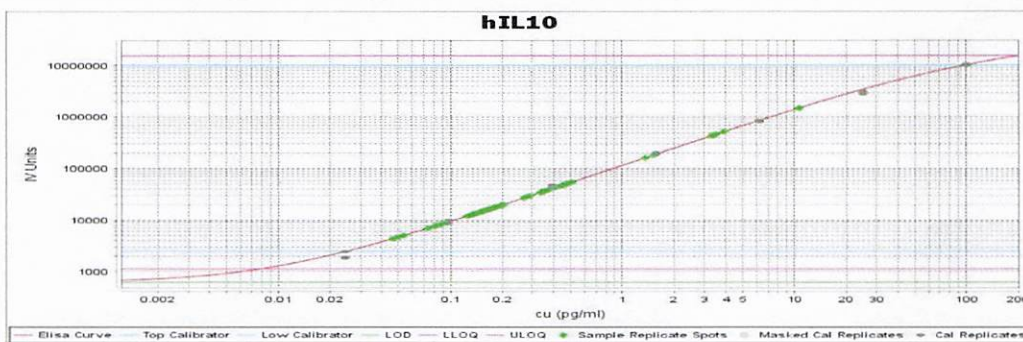
**IL-22 Calibrator Curve**



**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