

<b>Product Number:</b>	85-0401
<b>Lot Number:</b>	504104
<b>Expiration:</b>	26-Oct-2024
<b>Platform(s):</b>	SP-X



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

<b>CorPlex Cytokine Plate</b>	97-0401	319548P
<b>Biotinylated Antibody Reagent</b>	98-0401	319535B
<b>Sample Diluent 21</b>	1863329	319785
<b>IL-1β Calibrator</b>	1880100	319310
<b>IL-6 Calibrator</b>	1880101	319336
<b>IL-10 Calibrator</b>	1880102	319141
<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 Corplex Cytokine Panel 2 Kit Instructions 85-0401 (KI-0209).

Review/Approval

Kathleen Cornell  
 Name  
 Quality Control  
 Supervisor  
 Title

*Kathleen Cornell*  
 Signature/Date  
 29  
 10/2024

**Release Materials\***

**IL-1 $\beta$**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319012	4.86	3.88 - 9.76
<b>Panel 2</b>	319012	1.58	1.09 - 3.03
<b>Panel 3</b>	319012	0.510	0.303 - 1.09
<b>Panel 4</b>	319012	0.199	0.102 - 0.394
<b>Panel 5</b>	319012	0.083	0.030 - 0.172

**IL-6**

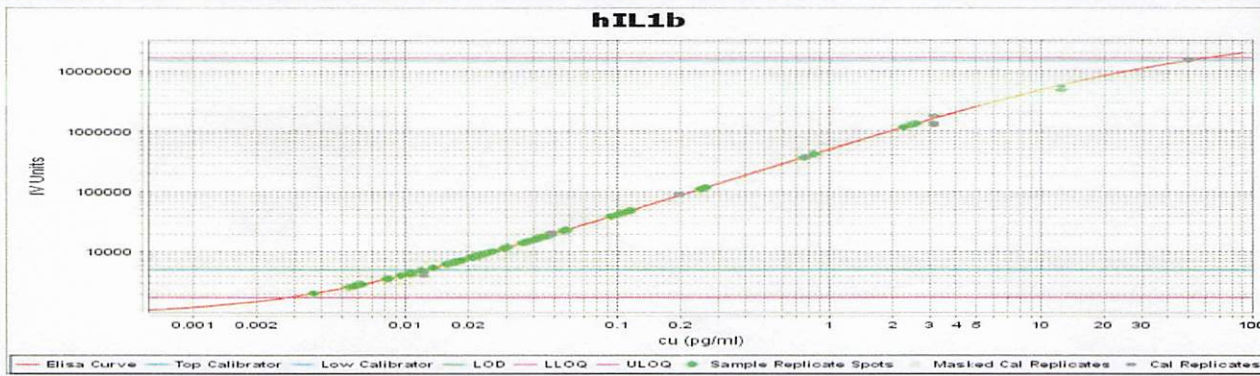
	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319012	23.6	17.0 - 37.8
<b>Panel 2</b>	319012	9.16	6.65 - 13.7
<b>Panel 3</b>	319012	3.66	1.84 - 5.81
<b>Panel 4</b>	319012	1.64	0.608 - 2.54
<b>Panel 5</b>	319012	0.825	0.336 - 1.16

**IL-10**

	<b>Lot</b>	<b>Result (pg/mL)</b>	<b>Spec. Range (pg/mL)</b>
<b>Panel 1</b>	319012	47.2	39.2 - 58.8
<b>Panel 2</b>	319012	14.8	13.1 - 19.6
<b>Panel 3</b>	319012	5.49	4.35 - 7.51
<b>Panel 4</b>	319012	2.07	1.48 - 2.82
<b>Panel 5</b>	319012	0.823	0.584 - 1.07

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

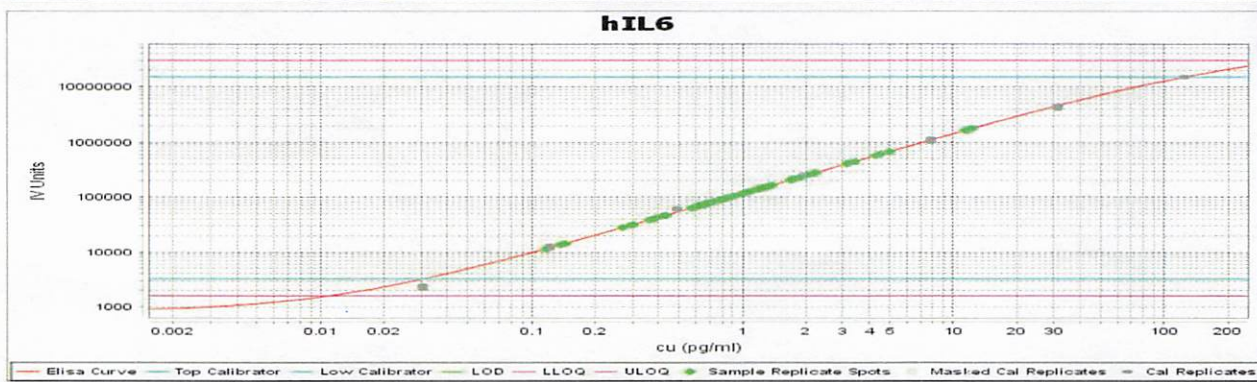
**IL-1 $\beta$  Calibrator Curve**



**Calibrator Levels  
(pg/mL)**

A	50.0
B	12.5
C	3.13
D	0.781
E	0.195
F	0.049
G	0.012
H	0.000

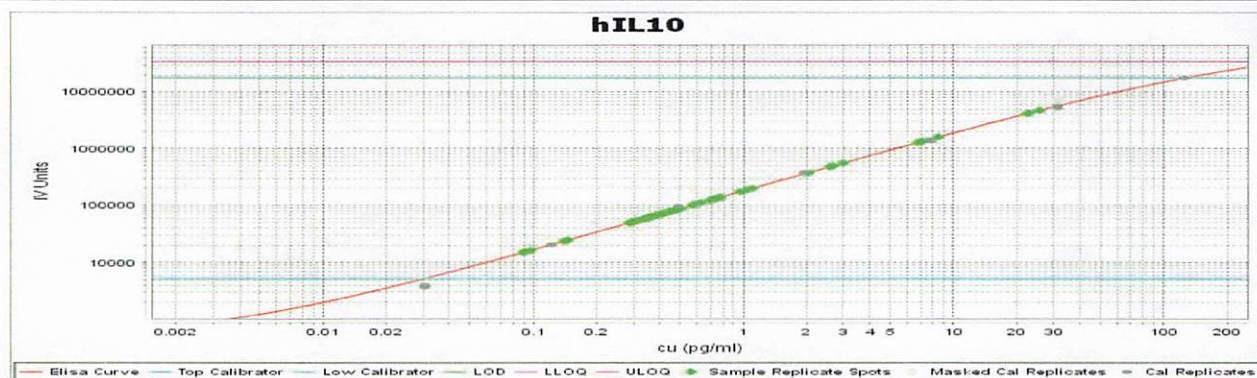
**IL-6 Calibrator Curve**



**Calibrator Levels  
(pg/mL)**

A	125
B	31.3
C	7.81
D	1.95
E	0.488
F	0.122
G	0.031
H	0.000

**IL-10 Calibrator Curve**



**Calibrator Levels  
(pg/mL)**

A	125
B	31.3
C	7.81
D	1.95
E	0.488
F	0.122
G	0.031
H	0.000