

Product Number:	85-0401
Lot Number:	319144
Expiration:	28-Jan-2022
Platform(s) :	SP-X



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

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0401	319144P
Biotinylated Antibody Rgt	98-0401	319220B
Sample Diluent 21	1863329	319213
IL-1 $\beta$ Calibrator	1880100	319197
IL-6 Calibrator	1880101	319198
IL-10 Calibrator	1880102	319199
Wash Buffer (25X) <sup>1</sup>	1860055	N/A
Streptavidin-HRP Reagent <sup>1</sup>	1858263	N/A
Super Signal <sup>®</sup> Luminol* <sup>1</sup>	1858089	N/A
Super Signal <sup>™</sup> Peroxide* <sup>1</sup>	1858090	N/A
Streptavidin-HRP Reagent <sup>1</sup>	1858263	N/A
Super Signal <sup>™</sup> Luminol* <sup>1</sup>	1858089	N/A
Super Signal <sup>™</sup> Peroxide* <sup>1</sup>	1858090	N/A

<sup>1</sup> Reagents are not Kit Lot Specific

\*SuperSignal<sup>®</sup> 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 the CorPlex Cytokine 2 Kit Instructions (85A-0401).

**Release Materials**

**IL-1 $\beta$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319012	6.557	5.207 - 9.540
Panel 2	319012	1.951	1.498 - 2.943
Panel 3	319012	0.621	0.438 - 1.088
Panel 4	319012	0.223	0.150 - 0.383
Panel 5	319012	0.085	0.051 - 0.161

**IL-6**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319012	23.54	17.88 - 37.07
Panel 2	319012	8.934	6.616 - 13.87
Panel 3	319012	3.278	2.317 - 5.420
Panel 4	319012	1.321	0.974 - 2.159
Panel 5	319012	0.654	0.460 - 1.120

**IL-10**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319012	49.89	40.40 - 60.60
Panel 2	319012	16.31	13.76 - 20.65
Panel 3	319012	5.770	4.964 - 7.447
Panel 4	319012	2.043	1.756 - 2.754
Panel 5	319012	0.777	0.705 - 1.063

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

Review/Approval

Shivani Goel

Quality Control  
Manager

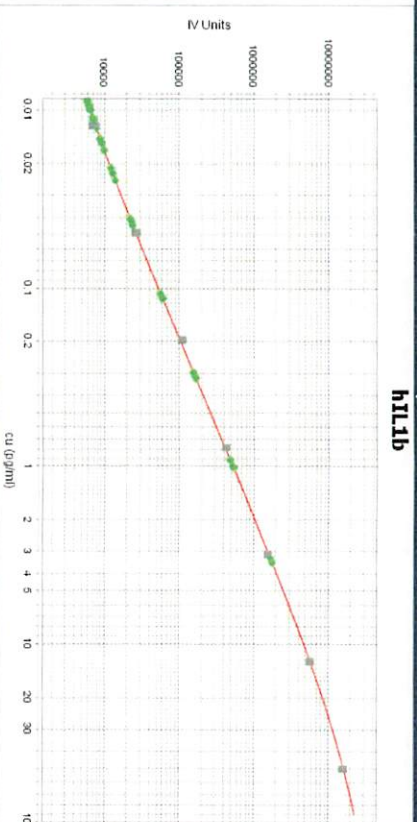
Shivani Goel 14 April 2021

Name

Title

Signature/Date

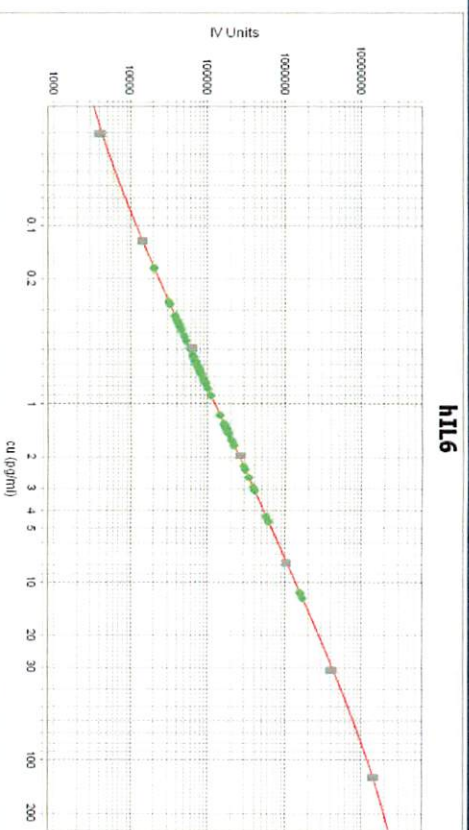
Ir-1 $\beta$  Calibration Curve  
**hIL1 $\beta$**



Calibrator Levels  
(pg/ml)

- 1 50
- 2 12.5
- 3 3.13
- 4 0.781
- 5 0.195
- 6 0.049
- 7 0.012
- 8 0.000

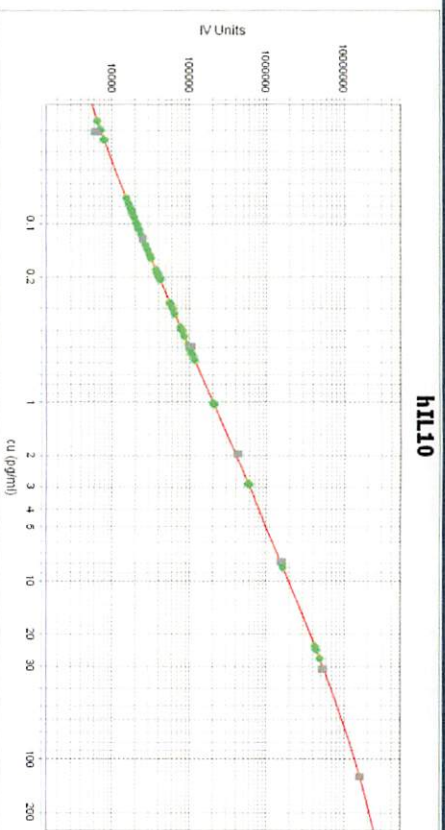
Ir-6 Calibration Curve  
**hIL6**



Calibrator Levels  
(pg/ml)

- 1 125
- 2 31.25
- 3 7.813
- 4 1.953
- 5 0.488
- 6 0.122
- 7 0.031
- 8 0.000

Ir-10 Calibration Curve  
**hIL10**



Calibrator Levels  
(pg/ml)

- 1 125
- 2 31.25
- 3 7.813
- 4 1.953
- 5 0.488
- 6 0.122
- 7 0.031
- 8 0.000