

Product Number:	85-0401
Lot Number:	319143
Expiration:	28-Jan-2022
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0401	319143P
Biotinylated Antibody Rgt	98-0401	319219B
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).



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

**Release Materials**

**IL-1 $\beta$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319012	7.311	5.207 - 9.540
Panel 2	319012	2.097	1.498 - 2.943
Panel 3	319012	0.679	0.438 - 1.088
Panel 4	319012	0.248	0.150 - 0.383
Panel 5	319012	0.088	0.051 - 0.161

**IL-6**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319012	25.98	17.88 - 37.07
Panel 2	319012	9.566	6.616 - 13.87
Panel 3	319012	3.468	2.317 - 5.420
Panel 4	319012	1.526	0.974 - 2.159
Panel 5	319012	0.767	0.460 - 1.120

**IL-10**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319012	51.09	40.40 - 60.60
Panel 2	319012	16.56	13.76 - 20.65
Panel 3	319012	6.051	4.964 - 7.447
Panel 4	319012	2.340	1.756 - 2.754
Panel 5	319012	0.878	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

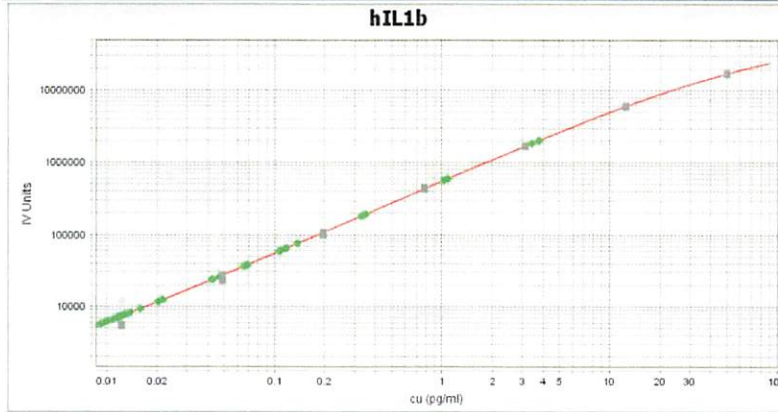
*Shivani Goel 14 Apr 2021*

Name

Title

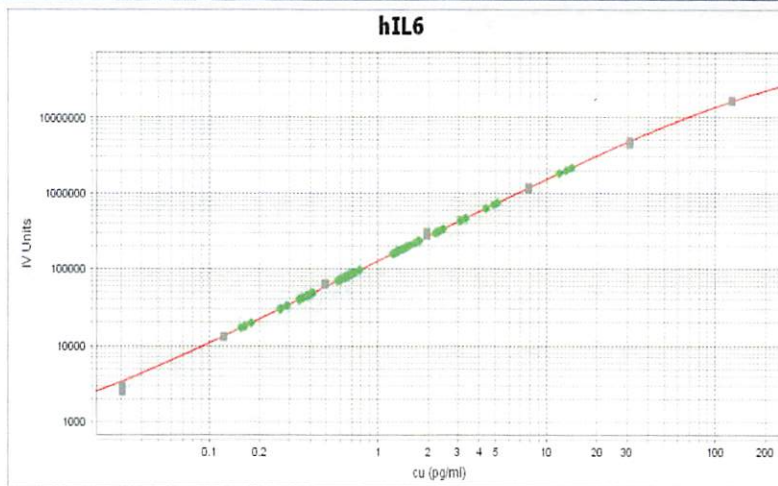
Signature/Date

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



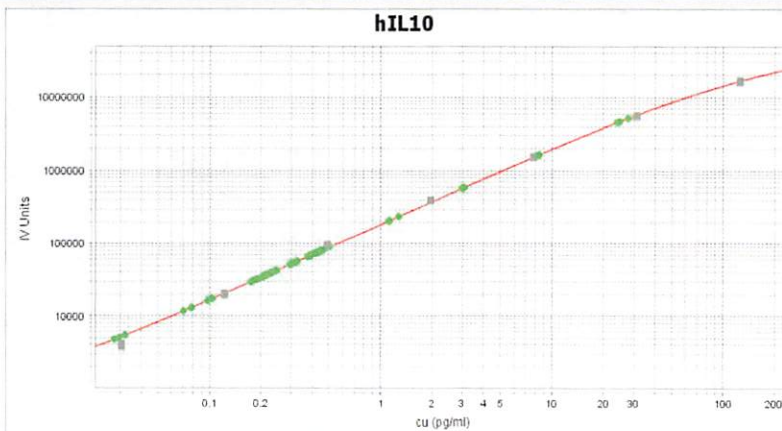
Calibrator Levels (pg/ml)	
1	50
2	12.5
3	3.13
4	0.781
5	0.195
6	0.049
7	0.012
8	0.000

**IL-6 Calibration Curve**



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

**IL-10 Calibration Curve**



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