

Product Number:	85-0401
Lot Number:	319312
Expiration:	12-Nov-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	319312P
Biotinylated Antibody Rgt	98-0401	319312B
Sample Diluent 21	1863329	319307
IL-1 β Calibrator	1880100	319197
IL-6 Calibrator	1880101	319198
IL-10 Calibrator	1880102	319199
Wash Buffer (25X) ¹	1860055	N/A
Streptavidin-HRP Reagent ¹	1858263	N/A
Super Signal [®] Luminol* ¹	1858089	N/A
Super Signal [™] Peroxide* ¹	1858090	N/A

¹ Reagents are not Kit Lot Specific

*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 the CorPlex Cytokine 2 Kit Instructions (85A-0401).

Release Materials

IL-1 β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319012	6.697	5.207 - 9.540
Panel 2	319012	1.880	1.498 - 2.943
Panel 3	319012	0.623	0.438 - 1.088
Panel 4	319012	0.214	0.150 - 0.383
Panel 5	319012	0.071	0.051 - 0.161

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319012	25.01	17.88 - 37.07
Panel 2	319012	9.23	6.616 - 13.87
Panel 3	319012	3.265	2.317 - 5.420
Panel 4	319012	1.410	0.974 - 2.159
Panel 5	319012	0.663	0.460 - 1.120

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319012	48.26	40.40 - 60.60
Panel 2	319012	16.01	13.76 - 20.65
Panel 3	319012	5.347	4.964 - 7.447
Panel 4	319012	2.000	1.756 - 2.754
Panel 5	319012	0.740	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

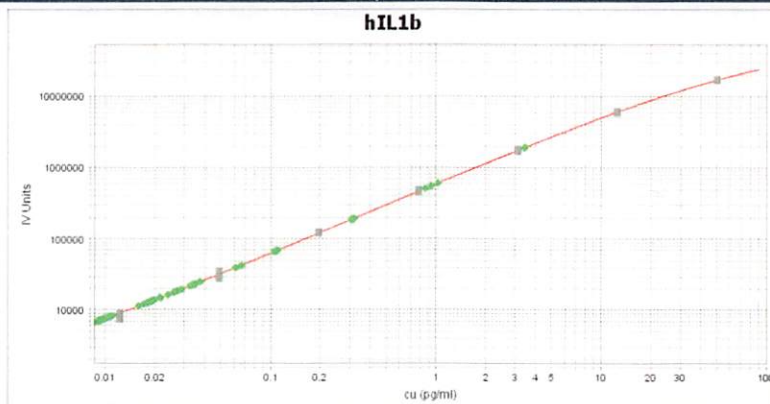
Shivani Goel 30 Jun 2021

Name

Title

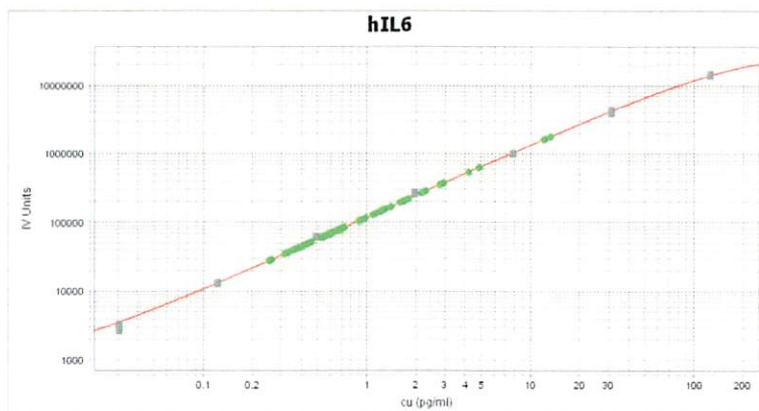
Signature/Date

IL-1 β Calibration Curve



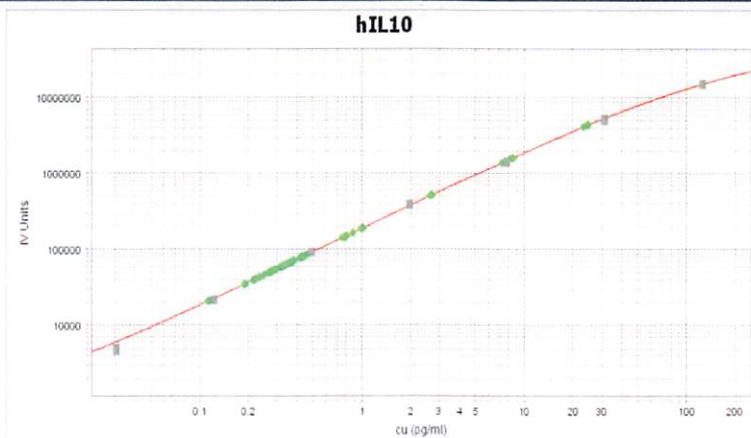
Calibrator Levels (pg/ml)	
1	50
2	12.5
3	3.13
4	0.781
5	0.195
6	0.049
7	0.012
8	0.000

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