

Product Number:	85-0426
Lot Number:	319331
Expiration:	12-Nov-2022
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0426	319331P
Biotinylated Antibody Rgt	98-0426	319331B
Sample Diluent 21	1863329	319307
IL-8 Calibrator	1858454	319011
sTNF-RII Calibrator	1880104	319337
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
Streptavidin-HRP Reagent ¹	1858263	N/A



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

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 3 Kit Instructions (85A-0426).

Release Materials

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319269	549.5	379.5 - 801.2
Panel 2	319269	199.8	141.3 - 278.6
Panel 3	319269	62.89	48.87 - 87.06
Panel 4	319269	21.22	17.02 - 29.78
Panel 5	319269	6.783	5.062 - 11.80

sTNFR-II

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319269	5428.0	3622.3 - 6105.0
Panel 2	319269	1782.0	1118.0 - 2215.2
Panel 3	319269	682.0	448.79 - 914.8
Panel 4	319269	289.8	218.35 - 454.9
Panel 5	319269	179.1	116.38 - 294.7

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

Review/Approval

Shivani Goel

Quality Control
Manager

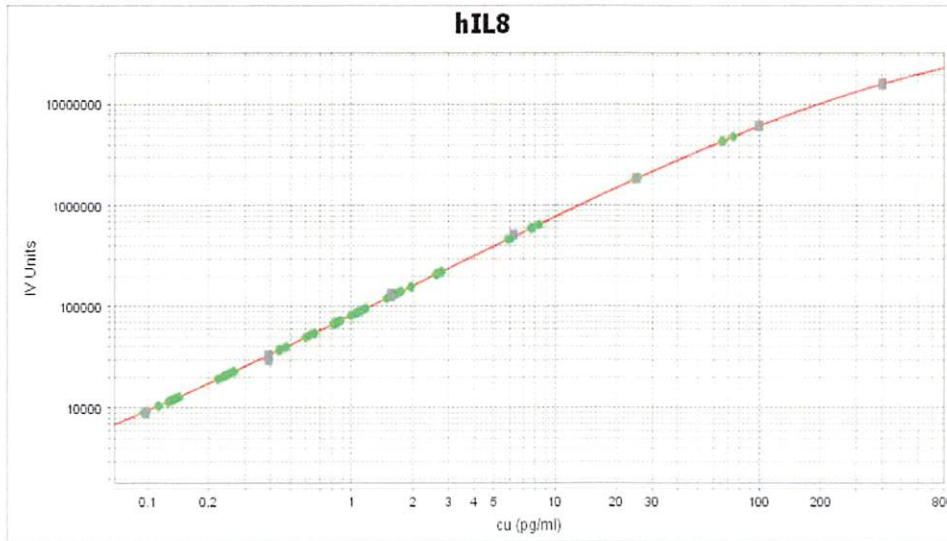
Shivani Goel 16 Jun 2021

Name

Title

Signature/Date

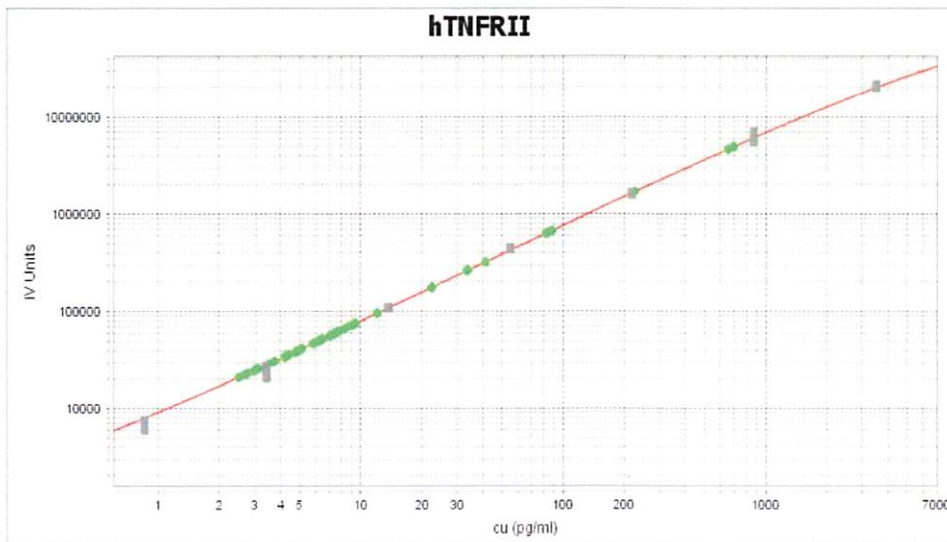
IL-8 Calibration Curve



Calibrator Levels (pg/ml)

1	400.0
2	100.0
3	25.00
4	6.250
5	1.563
6	0.391
7	0.098
8	0.000

sTNF-RII Calibration Curve



Calibrator Levels (pg/mL)

1	3500
2	875.0
3	218.8
4	54.69
5	13.67
6	3.418
7	0.854
8	0.000