

Product Number:	85-0329
Lot Number:	318404
Expiration:	9-Jun-2021
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	318404P
Biotinylated Antibody Rgt	98-0329	318404B
Sample Diluent 21 <sup>1</sup>	1863329	N/A
IL-12p70 Calibrator	18910034	317901
IL-1 $\beta$ Calibrator	18910040	318067
IL-4 Calibrator	1858452	317999
IL-5 Calibrator	1858462	317984
IFN $\gamma$ Calibrator	1858260	318097
IL-6 Calibrator	18910044	318089
IL-8 Calibrator	1858454	318081
IL-22 Calibrator	18910036	318408
TNF $\alpha$ Calibrator	1880012	318436
IL-10 Calibrator	18910043	318068
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

<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 Kit Instructions (85A-0329).



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

**Release Materials**

**IL-12p70**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	35.5	29.4 - 41.6
QC3 Mid	318591	12.8	9.62 - 13.5
QC3 Low	318591	1.33	1.02 - 1.43

**IL-1 $\beta$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	12.3	10.1 - 13.3
QC3 Mid	318591	4.53	3.50 - 4.55
QC3 Low	318591	0.624	0.552 - 0.759

**IL-4**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	28.0	21.5 - 32.0
QC3 Mid	318591	10.9	8.53 - 13.3
QC3 Low	318591	2.79	2.57 - 5.24

**IL-5**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	27.1	23.5 - 35.3
QC3 Mid	318591	10.5	7.79 - 10.8
QC3 Low	318591	0.992	1.02 - 1.51

**IFN $\gamma$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	0.951	0.856 - 1.20
QC3 Mid	318591	0.430	0.327 - 0.541
QC3 Low	318591	0.128	0.108 - 0.282

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

Release Materials

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	34.5	29.2 - 39.2
QC3 Mid	318591	12.2	10.3 - 13.7
QC3 Low	318591	2.17	1.99 - 2.62

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	66.2	57.8 - 83.4
QC3 Mid	318591	36.3	28.7 - 41.8
QC3 Low	318591	20.4	16.7 - 23.4

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	8.65	9.11 - 12.8
QC3 Mid	318591	3.49	3.40 - 4.86
QC3 Low	318591	0.764	0.632 - 0.904

TNF $\alpha$

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	58.2	50.6 - 70.6
QC3 Mid	318591	24.5	20.5 - 28.3
QC3 Low	318591	7.00	6.82 - 10.5

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	14.8	13.0 - 17.7
QC3 Mid	318591	5.34	4.66 - 6.48
QC3 Low	318591	0.932	0.800 - 1.19

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

Review/Approval

Patrizia Stadler

Associate  
Director of QC

*P. Stadler*

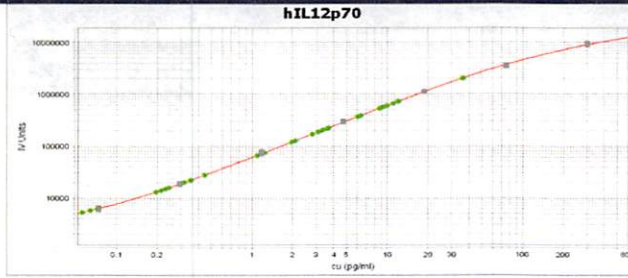
24 MAR 2020

Name

Title

Signature/Date

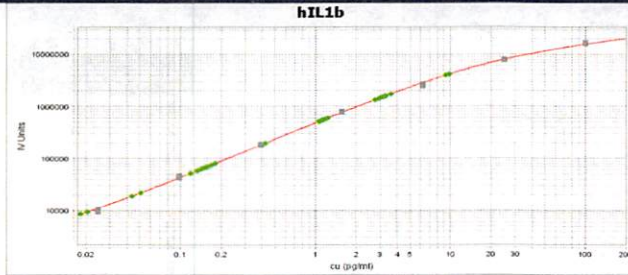
**IL-12p70 Calibration Curve**



**Calibrator Levels (pg/mL)**

1	300
2	75.0
3	18.8
4	4.69
5	1.17
6	0.293
7	0.073
8	0.000

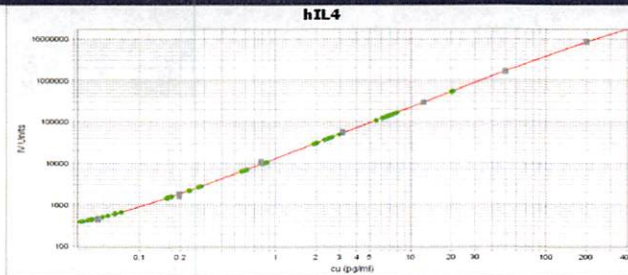
**IL-1 $\beta$  Calibration Curve**



**Calibrator Levels (pg/mL)**

1	100
2	25.0
3	6.25
4	1.56
5	0.391
6	0.098
7	0.024
8	0.000

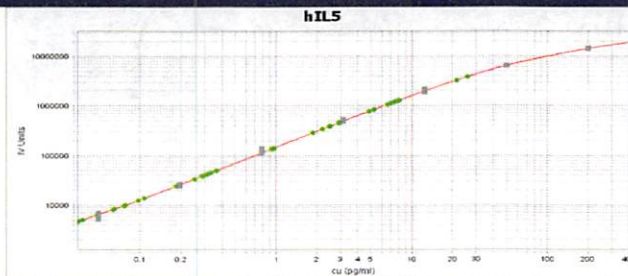
**IL-4 Calibration Curve**



**Calibrator Levels (pg/mL)**

1	200
2	50.0
3	12.5
4	3.13
5	0.781
6	0.195
7	0.049
8	0.000

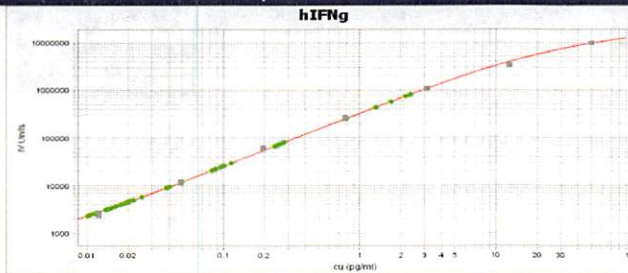
**IL-5 Calibration Curve**



**Calibrator Levels (pg/mL)**

1	200
2	50.0
3	12.5
4	3.13
5	0.781
6	0.195
7	0.049
8	0.000

**IFN $\gamma$  Calibration Curve**



**Calibrator Levels (pg/mL)**

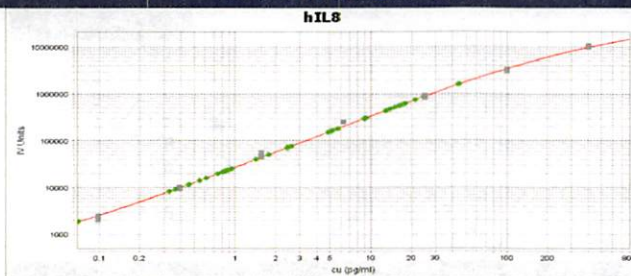
1	200
2	50.0
3	12.5
4	3.13
5	0.781
6	0.195
7	0.049
8	0.000

**IL-6 Calibration Curve**



Calibrator Levels (pg/mL)	
1	300
2	75.0
3	18.8
4	4.69
5	1.17
6	0.293
7	0.073
8	0.000

**IL-8 Calibration Curve**



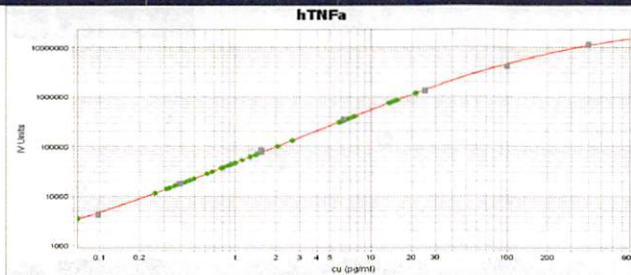
Calibrator Levels (pg/mL)	
1	400
2	100
3	25.0
4	6.25
5	1.56
6	0.391
7	0.098
8	0.000

**IL-22 Calibration Curve**



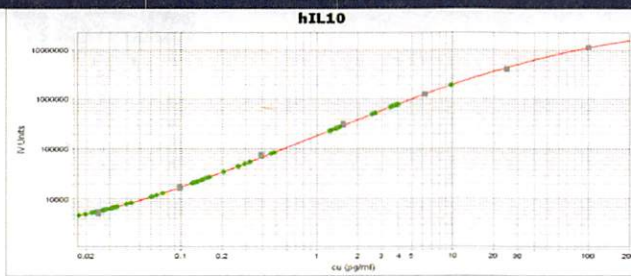
Calibrator Levels (pg/mL)	
1	100
2	25.0
3	6.25
4	1.56
5	0.391
6	0.098
7	0.024
8	0.000

**TNF $\alpha$  Calibration Curve**



Calibrator Levels (pg/mL)	
1	400
2	100
3	25.0
4	6.25
5	1.56
6	0.391
7	0.098
8	0.000

**IL-10 Calibration Curve**



Calibrator Levels (pg/mL)	
1	100
2	25.0
3	6.25
4	1.56
5	0.391
6	0.098
7	0.024
8	0.000