

Product Number:	85-0329
Lot Number:	318459
Expiration:	12-Aug-2021
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	318459P
Biotinylated Antibody Rgt	98-0329	318744B
Sample Diluent 21 ¹	1863329	N/A
IL-12p70 Calibrator	18910034	318629
IL-1b Calibrator	18910040	318673
IL-4 Calibrator	1858452	318480
IL-5 Calibrator	1858462	318465
IFNg Calibrator	1858260	318538
IL-6 Calibrator	18910044	318583
IL-8 Calibrator	1858454	318342
IL-22 Calibrator	18910036	318408
TNFa Calibrator	1880012	318436
IL-10 Calibrator	18910043	318630
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



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 Cytokine1 Kit Instructions (85A-0329).

Release Materials

IL-12p70

	Lot	Result (pg/mL)	Mean Range (pg/mL)		
Panel 1	318641	165.50	164	-	246
Panel 2	318641	56.30	54.9	-	83.5
Panel 3	318641	20.79	17.7	-	28.3
Panel 4	318641	7.77	5.19	-	10.40
Panel 5	318641	2.53	1.75	-	3.55

IL-1b

	Lot	Result (pg/mL)	Mean Range (pg/mL)		
Panel 1	318641	23.51	19.6	-	29.4
Panel 2	318641	8.01	6.56	-	9.85
Panel 3	318641	2.69	2.02	-	3.42
Panel 4	318641	0.96	0.611	-	1.28
Panel 5	318641	0.31	0.140	-	0.511

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)		
Panel 1	318641	64.36	69.3	-	111
Panel 2	318641	20.45	22.7	-	36.95
Panel 3	318641	7.37	6.82	-	14.1
Panel 4	318641	2.48	2.27	-	5.302
Panel 5	318641	0.77	0.885	-	1.69

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)		
Panel 1	318641	60.04	50.98	-	76.5
Panel 2	318641	19.44	16.7	-	26.3
Panel 3	318641	7.27	5.85	-	9.33
Panel 4	318641	2.50	1.86	-	3.47
Panel 5	318641	0.94	0.651	-	1.293

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

Release Materials

IFN γ

	Lot	Result (pg/mL)	Mean Range (pg/mL)		
Panel 1	318641	9.30	8.62	-	12.93
Panel 2	318641	3.55	3.05	-	4.57
Panel 3	318641	1.30	1.04	-	1.63
Panel 4	318641	0.55	0.340	-	0.664
Panel 5	318641	0.22	0.137	-	0.284

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)		
Panel 1	318641	109.40	118	-	187
Panel 2	318641	3.29	3.67	-	7.78
Panel 3	318641	1.19	1.19	-	2.68
Panel 4	318641	0.44	0.449	-	1.09
Panel 5	318641	0.29	0.214	-	0.437

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)		
Panel 1	318641	453.10	382	-	573
Panel 2	318641	135.10	116	-	194
Panel 3	318641	49.25	39.3	-	68.4
Panel 4	318641	17.32	12.1	-	24.91
Panel 5	318641	5.32	2.49	-	9.304

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)		
Panel 1	318641	19.47	16.8	-	34.4
Panel 2	318641	5.85	4.69	-	10.2
Panel 3	318641	2.11	1.604	-	3.42
Panel 4	318641	0.76	0.538	-	1.24
Panel 5	318641	0.24	0.168	-	0.434

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)		
Panel 1	318641	63.50	62.9	-	103
Panel 2	318641	21.98	22.4	-	36.04
Panel 3	318641	7.52	7.40	-	12.4
Panel 4	318641	2.60	2.43	-	4.48
Panel 5	318641	0.92	0.877	-	1.70

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)		
Panel 1	318641	54.04	46.4	-	69.6
Panel 2	318641	18.89	14.4	-	25.8
Panel 3	318641	7.23	5.38	-	8.82
Panel 4	318641	2.77	2.03	-	3.20
Panel 5	318641	1.01	0.714	-	1.23

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

Review/Approval

Patrizia Stadler

Associate Director of
QC

P. Stadler

02 Jun 2020

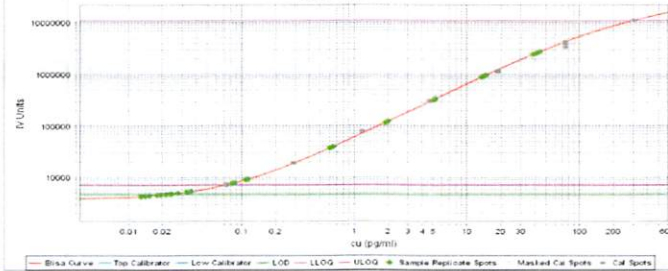
Name

Title

Signature/Date

IL-12p70 Calibration Curve

hIL12p70

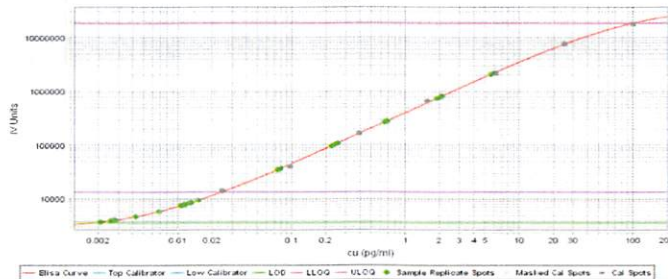


Calibrator Levels (pg/ml)

1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

IL-1 β Calibration Curve

hIL1 β

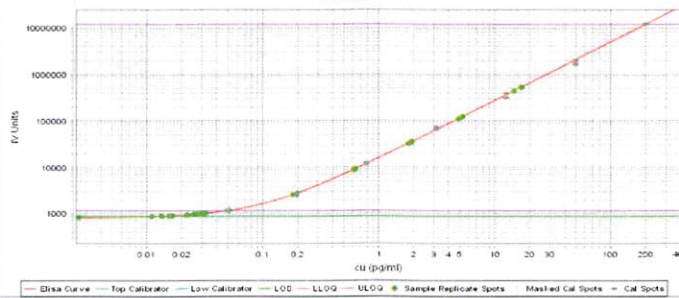


Calibrator Levels (pg/mL)

1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

IL-4 Calibration Curve

hIL4

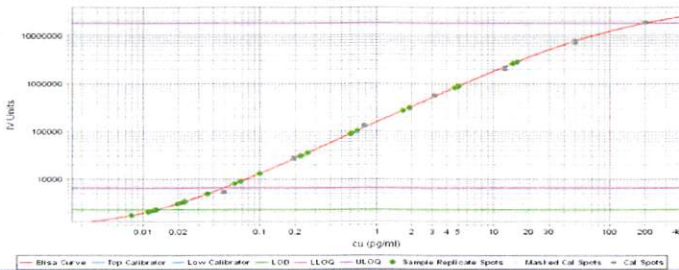


Calibrator Levels (pg/mL)

1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

IL-5 Calibration Curve

hIL5

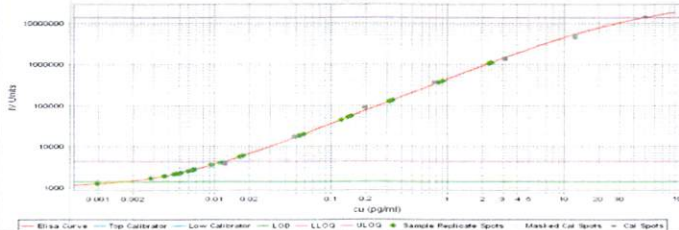


Calibrator Levels (pg/mL)

1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

IFN γ Calibration Curve

hIFN γ

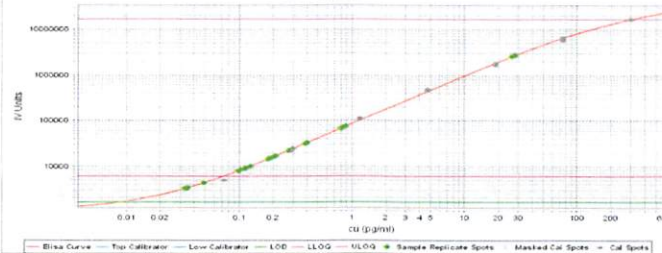


Calibrator Levels (pg/mL)

1	200
2	50
3	12.50
4	3.125
5	0.781
6	0.195
7	0.049
8	0.000

IL-6 Calibration Curve

hIL6

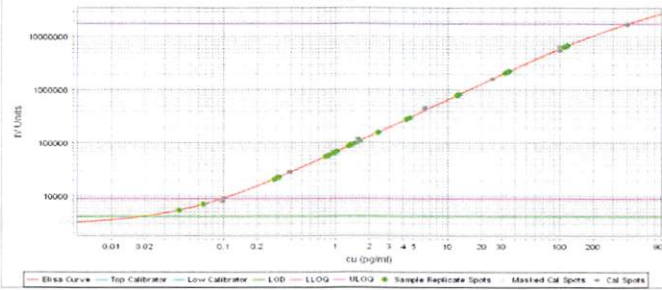


Calibrator Levels (pg/mL)

1	300
2	75
3	18.75
4	4.688
5	1.172
6	0.293
7	0.073
8	0.000

IL-8 Calibration Curve

hIL8

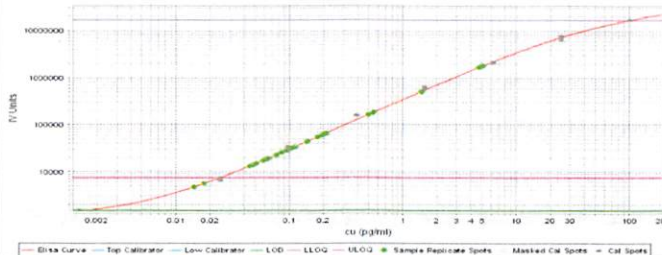


Calibrator Levels (pg/mL)

1	400
2	100
3	25.00
4	6.250
5	1.563
6	0.391
7	0.098
8	0.000

IL-22 Calibration Curve

hIL22

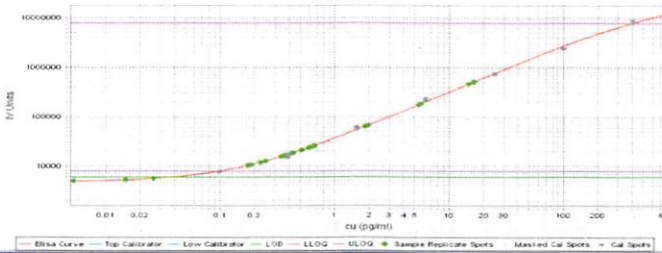


Calibrator Levels (pg/mL)

1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000

TNFα Calibration Curve

hTNFα

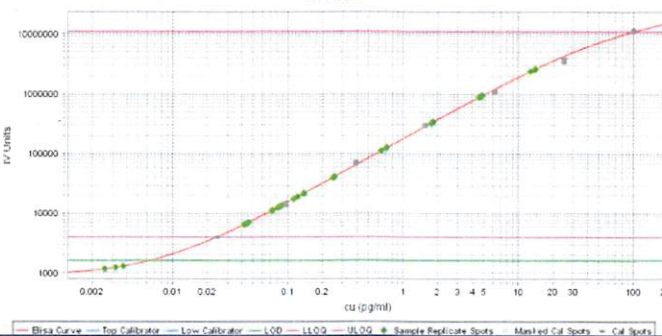


Calibrator Levels (pg/mL)

1	400
2	100
3	25.00
4	6.250
5	1.563
6	0.391
7	0.098
8	0.000

IL-10 Calibration Curve

hIL10



Calibrator Levels (pg/mL)

1	100
2	25
3	6.25
4	1.563
5	0.391
6	0.098
7	0.024
8	0.000