

Product Number:	85-0329
Lot Number:	319824
Expiration:	10-May-2023
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	319823P
Biotinylated Antibody Rgt	98-0329	319494B
Sample Diluent 21	1863329	319709
IL-12p70 Calibrator	18910034	319156
IL-1 β Calibrator	18910040	319564
IL-4 Calibrator	1858452	319458
IL-5 Calibrator	1858462	318837
IFN γ Calibrator	1858260	319711
IL-6 Calibrator	18910044	319484
IL-8 Calibrator	1858454	319575
IL-22 Calibrator	18910036	319459
TNF α Calibrator	1880012	319247
IL-10 Calibrator	18910043	319215
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 Cytokine Kit Instructions (85A-0329).

Release Materials

IL-12p70

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	149.0	133.1 - 204.3
Panel 2	318935	46.19	37.09 - 66.72
Panel 3	318935	15.37	13.48 - 21.97
Panel 4	318935	4.433	3.812 - 6.073
Panel 5	318935	1.303	1.188 - 2.091

IL-1 β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	24.95	12.49 - 32.68
Panel 2	318935	7.350	4.122 - 9.719
Panel 3	318935	2.586	1.349 - 3.475
Panel 4	318935	0.759	0.418 - 0.968
Panel 5	318935	0.246	0.133 - 0.345

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	57.66	51.07 - 77.41
Panel 2	318935	16.80	14.17 - 24.05
Panel 3	318935	6.107	5.217 - 8.887
Panel 4	318935	2.002	1.455 - 2.760
Panel 5	318935	0.741	0.418 - 1.250

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	81.83	43.59 - 101.9
Panel 2	318935	26.09	13.43 - 32.50
Panel 3	318935	9.118	4.948 - 11.09
Panel 4	318935	2.763	1.317 - 3.402
Panel 5	318935	0.801	0.473 - 1.224

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	318935	12.25	8.190 - 14.53
Panel 2	318935	3.719	2.393 - 4.771
Panel 3	318935	1.312	0.855 - 1.720
Panel 4	318935	0.417	0.277 - 0.494
Panel 5	318935	0.167	0.114 - 0.221

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	200.7	142.9 - 214.3
Panel 2	318935	88.24	62.87 - 99.85
Panel 3	318935	17.72	13.37 - 20.05
Panel 4	318935	8.256	5.951 - 8.927
Panel 5	318935	2.625	1.992 - 3.084

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	440.6	365.0 - 604.4
Panel 2	318935	118.9	92.02 - 191.0
Panel 3	318935	41.52	32.36 - 67.44
Panel 4	318935	11.42	9.991 - 18.54
Panel 5	318935	3.579	3.355 - 6.599

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	77.9	50.55 - 160.4
Panel 2	318935	13.18	7.411 - 28.41
Panel 3	318935	4.207	2.237 - 9.205
Panel 4	318935	1.178	0.738 - 2.619
Panel 5	318935	0.336	0.226 - 0.888

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	389.4	250.8 - 453.3
Panel 2	318935	109.3	68.46 - 156.1
Panel 3	318935	37.14	25.97 - 51.08
Panel 4	318935	10.93	7.455 - 14.81
Panel 5	318935	3.171	2.339 - 4.887

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	106.6	76.56 - 126.0
Panel 2	318935	30.16	21.73 - 39.58
Panel 3	318935	10.14	7.682 - 13.72
Panel 4	318935	3.095	2.408 - 3.832
Panel 5	318935	1.072	0.837 - 1.442

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

Review/Approval

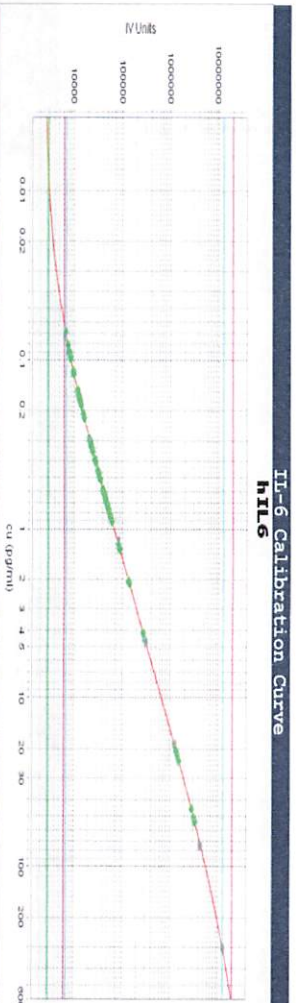
Matthew Smart

Quality Control
Director

Name

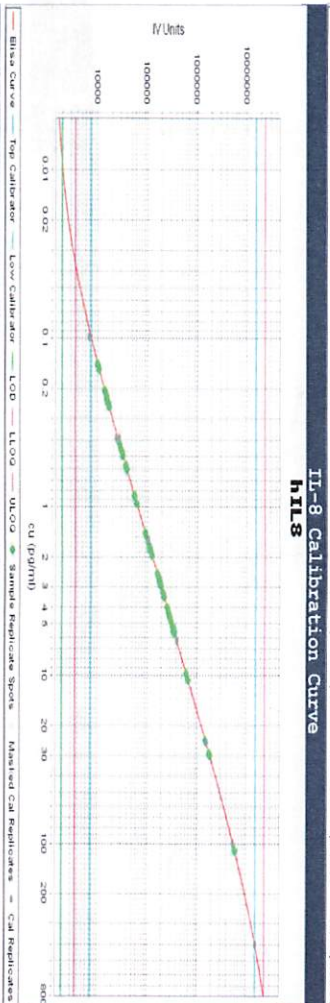
Title

M. Smart 17 Aug 2022
Signature/Date



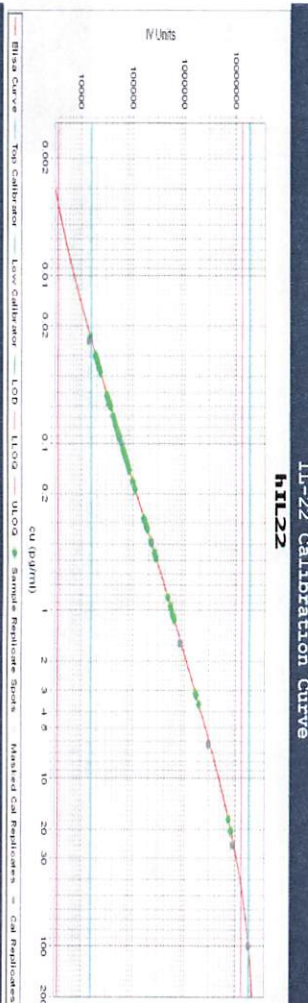
hIL-6 Calibration Curve

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



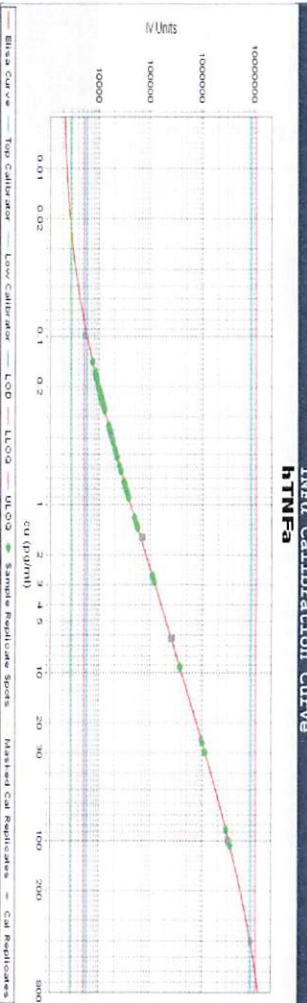
hIL-8 Calibration Curve

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



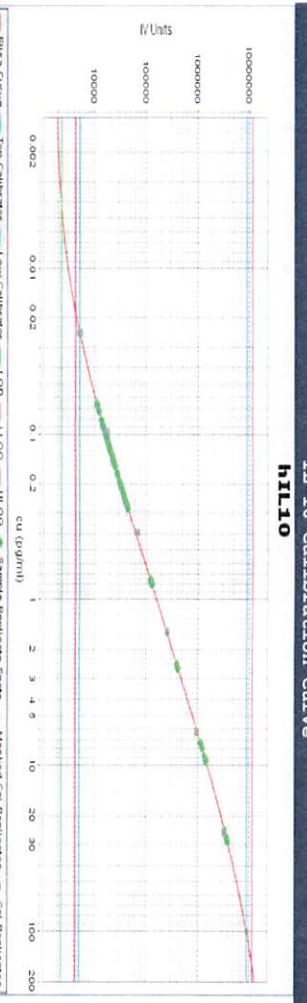
hIL-22 Calibration Curve

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



hTNF-α Calibration Curve

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



hIL-10 Calibration Curve

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

