

Product Number:	85-0329
Lot Number:	319779
Expiration:	10-May-2023
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	319779P
Biotinylated Antibody Rgt	98-0329	319779B
Sample Diluent 21	1863329	319709
IL-12p70 Calibrator	18910034	319156
IL-1β Calibrator	18910040	319564
IL-4 Calibrator	1858452	319458
IL-5 Calibrator	1858462	319223
IFNγ Calibrator	1858260	319711
IL-6 Calibrator	18910044	319484
IL-8 Calibrator	1858454	319575
IL-22 Calibrator	18910036	319176
TNFα 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

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).



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)
Panel 1	318935	139.5	133.1 - 204.3
Panel 2	318935	43.88	37.09 - 66.72
Panel 3	318935	15.08	13.48 - 21.97
Panel 4	318935	4.170	3.812 - 6.073
Panel 5	318935	1.330	1.188 - 2.091

IL-1β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	22.46	12.49 - 32.68
Panel 2	318935	7.183	4.122 - 9.719
Panel 3	318935	2.277	1.349 - 3.475
Panel 4	318935	0.631	0.418 - 0.968
Panel 5	318935	0.210	0.133 - 0.345

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	52.19	51.07 - 77.41
Panel 2	318935	16.94	14.17 - 24.05
Panel 3	318935	5.913	5.217 - 8.887
Panel 4	318935	1.824	1.455 - 2.760
Panel 5	318935	0.623	0.418 - 1.250

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	65.35	43.59 - 101.9
Panel 2	318935	20.42	13.43 - 32.50
Panel 3	318935	7.263	4.948 - 11.09
Panel 4	318935	1.932	1.317 - 3.402
Panel 5	318935	0.743	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	12.54	8.190 - 14.53
Panel 2	318935	3.957	2.393 - 4.771
Panel 3	318935	1.399	0.855 - 1.720
Panel 4	318935	0.415	0.277 - 0.494
Panel 5	318935	0.171	0.114 - 0.221

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	184.5	142.9 - 214.3
Panel 2	318935	86.55	62.87 - 99.85
Panel 3	318935	17.75	13.37 - 20.05
Panel 4	318935	7.845	5.951 - 8.927
Panel 5	318935	2.818	1.992 - 3.084

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	387.5	365.0 - 604.4
Panel 2	318935	119.1	92.02 - 191.0
Panel 3	318935	36.93	32.36 - 67.44
Panel 4	318935	11.39	9.991 - 18.54
Panel 5	318935	3.716	3.355 - 6.599

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	115.2	50.55 - 160.4
Panel 2	318935	17.52	7.411 - 28.41
Panel 3	318935	5.963	2.237 - 9.205
Panel 4	318935	1.471	0.738 - 2.619
Panel 5	318935	0.519	0.226 - 0.888

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	331.0	250.8 - 453.3
Panel 2	318935	109.6	68.46 - 156.1
Panel 3	318935	37.50	25.97 - 51.08
Panel 4	318935	10.26	7.455 - 14.81
Panel 5	318935	2.883	2.339 - 4.887

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	100.6	76.56 - 126.0
Panel 2	318935	29.61	21.73 - 39.58
Panel 3	318935	10.04	7.682 - 13.72
Panel 4	318935	3.025	2.408 - 3.832
Panel 5	318935	1.108	0.837 - 1.442

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

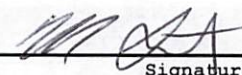
Review/Approval

Matthew Smart

Quality Control
Director

Name

Title

 2/15/2022
Signature/Date

