

Product Number:	85-0329
Lot Number:	319880
Expiration:	3-Feb-2024
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	319879P
Biotinylated Antibody Rgt	98-0329	319800B
Sample Diluent 21	1863329	319785
IL-12p70 Calibrator	18910034	318756
IL-1 β Calibrator	18910040	319564
IL-4 Calibrator	1858452	319458
IL-5 Calibrator	1858462	319838
IFN γ Calibrator	1858260	319843
IL-6 Calibrator	18910044	319484
IL-8 Calibrator	1858454	319436
IL-22 Calibrator	18910036	319176
TNF α Calibrator	1880012	319247
IL-10 Calibrator	18910043	319215
Wash Buffer (12.5X) ¹	1863332	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	319839	144.2	107.1 - 185.7
Panel 2	319839	45.48	33.39 - 50.13
Panel 3	319839	14.34	12.90 - 19.65
Panel 4	319839	5.174	2.913 - 5.970
Panel 5	319839	1.850	0.913 - 1.932

IL-1 β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	13.41	10.92 - 19.91
Panel 2	319839	4.175	3.366 - 5.402
Panel 3	319839	1.353	1.258 - 2.215
Panel 4	319839	0.535	0.312 - 0.698
Panel 5	319839	0.208	0.120 - 0.257

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	59.97	48.37 - 93.27
Panel 2	319839	18.16	15.53 - 23.53
Panel 3	319839	6.192	6.040 - 10.18
Panel 4	319839	2.543	1.688 - 3.269
Panel 5	319839	1.138	0.635 - 1.242

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	45.67	25.73 - 82.96
Panel 2	319839	14.55	8.338 - 21.52
Panel 3	319839	4.508	2.981 - 9.260
Panel 4	319839	2.167	1.095 - 2.573
Panel 5	319839	0.812	0.414 - 0.973

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	319839	9.421	5.142 - 12.43
Panel 2	319839	3.022	1.584 - 4.084
Panel 3	319839	1.061	0.609 - 1.848
Panel 4	319839	0.468	0.193 - 0.572
Panel 5	319839	0.248	0.069 - 0.296

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	227.5	132.5 - 318.7
Panel 2	319839	54.67	37.34 - 72.84
Panel 3	319839	17.63	13.54 - 30.82
Panel 4	319839	8.288	4.151 - 11.39
Panel 5	319839	3.461	1.669 - 4.008

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	207.4	146.6 - 414.1
Panel 2	319839	60.47	42.19 - 107.0
Panel 3	319839	18.66	17.48 - 41.71
Panel 4	319839	7.719	5.706 - 12.96
Panel 5	319839	3.199	2.437 - 4.581

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	80.02	40.01 - 100.4
Panel 2	319839	12.43	7.892 - 14.30
Panel 3	319839	5.910	3.903 - 7.620
Panel 4	319839	1.099	0.699 - 1.525
Panel 5	319839	0.365	0.176 - 0.454

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	60.54	33.25 - 108.2
Panel 2	319839	18.43	11.31 - 27.9
Panel 3	319839	6.090	4.165 - 11.41
Panel 4	319839	2.032	1.149 - 2.987
Panel 5	319839	0.845	0.525 - 1.051

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	37.56	26.13 - 64.76
Panel 2	319839	12.04	8.347 - 18.55
Panel 3	319839	4.038	3.175 - 8.174
Panel 4	319839	1.661	1.081 - 2.369
Panel 5	319839	0.780	0.435 - 1.031

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

Review/Approval

Tony Iem

Quality Control
Supervisor

 21NOV2021

Name

Title

Signature/Date



