

Product Number:	85-0329
Lot Number:	319791
Expiration:	10-May-2023
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	319790P
Biotinylated Antibody Rgt	98-0329	319779B
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	319223
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	177.6	133.1	204.3
Panel 2	318935	52.29	37.09	66.72
Panel 3	318935	16.03	13.48	21.97
Panel 4	318935	4.557	3.812	6.073
Panel 5	318935	1.152	1.188	2.091
IL-1β				
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Panel 1	318935	30.21	12.49	32.68
Panel 2	318935	8.229	4.122	9.719
Panel 3	318935	2.451	1.349	3.475
Panel 4	318935	0.691	0.418	0.968
Panel 5	318935	0.179	0.133	0.345
IL-4				
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Panel 1	318935	51.90	51.07	77.41
Panel 2	318935	15.95	14.17	24.05
Panel 3	318935	5.490	5.217	8.887
Panel 4	318935	1.638	1.455	2.760
Panel 5	318935	0.508	0.418	1.250
IL-5				
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Panel 1	318935	71.01	43.59	101.9
Panel 2	318935	20.76	13.43	32.50
Panel 3	318935	6.634	4.948	11.09
Panel 4	318935	2.066	1.317	3.402
Panel 5	318935	0.571	0.473	1.224

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