

Product Number:	85-0329
Lot Number:	319959
Expiration:	3-Feb-2024
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	319836P
Biotinylated Antibody Rgt	98-0329	319836B
Sample Diluent 21	1863329	319785
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	319843
IL-6 Calibrator	18910044	319484
IL-8 Calibrator	1858454	319575
IL-22 Calibrator	18910036	319459
TNF α Calibrator	1880012	319247
IL-10 Calibrator	18910043	319842
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 Cytokine1 Kit Instructions (85A-0329).

Release Materials

IL-12p70

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	159.4	101.2 - 193.3
Panel 2	319839	45.11	32.24 - 52.86
Panel 3	319839	17.50	12.45 - 20.78
Panel 4	319839	4.650	2.632 - 6.594
Panel 5	319839	1.512	0.764 - 2.272

IL-1 β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	17.31	10.37 - 21.04
Panel 2	319839	4.715	3.172 - 5.834
Panel 3	319839	1.723	1.201 - 2.331
Panel 4	319839	0.516	0.279 - 0.773
Panel 5	319839	0.206	0.096 - 0.301

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	75.66	45.70 - 99.67
Panel 2	319839	19.81	13.41 - 27.61
Panel 3	319839	7.580	5.612 - 11.10
Panel 4	319839	2.170	1.515 - 3.657
Panel 5	319839	0.877	0.559 - 1.400

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	48.73	27.65 - 80.96
Panel 2	319839	11.47	8.208 - 22.84
Panel 3	319839	5.248	2.848 - 9.877
Panel 4	319839	1.341	1.085 - 2.550
Panel 5	319839	0.513	0.320 - 1.170

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	11.46	5.622 - 14.47
Panel 2	319839	3.442	1.692 - 4.770
Panel 3	319839	1.435	0.731 - 2.000
Panel 4	319839	0.457	0.215 - 0.692
Panel 5	319839	0.220	0.071 - 0.348

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	239.4	141.9 - 312.1
Panel 2	319839	57.45	39.07 - 73.71
Panel 3	319839	21.25	14.61 - 30.61
Panel 4	319839	7.335	4.493 - 11.67
Panel 5	319839	2.914	1.407 - 4.765

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	289.6	164.0 - 399.7
Panel 2	319839	78.04	45.67 - 107.4
Panel 3	319839	27.44	18.96 - 41.09
Panel 4	319839	8.418	6.105 - 13.08
Panel 5	319839	3.288	2.260 - 5.125

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	64.30	36.22 - 101.9
Panel 2	319839	8.359	6.825 - 14.71
Panel 3	319839	4.463	3.227 - 7.790
Panel 4	319839	0.777	0.646 - 1.509
Panel 5	319839	0.224	0.180 - 0.453

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	86.88	37.50 - 105.1
Panel 2	319839	22.27	12.23 - 27.53
Panel 3	319839	8.240	4.430 - 11.09
Panel 4	319839	1.970	1.213 - 2.913
Panel 5	319839	0.744	0.543 - 1.056

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	52.68	28.83 - 63.34
Panel 2	319839	13.20	9.000 - 18.28
Panel 3	319839	5.192	3.397 - 7.891
Panel 4	319839	1.572	1.141 - 2.386
Panel 5	319839	0.671	0.457 - 1.051

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

Review/Approval

Tony Iem

Quality Control
Supervisor

 16 FEB 2013

Name

Title

Signature/Date



