

Product Number:	85-0329
Lot Number:	319930
Expiration:	3-Feb-2024
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-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	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	148.2	101.2 - 193.3
Panel 2	319839	46.71	32.24 - 52.86
Panel 3	319839	18.62	12.45 - 20.78
Panel 4	319839	5.279	2.632 - 6.594
Panel 5	319839	2.233	0.764 - 2.272

IL-1 β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	17.58	10.37 - 21.04
Panel 2	319839	5.449	3.172 - 5.834
Panel 3	319839	2.087	1.201 - 2.331
Panel 4	319839	0.619	0.279 - 0.773
Panel 5	319839	0.248	0.096 - 0.301

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	78.91	45.70 - 99.67
Panel 2	319839	26.46	13.41 - 27.61
Panel 3	319839	10.31	5.612 - 11.10
Panel 4	319839	3.353	1.515 - 3.657
Panel 5	319839	1.185	0.559 - 1.400

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	44.47	27.65 - 80.96
Panel 2	319839	15.03	8.208 - 22.84
Panel 3	319839	6.044	2.848 - 9.877
Panel 4	319839	1.829	1.085 - 2.550
Panel 5	319839	0.888	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	11.76	5.622 - 14.47
Panel 2	319839	3.956	1.692 - 4.770
Panel 3	319839	1.664	0.731 - 2.000
Panel 4	319839	0.600	0.215 - 0.692
Panel 5	319839	0.299	0.071 - 0.348

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	235.0	141.9 - 312.1
Panel 2	319839	63.92	39.07 - 73.71
Panel 3	319839	25.44	14.61 - 30.61
Panel 4	319839	9.883	4.493 - 11.67
Panel 5	319839	4.660	1.407 - 4.765

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	287.3	164.0 - 399.7
Panel 2	319839	88.12	45.67 - 107.4
Panel 3	319839	32.59	18.96 - 41.09
Panel 4	319839	10.69	6.105 - 13.08
Panel 5	319839	4.487	2.260 - 5.125

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	86.00	36.22 - 101.9
Panel 2	319839	11.83	6.825 - 14.71
Panel 3	319839	5.571	3.227 - 7.790
Panel 4	319839	1.109	0.646 - 1.509
Panel 5	319839	0.385	0.180 - 0.453

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	74.95	37.50 - 105.1
Panel 2	319839	21.20	12.23 - 27.53
Panel 3	319839	8.156	4.430 - 11.09
Panel 4	319839	2.038	1.213 - 2.913
Panel 5	319839	0.808	0.543 - 1.056

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	319839	47.49	28.83 - 63.34
Panel 2	319839	14.82	9.000 - 18.28
Panel 3	319839	5.379	3.397 - 7.891
Panel 4	319839	1.882	1.141 - 2.386
Panel 5	319839	0.849	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

 27 FEB 2021

Name

Title

Signature/Date



