

Product Number:	85-0329
Lot Number:	319797
Expiration:	10-May-2023
Platform(s):	SP-X



Data below represents results generated on the SP-X Imaging and Analysis System

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	319722P
Biotinylated Antibody Rgt	98-0329	319722B
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	318886
IL-6 Calibrator	18910044	319484
IL-8 Calibrator	1858454	319436
IL-22 Calibrator	18910036	319176
TNFα Calibrator	1880012	319460
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).

Release Materials

IL-12p70

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	164.7	133.1 - 204.3
Panel 2	318935	49.27	37.09 - 66.72
Panel 3	318935	16.24	13.48 - 21.97
Panel 4	318935	4.558	3.812 - 6.073
Panel 5	318935	1.549	1.188 - 2.091

IL-1β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	24.91	12.49 - 32.68
Panel 2	318935	7.266	4.122 - 9.719
Panel 3	318935	2.604	1.349 - 3.475
Panel 4	318935	0.771	0.418 - 0.968
Panel 5	318935	0.383	0.133 - 0.345

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	58.34	51.07 - 77.41
Panel 2	318935	16.92	14.17 - 24.05
Panel 3	318935	6.194	5.217 - 8.887
Panel 4	318935	1.837	1.455 - 2.760
Panel 5	318935	0.742	0.418 - 1.250

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	66.85	43.59 - 101.9
Panel 2	318935	19.72	13.43 - 32.50
Panel 3	318935	7.032	4.948 - 11.09
Panel 4	318935	2.116	1.317 - 3.402
Panel 5	318935	0.863	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	12.19	8.190 - 14.53
Panel 2	318935	3.943	2.393 - 4.771
Panel 3	318935	1.387	0.855 - 1.720
Panel 4	318935	0.437	0.277 - 0.494
Panel 5	318935	0.228	0.114 - 0.221

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	193.0	142.9 - 214.3
Panel 2	318935	85.09	62.87 - 99.85
Panel 3	318935	16.75	13.37 - 20.05
Panel 4	318935	7.290	5.951 - 8.927
Panel 5	318935	2.901	1.992 - 3.084

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	453.3	365.0 - 604.4
Panel 2	318935	132.2	92.02 - 191.0
Panel 3	318935	46.21	32.36 - 67.44
Panel 4	318935	12.20	9.991 - 18.54
Panel 5	318935	4.501	3.355 - 6.599

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	117.5	50.55 - 160.4
Panel 2	318935	17.96	7.411 - 28.41
Panel 3	318935	5.600	2.237 - 9.205
Panel 4	318935	1.682	0.738 - 2.619
Panel 5	318935	0.566	0.226 - 0.888

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	327.3	250.8 - 453.3
Panel 2	318935	106.4	68.46 - 156.1
Panel 3	318935	34.01	25.97 - 51.08
Panel 4	318935	9.38	7.455 - 14.81
Panel 5	318935	3.488	2.339 - 4.887

IL-10


	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	111.8	76.56 - 126.0
Panel 2	318935	31.46	21.73 - 39.58
Panel 3	318935	10.89	7.682 - 13.72
Panel 4	318935	3.080	2.408 - 3.832
Panel 5	318935	1.271	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

 11 Jun 2022

Name

Title

Signature/Date

