

Product Number:	85-0329
Lot Number:	319817
Expiration:	10-May-2023
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	319816P
Biotinylated Antibody Rgt	98-0329	319800B
Sample Diluent 21	1863329	319709
IL-12p70 Calibrator	18910034	319156
IL-1 $\beta$ Calibrator	18910040	319564
IL-4 Calibrator	1858452	319458
IL-5 Calibrator	1858462	318837
IFN $\gamma$ Calibrator	1858260	319711
IL-6 Calibrator	18910044	319484
IL-8 Calibrator	1858454	319575
IL-22 Calibrator	18910036	319459
TNF $\alpha$ Calibrator	1880012	319247
IL-10 Calibrator	18910043	319215
Wash Buffer (25X) <sup>1</sup>	1860055	N/A
Streptavidin-HRP Reagent <sup>1</sup>	1858263	N/A
Super Signal <sup>®</sup> Luminol* <sup>1</sup>	1858089	N/A
Super Signal <sup>®</sup> Peroxide* <sup>1</sup>	1858090	N/A



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

<sup>1</sup> Reagents are not Kit Lot Specific

\*SuperSignal<sup>®</sup> 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	148.7	133.1 - 204.3
Panel 2	318935	44.39	37.09 - 66.72
Panel 3	318935	16.20	13.48 - 21.97
Panel 4	318935	4.478	3.812 - 6.073
Panel 5	318935	1.409	1.188 - 2.091

**IL-1 $\beta$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	23.79	12.49 - 32.68
Panel 2	318935	7.254	4.122 - 9.719
Panel 3	318935	2.587	1.349 - 3.475
Panel 4	318935	0.747	0.418 - 0.968
Panel 5	318935	0.247	0.133 - 0.345

**IL-4**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	61.68	51.07 - 77.41
Panel 2	318935	18.63	14.17 - 24.05
Panel 3	318935	6.615	5.217 - 8.887
Panel 4	318935	1.964	1.455 - 2.760
Panel 5	318935	0.694	0.418 - 1.250

**IL-5**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	83.21	43.59 - 101.9
Panel 2	318935	28.12	13.43 - 32.50
Panel 3	318935	10.431	4.948 - 11.09
Panel 4	318935	3.145	1.317 - 3.402
Panel 5	318935	1.112	0.473 - 1.224

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

**Release Materials**

**IFN $\gamma$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	10.32	8.190 - 14.53
Panel 2	318935	3.302	2.393 - 4.771
Panel 3	318935	1.168	0.855 - 1.720
Panel 4	318935	0.386	0.277 - 0.494
Panel 5	318935	0.158	0.114 - 0.221

**IL-6**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	176.8	142.9 - 214.3
Panel 2	318935	80.84	62.87 - 99.85
Panel 3	318935	17.76	13.37 - 20.05
Panel 4	318935	8.029	5.951 - 8.927
Panel 5	318935	2.640	1.992 - 3.084

**IL-8**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	419.0	365.0 - 604.4
Panel 2	318935	115.5	92.02 - 191.0
Panel 3	318935	41.75	32.36 - 67.44
Panel 4	318935	12.39	9.991 - 18.54
Panel 5	318935	4.125	3.355 - 6.599

**IL-22**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	75.7	50.55 - 160.4
Panel 2	318935	12.08	7.411 - 28.41
Panel 3	318935	4.237	2.237 - 9.205
Panel 4	318935	1.163	0.738 - 2.619
Panel 5	318935	0.396	0.226 - 0.888

**TNF $\alpha$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	326.7	250.8 - 453.3
Panel 2	318935	103.8	68.46 - 156.1
Panel 3	318935	36.71	25.97 - 51.08
Panel 4	318935	11.15	7.455 - 14.81
Panel 5	318935	3.238	2.339 - 4.887

**IL-10**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	90.1	76.56 - 126.0
Panel 2	318935	27.69	21.73 - 39.58
Panel 3	318935	10.04	7.682 - 13.72
Panel 4	318935	3.016	2.408 - 3.832
Panel 5	318935	1.069	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

Name

Title

*M. Smart* 10 Aug 2022  
Signature/Date



