

Product Number:	85-0329
Lot Number:	319276
Expiration:	28-Jan-2022
Platform(s) :	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	319276P
Biotinylated Antibody Rgt	98-0329	319276B
Sample Diluent 21	1863329	319289
IL-12p70 Calibrator	18910034	318756
IL-1 $\beta$ Calibrator	18910040	318728
IL-4 Calibrator	1858452	318580
IL-5 Calibrator	1858462	318837
IFN $\gamma$ Calibrator	1858260	318886
IL-6 Calibrator	18910044	318724
IL-8 Calibrator	1858454	318500
IL-22 Calibrator	18910036	318813
TNF $\alpha$ Calibrator	1880012	318883
IL-10 Calibrator	18910043	318755
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	206.5	140.6 - 239.0
Panel 2	318935	56.76	44.36 - 73.70
Panel 3	318935	17.70	15.74 - 23.61
Panel 4	318935	5.106	4.400 - 6.601
Panel 5	318935	1.679	1.243 - 2.397

**IL-1 $\beta$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	23.79	16.60 - 33.00
Panel 2	318935	6.643	5.060 - 10.28
Panel 3	318935	2.114	1.780 - 3.430
Panel 4	318935	0.658	0.496 - 1.030
Panel 5	318935	0.239	0.151 - 0.367

**IL-4**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	62.99	53.27 - 82.46
Panel 2	318935	17.42	16.02 - 24.04
Panel 3	318935	6.060	6.011 - 9.016
Panel 4	318935	2.306	1.729 - 2.776
Panel 5	318935	1.136	0.434 - 1.312

**IL-5**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	100.60	54.07 - 121.1
Panel 2	318935	26.16	18.39 - 34.26
Panel 3	318935	8.47	6.909 - 11.42
Panel 4	318935	2.401	2.007 - 3.306
Panel 5	318935	1.012	0.615 - 1.271

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	13.05	9.152 - 15.15
Panel 2	318935	3.770	2.766 - 5.014
Panel 3	318935	1.252	1.021 - 1.670
Panel 4	318935	0.394	0.320 - 0.515
Panel 5	318935	0.169	0.142 - 0.223

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	196.1	140.8 - 249.6
Panel 2	318935	79.62	68.32 - 107.0
Panel 3	318935	14.90	14.11 - 21.96
Panel 4	318935	7.305	6.379 - 9.569
Panel 5	318935	2.657	1.951 - 3.380

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	554.8	287.6 - 812.3
Panel 2	318935	136.6	99.01 - 213.4
Panel 3	318935	46.73	33.87 - 79.01
Panel 4	318935	14.59	10.72 - 20.44
Panel 5	318935	5.123	3.706 - 7.504

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	150.5	73.64 - 264.5
Panel 2	318935	21.87	13.35 - 41.63
Panel 3	318935	7.096	6.048 - 11.67
Panel 4	318935	2.050	1.025 - 3.238
Panel 5	318935	0.700	0.478 - 1.052

TNF $\alpha$

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	378.7	213.6 - 599.8
Panel 2	318935	116.1	81.39 - 179.7
Panel 3	318935	37.72	30.07 - 60.48
Panel 4	318935	10.87	9.413 - 16.87
Panel 5	318935	3.671	2.187 - 6.149

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318935	117.7	86.85 - 130.3
Panel 2	318935	34.24	25.35 - 40.54
Panel 3	318935	10.42	9.246 - 13.87
Panel 4	318935	3.137	2.649 - 4.032
Panel 5	318935	1.141	0.890 - 1.546

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

Review/Approval

Shivani Goel

Quality Control  
Manager

*Shivani Goel 01 Jun 2021*

Name

Title

Signature/Date