

Product Number:	85-0329
Lot Number:	318855
Expiration:	10/25/2021 ²
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	318797P
Biotinylated Antibody Rgt	98-0329	318797B
Sample Diluent 21 ¹	1863329	N/A
IL-12p70 Calibrator	18910034	318629
IL-1β Calibrator	18910040	318728
IL-4 Calibrator	1858452	318480
IL-5 Calibrator	1858462	318465
IFNγ Calibrator	1858260	318631
IL-6 Calibrator	18910044	318724
IL-8 Calibrator	1858454	318342
IL-22 Calibrator	18910036	318408
TNFα Calibrator	1880012	318436
IL-10 Calibrator	18910043	318630
Wash Buffer (25X) ¹	18242300	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.

² If Wash Buffer lot # 1860055 is used kit expiration is 18Sep2021. If Wash buffer lot # 18242300 is used kit expiration is 25Oct 2021

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	318641	174.4	148.9	- 240.7
Panel 2	318641	57.96	51.04	- 81.68
Panel 3	318641	20.50	17.95	- 26.93
Panel 4	318641	6.972	5.525	- 9.670
Panel 5	318641	2.215	1.672	- 3.404

IL-1β

	Lot	Result (pg/mL)	Mean	Range (pg/mL)
Panel 1	318641	19.69	19.65	- 29.48
Panel 2	318641	6.343	6.475	- 10.24
Panel 3	318641	2.358	2.123	- 3.489
Panel 4	318641	0.747	0.649	- 1.279
Panel 5	318641	0.246	0.223	- 0.460

IL-4

	Lot	Result (pg/mL)	Mean	Range (pg/mL)
Panel 1	318641	78.91	66.79	- 105.1
Panel 2	318641	24.95	20.41	- 35.27
Panel 3	318641	7.794	6.722	- 12.50
Panel 4	318641	2.522	2.180	- 4.624
Panel 5	318641	0.815	0.815	- 1.597

IL-5

	Lot	Result (pg/mL)	Mean	Range (pg/mL)
Panel 1	318641	68.00	50.55	- 76.98
Panel 2	318641	22.38	16.93	- 25.93
Panel 3	318641	7.808	5.920	- 9.298
Panel 4	318641	2.692	1.943	- 3.293
Panel 5	318641	0.896	0.700	- 1.238

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	318641	12.41	8.541	- 12.86
Panel 2	318641	4.363	2.940	- 4.681
Panel 3	318641	1.536	0.994	- 1.686
Panel 4	318641	0.565	0.348	- 0.647
Panel 5	318641	0.219	0.138	- 0.270

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Panel 1	318641	121.7	97.18	- 184.2
Panel 2	318641	3.616	2.780	- 7.433
Panel 3	318641	1.277	0.928	- 2.564
Panel 4	318641	0.476	0.330	- 1.033

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Panel 1	318641	395.9	372.9	- 559.4
Panel 2	318641	135.8	109.4	- 188.6
Panel 3	318641	44.19	36.67	- 64.61
Panel 4	318641	15.27	11.28	- 24.27
Panel 5	318641	5.169	3.867	- 7.799

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Panel 1	318641	22.71	14.24	- 33.29
Panel 2	318641	5.999	4.355	- 9.706
Panel 3	318641	2.047	1.528	- 3.283
Panel 4	318641	0.701	0.462	- 1.223
Panel 5	318641	0.208	0.156	- 0.428

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Panel 1	318641	67.11	59.79	- 100.9
Panel 2	318641	22.99	20.44	- 35.53
Panel 3	318641	7.447	6.956	- 12.29
Panel 4	318641	2.605	2.303	- 4.312
Panel 5	318641	0.888	0.547	- 1.821

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Panel 1	318641	69.33	48.59	- 73.50
Panel 2	318641	22.96	15.68	- 27.33
Panel 3	318641	8.042	5.500	- 9.707
Panel 4	318641	2.954	2.011	- 3.603
Panel 5	318641	1.022	0.706	- 1.339

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

Review/Approval

Patrizia Stadler Associate
Director of QC

P. Stadler

05 Aug 2020

Name

Title

Signature/Date