

Product Number:	85-0329
Lot Number:	318733
Expiration:	18-Sep-2021
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	318627P
Biotinylated Antibody Rgt	98-0329	318718B
Sample Diluent 2l ¹	1863329	N/A
IL-12p70 Calibrator	18910034	318571
IL-1 β Calibrator	18910040	318673
IL-4 Calibrator	1858452	318480
IL-5 Calibrator	1858462	318465
IFN γ Calibrator	1858260	318538
IL-6 Calibrator	18910044	318583
IL-8 Calibrator	1858454	318342
IL-22 Calibrator	18910036	318408
TNF α Calibrator	1880012	318436
IL-10 Calibrator	18910043	318630
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



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	318641	181.9	74.90 - 312.8
Panel 2	318641	64.55	25.27 - 105.7
Panel 3	318641	22.53	8.118 - 35.35
Panel 4	318641	7.460	2.411 - 12.41
Panel 5	318641	2.440	0.828 - 4.244

IL-1 β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	27.41	9.353 - 36.93
Panel 2	318641	9.540	3.051 - 12.44
Panel 3	318641	3.100	0.942 - 4.185
Panel 4	318641	1.100	0.288 - 1.506
Panel 5	318641	0.400	0.053 - 0.563

IL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	80.07	31.71 - 139.3
Panel 2	318641	27.34	10.51 - 46.13
Panel 3	318641	9.118	3.155 - 16.82
Panel 4	318641	3.398	1.070 - 6.178
Panel 5	318641	1.271	0.416 - 2.046

IL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	318641	58.93	23.48 - 97.00
Panel 2	318641	20.04	7.797 - 32.93
Panel 3	318641	6.634	2.673 - 11.72
Panel 4	318641	2.296	0.879 - 4.207
Panel 5	318641	0.770	0.306 - 1.549

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	10.67	4.077	-	16.34
Panel 2	318641	3.753	1.401	-	5.834
Panel 3	318641	1.364	0.475	-	2.058
Panel 4	318641	0.510	0.163	-	0.798
Panel 5	318641	0.200	0.067	-	0.338

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)		
Panel 1	318641	138.8	53.84	-	235.8
Panel 2	318641	5.098	1.795	-	9.227
Panel 3	318641	1.700	0.585	-	3.138
Panel 4	318641	0.643	0.219	-	1.262
Panel 5	318641	0.222	0.214	-	0.437

IL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)		
Panel 1	318641	476.3	175.2	-	729.0
Panel 2	318641	146.5	53.79	-	240.7
Panel 3	318641	49.12	18.28	-	84.37
Panel 4	318641	16.55	5.861	-	29.62
Panel 5	318641	5.400	0.997	-	10.30

IL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)		
Panel 1	318641	23.35	8.132	-	41.07
Panel 2	318641	7.093	2.348	-	11.98
Panel 3	318641	2.257	0.776	-	4.049
Panel 4	318641	0.795	0.274	-	1.445
Panel 5	318641	0.244	0.080	-	0.499

TNF α

	Lot	Result (pg/mL)	Mean Range (pg/mL)		
Panel 1	318641	87.93	28.83	-	127.9
Panel 2	318641	31.86	10.37	-	45.07
Panel 3	318641	10.69	3.411	-	15.41
Panel 4	318641	3.347	1.115	-	5.442
Panel 5	318641	1.381	0.402	-	2.035

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)		
Panel 1	318641	65.36	22.96	-	85.32
Panel 2	318641	23.64	7.491	-	30.02
Panel 3	318641	8.435	2.644	-	10.62
Panel 4	318641	2.950	0.964	-	3.944
Panel 5	318641	1.123	0.339	-	1.486

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

21 May 2020

Name

Title

Signature/Date