

Product Number:	100-0487
Lot Number:	503898
Expiration:	1-May-2024
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	197-0487	319567P
Biotinylated Antibody Rgt	198-0487	319567B
Sample Diluent 22	1880098	312104A
IL-2 Calibrator	1880099	319311
Wash Buffer (12.5X) ¹	1863332	N/A
Streptavidin-HRP Reagent ¹	1858263	N/A
Super Signal [®] Luminol ^{1,2}	1858089	N/A
Super Signal [®] Peroxide ^{1,2}	1858090	N/A

¹ Reagents are not Kit Lot Specific

² 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 IL-2 Kit Instructions (KI-0223).

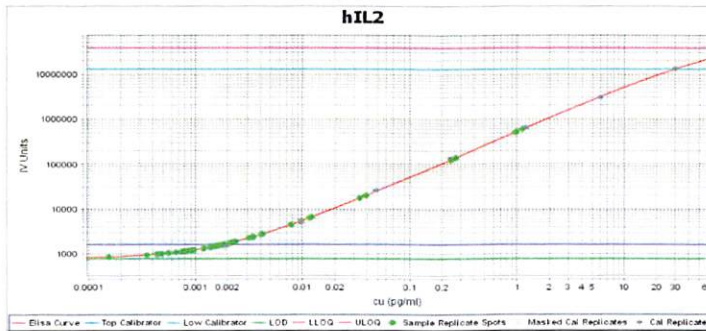
Release Materials

IL-2

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	029701	4.083	2.455 - 4.880
Panel 2	029701	1.010	0.568 - 1.349
Panel 3	029701	0.149	0.085 - 0.312
Panel 4	029701	0.048	0.019 - 0.067

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

IL-2 Calibration Curve



Calibrator Levels (pg/ml)

1	30
2	6.0
3	1.20
4	0.240
5	0.048
6	0.010
7	0.002
8	0.000

Review/Approval

Robert Somogyi

Quality Control
Manager

Name

Title

Signature/Date

[Handwritten Signature] 29 Jun 2023