

Product Number:	100-0487
Lot Number:	503998
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	327707P
Biotinylated Antibody Reagent	198-0487	327604B
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

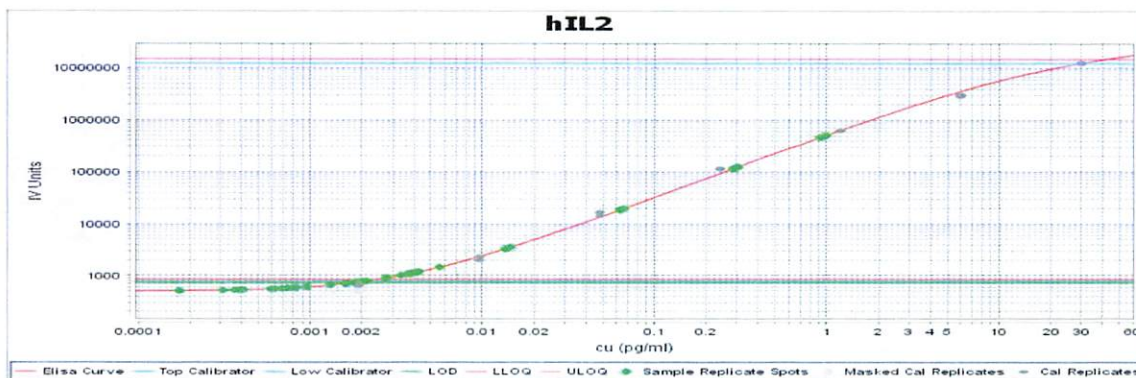
² SuperSignal® technology is protected by US Patent #6,432,662 and is a registered trademark of ThermoFisher Scientific.

All reagents are specific for use on the SP-X platform - refer to KI-0223 Corplex IL-2 Kit Instructions 100-0487.

Release Materials*			
IL-2			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	029701	3.768	2.455 - 4.880
Panel 2	029701	1.106	0.568 - 1.349
Panel 3	029701	0.258	0.085 - 0.312
Panel 4	029701	0.057	0.019 - 0.067

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

IL-2 Calibrator Curve



Calibrator Levels (pg/mL)

- 1 30.00
- 2 6.000
- 3 1.200
- 4 0.240
- 5 0.048
- 6 0.010
- 7 0.002
- 8 0.000

Review/Approval

Kathleen Cornell

Quality Control Supervisor

[Signature] 17 Nov 2023

Name

Title

Signature/Date