

Product Number:	100-0487
Lot Number:	504323
Expiration:	21-Mar-2025
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 Reagent	198-0487	420603B
Sample Diluent 22	1880098	408105
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

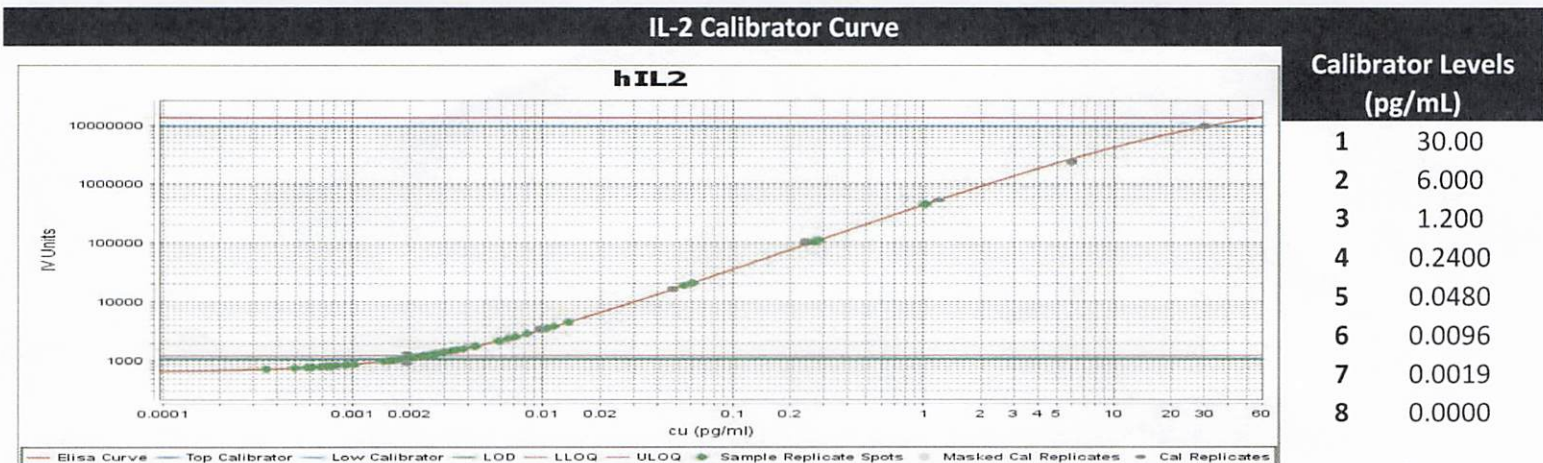
¹ Reagents are not Kit Lot Specific

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	4.089	2.660 - 5.086
Panel 2	029701	1.096	0.6062 - 1.496
Panel 3	029701	0.2357	0.1102 - 0.3334
Panel 4	029701	0.0472	0.0197 - 0.0801

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



Review/Approval	Kathleen Cornell Name	Quality Control Supervisor Title	 05 Aug 2024 Signature/Date
-----------------	--------------------------	--	-----------------------------------