

Product Number:	100-0487
Lot Number:	504203
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	327707P
Biotinylated Antibody Reagent	198-0487	327604B
Sample Diluent 22	1880098	408105
IL-2 Calibrator	1880099	319311
Wash Buffer (12.5X) <sup>1</sup>	1863332	N/A
Streptavidin-HRP Reagent <sup>1</sup>	1858263	N/A
Super Signal® Luminol <sup>1,2</sup>	1858089	N/A
Super Signal® Peroxide <sup>1,2</sup>	1858090	N/A

<sup>1</sup> Reagents are not Kit Lot Specific

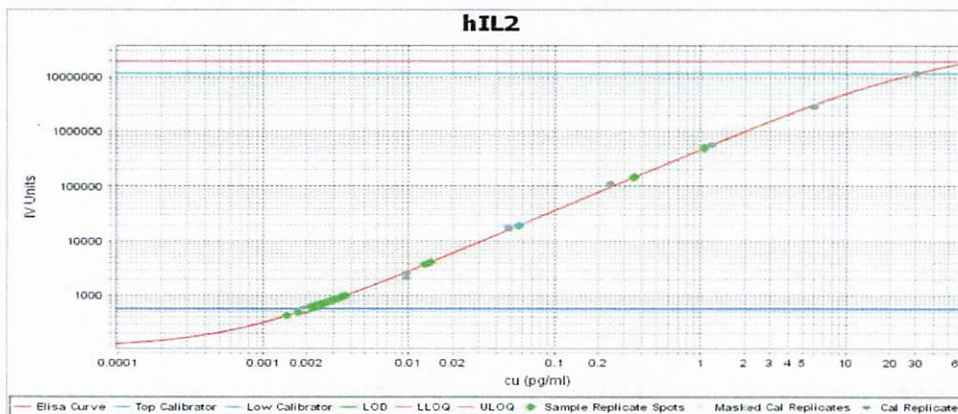
<sup>2</sup> SuperSignal™ technology is protected by US Patent #6,432,662 and is a registered trademark of ThermoFisher Scientific.

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	4.236	2.660 - 5.086
Panel 2	029701	1.403	0.6062 - 1.496
Panel 3	029701	0.2262	0.1102 - 0.3334
Panel 4	029701	0.0539	0.0197 - 0.0801

\* 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	0.2400
5	0.0480
6	0.0096
7	0.0019
8	0.0000

Review/Approval

Kathleen Cornell

Quality Control  
Supervisor

*Kathleen Cornell* 16 May 2024

Name

Title

Signature/Date