

Product Number:	100-0487
Lot Number:	536610
Expiration:	11-May-2026
Platform(s):	SP-X



Data below represents results generated on the SP-X Imaging and Analysis System

Component	Part Number	Lot Number	Control Lot Number
CorPlex Cytokine Plate	197-0487	504402P	504402AP
Biotinylated Antibody Reagent	198-0487	503402B	N/A
Sample Diluent 22	1880098	527902	N/A
IL-2 Calibrator	1880099	319311	N/A
Wash Buffer (12.5X) ¹	1863332	N/A	N/A
Streptavidin-HRP Reagent ¹	1858263	N/A	N/A
Super Signal® Luminol ^{1,2}	1858089	N/A	N/A
Super Signal® Peroxide ^{1,2}	1858090	N/A	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.

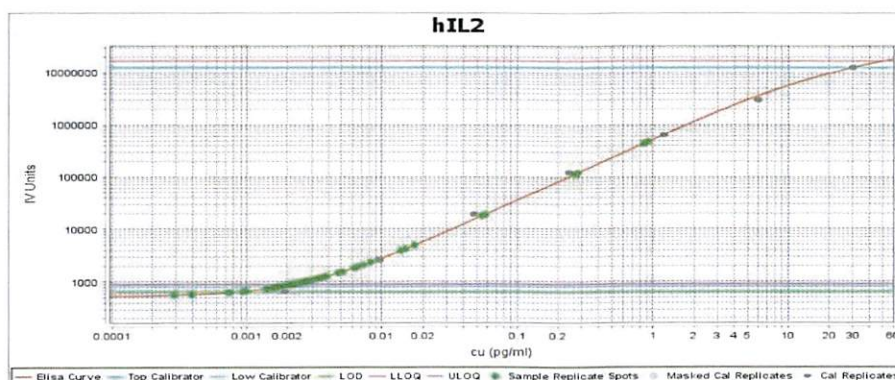
Release Materials*

IL-2

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	029701	3.550	2.660 - 5.086
Panel 2	029701	1.081	0.6062 - 1.496
Panel 3	029701	0.2243	0.1102 - 0.3334
Panel 4	029701	0.0612	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

Dhruvi Patel *QA Specialist II* *SDP* / 11 Dec 2025

Name

Title

Signature/Date

Quanterix Corporation
900 Middlesex Turnpike, Building 1, Billerica, 01821
techsupport@quantexix.com

©2025 Quanterix Corp. Simoa® is a registered trademark of Quanterix Corp.

For research use only. Not for use in diagnostic procedures.

Release

COA-0067 02