

Product Number:	100-0487
Lot Number:	319301
Expiration:	20-Jul-2022
Platform(s):	SP-X



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

Component	Part Number	Lot Number
CorPlex Cytokine Plate	197-0487	319301P
Biotinylated Antibody Rgt	198-0487	319301B
Sample Diluent 22	1880098	319304
IL-2 Calibrator	1880099	319311
Wash Buffer (25X) ¹	1860055	N/A
Streptavidin-HRP Reagent ¹	1858263	N/A
Super Signal [®] Luminol* ¹	1858089	N/A
Super Signal [™] Peroxide* ¹	1858090	N/A
Streptavidin-HRP Reagent ¹	1858263	N/A

¹ Reagents are not Kit Lot Specific

*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 Cytokine 2 Kit Instructions (85A-0401).

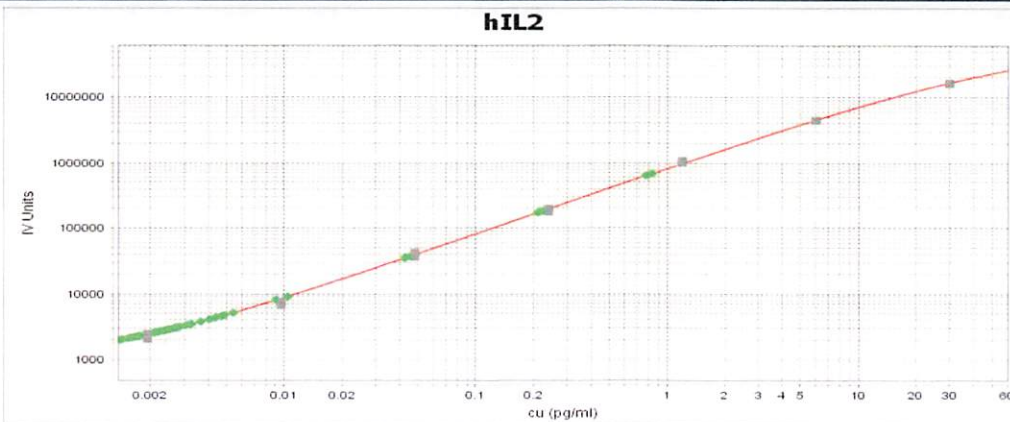
Release Materials

IL-2

Panel	Lot	Result (pg/mL)	Mean Range (pg/mL)
Panel 1	029701	3.220	2.557 - 5.086
Panel 2	029701	0.864	0.644 - 1.390
Panel 3	029701	0.175	0.119 - 0.317
Panel 4	029701	0.038	0.027 - 0.068

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

Shivani Goel

Quality Control
Manager

Shivani Goel 09 Jun 2021

Name

Title

Signature/Date