

Product Number:	85-0348
Lot Number:	504337
Expiration:	5-Jul-2025
Platform(s):	SP-X



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

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	419801P
Biotinylated Antibody Reagent	98-0329	403307B
Sample Diluent 21	1863329	403020
IL-1β Calibrator	18910040	319564
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

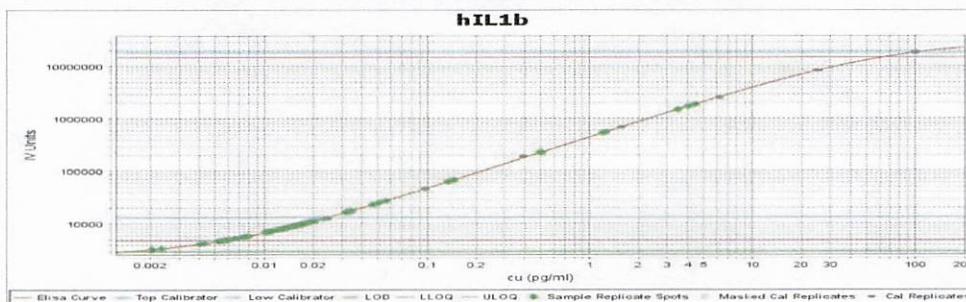
² SuperSignal™ technology is protected by US Patent #6,432,662 and is a registered trademark of ThermoFisher Scientific.

All reagents are specific for use on the SP-X platform - refer to KI-0286 Simoa Complex Cytokine Panel 1 IL-1b Instruction Book 85-0348.

Release Materials*			
IL-1β			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	15.9	9.79 - 20.4
Panel 2	319839	4.90	2.85 - 5.89
Panel 3	319839	2.00	1.06 - 2.31
Panel 4	319839	0.554	0.238 - 0.745
Panel 5	319839	0.204	0.084 - 0.274

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

IL-1β Calibrator Curve



Calibrator Levels (pg/mL)

A	100
B	25.0
C	6.25
D	1.56
E	0.391
F	0.098
G	0.024
H	0.000

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title

Signature/Date

[Signature] 13 Aug 2024