

Product Number:	100-0217
Lot Number:	319015
Expiration:	27-Nov-2021
Platform(s):	SP-X



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

Component	Part Number	Lot Number
IL-6 Plate	197-0217	319015P
Biotinylated Antibody Reagent	198-0217	319015B
IL-6 Calibrator	1861002	318801
Sample Diluent 16	1863325	N/A
Wash Buffer (25X) ¹	1860055	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

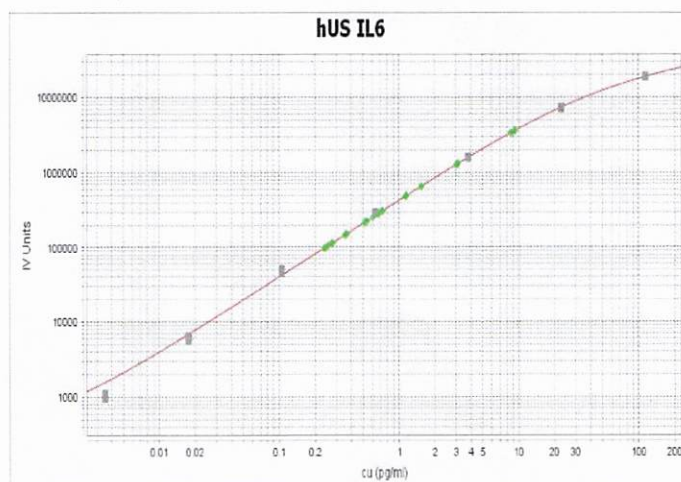
¹ Reagents are not Kit Lot Specific

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 the Simoa® IL-6 Developer Kit Instructions (KI-0196).

Release Materials		
IL-6		
	Result (pg/mL)	Spec. Range (pg/mL)
Control 2	9.01	4.53 - 21.5
Control 3	3.07	1.59 - 5.51

IL-6 Calibrator Curve



Calibrator Levels (pg/mL)	
1	114
2	22.8
3	3.80
4	0.633
5	0.106
6	0.0176
7	0.00350
0	0.000

Review/Approval

Kyle Giroux

Name

Senior Document Control Specialist, QA

Title

Kyle Giroux 01 Apr 2025

Signature/Date