

Product Number:	100-0217
Lot Number:	504722
Expiration:	See individual components below.
Platform(s):	SP-X

Component	Part Number	Lot Number	Expiration Date
IL-6 Plate	197-0217	524713P	4-Sep-30
Biotinylated Antibody Reagent	198-0217	525115B	8-Sep-28
IL-6 Calibrator	1861002	426401	24-Sep-29
Sample Diluent 16	1863325	N/A	
Wash Buffer (12.5X)	1863332	N/A ¹	
Streptavidin-HRP Reagent	1858263	N/A ¹	
Super Signal® Luminol	1858089	N/A ^{1,2}	
Super Signal® Peroxide	1858090	N/A ^{1,2}	

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.

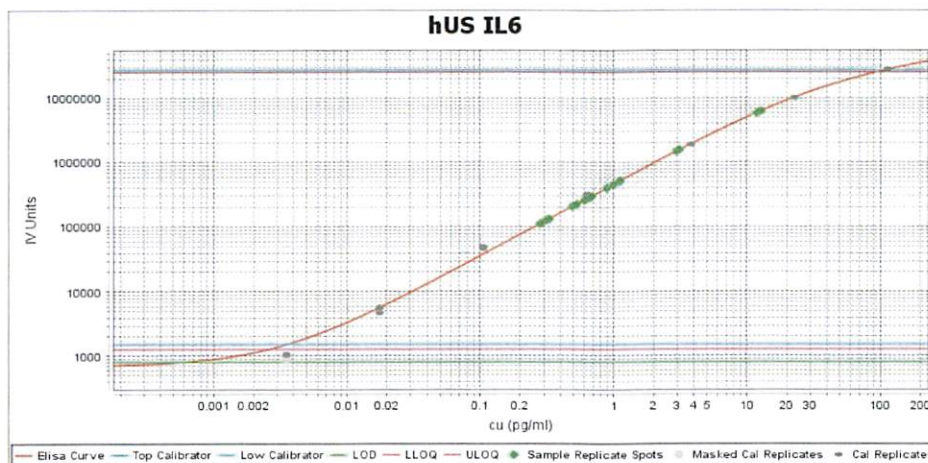


Data below represents results generated on the SP-X™ Imaging and Analysis System

All reagents are specific for use on the SP-X™ platform - refer to the Simoa® IL-6 Developer Kit Instructions (KI-0196).

Release Materials			* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
IL-6			
	Result (pg/mL)	Spec. Range (pg/mL)	
Control 2	12.3	4.53 - 21.5	
Control 3	3.05	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

Dhruvi Patel *QA Specialist II, QA* *DP* *18Sep2025*
Name Title Signature/Date