

Product Number:	100-0217
Lot Number:	319916
Expiration:	3-Feb-2024
Platform(s):	SP-X

IL-6 Plate	197-0217	319916P
Biotinylated Antibody Reagent	198-0217	319916B
IL-6 Calibrator	1861002	319912
Sample Diluent 16	1863325	N/A
Wash Buffer (12.5X)¹	1863332	N/A
Streptavidin-HRP Reagent¹	1858263	N/A
Super Signal® Luminol^{1,2}	1858089	N/A
Super Signal® Peroxide^{1,2}	1858090	N/A

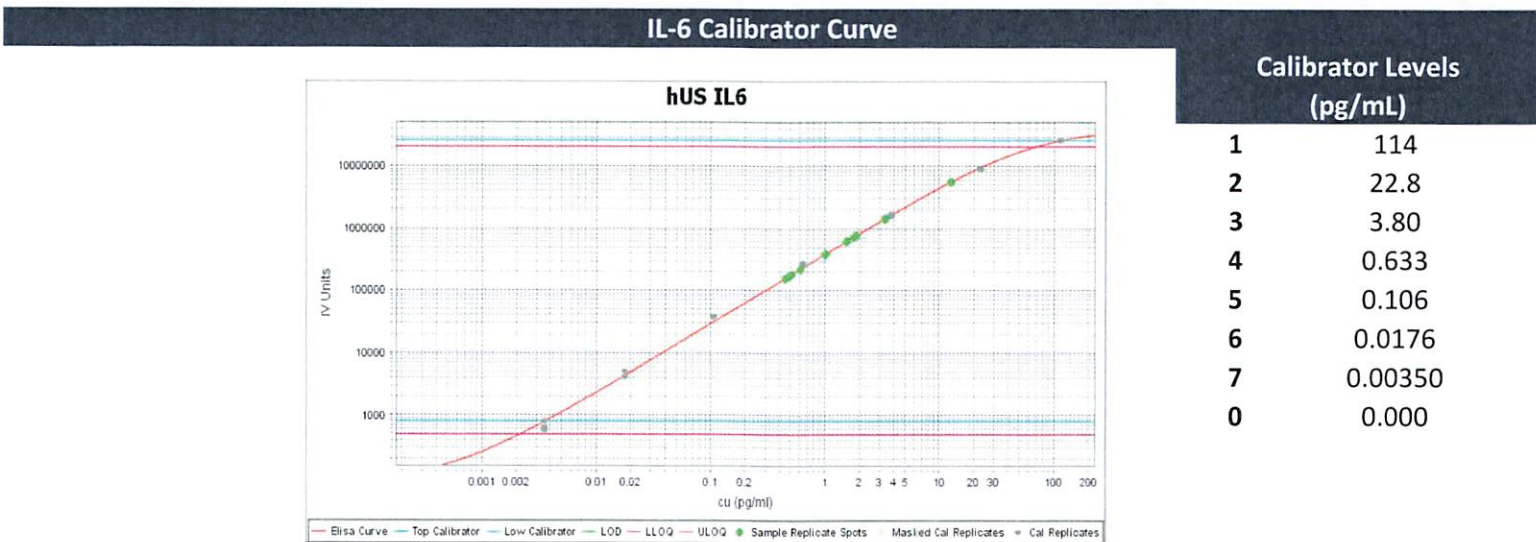


¹ Reagents are not Kit Lot Specific

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	12.6	4.53 - 21.5
Control 3	3.40	1.59 - 5.51



Review/Approval

Sonia Khillan

Quality Control Manager

Sonia Khillan 28 MAR 2023

Name

Title

Signature/Date

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

©Quanterix, Inc. Simoa® is a registered trademark of Quanterix, Inc.
 For research use only. Not for use in diagnostic procedures.