

Product Number:	100-0461
Lot Number:	504470C
Expiration:	15-Oct-2025
Platform(s):	SP-X



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

Component	Part Number	Lot Number
Homebrew Plate	197-0461	431304P
Diluent A ¹	1880090	419006
Lyophilized Homebrew Tag 1 ¹	18910024	513654000002540
EZ-Link NHS-PEG4-Biotin ¹	100348	RM2024100000053
Sulfo-SMCC ¹	1870039	513654000002781
PBS pH 7.2 Pack ¹	1870037	513654000002815
1M Tris HCl pH 7.4 ¹	1870036	429710
Microclimate Plate Lid ¹	1825107	N/A
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.

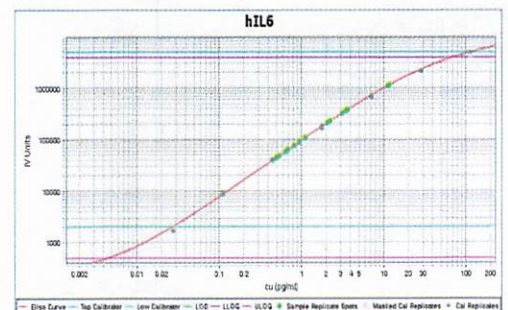
All reagents are specific for use on the SP-X platform - refer to Simoa Planar Array Kit Instructions (KI-0029).

The results below represent data from the functional testing of the homebrew plate using hIL-6:

Release Materials*		
	Result (pg/mL)	Spec. Range (pg/mL)
Control 2	11.6	4.53 - 21.5
Control 3	3.57	1.59 - 5.51

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

Curve Fit Analysis		
Concentration (pg/ml)	Backfit	% Recovery
114	115	100.8%
28.5	26.1	91.7%
7.13	6.79	95.3%
1.78	1.77	99.2%
0.445	0.465	104.4%
0.111	0.115	103.2%
0.028	0.023	82.7%
0	0.000	N/A



Review/Approval

Kyle Giroux

Senior Document Control Specialist, QA

[Signature] 25 Feb 2025

Name

Title

Signature/Date