

Product Number:	101622
Lot Number:	502687
Expiration:	7-Oct-2021
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	101624	034502
Detector Reagent	101623	034607
SBG Reagent	101625	034610
Sample Diluent	103732	028120
Calibrator Diluent	101627	034914
Calibrator Concentrate	100652	006408
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



Data below represents results generated on the Simoa™ HD-1 Analyzer.

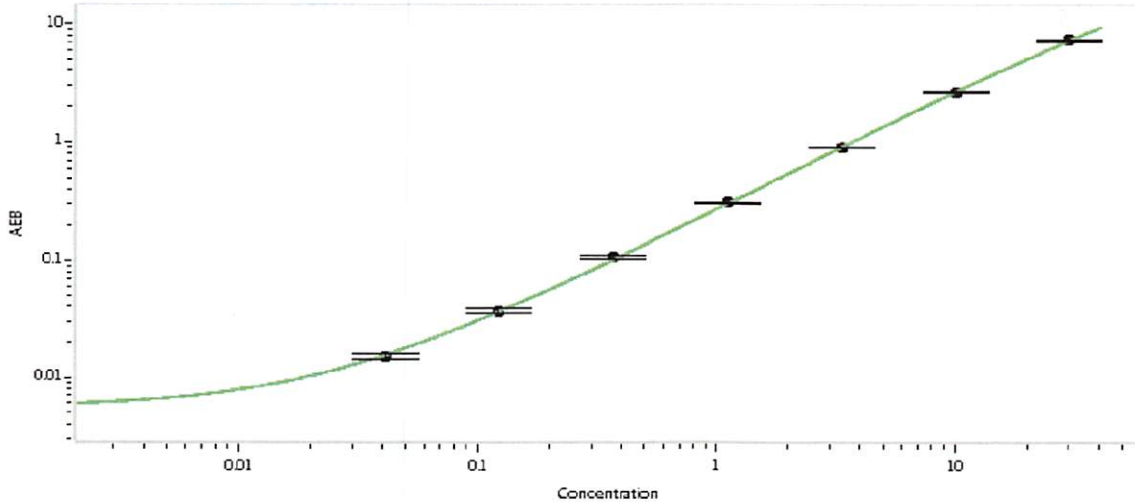
Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	901702	1.04	0.944-1.42
Control 2	901702	15.6	15.4-23.1
Panel 1	029301	44.8	39.8-59.7
Panel 2	029401	4.94	4.35-6.53
Panel 3	029301	1.37	1.22-1.82

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

Calibrator Concentrate Value Assignment (pg/mL)	675
---	-----

Calibration Curve

2020-12-21 IL-6 Kit (Lot 502657) Release QC (CU-122) Curve 1



Calibrator Levels (pg/mL)

A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 13 Jan 2021
Signature/Date