

Product Number:	101622
Lot Number:	502657
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	021203
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific



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

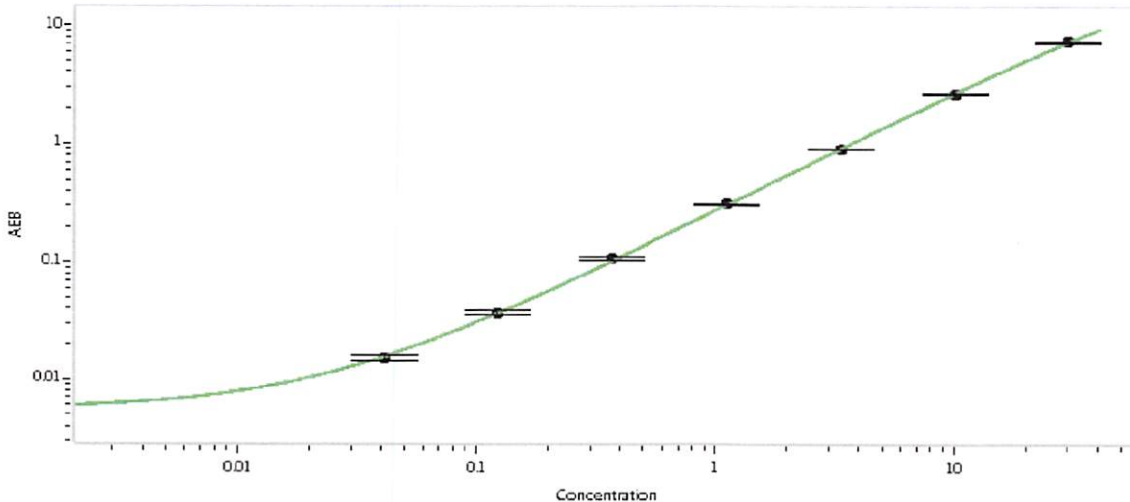
Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	901702	1.16	0.944-1.42
Control 2	901702	16.0	15.4-23.1
Panel 1	029301	47.4	39.8-59.7
Panel 2	029401	5.58	4.35-6.53
Panel 3	029301	1.58	1.22-1.82

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

Calibrator Concentrate Value Assignment (pg/mL)	606
---	-----

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

QC Supervisor

*[Signature]* 31 Dec 2020

Name

Title

Signature/Date