

Product Number:	102732
Lot Number:	502809
Expiration:	5-Mar-2022
Platform(s):	HD-1/HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	102722	107711
Detector Reagent	102724	107902
SBG Reagent	102726	108216
Sample Diluent	102728	106609
Calibrator Diluent	102731	106421
Calibrator Concentrate	100658	106608
RGP Reagent	103159	N/A <sup>1</sup>

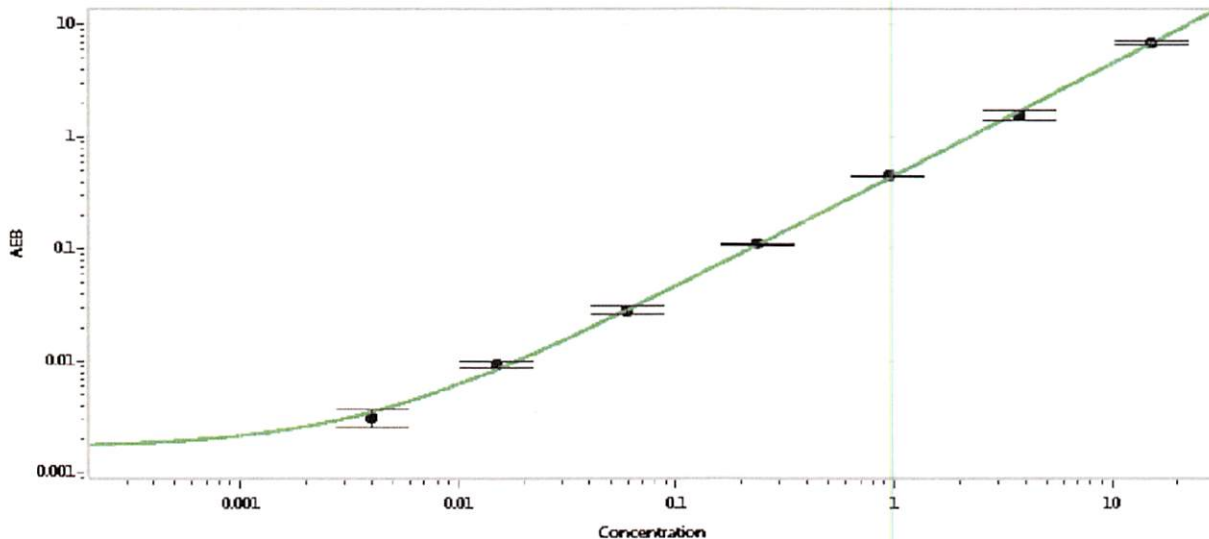
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	909502	0.133	0.093-0.167
Control 2	909502	3.68	2.83-4.48
Panel 1	911401	0.028	0.017-0.050
Panel 3	911403	1.83	1.18-2.02
Panel 4	911404	6.15	3.80-7.25

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

Calibrator Concentrate Value Assignment (pg/mL)	444
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.004
C	0.015
D	0.059
E	0.234
F	0.938
G	3.75
H	15.0

Review/Approval

Junpil Park

QC Supervisor

*[Signature]* 16 Apr 2021

Name

Title

Signature/Date