

Product Number:	102860
Lot Number:	503025
Expiration:	30-Dec-2021
Platform(s):	HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	102852	118106
Detector Reagent	102854	121036
SBG Reagent	102855	118604
Sample Diluent	102856	118507
Calibrator Diluent	102859	118509
Calibrator Concentrate	102857	121037
RGP Reagent	103159	N/A <sup>1</sup>

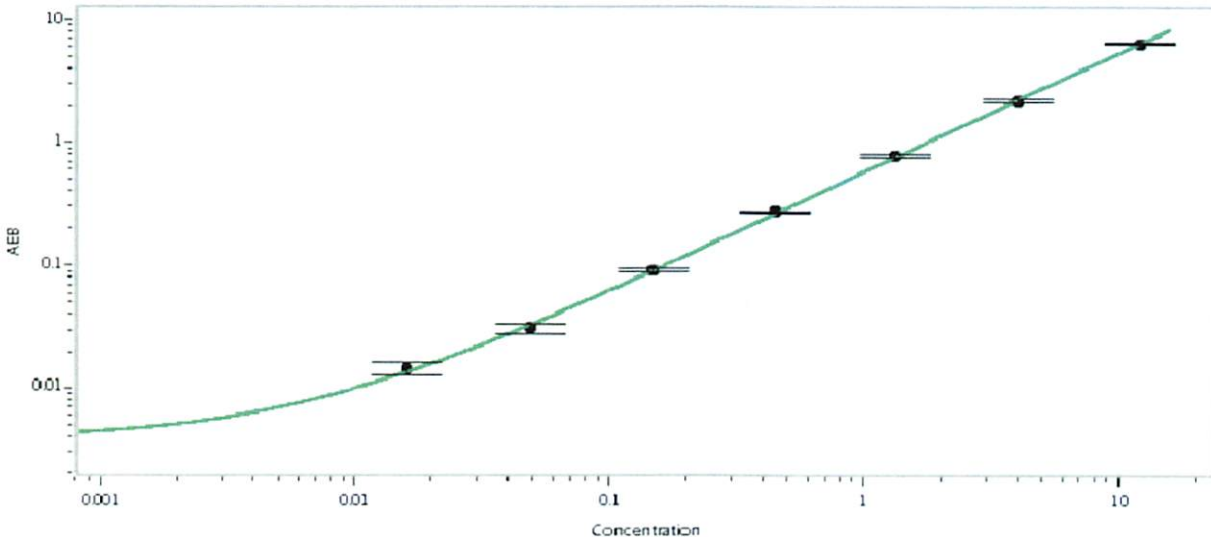
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	121506	0.310	0.229-0.344
Control 2	121506	5.00	3.87-5.81
Panel 1	033802	0.180	0.123-0.241
Panel 2	033802	2.27	1.70-2.65
Panel 3	033802	10.2	8.00-12.0

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

Calibrator Concentrate Value Assignment (pg/mL)	396
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.016
C	0.049
D	0.148
E	0.444
F	1.33
G	4.00
H	12.0

Review/Approval

Junpil Park

QC Supervisor

30 Aug 2021

Name

Title

Signature/Date