

Product Number:	102732
Lot Number:	502936
Expiration:	20-Apr-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	112505
Detector Reagent	102724	112610
SBG Reagent	102726	112611
Sample Diluent	102728	111021
Calibrator Diluent	102731	111022
Calibrator Concentrate	100658	112612
RGP Reagent	103159	N/A <sup>1</sup>

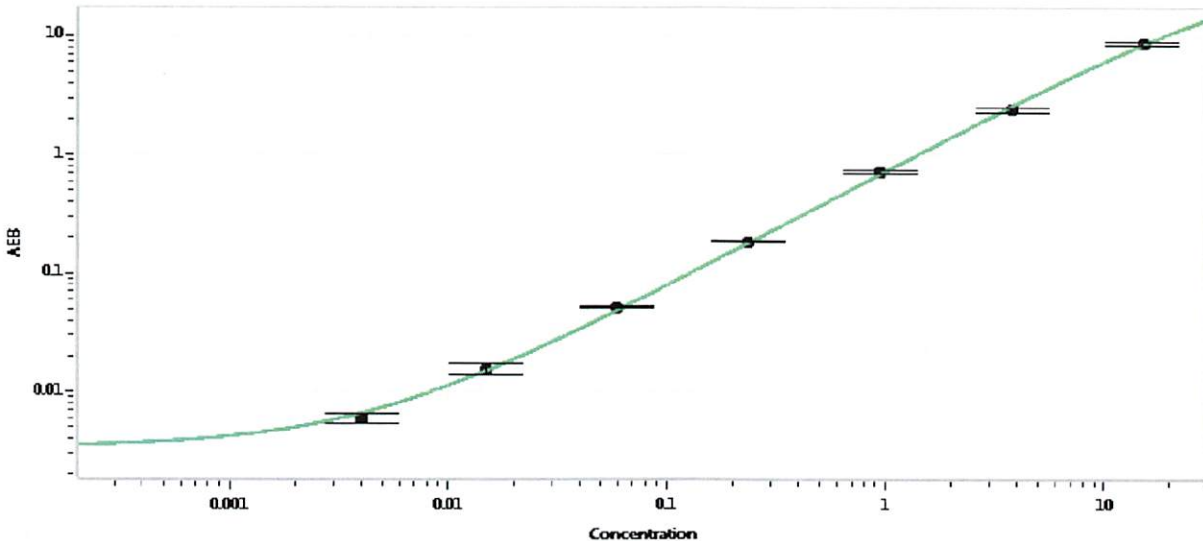
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	114108	0.076	0.067-0.159
Control 2	114108	2.57	1.91-3.81
Panel 2	112714	0.028	0.010-0.049
Panel 3	112714	1.27	1.02-1.87
Panel 4	112714	4.59	2.79-7.70

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

Calibrator Concentrate Value Assignment (pg/mL)	470
---	-----

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 Jun 2021

Name

Title

Signature/Date