

Product Number:	102860
Lot Number:	502926
Expiration:	12-Nov-2021
Platform(s):	HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	102852	113202
Detector Reagent	102854	113308
SBG Reagent	102855	113309
Sample Diluent	102856	113310
Calibrator Diluent	102859	113311
Calibrator Concentrate	102857	008305
RGP Reagent	103159	N/A ¹

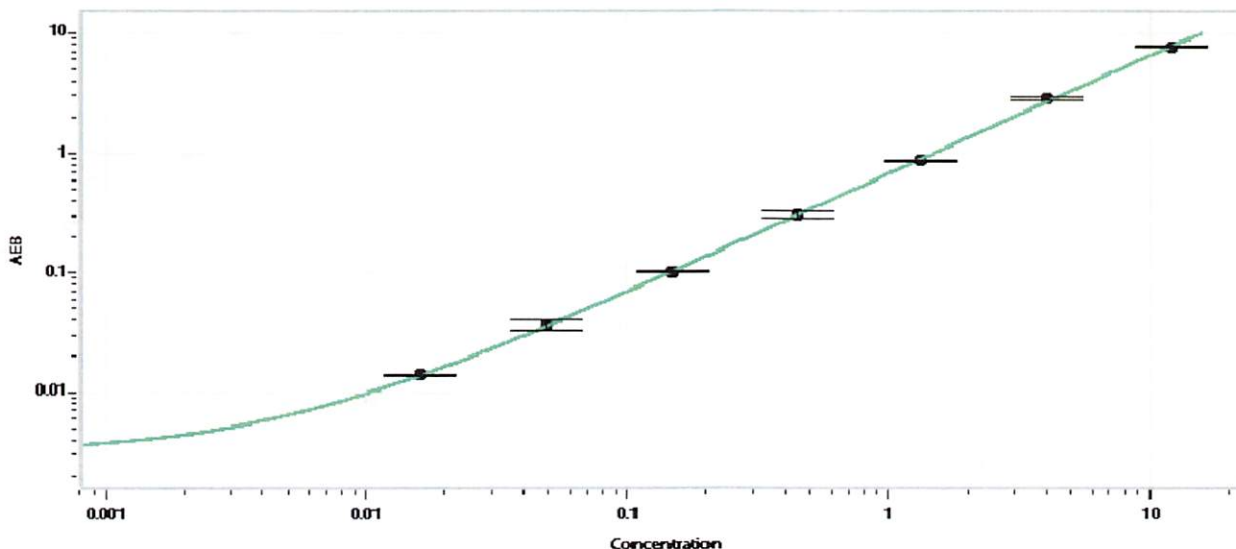
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	008305	0.241	0.180-0.300
Control 2	008305	4.86	3.60-6.00
Panel 1	033802	0.190	0.118-0.249
Panel 2	033802	2.00	1.64-2.65
Panel 3	033802	10.4	7.98-12.0

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

Calibrator Concentrate Value Assignment (pg/mL)	386
---	-----

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

[Signature] 03 Jun 2021

Name

Title

Signature/Date