

Product Number:	101599
Lot Number:	502852
Expiration:	22-Mar-2022
Platform(s):	HD-1/HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101601	109101
Detector Reagent	102012	109113
SBG Reagent	101602	109206
Sample Diluent	103886	109207
Calibrator Diluent	102014	108130
Calibrator Concentrate	100778	108911
RGP Reagent	103159	N/A ¹

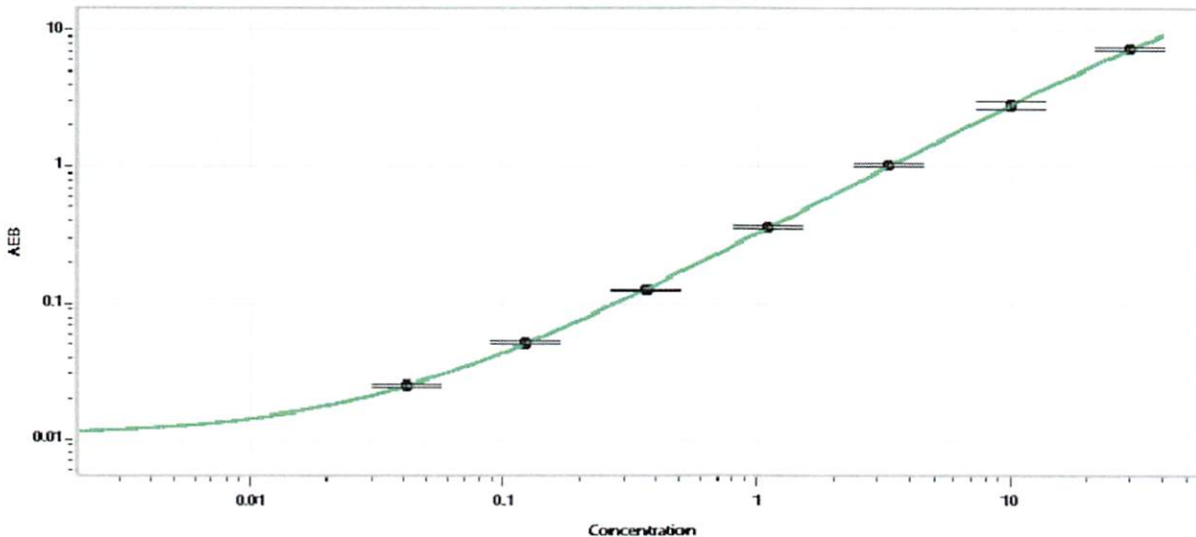
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108214	0.250	0.223-0.374
Control 2	109503	40.4	29.4-46.0
Panel 3	034908	10.2	7.76-12.7
Panel 4	034908	32.5	20.9-45.7

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

Calibrator Concentrate Value Assignment (pg/mL)	617
---	-----

Calibration Curve



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
Name

QC Supervisor
Title

[Signature]
Signature/Date

19 Apr 2021