

Product Number:	101580
Lot Number:	503011
Expiration:	26-May-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	101582	117108
Detector Reagent	101581	117111
SBG Reagent	101583	117219
Sample Diluent	103735	117114
Calibrator Diluent	101585	117209
Calibrator Concentrate	100655	117208
RGP Reagent	103159	N/A ¹

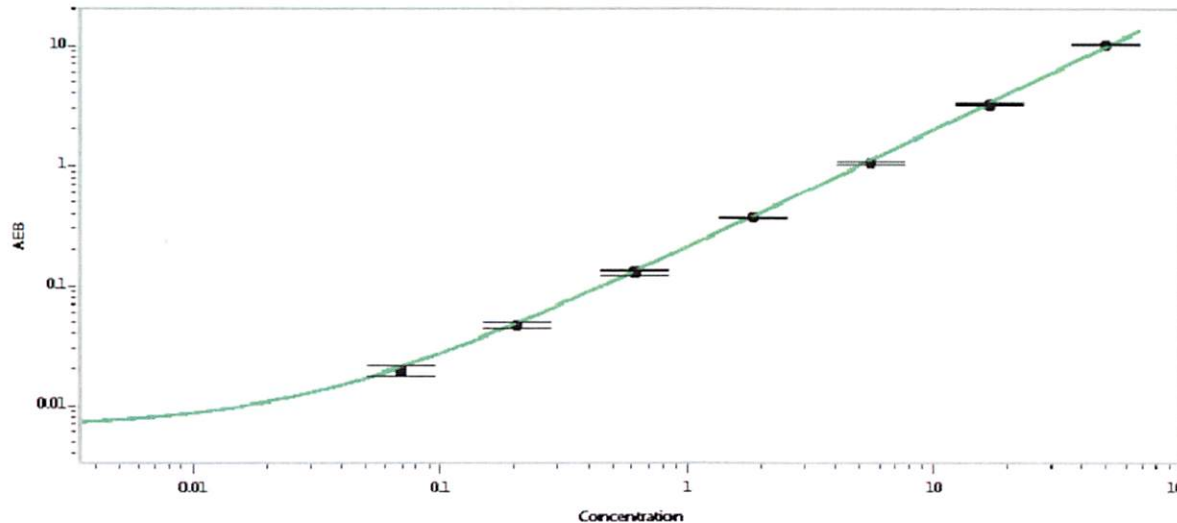
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	024603	1.35	0.762-2.33
Control 2	024603	24.2	19.1-34.9
Panel 2	115317	1.98	0.908-4.22
Panel 3	116301	27.8	18.4-39.3

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

Calibrator Concentrate Value Assignment (pg/mL)	1240
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.069
C	0.206
D	0.617
E	1.85
F	5.56
G	16.7
H	50.0

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 20 Jun 2021
Signature/Date