

Product Number:	103337
Lot Number:	502601
Expiration:	21-Oct-2021
Platform(s):	HD-1/HD-X & SR-X



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

Component	Item Number	Lot Number
Bead Reagent	103334	029504
Detector Reagent	101091	030701
SBG Reagent	101101	030710
Sample Diluent	103336	030709
Calibrator Diluent	101095	029716
Calibrator Concentrate	101093	032802
RGP Reagent	103159	N/A ¹

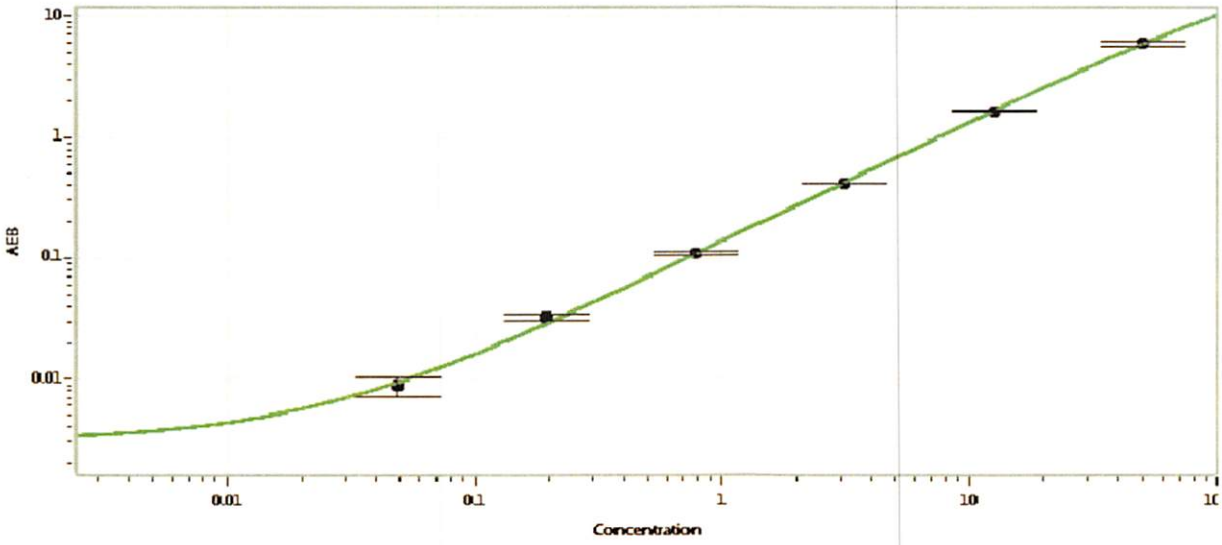
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	911603	1.92	1.19-2.10
Control 2	911603	38.4	20.2-43.8
Panel 1	024004	3.78	1.22-4.72
Panel 2	024004	4.21	3.07-4.61
Panel 3	024004	52.9	37.4-56.1

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

Calibrator Concentrate Value Assignment (pg/mL)	1,750
---	-------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.049
C	0.195
D	0.781
E	3.13
F	12.5
G	50.0

Review/Approval

Junpil Park
Name
QC Supervisor
Title

[Signature] 04 Dec 2020
Signature/Date