

Product Number:	103337
Lot Number:	502905
Expiration:	2-Apr-2022
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	111209
Detector Reagent	101091	111211
SBG Reagent	101101	111212
Sample Diluent	103336	111213
Calibrator Diluent	101095	109216
Calibrator Concentrate	101093	112008
RGP Reagent	103159	N/A ¹

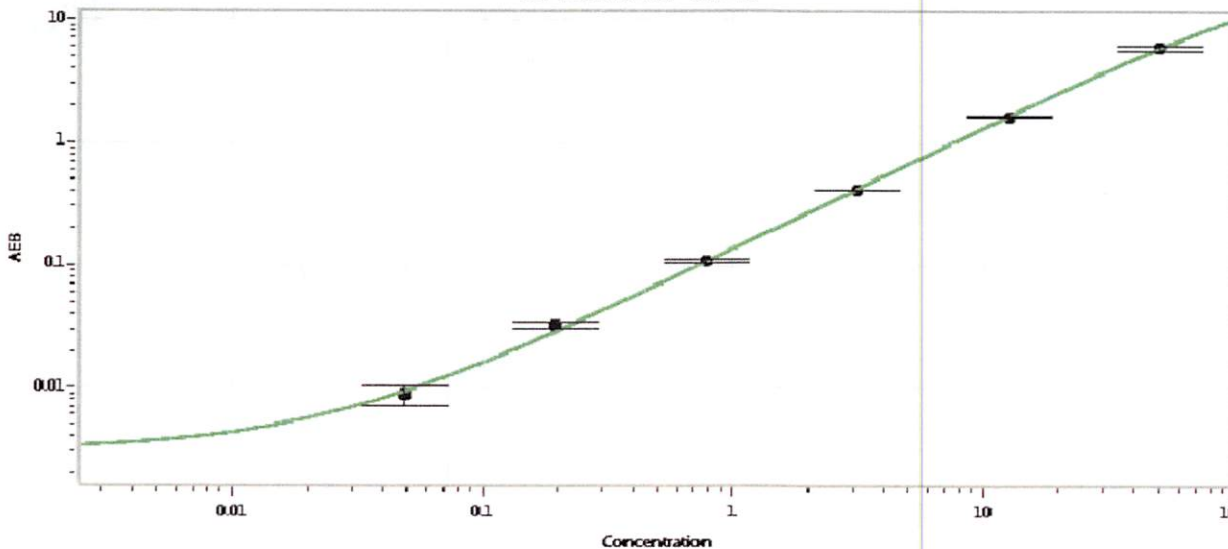
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	112008	0.959	0.750-1.25
Control 2	112008	18.2	15.0-25.0
Panel 2	024004	2.68	2.26-4.95
Panel 3	024004	37.6	32.5-57.3

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

Calibrator Concentrate Value Assignment (pg/mL)	1,277
---	-------

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/Date
24 Jun 2021