

Product Number:	101154
Lot Number:	502751
Expiration:	28-Dec-2021
Platform(s):	HD-1/HDX



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

Component	Part Number	Lot Number
Bead Reagent	101141	104317
Detector Reagent	101143	104315
SBG Reagent	101153	104331
Sample Diluent	103320	036319
Calibrator Diluent	101147	102711
Calibrator Concentrate	101145	105003
RGP Reagent	103159	N/A ¹

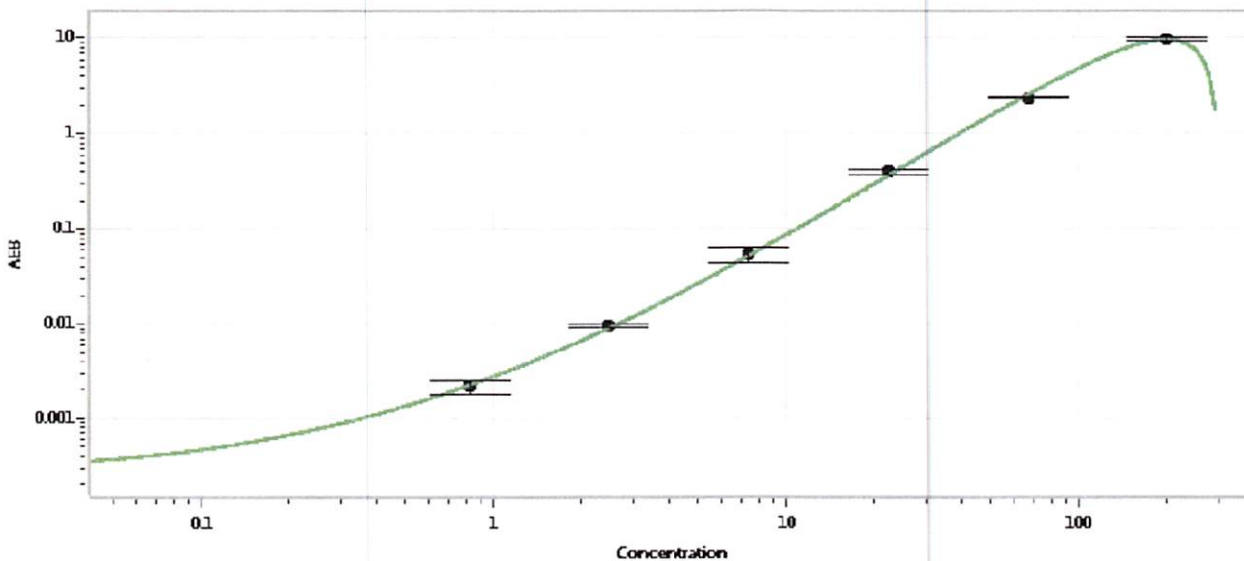
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	918402	40.4	31.7-50.4
Control 2	918402	303	251-461
Panel 1	909901	81.2	66.3-112
Panel 2	909902	70.4	58.0-102
Panel 3	909903	235	161-367
Panel 4	909904	220	166-314

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

Calibrator Concentrate Value Assignment (pg/mL)	3841
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.823
C	2.47
D	7.41
E	22.2
F	66.7
G	200

Review/Approval

Junpil Park

QC Supervisor

[Signature] 16 Mar 2021

Name

Title

Signature/Date