

Product Number:	103337
Lot Number:	502523
Expiration:	30-May-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	017010
Detector Reagent	101091	017001
SBG Reagent	101101	017014
Sample Diluent	103336	017003
Calibrator Diluent	101095	015114
Calibrator Concentrate	101093	906801
RGP Reagent	103159	N/A <sup>1</sup>

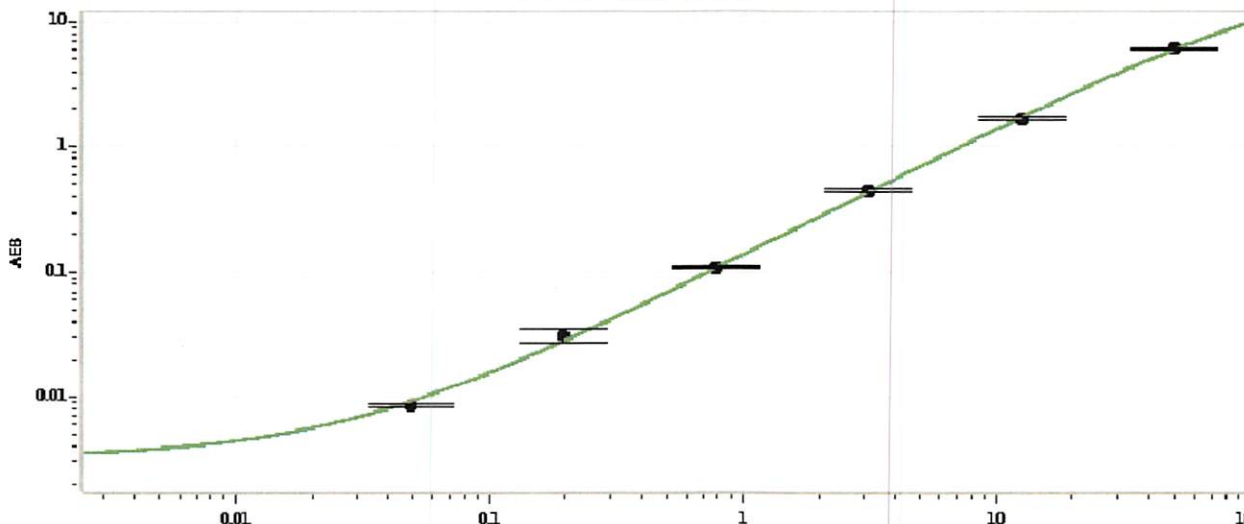
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	911603	1.82	1.25-1.88
Control 2	911603	36.6	21.4-38.9
Panel 1	024004	1.72	1.09-3.93
Panel 2	024004	4.28	2.97-4.45
Panel 3	024004	54.1	36.4-54.6

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

Calibrator Concentrate Value Assignment (pg/mL)	1,325
---	-------

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

Shivani Goel

QC Manager

Shivani Goel 14 Sep 2020

Name

Title

Signature/Date