

Product Number:	100196
Lot Number:	501958
Expiration:	30-Apr-2020
Platform(s):	HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	100391	921011
Detector Reagent	100315	921204
SBG Reagent	100316	912005
Sample Diluent	100317	912022
Calibrator Diluent	100665	921111
Calibrator Concentrate	100664	907102
RGP Reagent	103159	N/A <sup>1</sup>

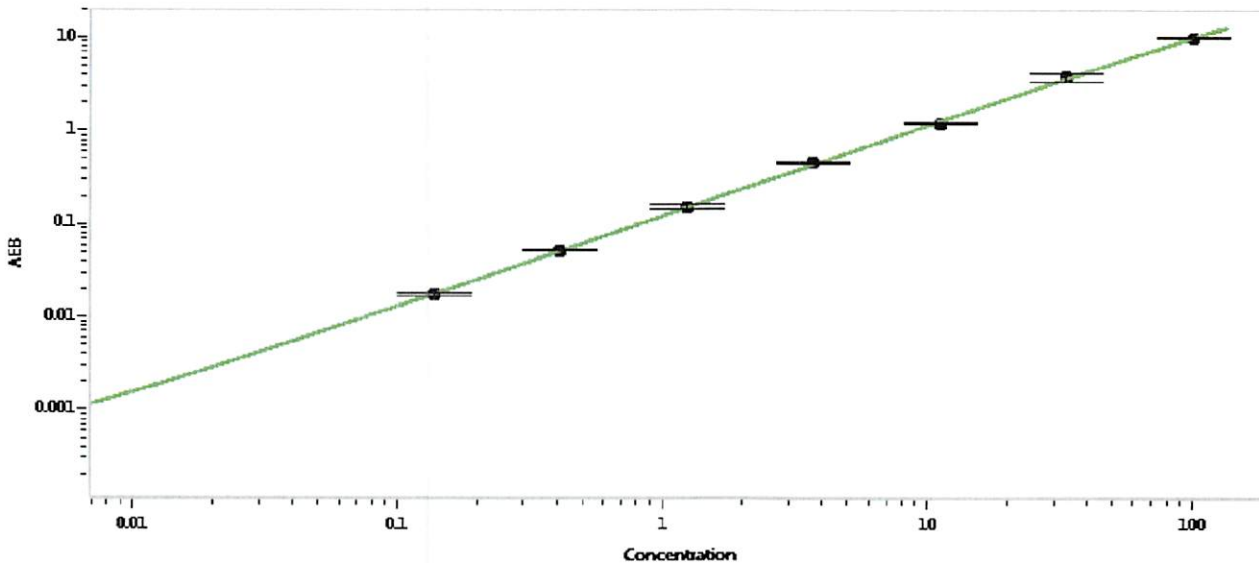
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	52521	118	73.1-125
Control 2	52521	11.4	7.56-12.1
Panel 1a	MF-062018-01	0.529	0.324-1.18
Panel 2	MF-062018-02	9.89	7.29-10.9
Panel 3	MF-062018-03	131	91.2-141

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

Calibrator Concentrate Value Assignment (pg/mL)	2290
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval

Daniel Ritchie

QC Senior Analyst

*[Signature]* / 18 Aug 2019

Name

Title

Signature/Date