

Product Number:	101622
Lot Number:	502251
Expiration:	4-Dec-2020
Platform(s):	HD-1/HD-x & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101624	005110
Detector Reagent	101623	005102
SBG Reagent	101625	004908
Sample Diluent	103203	933819
Calibrator Diluent	101627	003513
Calibrator Concentrate	100652	005601
RGP Reagent	103159	N/A <sup>1</sup>

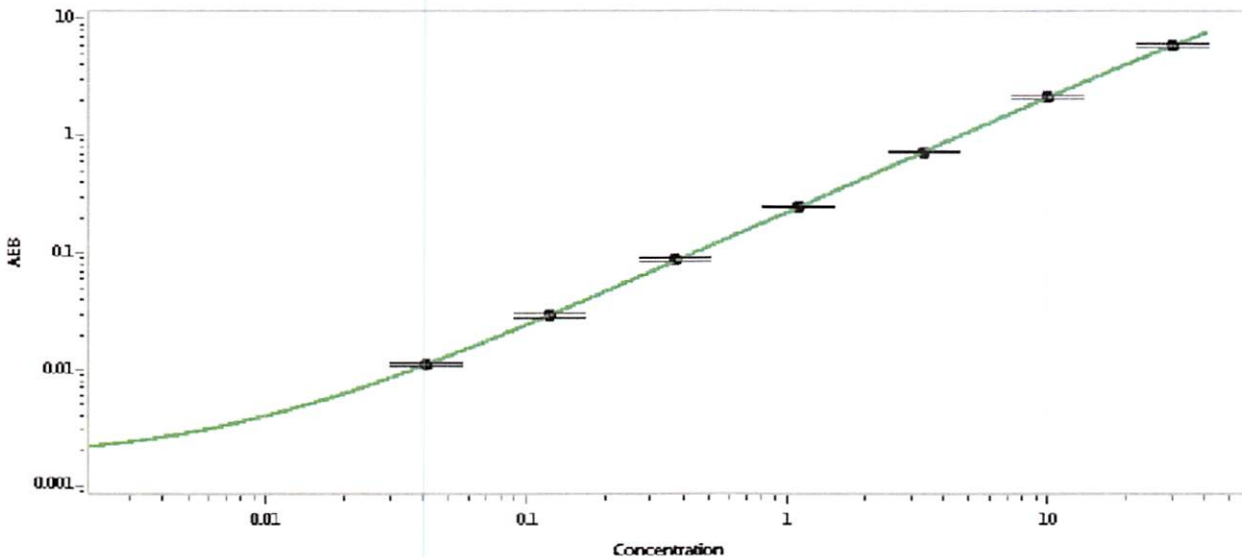
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	901702	1.10	0.944-1.42
Control 2	901702	17.3	15.4-23.1
Panel 1	MF-050918-01	51.5	41.6-67.5
Panel 2	MF-050918-02	7.01	5.03-8.84
Panel 3	MF-050918-03	0.940	0.644-1.24

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

Calibrator Concentrate Value Assignment (pg/mL)	645
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval: Patrizia Stadler Assoc. Director, QC *r. Stadler* 28 Feb 2020  
Name Title Signature/Date