

Product Number:	101605
Lot Number:	501575
Expiration:	25-Sep-2019
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	101607	900710
Detector Reagent	101606	900920
SBG Reagent	101608	833705
Sample Diluent	103200	833304
Calibrator Diluent	101610	826812
Calibrator Concentrate	100895	828001
RGP Reagent	103159	N/A <sup>1</sup>

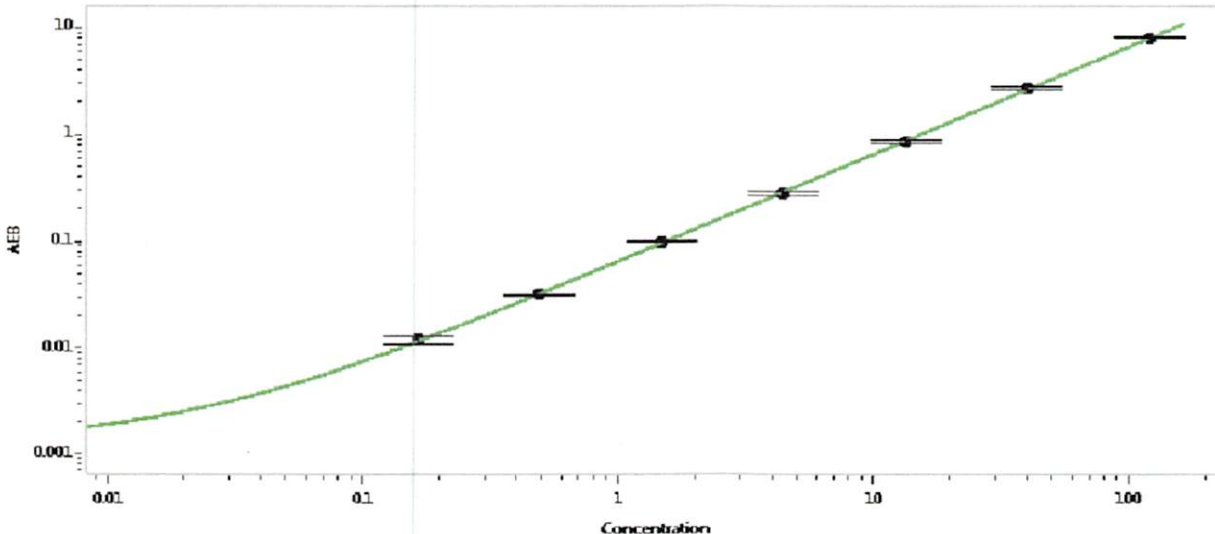
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-112515-02	5.77	3.13-7.30
Control 2	MF-112515-02	172	93.3-218
Panel 1	MF-112315-03	74.7	39.4-91.9
Panel 2	MF-112315-04	16.8	8.67-20.2
Panel 3	MF-112315-05	26.9	14.3-33.3

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

Calibrator Concentrate Value Assignment (pg/mL)	3,725
-------------------------------------------------	-------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.165
C	0.494
D	1.48
E	4.44
F	13.3
G	40.0
H	120

Review/Approval	Patrizia Stadler	Associate Director, QC	<i>P. Stadler</i>	23 Jan 2019
	Name	Title	Signature/Date	