

Product Number:	101664
Lot Number:	501559
Expiration:	8-Dec-2019
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101666	834810
Detector Reagent	101665	834202
SBG Reagent	101667	834203
Sample Diluent	101668	834204
Calibrator Diluent	101669	834205
Calibrator Concentrate	103217	834607
RGP Reagent	103159	N/A <sup>1</sup>

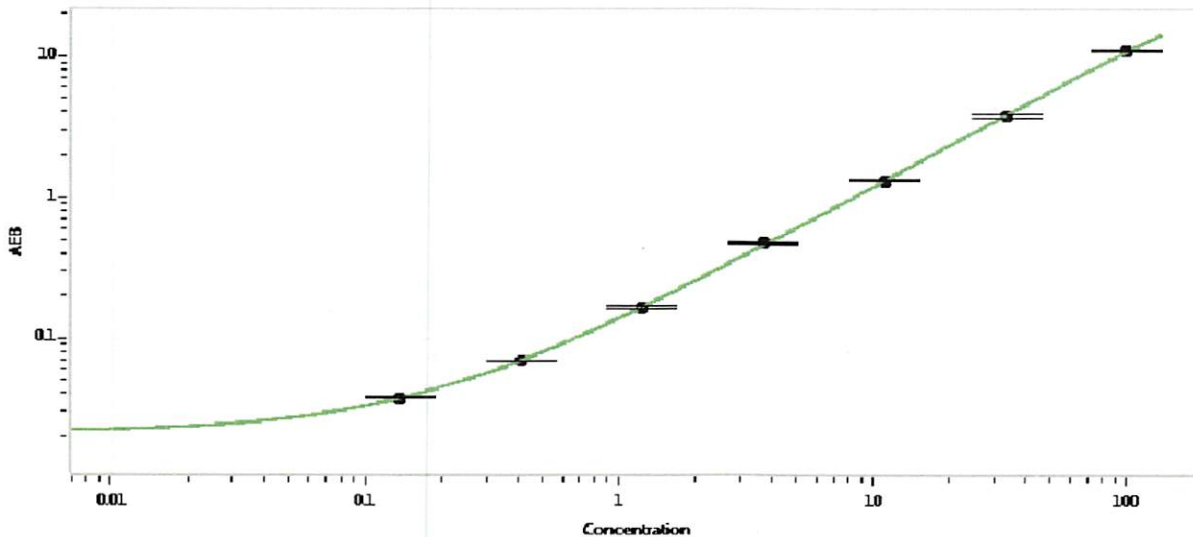
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	52520	1.01	0.740-1.33
Control 2	52520	33.4	27.1-40.6
Panel 1	MF-082015-01	1.20	0.942-1.53
Panel 2	MF-081415-02	10.3	7.46-11.2
Panel 3	MF-081415-03	11.1	9.20-15.0
Panel 4	MF-090315-09	41.5	27.6-66.8

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

Calibrator Concentrate Value Assignment (pg/mL)	1427
---	------

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

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]* 20/DEC/2018  
Signature/Date