

Product Number:	100196
Lot Number:	501174
Expiration:	1-Aug-2018
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	100391	80743
Detector Reagent	100315	80731
SBG Reagent	100316	80730
Sample Diluent	100317	80204
Calibrator Diluent	100665	80323
Calibrator Concentrate	100664	72701
RGP Reagent	103159	N/A <sup>1</sup>

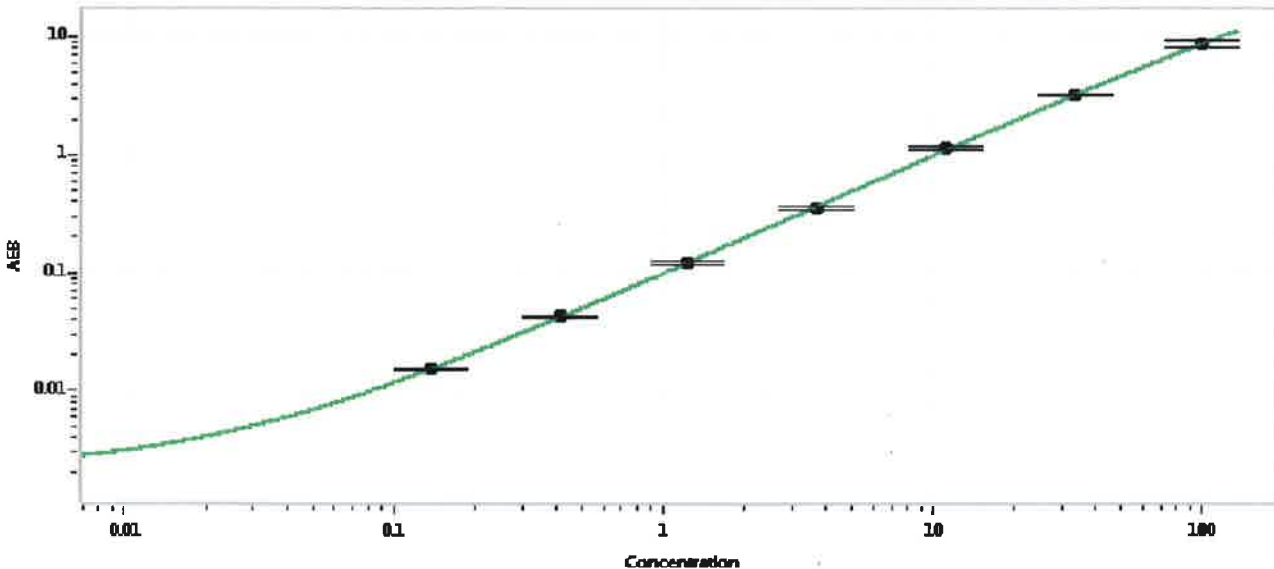
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	52521	105	75.6-118
Control 2	52521	10.2	7.72-11.6
Panel 1	MF-012116-05	1.56	0.871-1.73
Panel 2	MF-090815-18	9.00	4.45-9.13
Panel 3	MF-090815-19	90.5	53.9-94.9

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

Calibrator Concentrate Value Assignment (pg/mL)	2154
---	------

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      QC Supervisor      *[Signature]* 29 MAR 2018  
Name      Title      Signature/Date