

Product Number:	100199
Lot Number:	501325
Expiration:	3-Jan-2019
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	101045	819206
Detector Reagent	101047	819004
SBG Reagent	101057	819003
Sample Diluent	101055	818416
Calibrator Diluent	101051	816307
Calibrator Concentrate	101049	81302
RGP Reagent	103159	N/A ¹

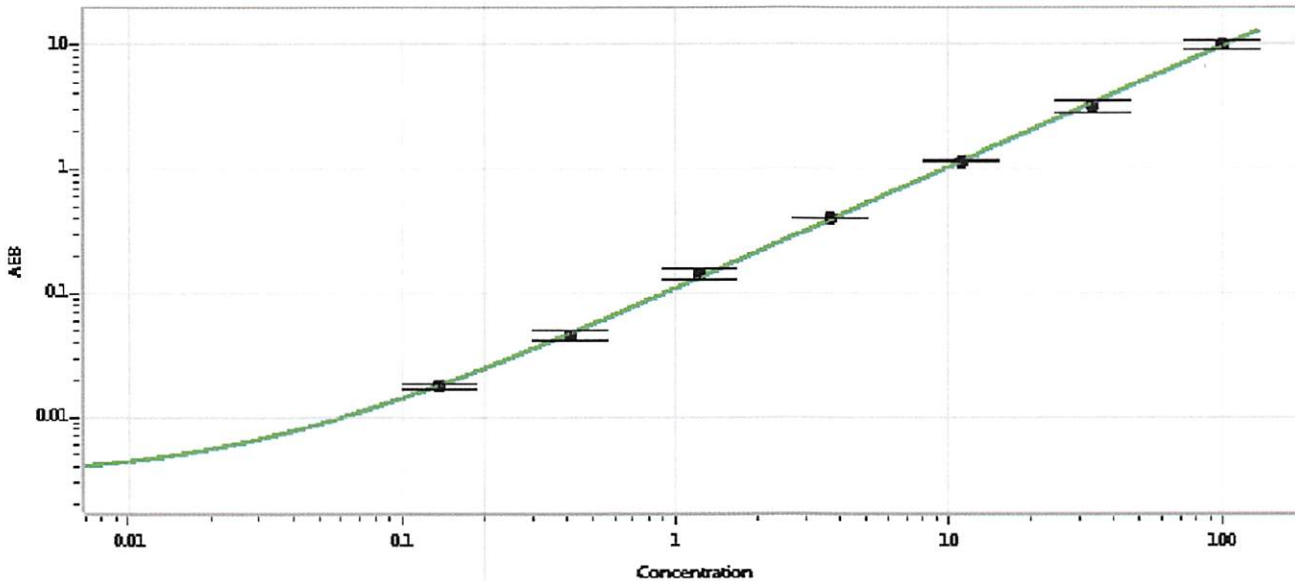
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-032017-04	1.61	1.30-1.95
Control 2	MF-032017-04	51.7	42.9-64.4
Panel 1	818701	0.513	0.337-1.33
Panel 2	MF-060118-02	2273	1539-2309
Panel 3	MF-060118-03	3337	2228-3341

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

Calibrator Concentrate Value Assignment (pg/mL)	1,832
---	-------

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] 22/1/2018
Signature/Date