

Product Number:	100199
Lot Number:	501359
Expiration:	9-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	821401
Detector Reagent	101047	821501
SBG Reagent	101057	821305
Sample Diluent	101055	819016
Calibrator Diluent	101051	819815
Calibrator Concentrate	101049	81302
RGP Reagent	103159	N/A ¹

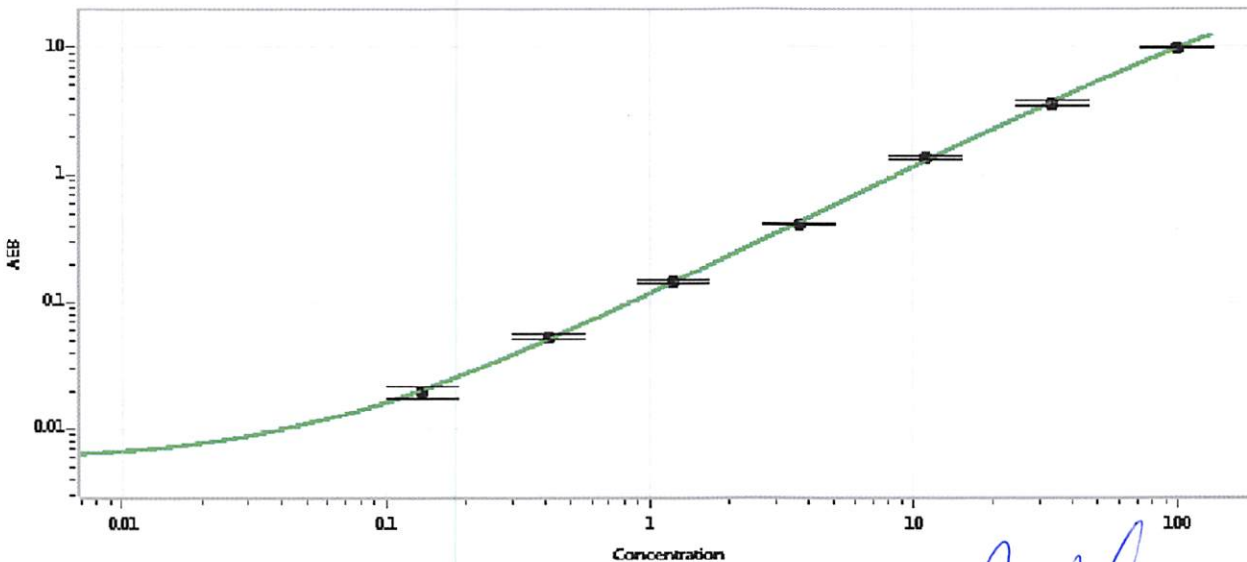
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-032017-04	1.60	1.30-1.95
Control 2	MF-032017-04	49.2	42.7-64.0
Panel 1	818701	0.867	0.354-1.32
Panel 2	MF-060118-02	1850	1534-2301
Panel 3	MF-060118-03	2824	2230-3345

* 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

James Drumme
Senior Director, Manufacturing Operations
On behalf of:

[Signature]
Aug 10, 2018

Review/Approval

Sivasankari Murugan QC Supervisor
Name Title

Signature/Date