

Product Number:	100199
Lot Number:	501250
Expiration:	9-Nov-2018
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	101045	81301
Detector Reagent	101047	81352
SBG Reagent	101057	81300
Sample Diluent	101055	81296
Calibrator Diluent	101051	81231
Calibrator Concentrate	101049	81302
RGP Reagent	103159	N/A ¹

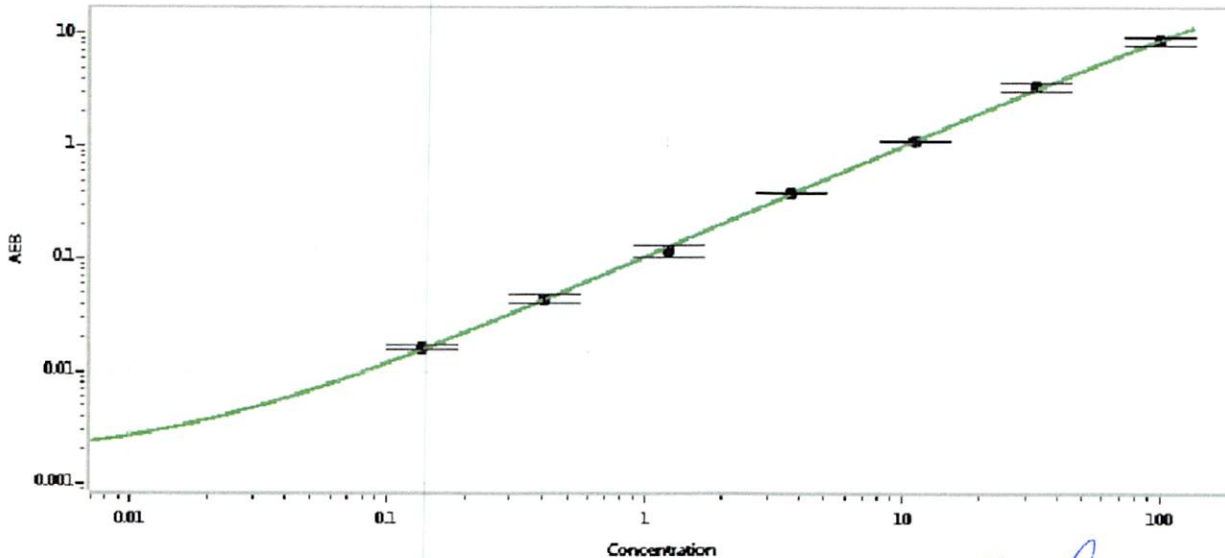
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-032017-04	1.57	1.30-1.95
Control 2	MF-032017-04	58.7	42.9-64.4
Panel 1	MF-032217-08	466	363-545
Panel 2	MF-032217-09	2257	2056-3084
Panel 3	MF-032217-10	504	343-515

* 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 Drummey
Senior Director, Manufacturing Operations
On behalf of:

Review/Approval: Sivasankari Murugan, QC Supervisor
Name: Sivasankari Murugan, Title: QC Supervisor, Signature/Date: [Signature], 22 May 2018