

Product Number:	100200
Lot Number:	501461
Expiration:	10-Jan-2019
Platform(s):	HD-1 & SR-X



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

Component	Item Number	Lot Number
Bead Reagent	101089	820601
Detector Reagent	101091	821308
SEB Reagent	101101	819104
Sample Diluent	101099	818005
Calibrator Diluent	101095	816308
Calibrator Concentrate	101093	828403
RGP Reagent	103159	N/A ¹

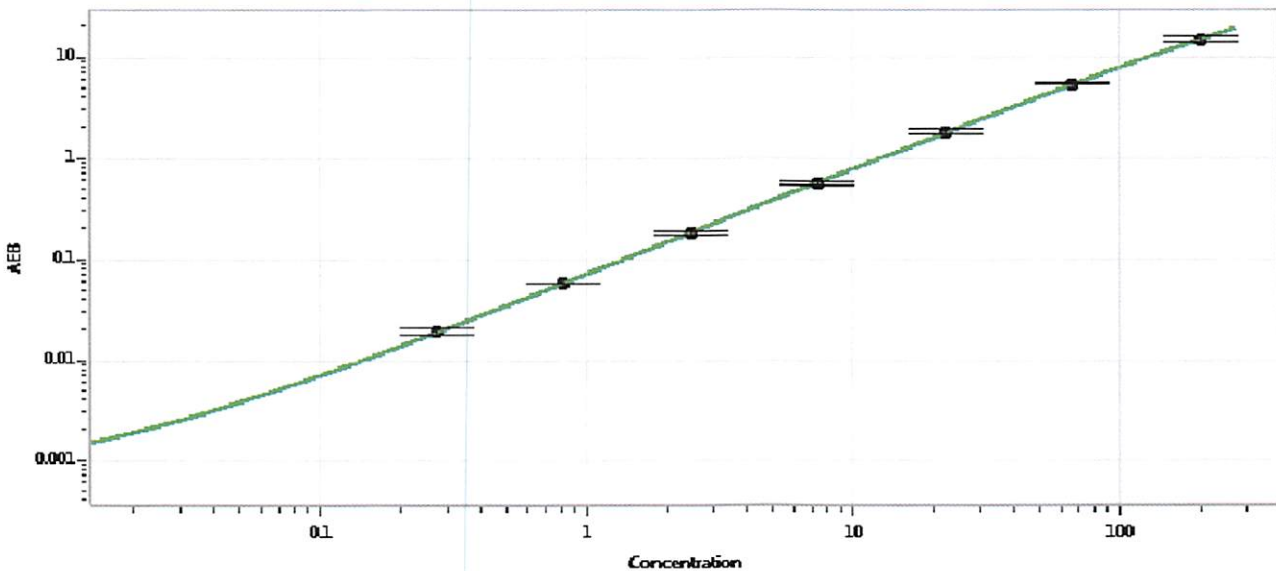
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-092215-03	9.72	6.99-11.4
Control 2	MF-092215-03	97.8	67.3-112
Panel 1	MF-012116-06	2.46	1.17-2.99
Panel 2	MF-093015-02	4.69	1.45-5.96
Panel 3	MF-093015-03	40.2	24.1-59.5

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

Calibrator Concentrate Value Assignment (pg/mL)	1,432
---	-------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Sivasankari Murugar QC Supervisor
Name Title

[Signature] 18 OCT 2018
Signature/Date