

Product Number:	101605
Lot Number:	501444
Expiration:	25-Sep-2019
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	101607	827610
Detector Reagent	101606	827105
SBG Reagent	101608	827106
Sample Diluent	103200	827002
Calibrator Diluent	101610	826809
Calibrator Concentrate	100895	828001
RGP Reagent	103159	N/A ¹

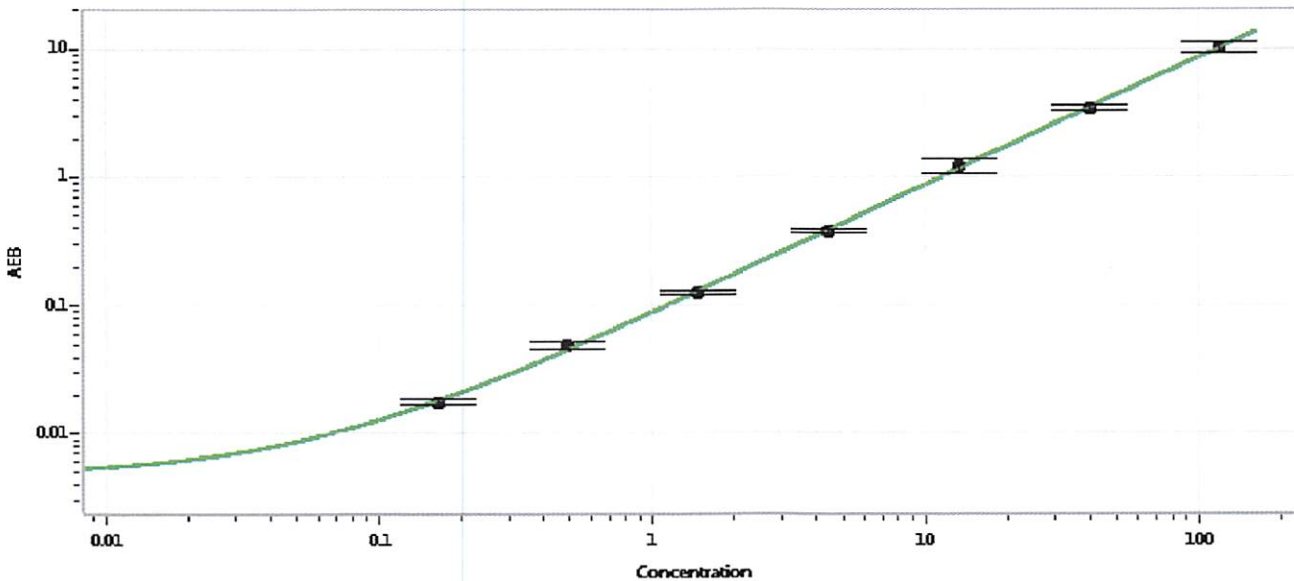
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-112515-02	6.16	3.13-7.30
Control 2	MF-112515-02	176	93.3-218
Panel 1	MF-112315-03	77.8	39.4-91.9
Panel 2	MF-112315-04	19.4	8.67-20.2
Panel 3	MF-112315-05	26.9	14.3-33.3

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

Calibrator Concentrate Value Assignment (pg/mL)	3,725
---	-------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.165
C	0.494
D	1.48
E	4.44
F	13.3
G	40.0
H	120

Review/Approval

Sivasankari Murugan

QC Supervisor

[Signature] 10/09/2018

Name

Title

Signature/Date