

Product Number:	101605
Lot Number:	501202
Expiration:	31-Jan-2019
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	101607	80333
Detector Reagent	101606	80322
SBG Reagent	101608	80311
Sample Diluent	101609	80321
Calibrator Diluent	101610	80944
Calibrator Concentrate	100895	80334
RGP Reagent	103159	N/A ¹

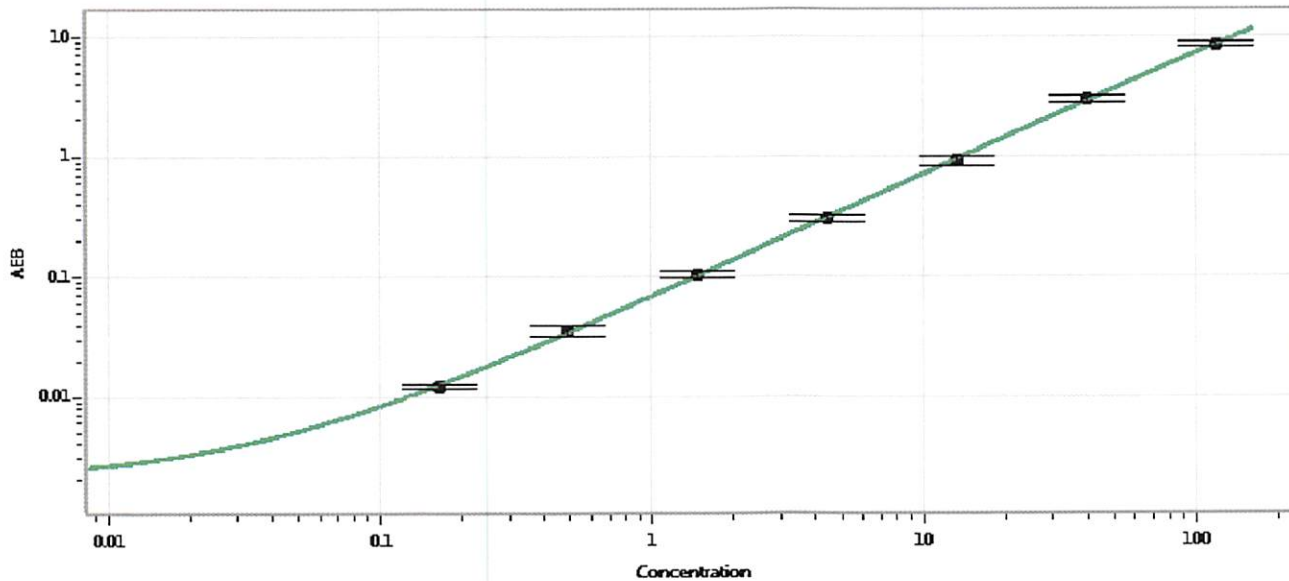
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-112515-02	6.38	3.13-7.30
Control 2	MF-112515-02	192	93.3-218
Panel 1	MF-112315-03	82.8	39.4-91.9
Panel 2	MF-112315-04	19.8	8.67-20.2
Panel 3	MF-112315-05	32.3	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)	5183
-------------------------------------------------	------

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
Name

QC Supervisor
Title

[Signature]
Signature/Date