

Product Number:	101605
Lot Number:	501239
Expiration:	12-Apr-2019
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	101607	81216
Detector Reagent	101606	81173
SBG Reagent	101608	81172
Sample Diluent	103200	81200
Calibrator Diluent	101610	80945
Calibrator Concentrate	100895	80334
RGP Reagent	103159	N/A ¹

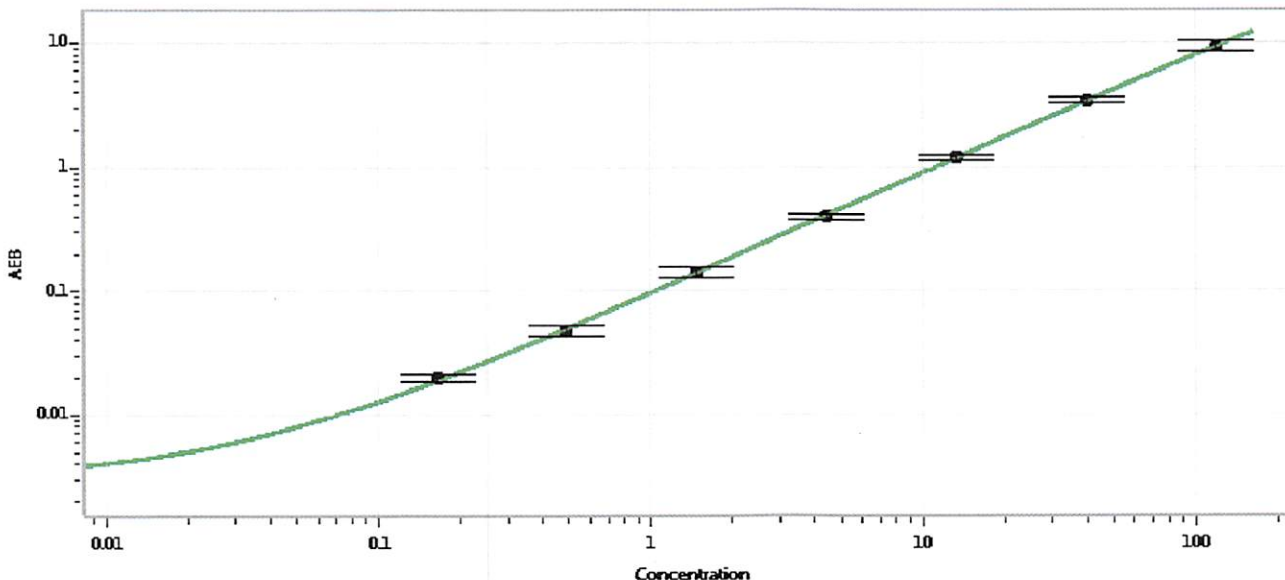
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-112515-02	6.01	3.13-7.30
Control 2	MF-112515-02	203	93.3-218
Panel 1	MF-112315-03	83.2	39.4-91.9
Panel 2	MF-112315-04	19.0	8.67-20.2
Panel 3	MF-112315-05	31.8	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
11 MAY 2018