

Product Number:	100198
Lot Number:	501100
Expiration:	26-Jul-2018
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	100815	80260
Detector Reagent	100816	80267
SBG Reagent	100817	80262
Sample Diluent	100818	80270
Calibrator Diluent	100829	80264
Calibrator Concentrate	100828	73380
RGP Reagent	103159	N/A <sup>1</sup>

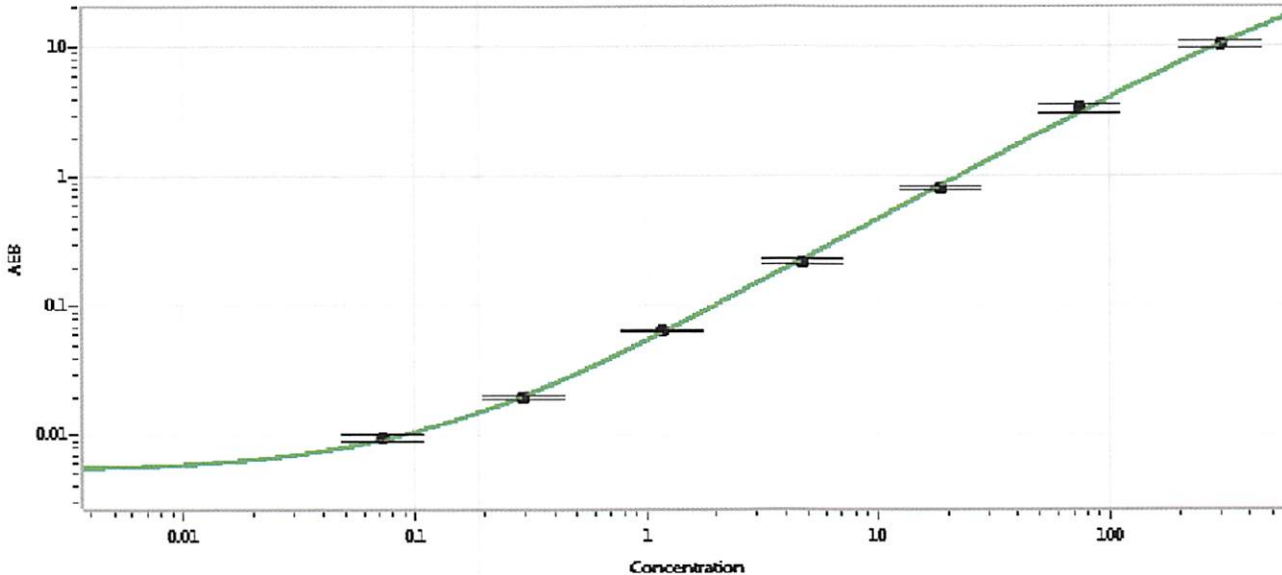
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-091217-02	7.96	6.51-9.77
Control 2	MF-091217-02	165	113-178
Panel 1	MF-061617-01	3.58	3.12-4.90
Panel 2	MF-061617-02	42.0	34.9-52.3
Panel 3	MF-061617-03	335	183-361

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

Calibrator Concentrate Value Assignment (pg/mL)	15328
---	-------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.073
C	0.293
D	1.17
E	4.69
F	18.8
G	75.0
H	300

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]* 20 FEB 2018  
Signature/Date