

Product Number:	102258
Lot Number:	501123
Expiration:	3-Nov-2018
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	102246	80390
Detector Reagent	102248	80340
SBG Reagent	102250	80341
Sample Diluent	102252	80342
Calibrator Diluent	102620	80343
Calibrator Concentrate	102619	73520
RGP Reagent	103159	N/A ¹

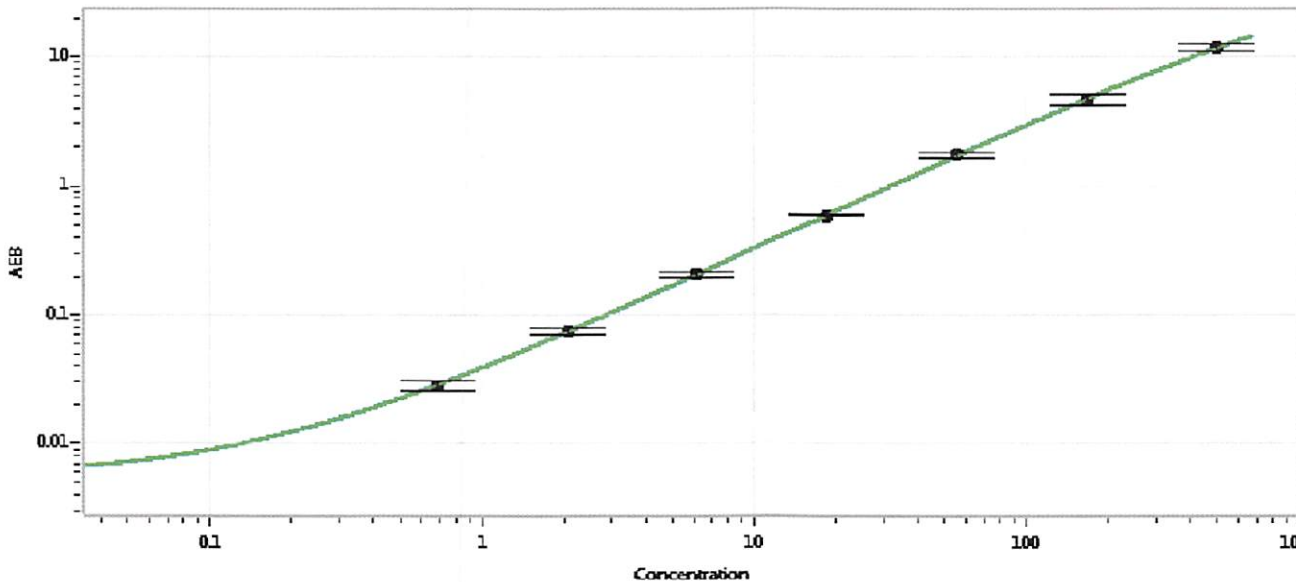
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	72720	4.18	3.70-5.82
Control 2	72720	179	117-224
Panel 1	MF-092717-01	4.69	4.27-7.42
Panel 2	MF-092717-02	53.0	46.8-78.1
Panel 3	MF-092717-03	197	124-291

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

Calibrator Concentrate Value Assignment (pg/mL)	7,585
---	-------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.686
C	2.06
D	6.17
E	18.5
F	55.6
G	167
H	500

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 26 FEB 2018
Signature/Date