

Product Number:	100200
Lot Number:	501148
Expiration:	9-May-2018
Platform(s):	HD-1 & SR-X



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

Component	Item Number	Lot Number
Bead Reagent	101089	73323
Detector Reagent	101091	80533
SBG Reagent	101101	73331
Sample Diluent	101099	73136
Calibrator Diluent	101095	73137
Calibrator Concentrate	101093	63560
RGP Reagent	103159	N/A ¹

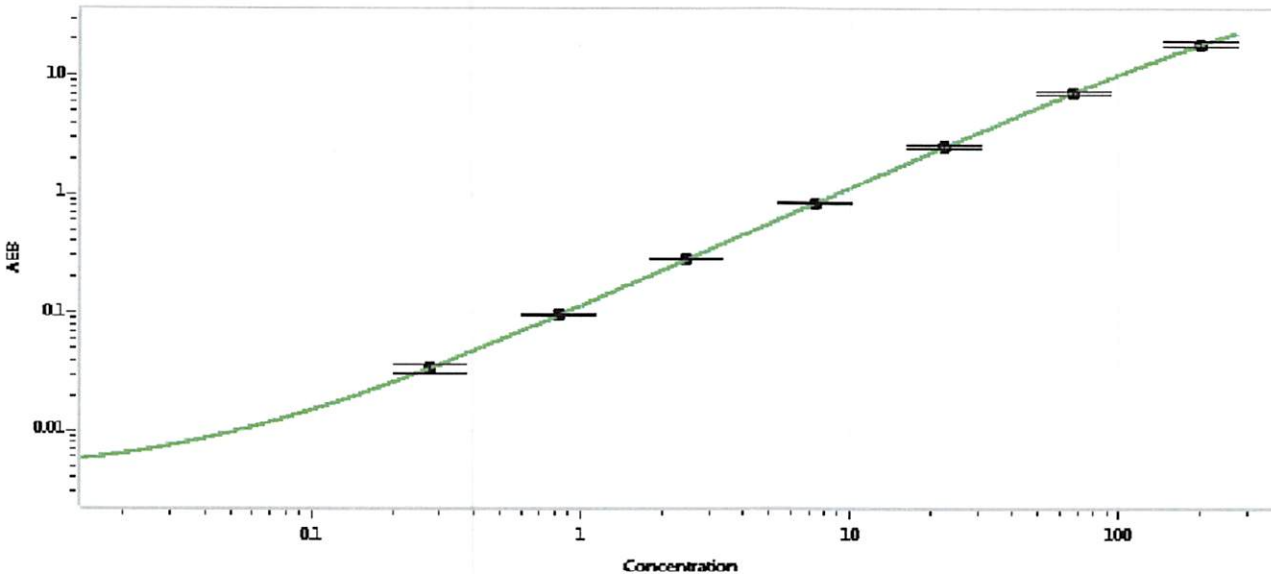
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-092215-03	10.2	6.99-11.4
Control 2	MF-092215-03	103	67.3-112
Panel 1	MF-012116-06	2.12	1.58-2.80
Panel 2	MF-093015-02	3.17	1.45-5.96
Panel 3	MF-093015-03	52.3	24.1-59.5

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

Calibrator Concentrate Value Assignment (pg/mL)	1,751
---	-------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date 22 MAR 2018