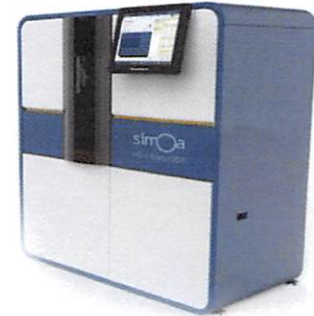


Product Number:	100200
Lot Number:	501261
Expiration:	5-Oct-2018
Platform(s):	HD-1 & SR-X



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

Component	Item Number	Lot Number
Bead Reagent	101089	80960
Detector Reagent	101091	81422
SBG Reagent	101101	80954
Sample Diluent	101099	80865
Calibrator Diluent	101095	80794
Calibrator Concentrate	101093	63560
RGP Reagent	103159	N/A ¹

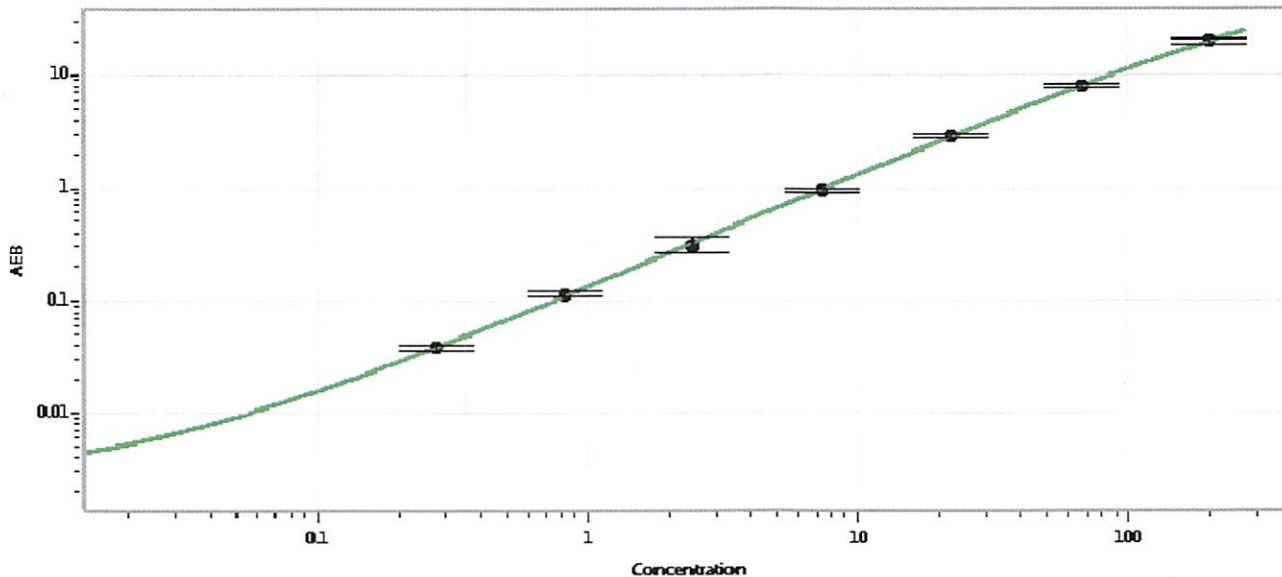
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-092215-03	7.73	6.99-11.4
Control 2	MF-092215-03	68.6	67.3-112
Panel 1	MF-012116-06	1.48	1.17-2.99
Panel 2	MF-093015-02	2.66	1.45-5.96
Panel 3	MF-093015-03	33.4	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

S. Murugan
Signature/Date 25 May 2018