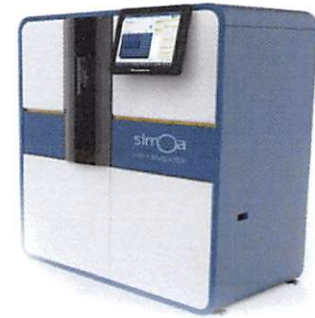


Product Number:	101154
Lot Number:	501246
Expiration:	27-Oct-2018
Platform(s):	HD-1



Component	Part Number	Lot Number
Bead Reagent	101141	81220
Detector Reagent	101143	81228
SBG Reagent	101153	81201
Sample Diluent	101151	81174
Calibrator Diluent	101147	81001
Calibrator Concentrate	101145	73524
RGP Reagent	103159	N/A ¹

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

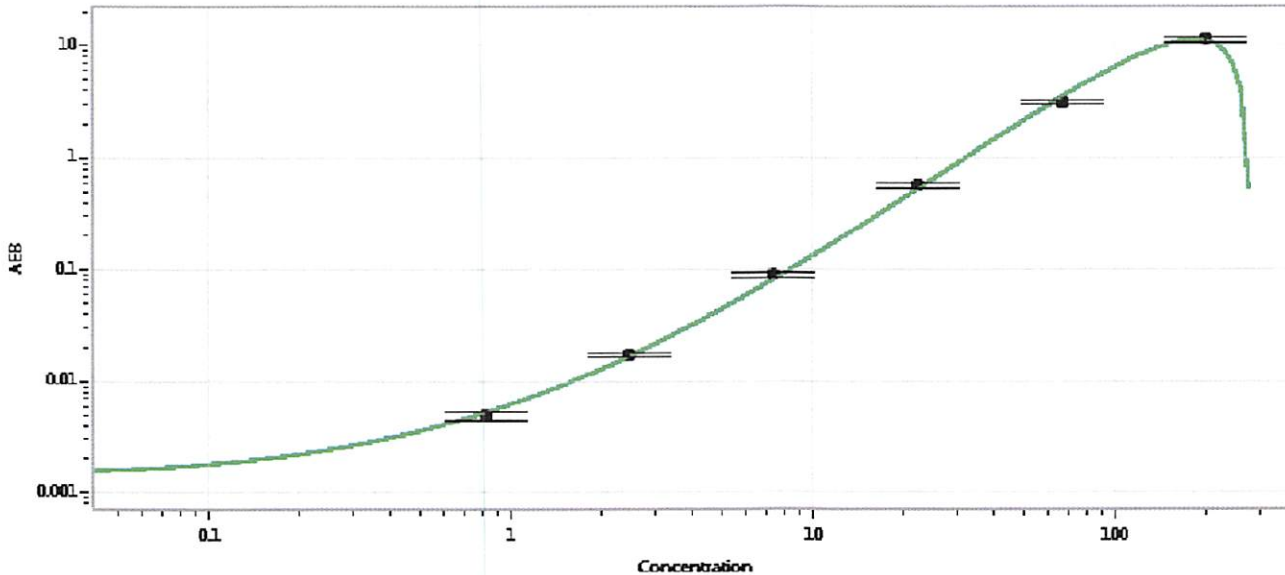
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	41952	69.4	54.1-81.2
Control 2	41952	374	277-415
Panel 1	1816	103	77.3-116
Panel 2	1712	93.1	68.1-103
Panel 3	2004	163	120-200

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

Calibrator Concentrate Value Assignment (pg/mL)	14,199
---	--------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.823
C	2.47
D	7.41
E	22.2
F	66.7
G	200

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 27 May 2018
Signature/Date