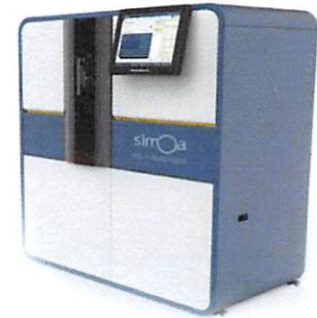


Product Number:	101132
Lot Number:	501301
Expiration:	9-Nov-2018
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	101119	81762
Detector Reagent	101121	81760
SBG Reagent	101131	81761
Sample Diluent	101129	81297
Calibrator Diluent	101125	81633
Calibrator Concentrate	101126	81771
RGP Reagent	103159	N/A <sup>1</sup>

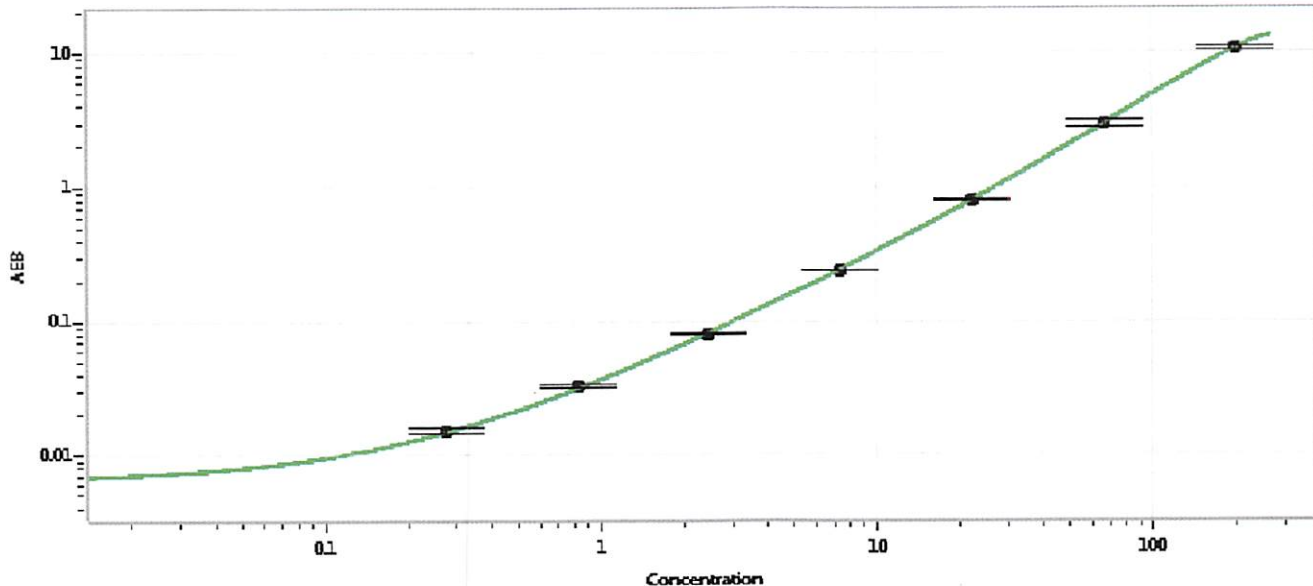
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-051716-04	2.18	1.35-3.54
Control 2	MF-051716-04	251	131-447
Panel 1	MF-051916-03	249	215-354
Panel 2	MF-051916-04	213	207-352

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

Calibrator Concentrate Value Assignment (pg/mL)	55,360
---	--------

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