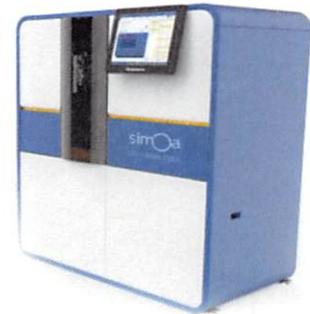


Product Number:	102732
Lot Number:	501130
Expiration:	9-May-2018
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	102722	80320
Detector Reagent	102724	73242
SBG Reagent	102726	73243
Sample Diluent	102728	73138
Calibrator Diluent	102731	73139
Calibrator Concentrate	100658	73217
RGP Reagent	103159	N/A ¹

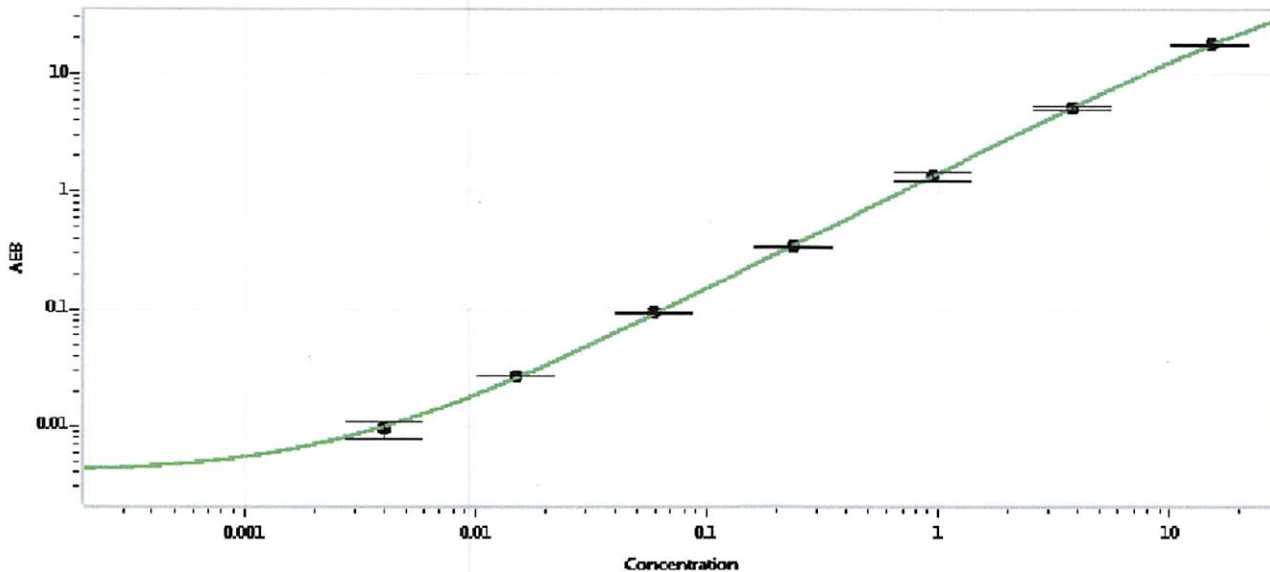
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-061217-01	0.160	0.132-0.293
Control 2	MF-061217-01	3.41	2.50-6.45
Panel 1	MF-061317-03	0.043	0.037-0.085
Panel 2	MF-061317-04	0.230	0.183-0.499
Panel 3	MF-061317-05	4.25	3.18-8.05

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

Calibrator Concentrate Value Assignment (pg/mL)	194
---	-----

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.004
C	0.015
D	0.059
E	0.234
F	0.938
G	3.75
H	15.0

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 16 FEB 2018
Signature/Date