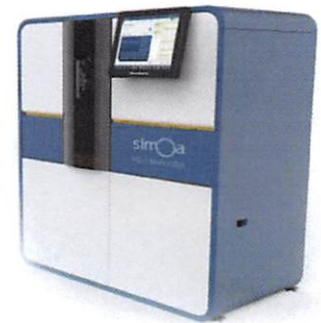


Product Number:	101599
Lot Number:	501404
Expiration:	31-May-2019
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101601	825402
Detector Reagent	102012	825501
SBG Reagent	101602	824307
Sample Diluent	101603	824303
Calibrator Diluent	102014	822014
Calibrator Concentrate	100778	81562
RGP Reagent	103159	N/A ¹

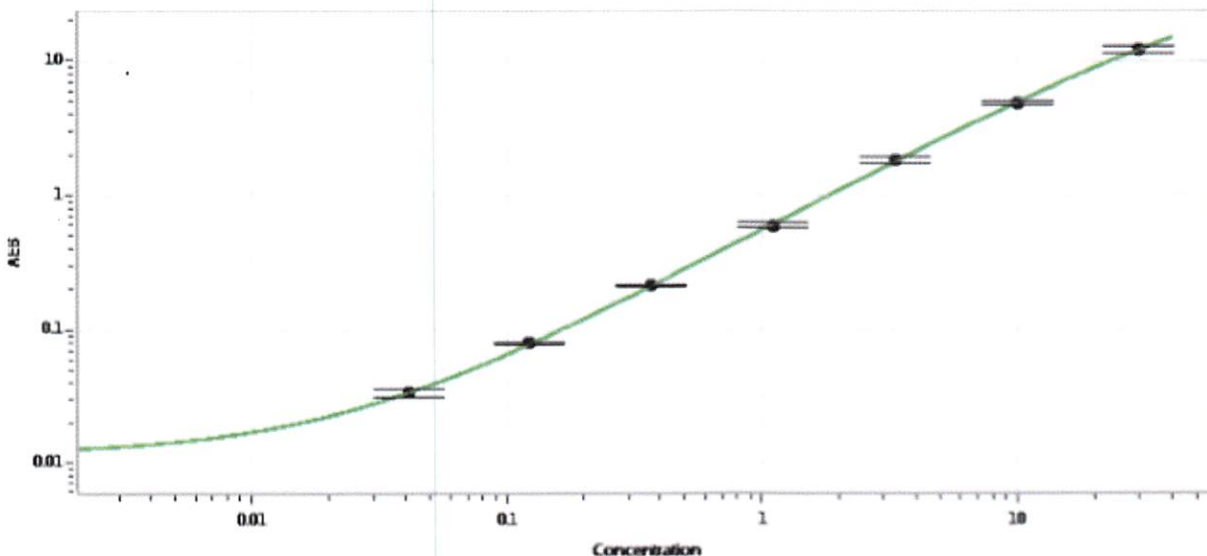
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-012518-02	0.431	0.370-0.555
Control 2	MF-012518-02	22.6	14.8-24.0
Panel 1	MF-071615-01	0.340	0.325-0.513
Panel 2	MF-071615-02	7.16	5.58-8.37
Panel 3	MF-071615-03	2.13	1.68-2.52

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

Calibrator Concentrate Value Assignment (pg/mL)	800
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Sivasankari Murugan QC Supervisor

Sivasankari Murugan 14 SEP 2018

Name

Title

Signature/Date