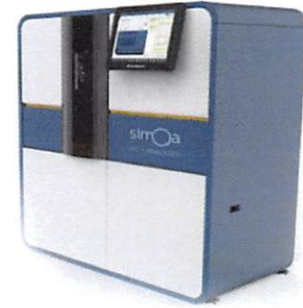


Product Number:	101672
Lot Number:	501393
Expiration:	9-Jan-2019
Platform(s):	HD-1 & SR-X



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

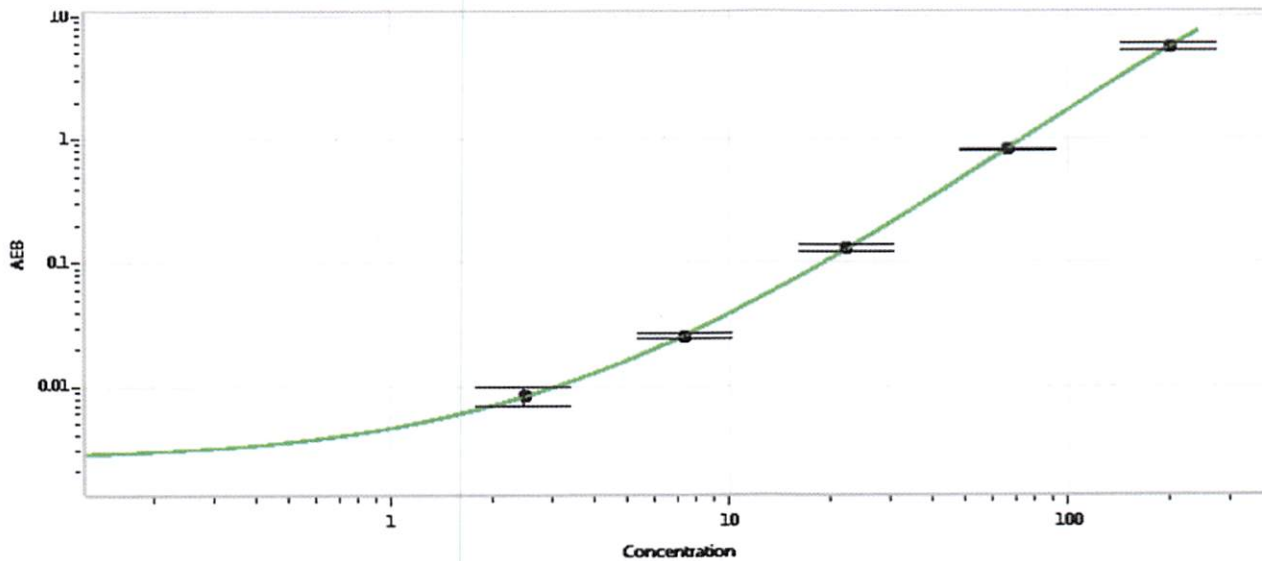
Component	Part Number	Lot Number
Bead Reagent	101674	819706
Detector Reagent	101673	824102
SBG Reagent	101675	819007
Sample Diluent	101676	818004
Calibrator Diluent	101677	816306
Calibrator Concentrate	101734	73205
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	MF-021816-02	19.2	17.1-28.3	
Control 2	MF-021816-02	316	263-434	
Panel 1	MF-040518-01	168	141-211	
Panel 2	MF-040518-02	43.0	35.0-58.0	

Calibrator Concentrate Value Assignment (pg/mL)	2711
---	------

Calibration Curve



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

Review/Approval

Sivasankari Murugar QC Supervisor
Name Title

[Signature] 06 SEP 2018
Signature/Date