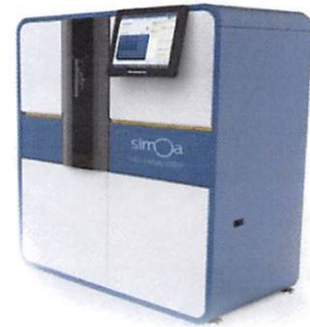


Product Number:	103186
Lot Number:	501386
Expiration:	20-Jul-2019
Platform(s):	HD-1 & SR-X



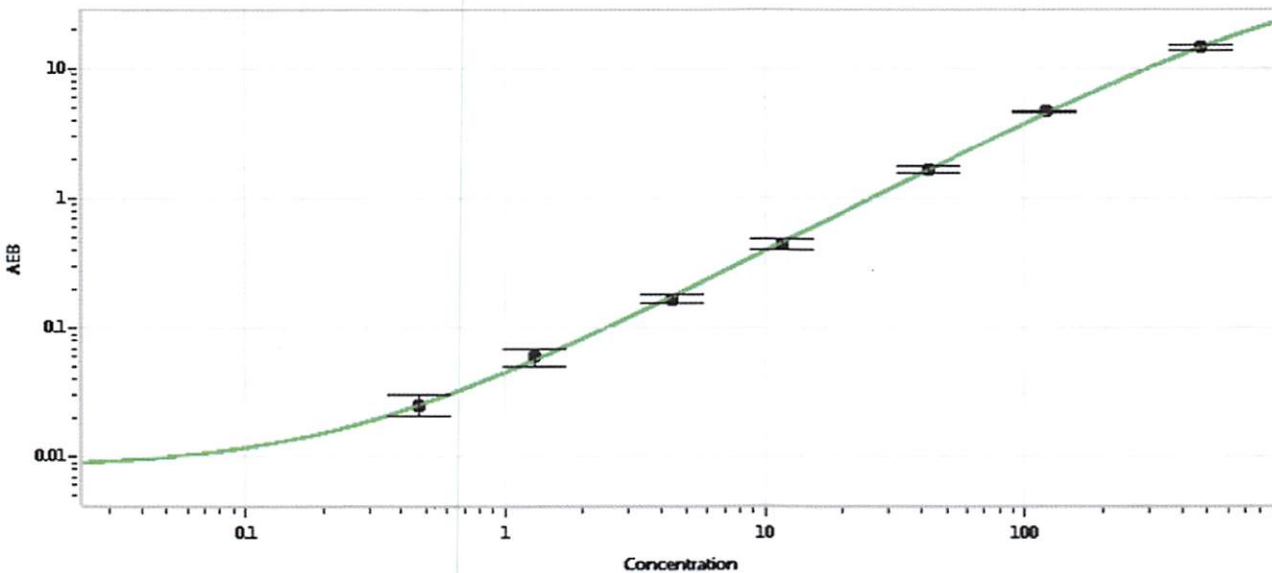
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102246	820703
Detector Reagent	102248	820406
SBG Reagent	102250	820103
Sample Diluent	102252	820104
Reference Calibrators	102255	822501
Controls	102256 - 102257	822806 823603
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	822806	4.49	3.57-5.62	
Control 2	823603	154	117-175	
Panel 1	823301	5.63	3.99-6.77	
Panel 2	823601	78.3	60.8-91.2	
Panel 3	823602	297	223-396	

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.467
C	1.31
D	4.38
E	11.6
F	42.5
G	120
H	472

Review/Approval

Sivasankari Murugar QC Supervisor
Name Title

[Signature] 29/06/2018
Signature/Date