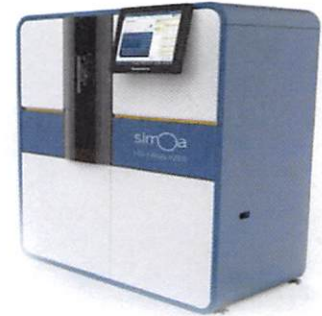


Product Number:	102215
Lot Number:	501364
Expiration:	1-Feb-2019
Platform(s):	HD-1 & SR-X



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

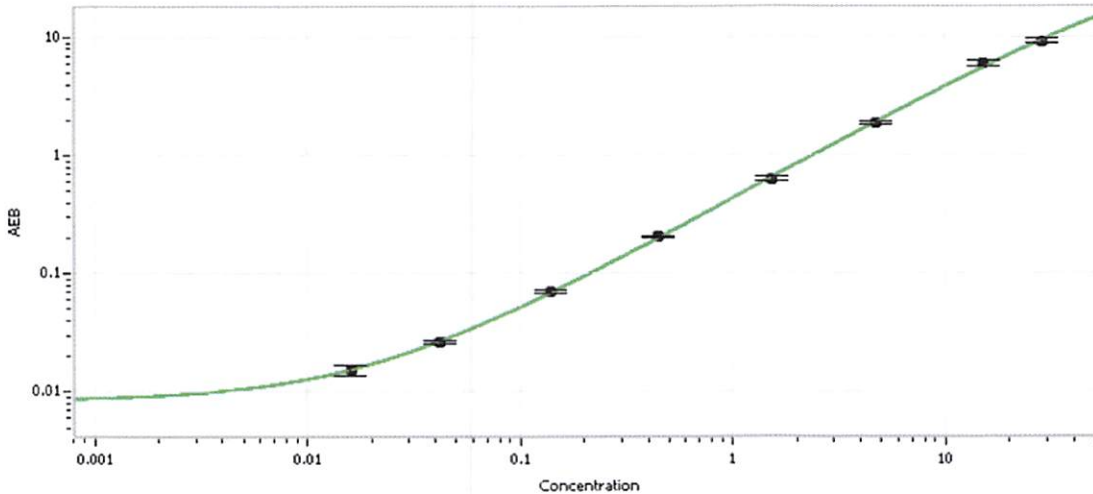
Component	Part Number	Lot Number
Bead Reagent	102211	822204
Detector Reagent	102212	821303
SBG Reagent	102213	821304
Sample Diluent	102214	821301
Calibrators	101310	81511/81783
Controls	100702-100703	81512
RGP Reagent	103159	N/A <sup>1</sup>

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

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	81512	0.152	0.117-0.175
Control 2	81512	16.3	12.6-19.3
Panel 1	MF-092217-01	0.282	0.177-0.353
Panel 2	MF-092217-02	2.85	1.99-3.51
Panel 3	MF-101117-01	17.7	12.8-20.8

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.016
C	0.042
D	0.140
E	0.446
F	1.54
G	4.77
H	15.3
I	28.6

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date 22 AUG 2018