

Product Number:	103186
Lot Number:	501496
Expiration:	5-Dec-2019
Platform(s):	HD-1 & SR-X



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

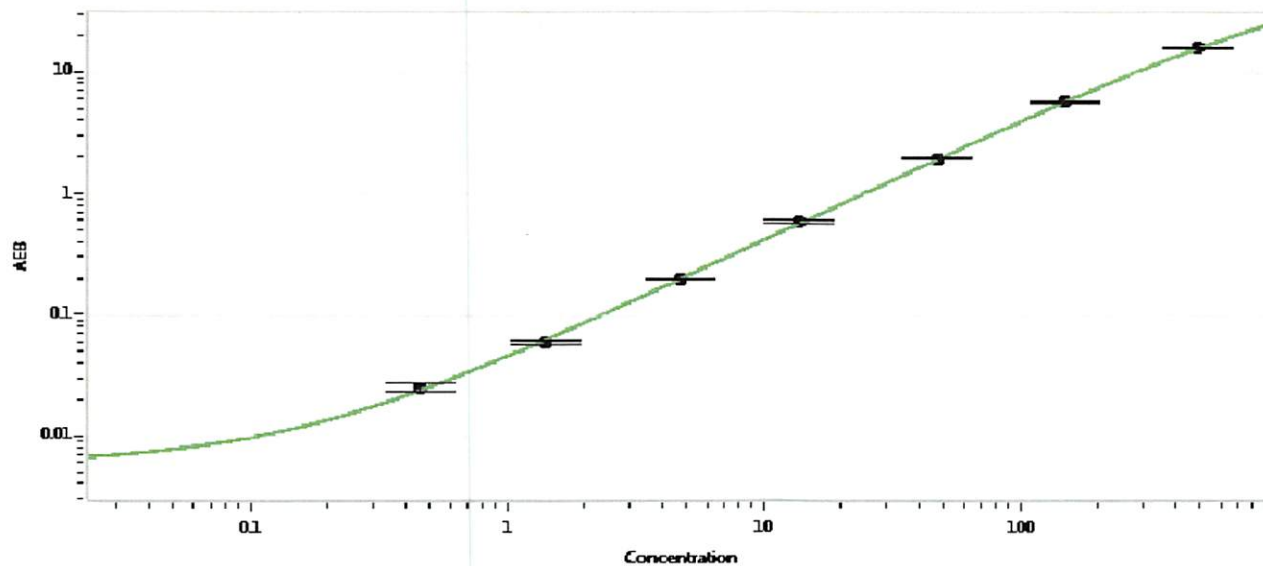
Component	Part Number	Lot Number
Bead Reagent	102246	834708
Detector Reagent	102248	834705
SBG Reagent	102250	833911
Sample Diluent	102252	834507
Reference Calibrators	102255	833201
Controls	102256 - 102257	830201
RGP Reagent	103159	N/A <sup>1</sup>

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

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	830201	4.21	3.16-5.11
Control 2	830201	160	125-187
Panel 1-A	823301	5.36	4.08-6.68
Panel 1-B	823302	6.39	4.90-7.61
Panel 2	823601	89.8	59.3-101
Panel 3	823602	336	236-402

\* 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	0.458
C	1.42
D	4.74
E	13.7
F	47.4
G	148
H	488

Review/Approval: Sivasankari Murugan QC Supervisor  
 Name: Sivasankari Murugan Title: QC Supervisor Signature/Date: [Signature] 11DEC2018