

Product Number:	103186
Lot Number:	501555
Expiration:	5-Dec-2019
Platform(s):	HD-1* & SR-X



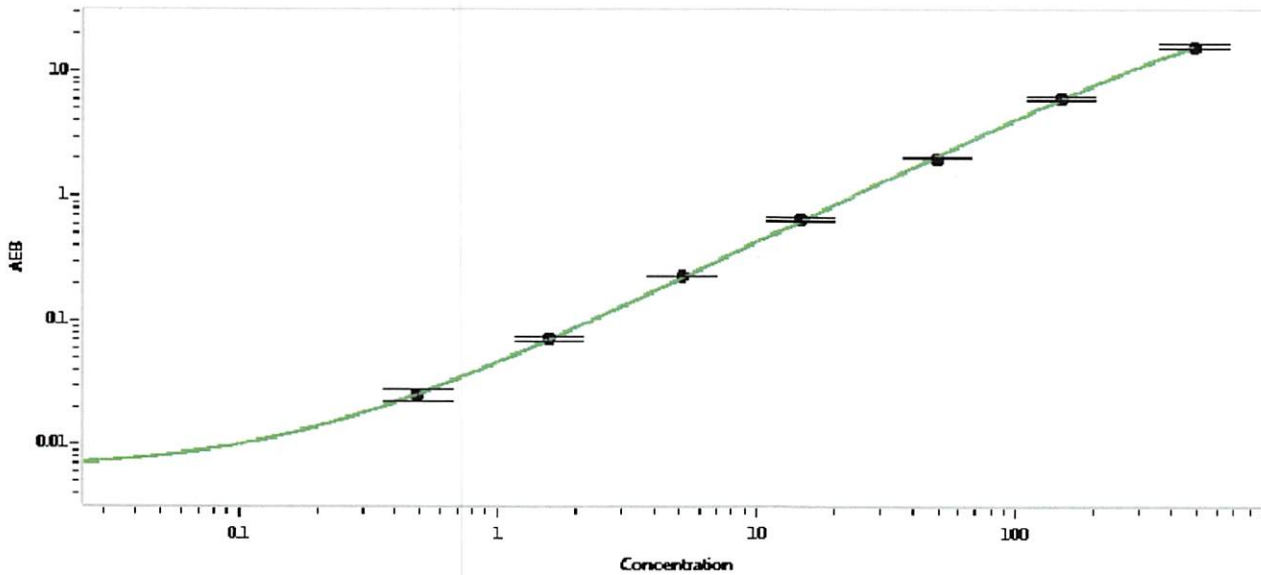
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102246	834708
Detector Reagent	102248	834705
SBG Reagent	102250	833911
Sample Diluent	102252	834507
Reference Calibrators	102255	831901
Controls	102256 - 102257	830201
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> 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	830201	4.39	3.16-5.11	
Control 2	830201	161	125-187	
Panel 1-A	823301	6.01	4.08-6.68	
Panel 1-B	823302	6.66	4.90-7.61	
Panel 2	823601	90.0	59.3-101	
Panel 3	823602	321	236-402	

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.494
C	1.58
D	5.10
E	14.8
F	49.3
G	150
H	484

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]* 24 DEC 2018  
Signature/Date