

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.

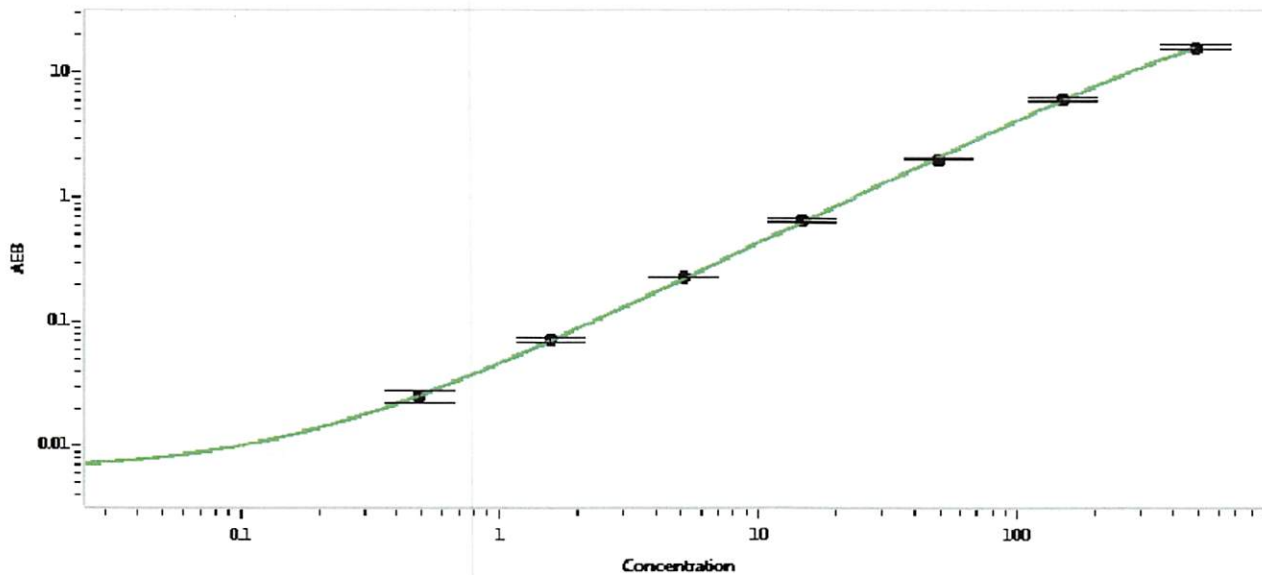
Component	Part Number	Lot Number
Bead Reagent	102246	834708
Detector Reagent	102248	834705
SBG Reagent	102250	833911
Sample Diluent	102252	834507
Reference Calibrators	102255	831901
Controls	102256 - 102257	830201
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	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

* 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	0.494
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