

Product Number:	103186
Lot Number:	501495
Expiration:	29-Aug-2019
Platform(s):	HD-1 & SR-X



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

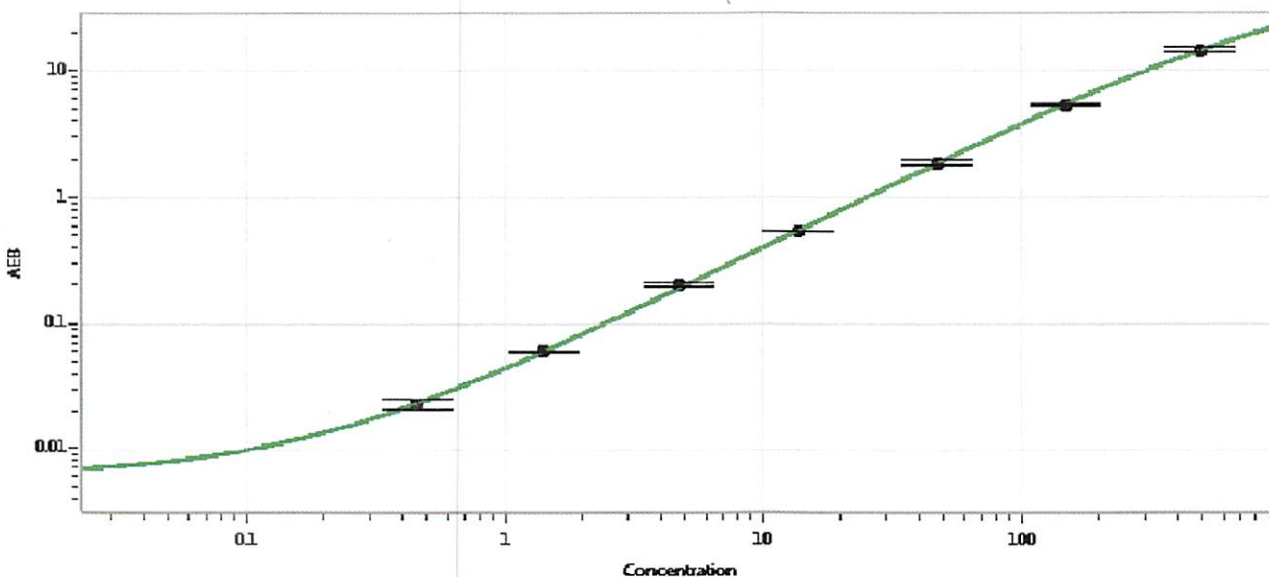
Component	Part Number	Lot Number
Bead Reagent	102246	831710
Detector Reagent	102248	831703
SBG Reagent	102250	824101
Sample Diluent	102252	826217
Reference Calibrators	102255	833201
Controls	102256 - 102257	830201
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	830201	4.23	3.16-5.11
Control 2	830201	167	125-187
Panel 1-A	823301	5.95	4.08-6.68
Panel 1-B	823302	7.40	4.90-7.61
Panel 2	823601	90.9	59.3-101
Panel 3	823602	354	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
Name

QC Supervisor
Title

S. Murugan 18 DEC 2018
Signature/Date