

Product Number:	101599
Lot Number:	501498
Expiration:	9-Aug-2019
Platform(s):	HD-1 & SR-X



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

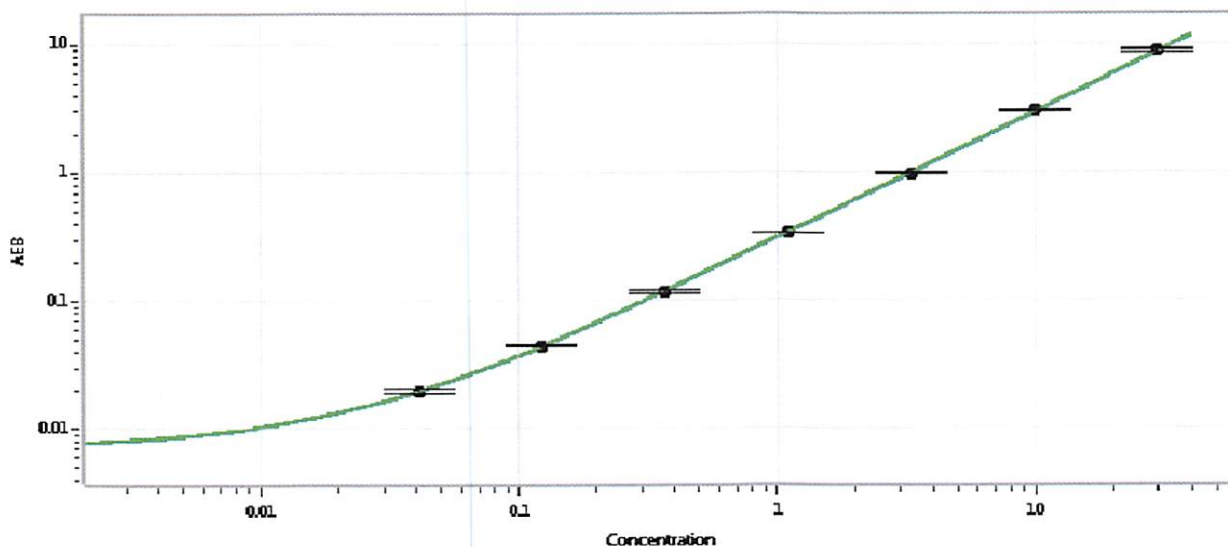
Component	Part Number	Lot Number
Bead Reagent	101601	831607
Detector Reagent	102012	831702
SBG Reagent	101602	831307
Sample Diluent	101603	831306
Calibrator Diluent	102014	829711
Calibrator Concentrate	100778	81562
RGP Reagent	103159	N/A ¹

¹ 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	MF-012518-02	0.393	0.363-0.545	
Control 2	MF-012518-02	22.1	14.7-24.7	
Panel 1	MF-120115-02	0.826	0.619-1.29	
Panel 2	MF-071615-02	7.70	5.58-8.37	
Panel 3	MF-071615-03	2.25	1.68-2.52	

Calibrator Concentrate Value Assignment (pg/mL)	800
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Sivasankari Murugan QC Supervisor

Name

Title

Sivasankari Murugan
Signature/Date