

Product Number:	101478
Lot Number:	501488
Expiration:	16-Feb-2019
Platform(s):	HD-1 & SR-X



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

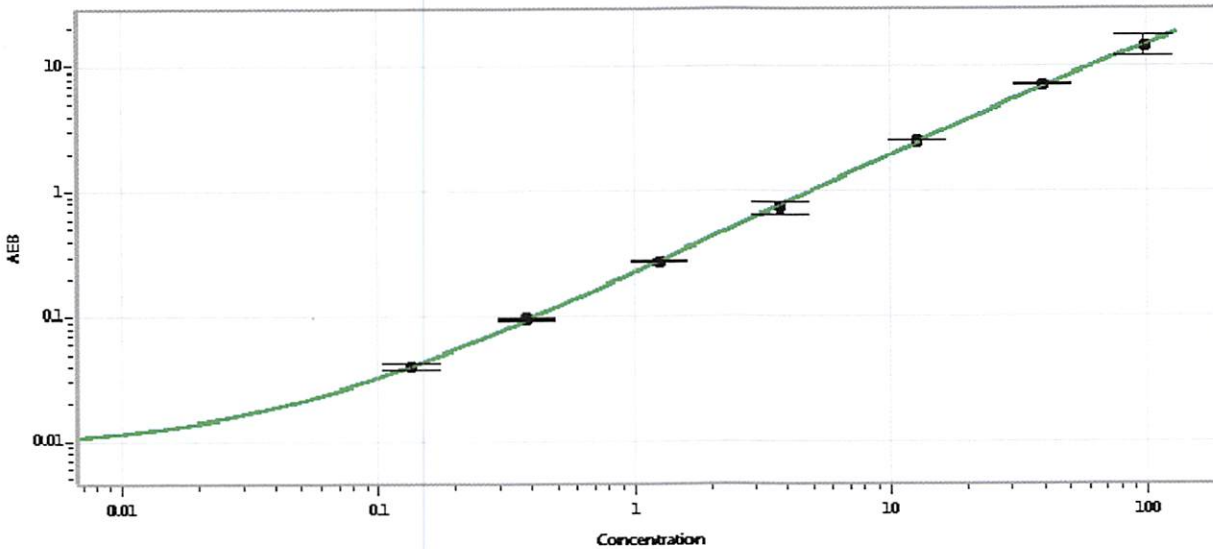
Component	Part Number	Lot Number
Bead Reagent	101480	826705
Detector Reagent	101481	826212
SBG Reagent	101482	822810
Sample Diluent	101483	827001
Calibrators	100758-100765	73180
Controls	100768 & 100769	73181
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	73181	4.18	3.43-5.14
Control 2	73181	95.2	74.2-111
Panel 1	MF-051017-04	0.680	0.508-0.855
Panel 2	MF-042117-06	4.92	3.89-6.07
Panel 3	MF-042117-07	58.9	45.5-73.7

* 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	0.133
C	0.381
D	1.25
E	3.70
F	12.8
G	39.5
H	98.1

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date
15 NOV 2018