

Product Number:	101212
Lot Number:	501557
Expiration:	25-Sep-2019
Platform(s):	HD-1



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

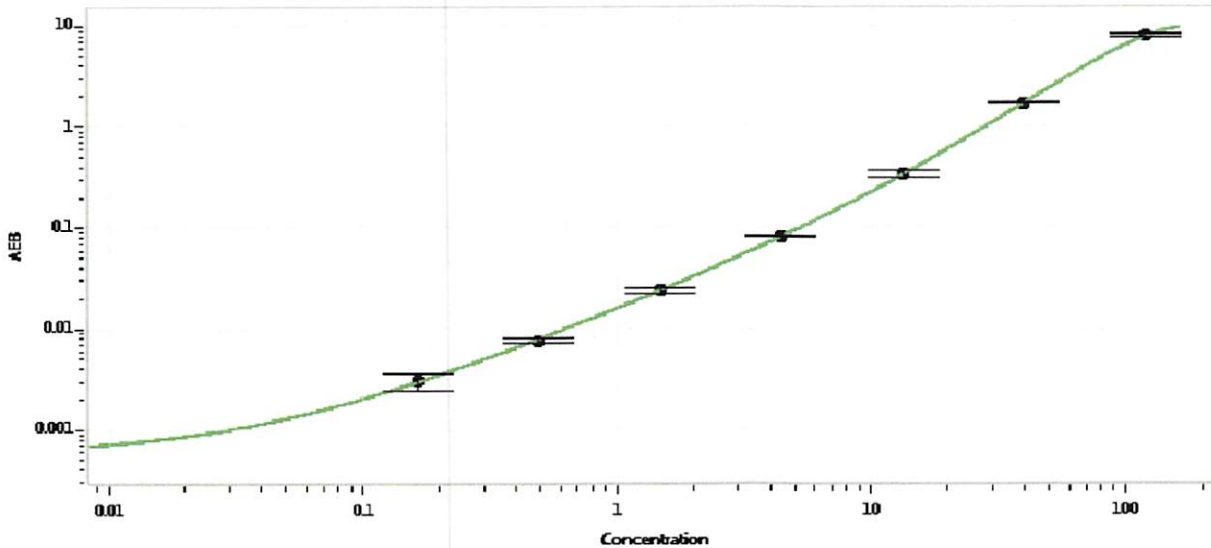
Component	Part Number	Lot Number
Bead Reagent	101198	834807
Detector Reagent	101200	834402
SBG Reagent	101211	834504
Sample Diluent	101209	834602
Calibrator Diluent	101204	826813
Calibrator Concentrate	101202	72780
RGP Reagent	103159	N/A ¹

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	828802	65.1	55.1-82.6	
Panel 1	MF-052318-05	192	141-211	
Panel 2	MF-052318-06	43.5	35.9-53.8	
Panel 3	MF-052318-07	97.8	72.7-109	

Calibrator Concentrate Value Assignment (pg/mL)	2,032
---	-------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.165
C	0.494
D	1.48
E	4.44
F	13.3
G	40.0
H	120

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date
27 DEC 2018