

Product Number:	100860
Lot Number:	501326
Expiration:	9-Jan-2019
Platform(s):	HD-1



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

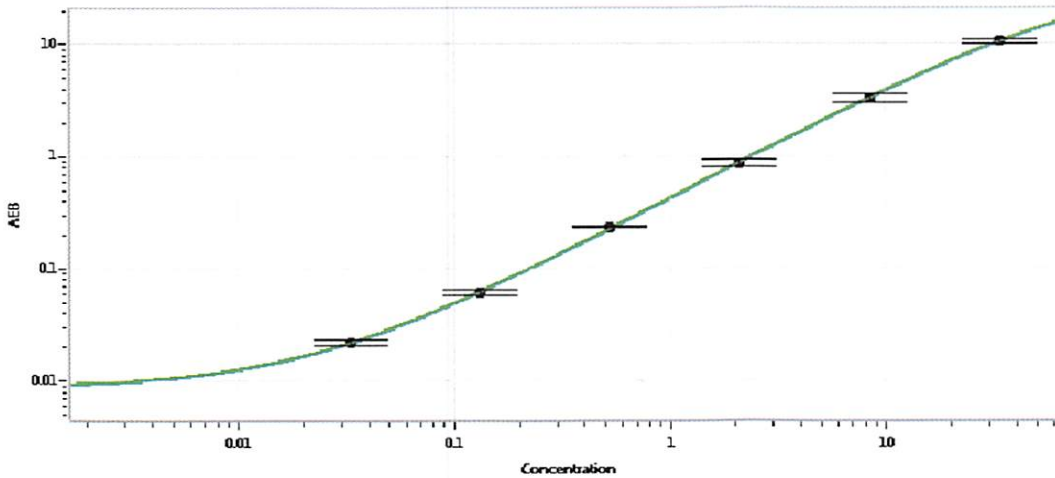
Component	Part Number	Lot Number
Bead Reagent	100865	819203
Detector Reagent	100862	819105
SBG Reagent	100868	819001
Sample Diluent	100870	819002
Calibrators	101111	81421
Controls	100875 & 100876	73110
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	73110	18.6	8.05-24.2
Control 2	73110	0.343	0.137-0.411
Panel 1	MF-011916-05	2.22	0.919-2.76
Panel 2	MF-011916-06	16.8	5.84-17.5

* 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	0.030
C	0.106
D	0.416
E	1.73
F	6.86
G	26.7

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Handwritten Signature]
Signature/Date
9/10/2018