

Product Number:	100860
Lot Number:	501219
Expiration:	14-Oct-2018
Platform(s):	HD-1



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

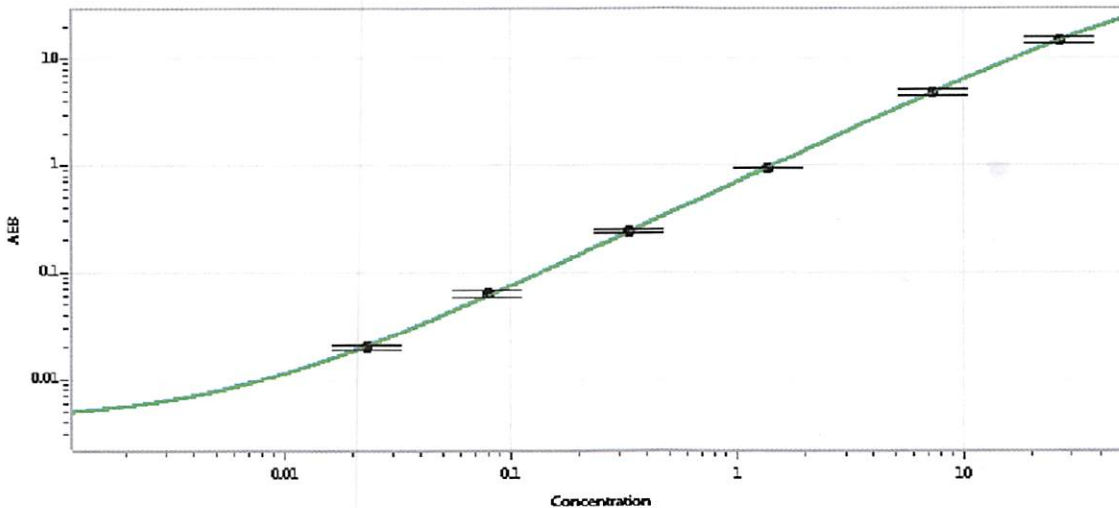
Component	Part Number	Lot Number
Bead Reagent	100865	81060
Detector Reagent	100862	81040
SBG Reagent	100868	81041
Sample Diluent	100870	81042
Calibrators	101111	72500/72790
Controls	100875 & 100876	73110
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	73110	17.3	12.7-19.1
Control 2	73110	0.288	0.217-0.325
Panel 1	MF-011916-05	2.18	1.25-2.42

* 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 0.023
C 0.079
D 0.335
E 1.38
F 7.37
G 26.7

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

S. Murugan
Signature/Date 01 MAY 2018