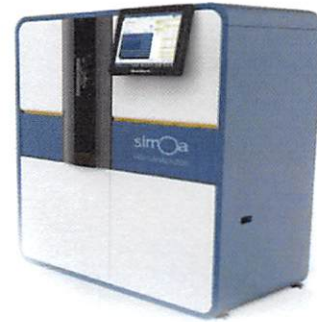


Product Number:	100860
Lot Number:	501272
Expiration:	23-Nov-2018
Platform(s):	HD-1



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

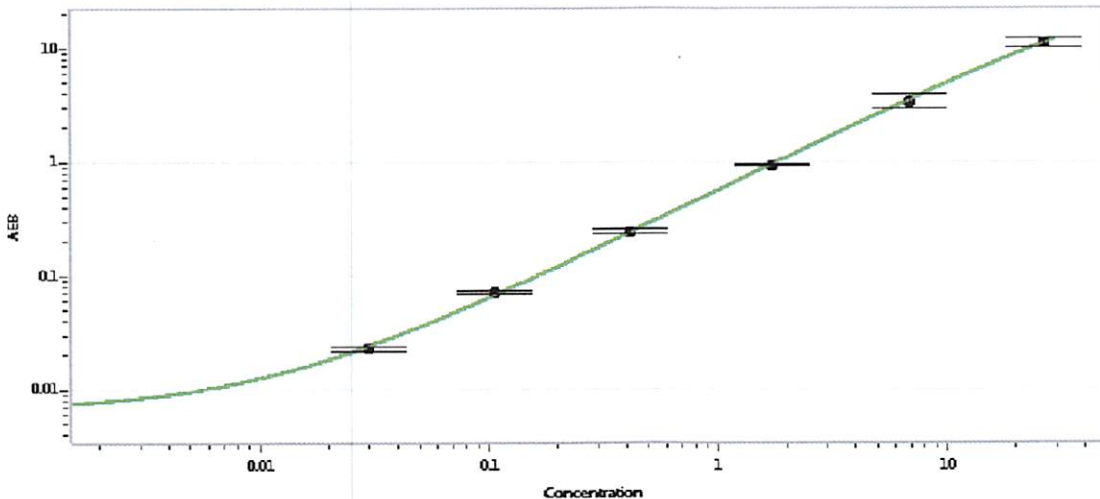
Component	Part Number	Lot Number
Bead Reagent	100865	81454
Detector Reagent	100862	81451
SBG Reagent	100868	81452
Sample Diluent	100870	81430
Calibrators	101111	81421
Controls	100875 & 100876	73110
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	73110	17.5	12.7-19.1
Control 2	73110	0.303	0.217-0.325
Panel 1	MF-011916-05	2.15	1.25-2.42
Panel 2	MF-011916-06	12.9	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	0.030
C	0.106
D	0.416
E	1.73
F	6.86
G	26.7

Review/Approval Sivasankari Murugan QC Supervisor [Signature] 05 Jun 2018
Name Title Signature/Date