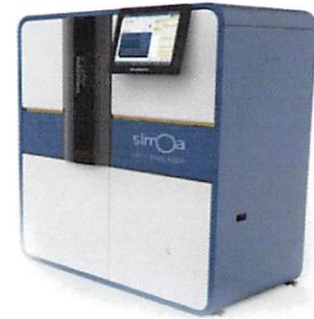


Product Number:	103186
Lot Number:	501213
Expiration:	27-Dec-2018
Platform(s):	HD-1 & SR-X



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

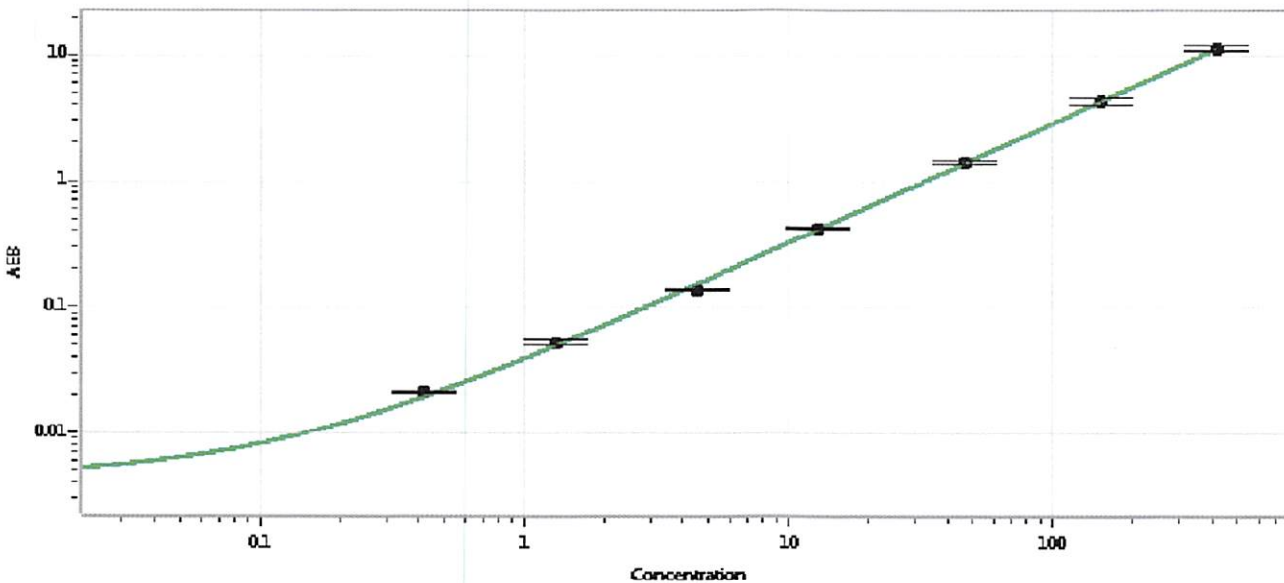
Component	Part Number	Lot Number
Bead Reagent	102246	81000
Detector Reagent	102248	80862
SBG Reagent	102250	80861
Sample Diluent	102252	80860
Reference Calibrators	102255	80250
Controls	102256 - 102257	80251
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80251	3.44	2.49-4.60
Control 2	80251	164	120-199
Panel 1	MF-092717-01	5.48	4.27-7.42
Panel 2	MF-092717-02	63.4	46.8-78.1
Panel 3	MF-092717-03	257	124-291

* 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	0.420
C	1.32
D	4.55
E	13.0
F	46.4
G	153
H	413

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date
18 APR 2018