

Product Number:	103186
Lot Number:	501350
Expiration:	20-Jul-2019
Platform(s):	HD-1 & SR-X



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

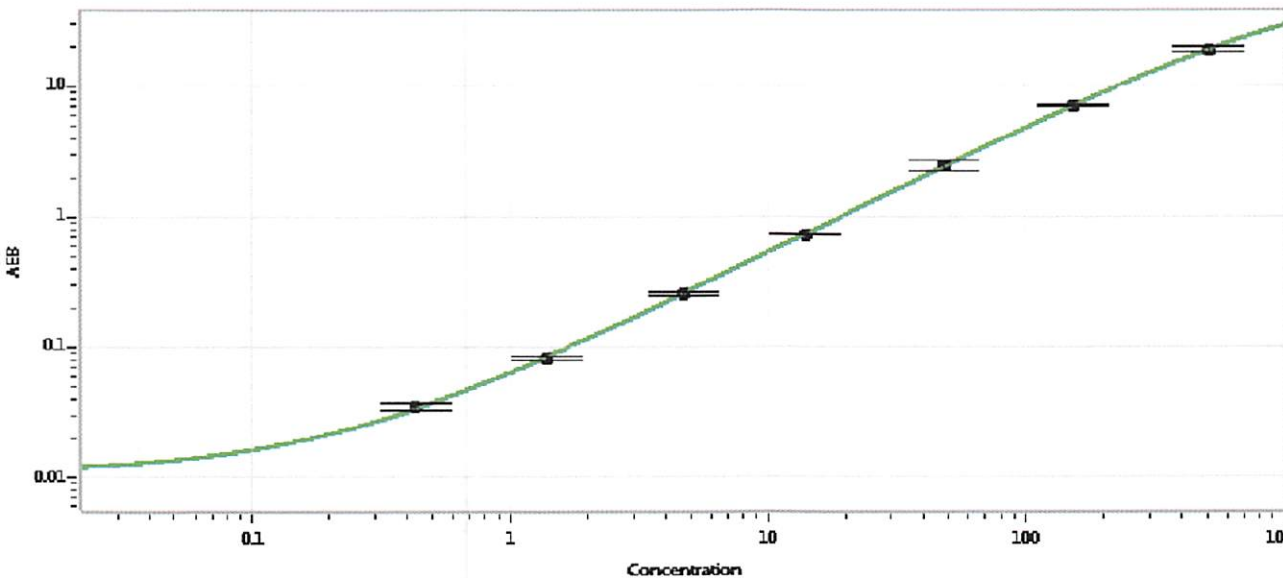
Component	Part Number	Lot Number
Bead Reagent	102246	820703
Detector Reagent	102248	820406
SBG Reagent	102250	820103
Sample Diluent	102252	820104
Reference Calibrators	102255	81570
Controls	102256 - 102257	81090
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	81090	4.47	3.26-6.87
Control 2	81090	176	128-257
Panel 1	MF-092717-01	6.14	4.27-7.42
Panel 2	MF-092717-02	66.3	46.8-78.1
Panel 3	MF-092717-03	249	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	0.431
C	1.40
D	4.75
E	14.0
F	48.5
G	153
H	516

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 01 Aug 2018
Signature/Date