

Product Number:	101580
Lot Number:	501307
Expiration:	10-Apr-2019
Platform(s):	HD-1 & SR-X



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

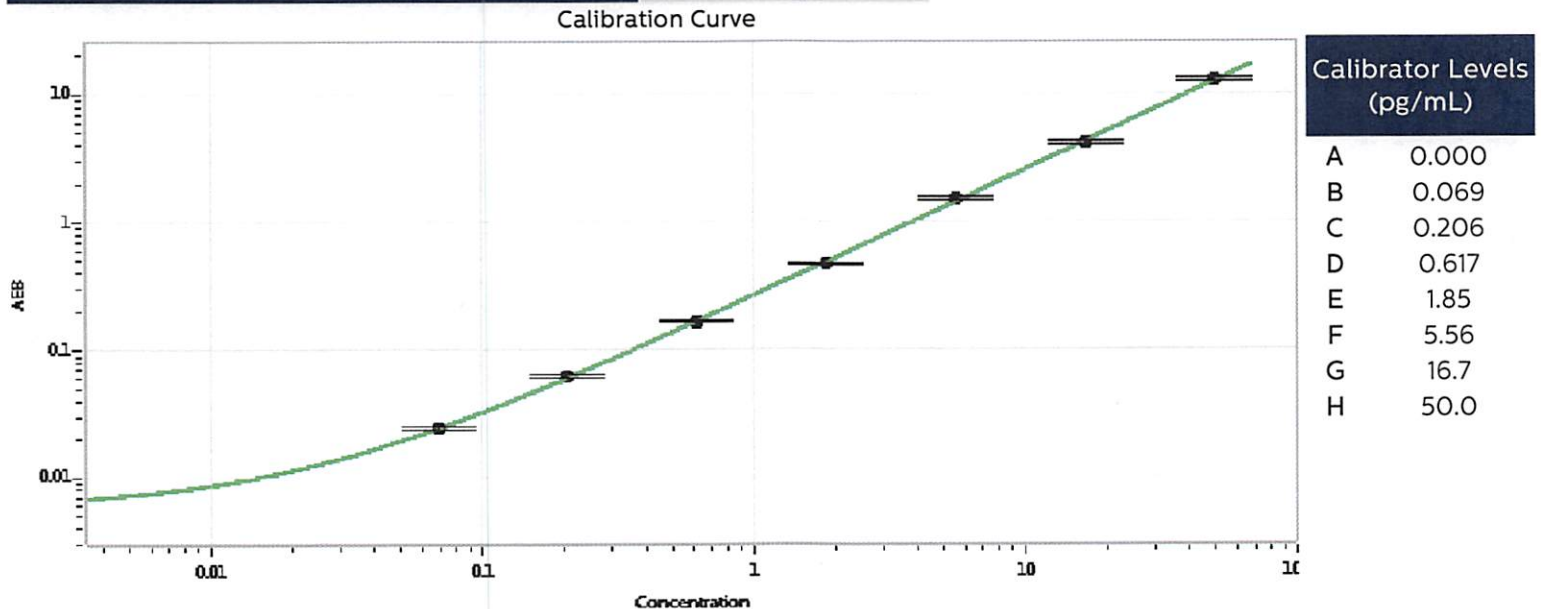
Component	Part Number	Lot Number
Bead Reagent	101582	818407
Detector Reagent	101581	81786
SBG Reagent	101583	81782
Sample Diluent	103201	81006
Calibrator Diluent	101585	81781
Calibrator Concentrate	100655	71510
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-042817-01	2.21	2.05-3.08
Control 2	MF-051617-02	44.0	35.8-53.7
Panel 1	MF-051017-03	25.3	20.9-32.4
Panel 2	MF-032217-11	1.31	1.17-2.07
Panel 3	MF-050117-05	1.98	1.19-2.11

\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Value Assignment (pg/mL)	1205
---	------



Review/Approval: Sivasankari Murugan QC Supervisor [Signature] 06/09/2018  
Name Title Signature/Date