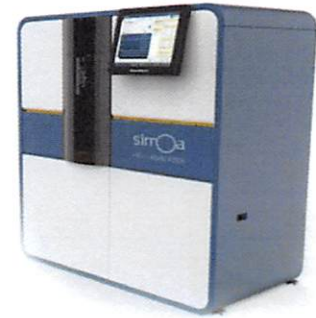


Product Number:	103186
Lot Number:	501279
Expiration:	2-Feb-2019
Platform(s):	HD-1 & SR-X



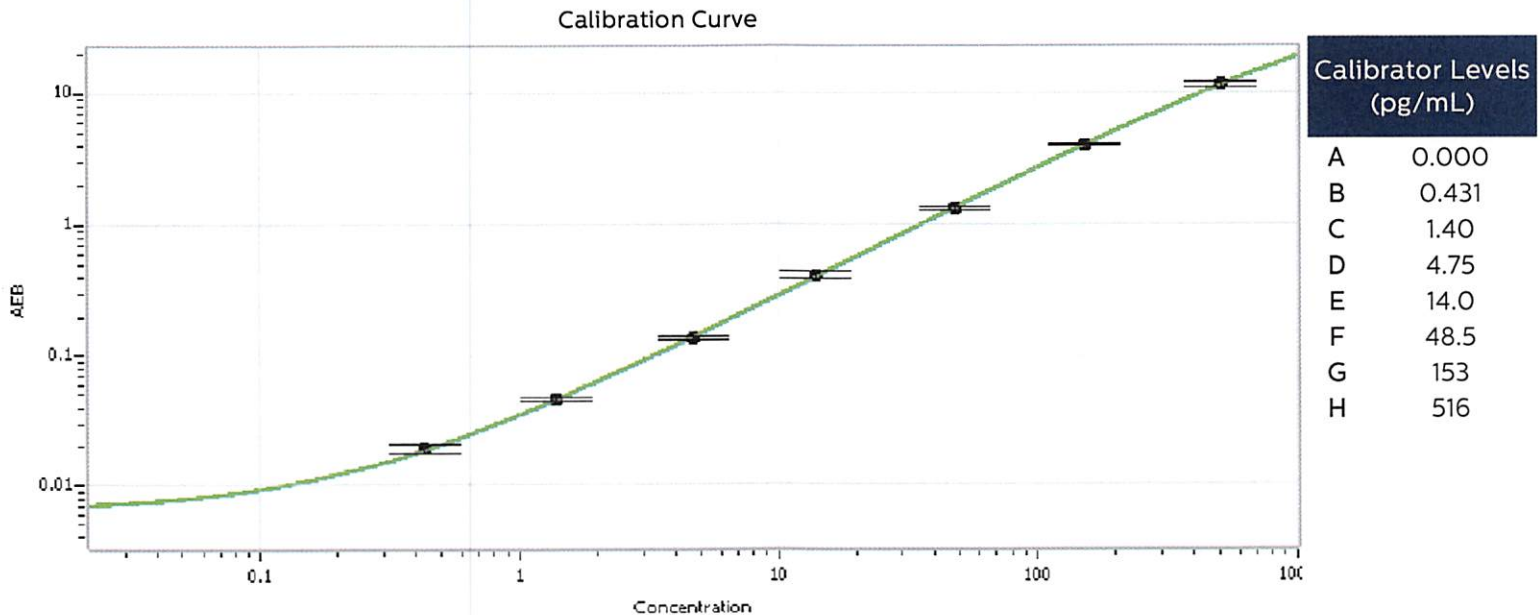
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102246	81552
Detector Reagent	102248	81450
SBG Reagent	102250	81227
Sample Diluent	102252	81223
Reference Calibrators	102255	81570
Controls	102256 - 102257	81090
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	81090	5.19	3.34-6.97
Control 2	81090	196	128-262
Panel 1	MF-092717-01	6.40	4.27-7.42
Panel 2	MF-092717-02	68.2	46.8-78.1
Panel 3	MF-092717-03	255	124-291

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



Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date  
14/Jan/2018