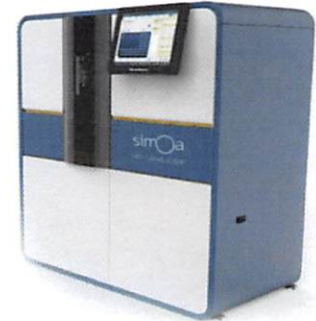


Product Number:	103186
Lot Number:	501419
Expiration:	29-Aug-2019
Platform(s):	HD-1 & SR-X



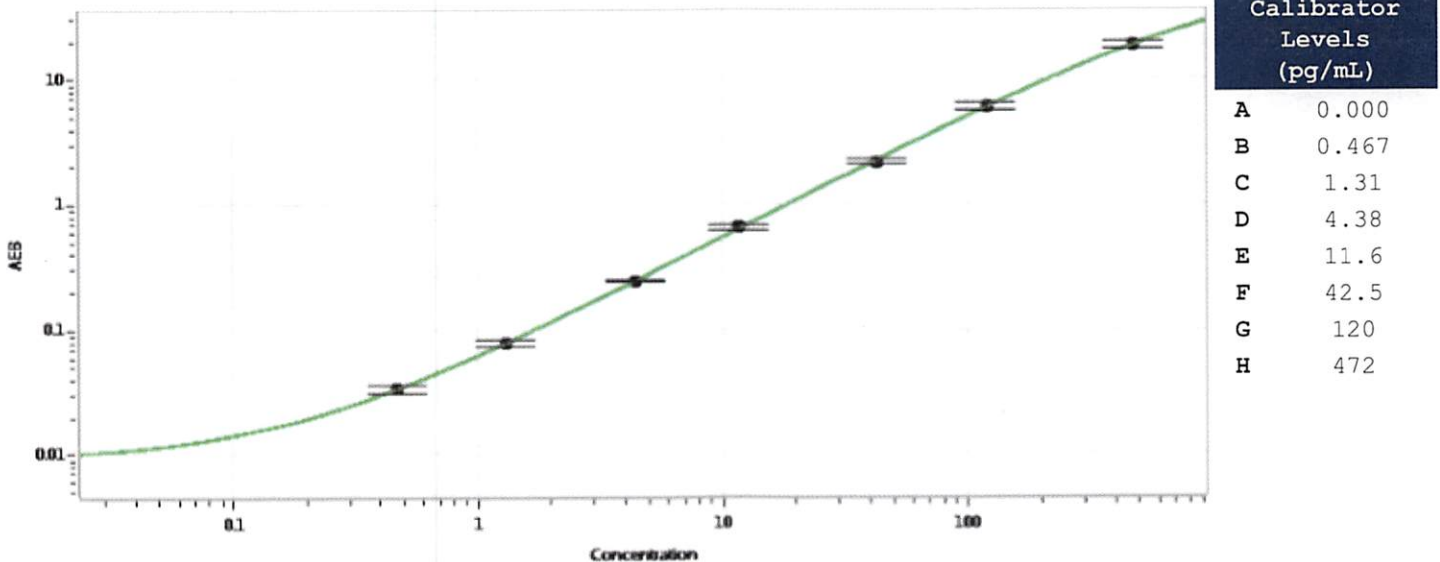
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102246	824903
Detector Reagent	102248	824203
SBG Reagent	102250	824101
Sample Diluent	102252	826205
Reference Calibrators	102255	822501
Controls	102256 - 102257	822806   823603
RGP Reagent	103159	N/A <sup>1</sup>

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

	Release Materials*			* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	822806	4.78	3.62-5.62	
Control 2	823603	154	118-176	
Panel 1-A	823301	5.07	3.99-6.68	
Panel 1-B	823302	6.47	4.74-7.59	
Panel 2	823601	93.7	56.9-100	
Panel 3	823602	341	226-402	

Calibration Curve



Review/Approval

Sivasankari Murugan QC Supervisor  
Name Title

*[Signature]* 01 OCT 2018  
Signature/Date