

Product Number:	102215
Lot Number:	501099
Expiration:	18-Jul-2018
Platform(s):	HD-1 & SR-X



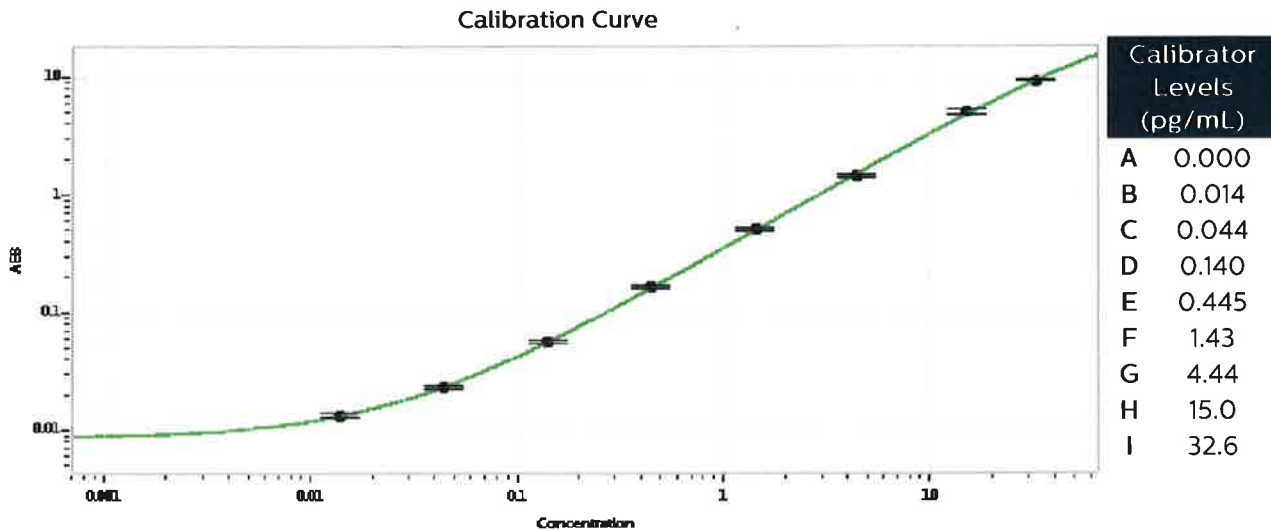
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102211	80181
Detector Reagent	102212	80240
SBG Reagent	102213	80180
Sample Diluent	102214	80190
Calibrators	101310	73250/ 73440
Controls	100702-100703	73260
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	73260	0.157	0.120-0.190
Control 2	73260	16.6	12.9-22.5
Panel 1	MF-092217-01	0.269	0.177-0.353
Panel 2	MF-092217-02	2.69	1.99-3.51
Panel 3	MF-101117-01	16.7	12.8-20.8

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



Review/Approval Sivasankari Murugan QC Supervisor [Signature] 05 FEB 2018  
 Name Title Signature/Date