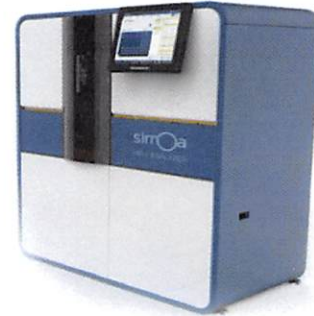


Product Number:	102215
Lot Number:	501230
Expiration:	20-Oct-2018
Platform(s):	HD-1 & SR-X



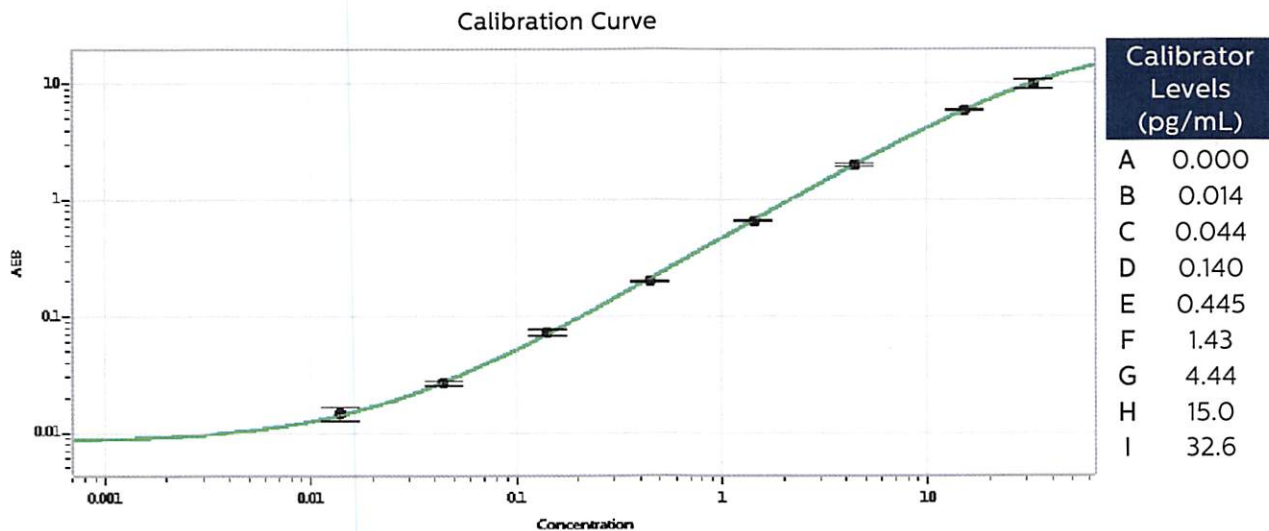
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102211	81101
Detector Reagent	102212	81135
SBG Reagent	102213	81103
Sample Diluent	102214	81102
Calibrators	101310	73250/ 73440
Controls	100702-100703	73260
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	73260	0.142	0.118-0.188
Control 2	73260	16.3	12.8-22.3
Panel 1	MF-092217-01	0.265	0.177-0.353
Panel 2	MF-092217-02	2.53	1.99-3.51
Panel 3	MF-101117-01	15.7	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
Name

QC Supervisor
Title

S. Murugan
Signature/Date 01MAY2018