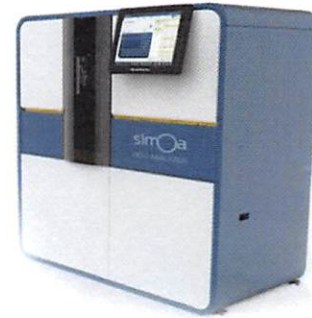


Product Number:	102215
Lot Number:	501270
Expiration:	20-Oct-2018
Platform(s):	HD-1 & SR-X



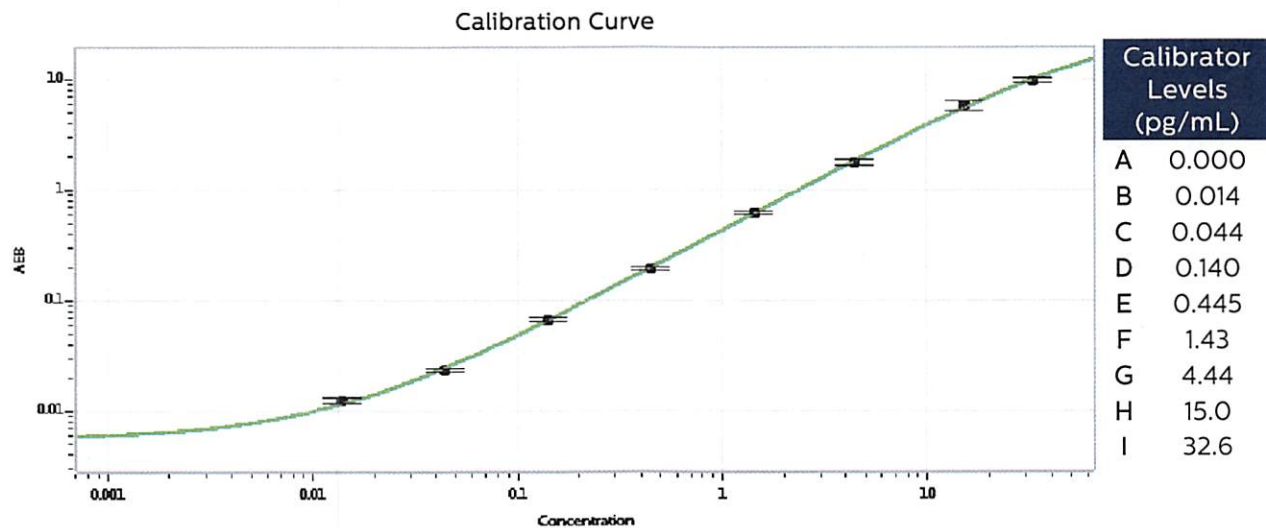
Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	73260/71301
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	0.152	0.118-0.188
Control 2	71301	14.6	11.8-18.3
Panel 1	MF-092217-01	0.250	0.177-0.353
Panel 2	MF-092217-02	2.58	1.99-3.51
Panel 3	MF-101117-01	17.0	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] 04/09/2018
 Name Title Signature/Date