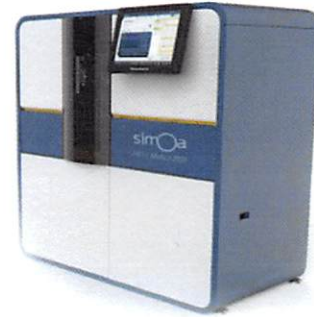


Product Number:	101478
Lot Number:	501286
Expiration:	10-Nov-2018
Platform(s):	HD-1 & SR-X



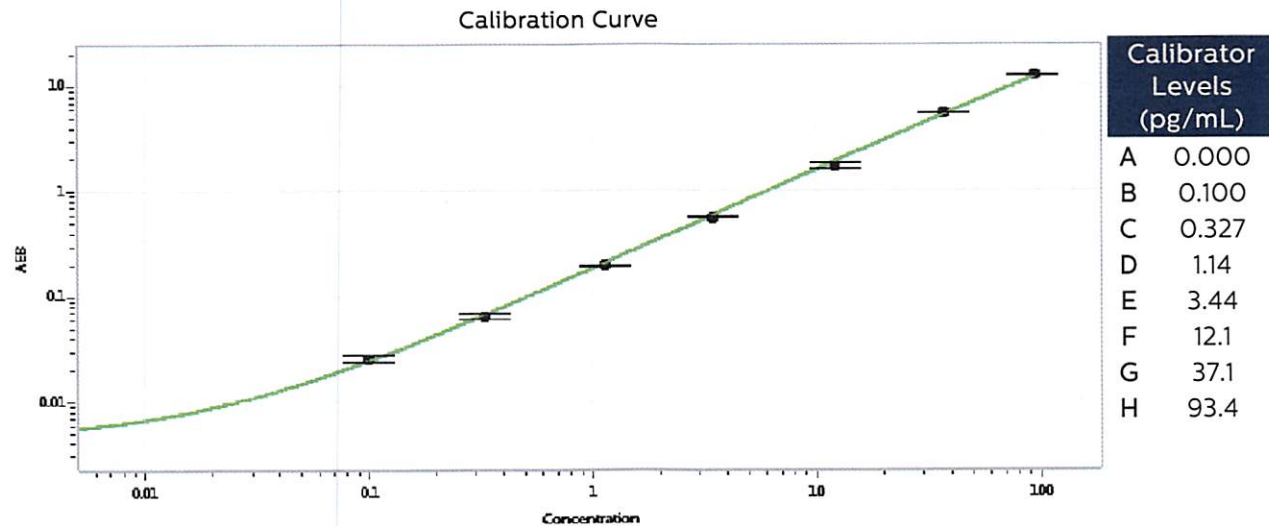
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101480	81620
Detector Reagent	101481	81303
SBG Reagent	101482	81304
Sample Diluent	101483	81295
Calibrators	100758-100765	80230
Controls	100768 & 100769	80231
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	80231	3.30	2.55-3.83
Control 2	80231	78.8	57.9-88.1
Panel 1	MF-051017-04	0.607	0.508-0.855
Panel 2	MF-042117-06	4.49	3.89-6.07
Panel 3	MF-042117-07	62.4	45.5-73.7

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



Review/Approval Sivasankari Murugan QC Supervisor [Signature] 14 Nov 2018
Name Title Signature/Date