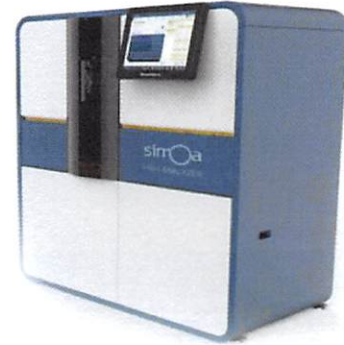


Product Number:	102215
Lot Number:	501296
Expiration:	20-Oct-2018
Platform(s):	HD-1 & SR-X



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

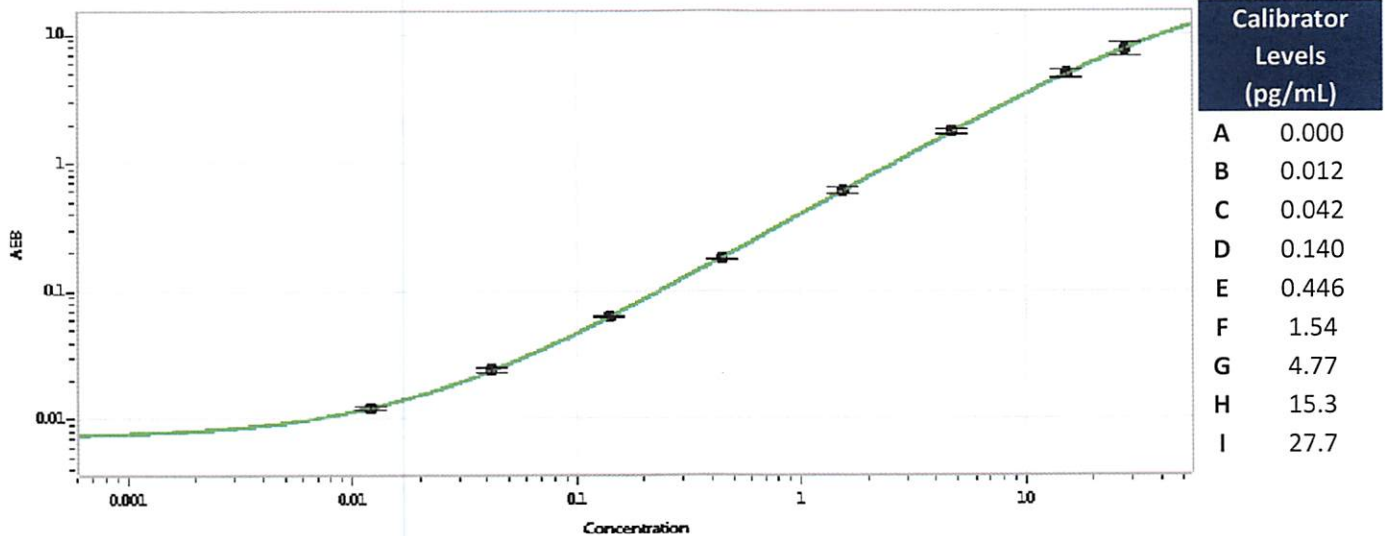
Component	Part Number	Lot Number
Bead Reagent	102211	81101
Detector Reagent	102212	81135
SBG Reagent	102213	81103
Sample Diluent	102214	81102
Calibrators	101310	81511/71300
Controls	100702-100703	81512
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	81512	0.142	0.115-0.173
Control 2	81512	17.9	12.1-19.3
Panel 1	MF-092217-01	0.263	0.177-0.353
Panel 2	MF-092217-02	2.85	1.99-3.51
Panel 3	MF-101117-01	20.8	12.8-20.8

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

Calibration Curve



Review/Approval: Sivasankari Murugan, QC Supervisor. Signature: [Signature], Date: 02/11/2018.