

Product Number:	102215
Lot Number:	501271
Expiration:	20-Oct-2018
Platform(s):	HD-1 & SR-X



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

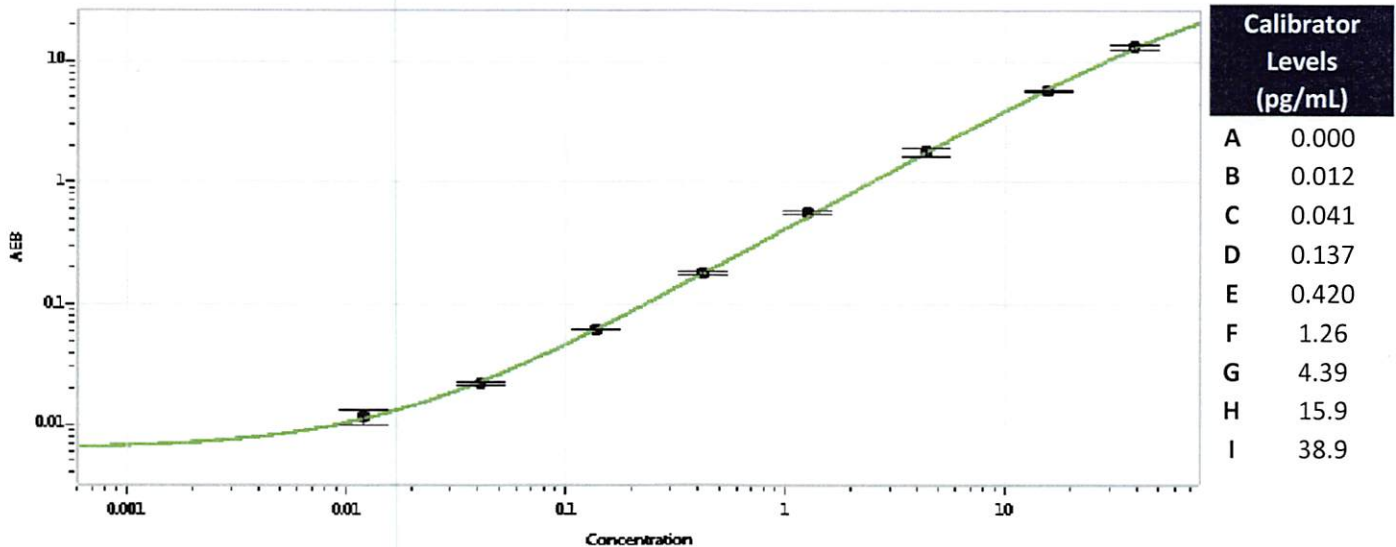
Component	Part Number	Lot Number
Bead Reagent	102211	81101
Detector Reagent	102212	81135
SBG Reagent	102213	81103
Sample Diluent	102214	81102
Calibrators	101310	70330/71300
Controls	100702-100703	70331
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	70331	0.142	0.105-0.157
Control 2	70331	13.9	11.5-17.4
Panel 1	MF-092217-01	0.265	0.177-0.353
Panel 2	MF-092217-02	2.74	1.99-3.51
Panel 3	MF-101117-01	15.2	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] 09 Jun 2018
 Name Title Signature/Date