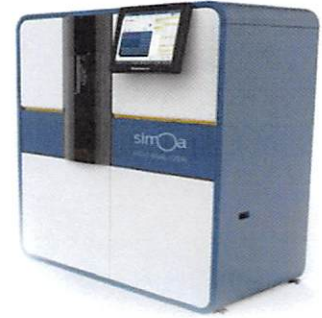


Product Number:	102215
Lot Number:	501424
Expiration:	10-Mar-2019
Platform(s):	HD-1 & SR-X



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

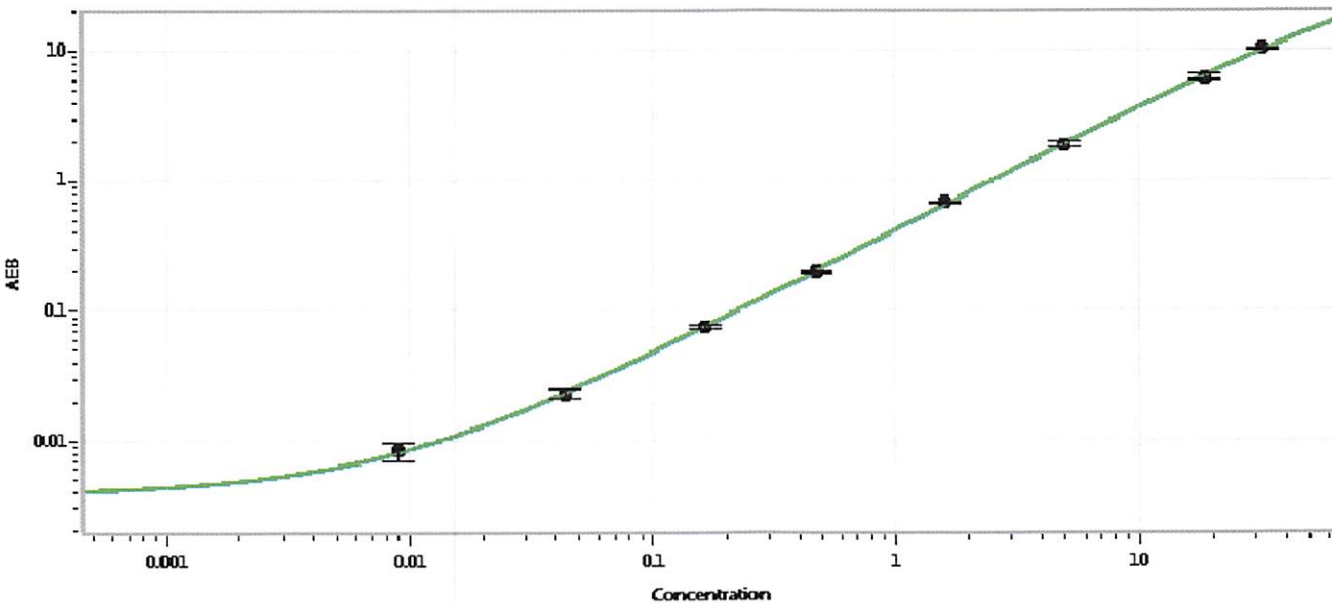
Component	Part Number	Lot Number
Bead Reagent	102211	826305
Detector Reagent	102212	826104
SBG Reagent	102213	825307
Sample Diluent	102214	825403
Calibrators	101310	829501
Controls	100702-100703	829502
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	829502	0.151	0.115-0.209
Control 2	829502	18.9	14.0-22.2
Panel 1	MF-092217-01	0.278	0.177-0.353
Panel 2	MF-092217-02	2.80	1.99-3.51
Panel 3	MF-101117-01	18.9	12.8-20.8

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.009
C	0.044
D	0.165
E	0.472
F	1.60
G	4.97
H	18.8
I	32.5

Review/Approval: Sivasankari Murugan (Name), QC Supervisor (Title), [Signature] (Signature), 29 OCT 2018 (Date)