

Product Number:	102530
Lot Number:	501199
Expiration:	9-Sep-2018
Platform(s):	HD-1



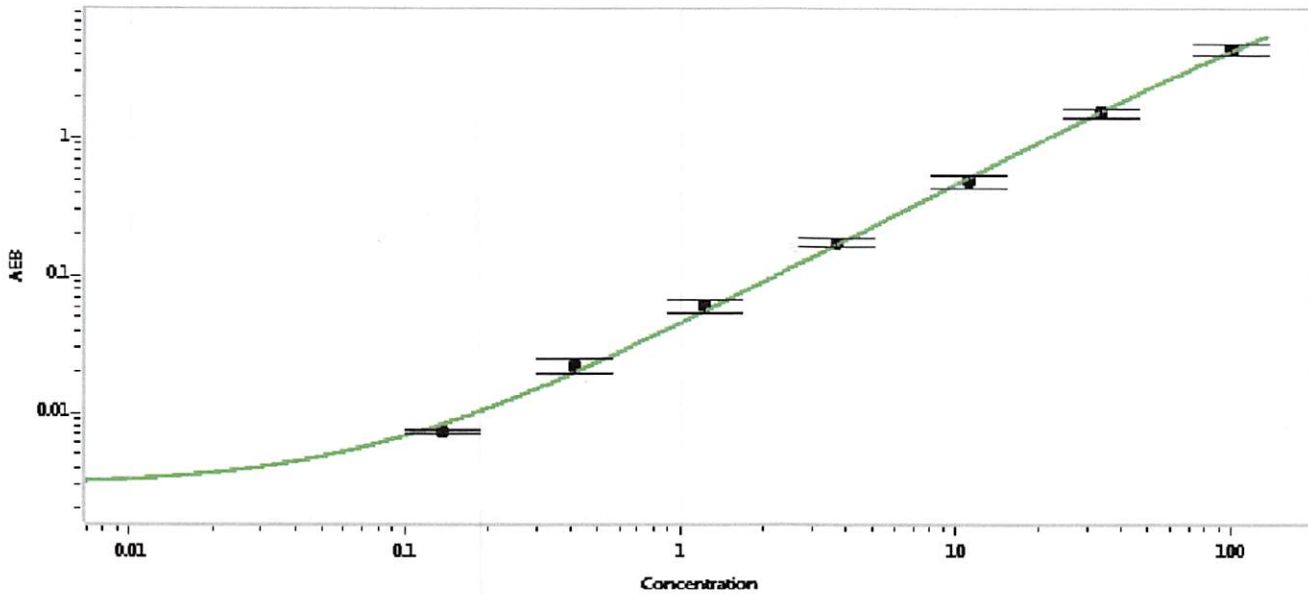
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102523	MF-032218-06
Detector Stock	102524	MF-032218-05
Calibrator Concentrate	102527	MF-040318-03
Calibrator Diluent	102525	MF-030918-02
Sample Diluent	102526	MF-032718-08

Generic Components - Included With Kit But Not Lot Number Specific		
Discovery Bead Diluent		101954
Discovery Detector Diluent		101955
Discovery SBG Diluent		101956
Discovery SBG Concentrate		101957
RGP Reagent		103159

Calibrator Concentrate Target Value	2 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval: Sivasankari Murugan (Name), QC Supervisor (Title), [Signature] 09 APR 2018 (Signature/Date)