

Product Number:	102669
Lot Number:	501378
Expiration:	13-Jan-2019
Platform(s):	HD-1



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

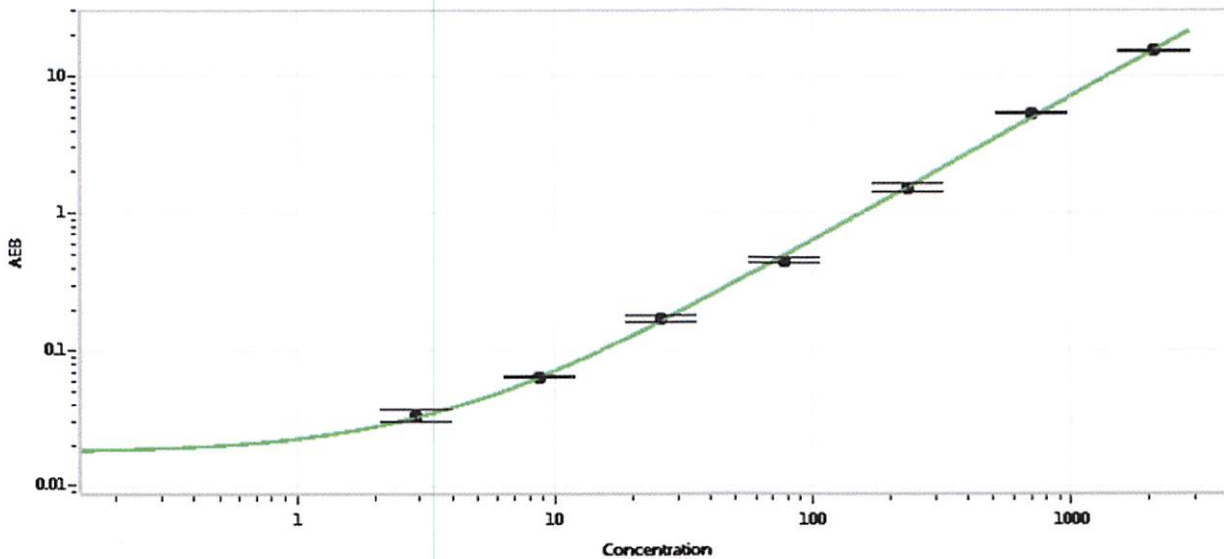
Component	Part Number	Lot Number
Bead Stock	102662	822007
Detector Stock	102663	821908
Calibrator Concentrate	102666	MF-020718-03
Calibrator Diluent	102664	819405
Sample Diluent	102665	821901

**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	60 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.88
C	8.64
D	25.9
E	77.8
F	233
G	700
H	2100

Review/Approval

Sivasankari Murugan QC Supervisor  
Name Title

*[Signature]* 10/11/2018  
Signature/Date