

Product Number:	102209
Lot Number:	501303
Expiration:	2-Dec-2018
Platform(s):	HD-1 and SR-X



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

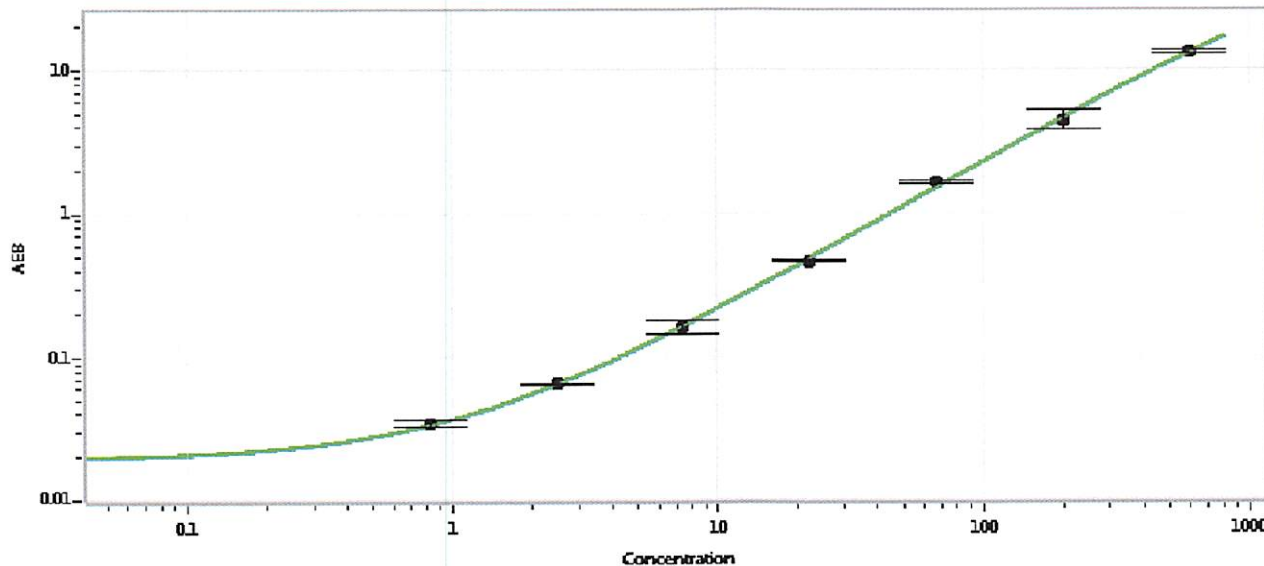
Component	Part Number	Lot Number
Bead Stock	102201	MF-062518-02
Detector Stock	102202	MF-062718-01
Calibrator Concentrate	102205	MF-120216-10
Calibrator Diluent	102203	MF-061918-04
Sample Diluent	102204	MF-061918-03

NOTE: For use on the HD-1 platform - refer to the Package Insert (PI-0077).  
For use on the SR-X platform - refer to the SR-X Kit Instructions (KI-0083).

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	12 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.823
C	2.47
D	7.41
E	22.2
F	66.7
G	200
H	600

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]* 28 Oct 2018  
Signature/Date