

Product Number:	102517
Lot Number:	503043
Expiration:	23-Dec-2021
Platform(s):	HD-1/HD-X & SR-X

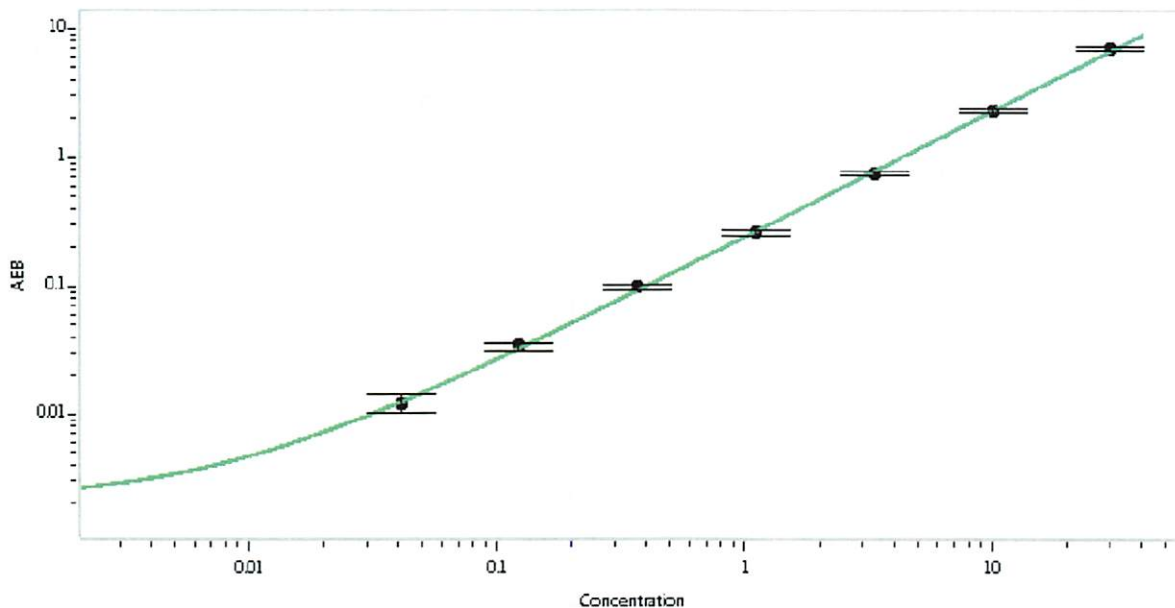


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102510	119201
Detector Stock	102511	119312
Calibrator Concentrate	102514	936010
Calibrator Diluent	102512	118311
Sample Diluent	102513	118310

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159

Calibrator Concentrate Target Value	600 pg/mL
-------------------------------------	-----------



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Junpil Park
Name

QC Supervisor
Title

Signature/Date 28 Jul 2021