

Product Number:	102794
Lot Number:	503308
Expiration:	15-Aug-2022
Platform(s):	HD-1/HD-X



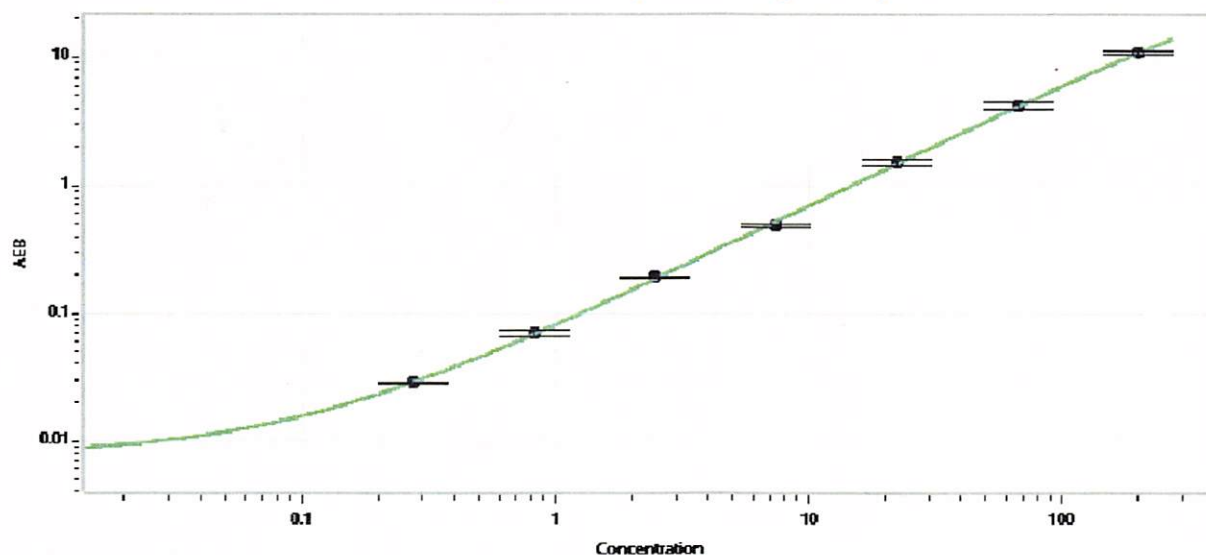
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102787	202205
Detector Stock	102788	203109
Calibrator Concentrate	102791	202306
Calibrator Diluent	102789	133918
Sample Diluent	102790	202304

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	4 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Shivani Goel

QC Manager

*Shivani Goel*

Name

Title

Signature/Date