

Product Number:	102318
Lot Number:	503075
Expiration:	20-May-2022
Platform(s):	HD-1/HDX

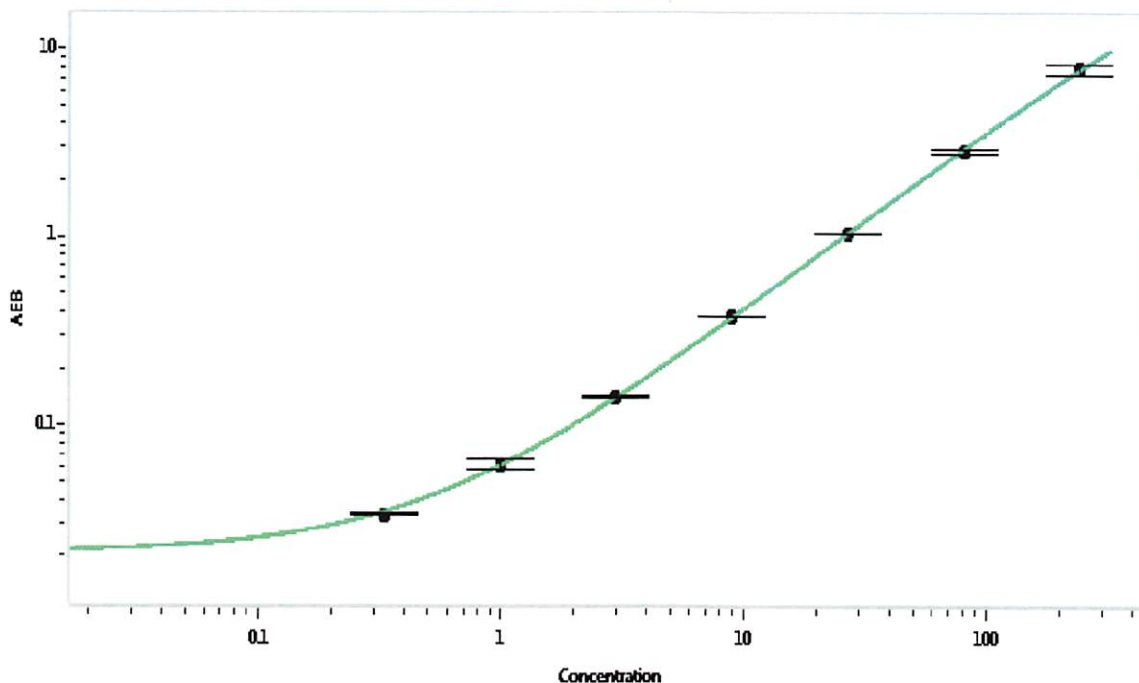
Component	Part Number	Lot Number
Bead Stock	102309	120905
Detector Stock	102310	121025
Calibrator Concentrate	102314	119507
Calibrator Diluent	102311	121027
Sample Diluent	102312	121026

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



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

Calibrator Concentrate Target Value	4,800 pg/mL
-------------------------------------	-------------



Calibrator Levels (pg/mL)	
A	0.000
B	0.329
C	0.988
D	2.96
E	8.89
F	26.7
G	80.0
H	240

Review/Approval

Junpil Park
Name

Supervisor, QC
Title

[Signature] 26 Aug 2021
Signature/Date