

Product Number:	102318
Lot Number:	503027
Expiration:	30-Apr-2022
Platform(s):	HD-1/HDX

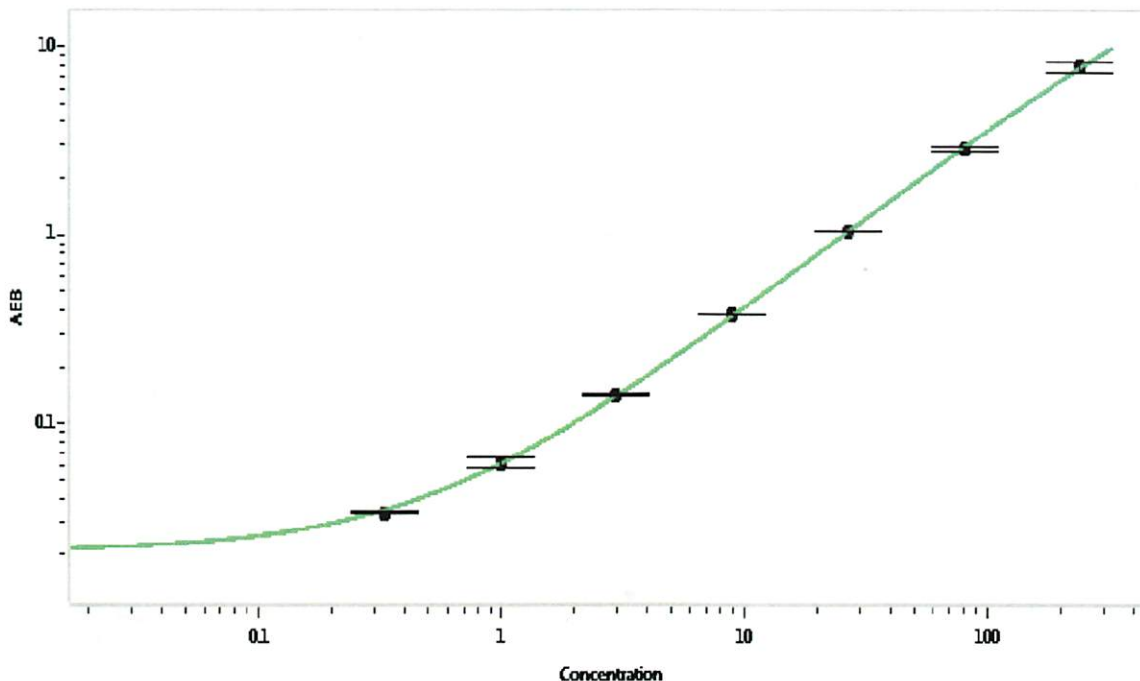


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102309	118401
Detector Stock	102310	118709
Calibrator Concentrate	102314	119507
Calibrator Diluent	102311	118407
Sample Diluent	102312	118406

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	4,800 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

Supervisor, QC

[Signature] 30 Jul 2021

Name

Title

Signature/Date