

Product Number:	102318
Lot Number:	503173
Expiration:	5-Aug-2022
Platform(s):	HD-1/HDX

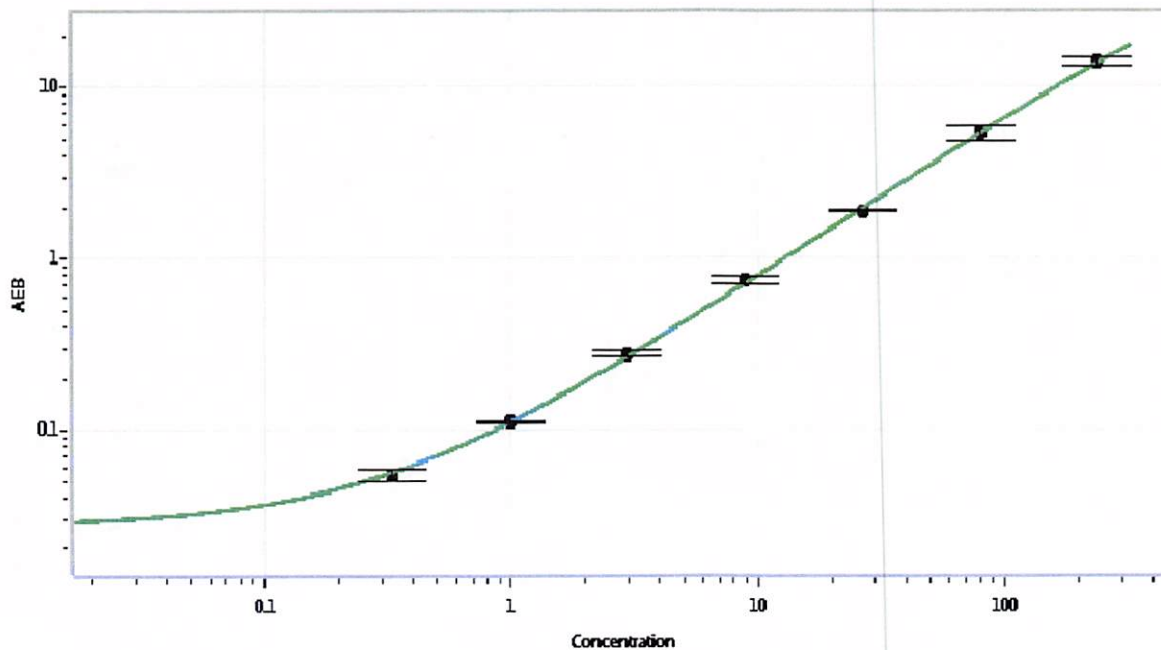
Component	Part Number	Lot Number
Bead Stock	102309	128013
Detector Stock	102310	128101
Calibrator Concentrate	102314	119507
Calibrator Diluent	102311	128103
Sample Diluent	102312	128102



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

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

Shivani Goel

Manager, QC

*Shivani Goel 26oct2021*

Name

Title

Signature/Date