

Product Number:	102343
Lot Number:	503326
Expiration:	3-Sep-2022
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	102337	203106
Detector Stock	102338	203602
Calibrator Concentrate	102342	132307
Calibrator Diluent	102339	133921
Sample Diluent	102340	203502

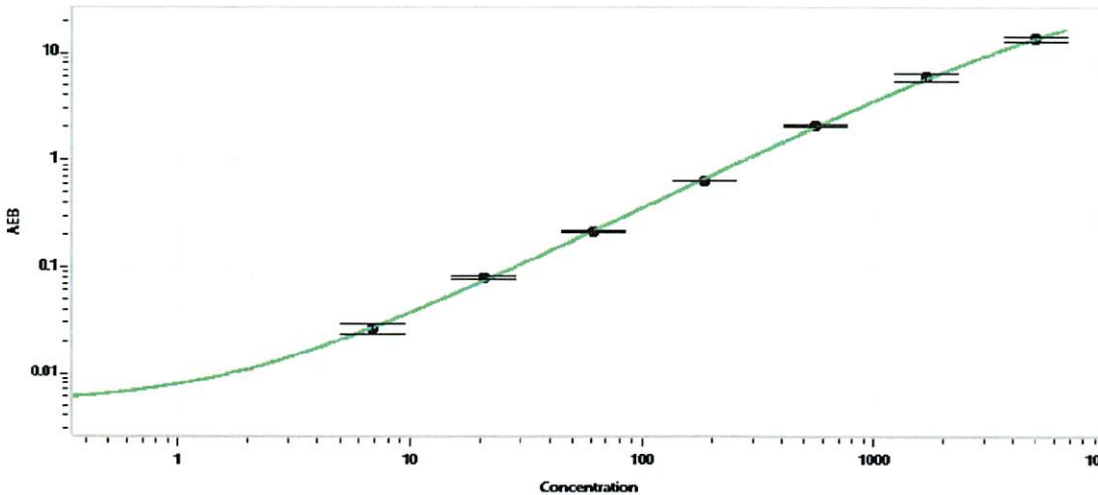
Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159



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

Calibrator Concentrate Target Value	100 ng/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.00
B	6.86
C	20.6
D	61.7
E	185
F	556
G	1667
H	5000

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

*Shivani Goel 18Feb2022*  
Signature/Date