

Product Number:	101984
Lot Number:	503351
Expiration:	8-Nov-2022
Platform(s):	HD-1/HD-X



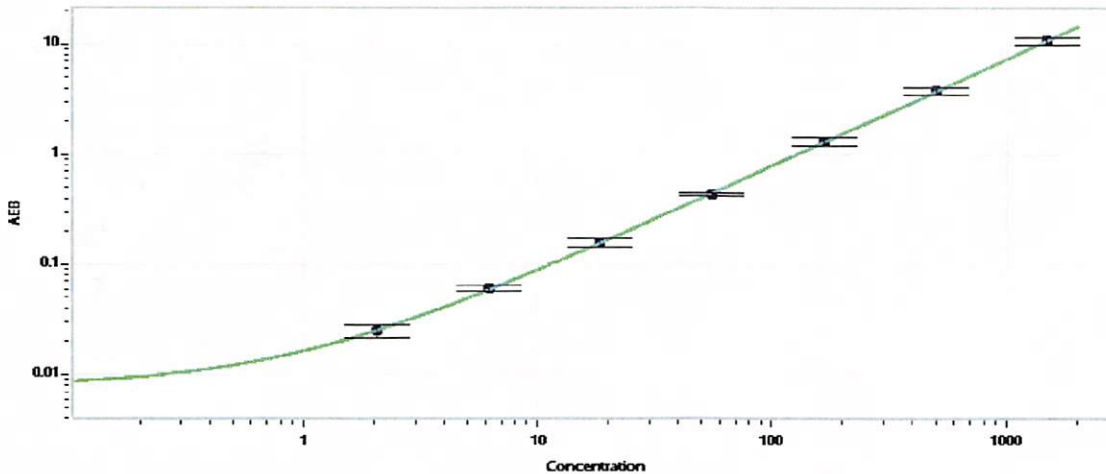
Component	Part Number	Lot Number
Bead Stock	101910	135504
Detector Stock	101911	204501
Calibrator Concentrate	101983	108306
Calibrator Diluent	101908	135604
Sample Diluent	101909	135603

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 / HD-X Analyzer.

Calibrator Concentrate Target Value	30 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	2.06
C	6.17
D	18.5
E	55.6
F	167
G	500
H	1500

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 31 Mar 2022  
Signature/Date