

Product Number:	102336
Lot Number:	504298
Expiration:	2-Mar-25
Platform(s):	SR-X and HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	102330	326406
Detector Stock	102331	326905
Calibrator Concentrate	102335	326601
Calibrator Diluent	102332	324511
Sample Diluent	102333	326809

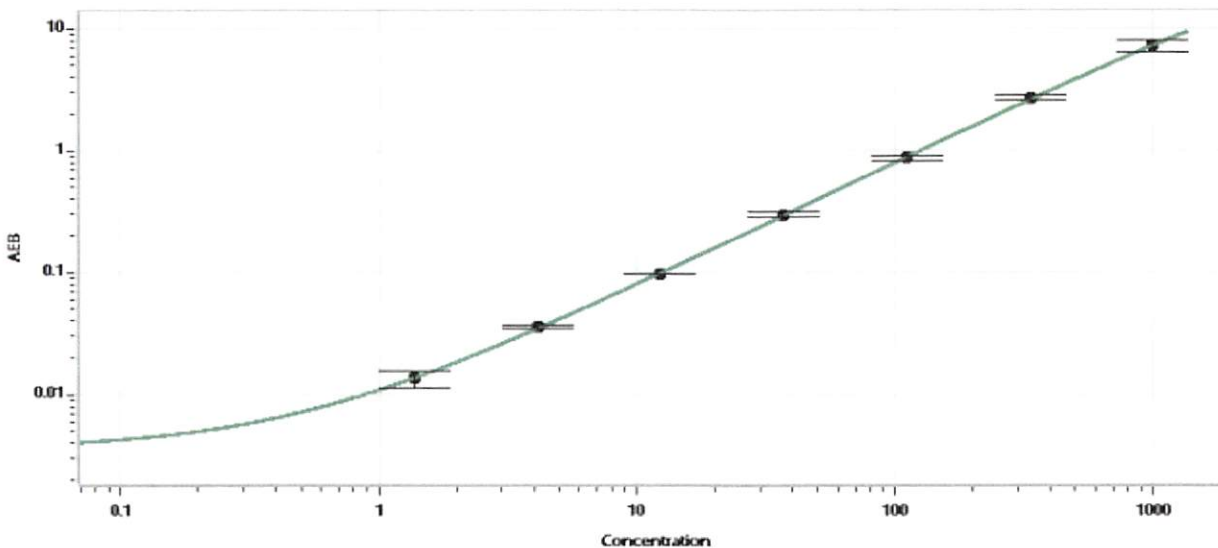
Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159



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

Calibrator Concentrate Target Value	20
--	----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.12
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval Cody Carlson QC Manager [Signature] 26 JUN 2024
 Name Title Signature/Date