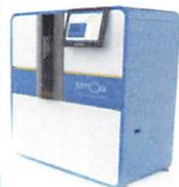


Product Number:	102343
Lot Number:	503531
Expiration:	20-Apr-2023
Platform(s):	HD-1/HD-X



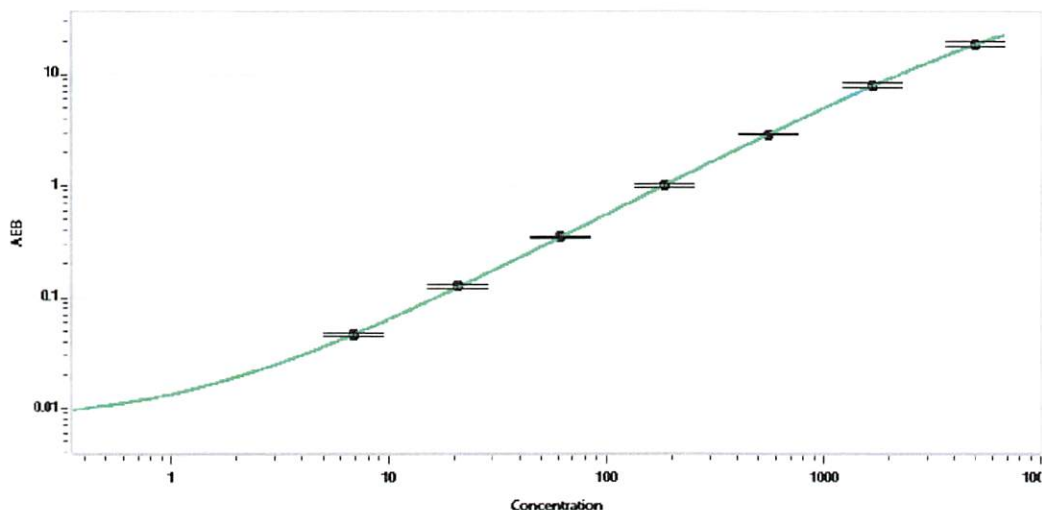
Component	Part Number	Lot Number
Bead Stock	102337	217805
Detector Stock	102338	217907
Calibrator Concentrate	102342	132307
Calibrator Diluent	102339	212417
Sample Diluent	102340	217908

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

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

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan / 14 JUL 2022
Signature/Date