

Product Number:	102648
Lot Number:	503556
Expiration:	10-Aug-2023
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Stock	102641	234310
Detector Stock	102642	225104
Calibrator Concentrate	102645	204607
Calibrator Diluent	102643	223013
Sample Diluent	102644	227216
Detector Diluent	102734	225103

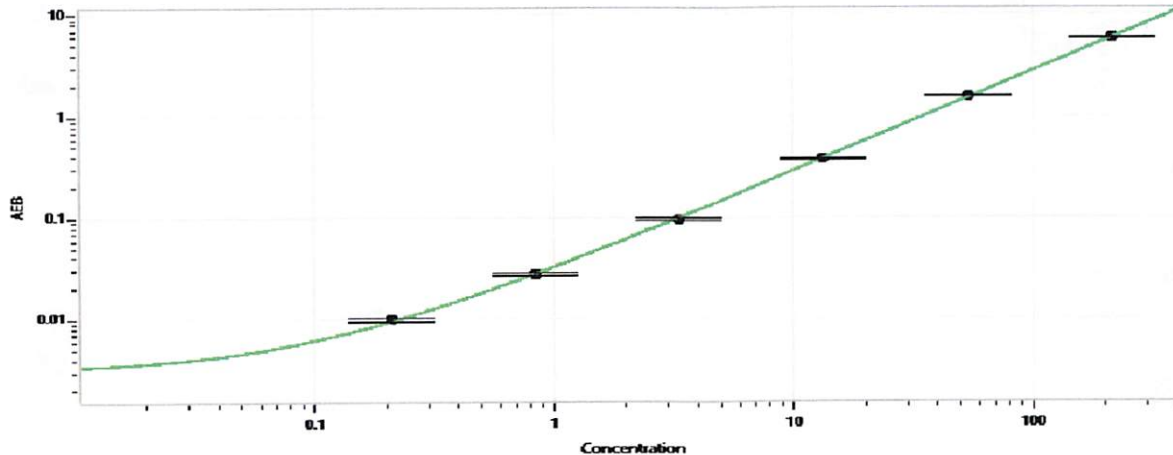
Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159

Calibrator Concentrate Target Value (pg/mL)	3,000
--	-------



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

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.210
C	0.840
D	3.36
E	13.4
F	53.8
G	215

Review/Approval

Sonia Khillan
Name

QC
Title

Sonia Khillan 12/1/2022
Signature/Date