

Product Number:	102209
Lot Number:	503370
Expiration:	24-Feb-2023
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Stock	102201	205714
Detector Stock	102202	205514
Calibrator Concentrate	102205	222709
Calibrator Diluent	102203	205518
Sample Diluent	102204	205517

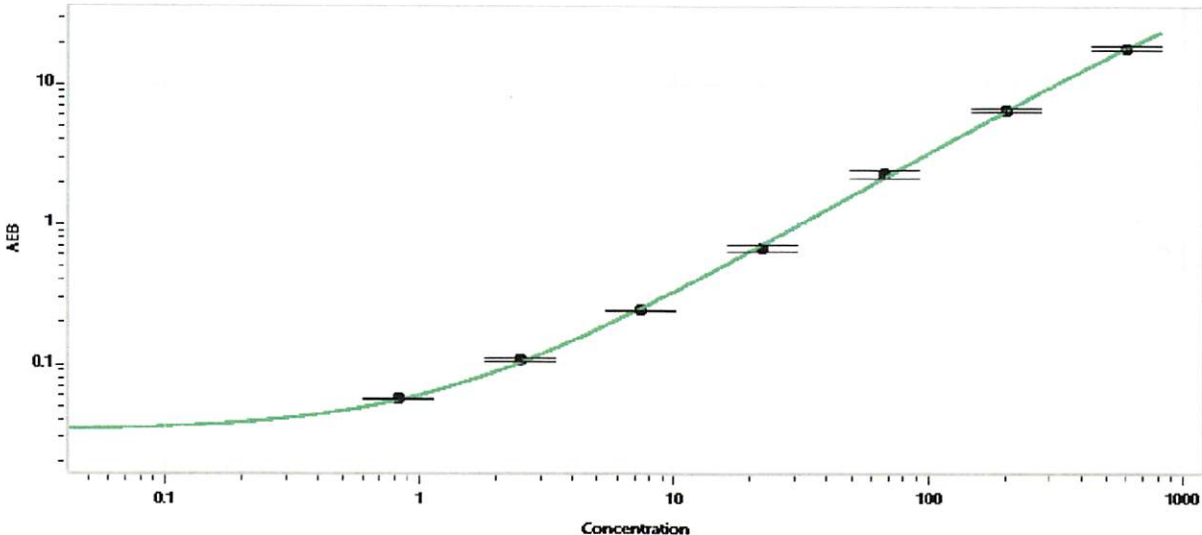
Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159



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

Calibrator Concentrate Target Value (ng/mL)	12 ng/mL
---	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.823
C	2.47
D	7.41
E	22.2
F	66.7
G	200
H	600

Review/Approval

Sonia Khillan
Name

QC Manager Sonia Khillan
Title

Sonia Khillan / 26 OCT 2022
Signature/Date