

Product Number:	102794
Lot Number:	503442
Expiration:	23-Apr-2023
Platform(s):	HD-1/HD-X



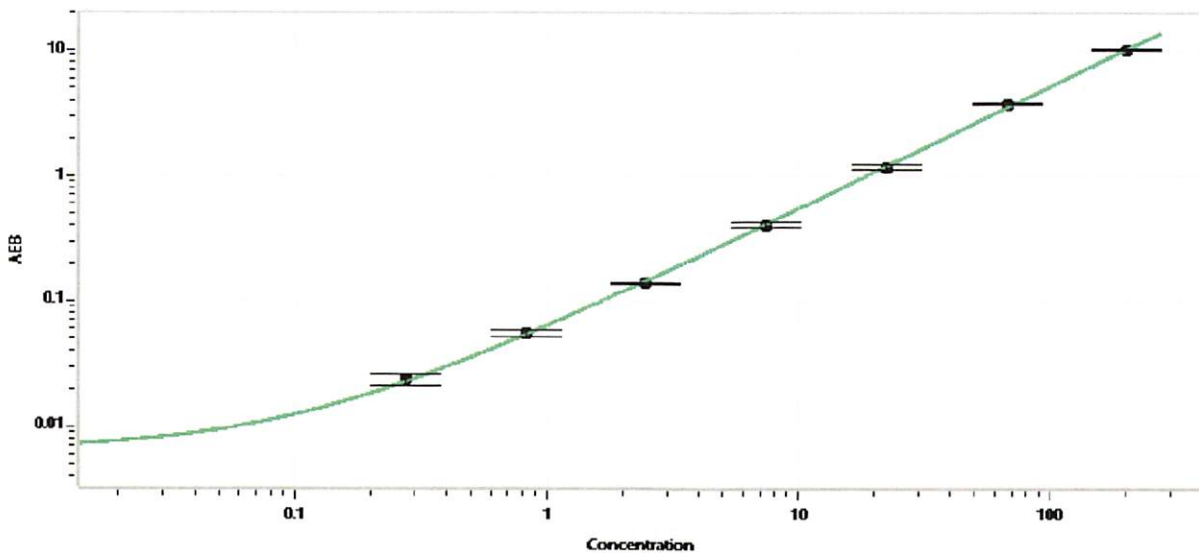
Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number
Bead Stock	102787	211301
Detector Stock	102788	211309
Calibrator Concentrate	102791	211006
Calibrator Diluent	102789	230515
Sample Diluent	102790	211310

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

Calibrator Concentrate Target Value (ng/mL)	4.0
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Sonia Khillan
Name

QC
Title

Sonia Khillan 1/10NOV2022
Signature/Date