

Product Number:	102635
Lot Number:	503765
Expiration:	10-Aug-2023
Platform(s):	HD-1/ HD-X



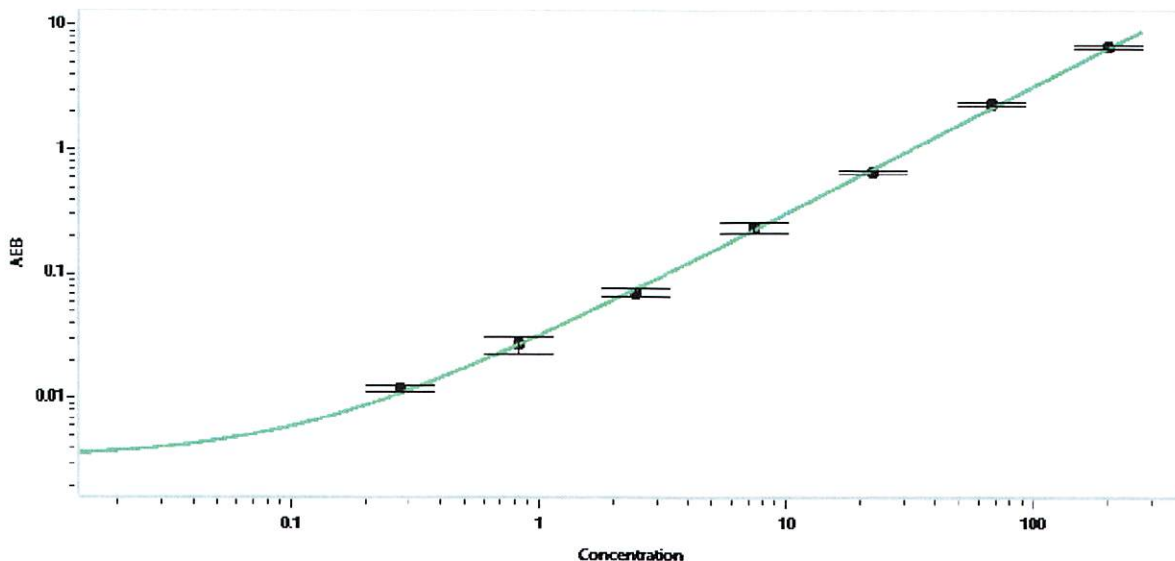
Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number
Bead Stock	102628	234805
Detector Stock	102629	235002
Calibrator Concentrate	102632	205503
Calibrator Diluent	102630	234017
Sample Diluent	102631	300602

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 ng/mL
--	---------

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 Manager
Title

Sonia Khillan 23 JAN 2023
Signature/Date