

Product Number:	101824
Lot Number:	503712
Expiration:	26-Jul-2023
Platform(s):	HD-1/ HD-X



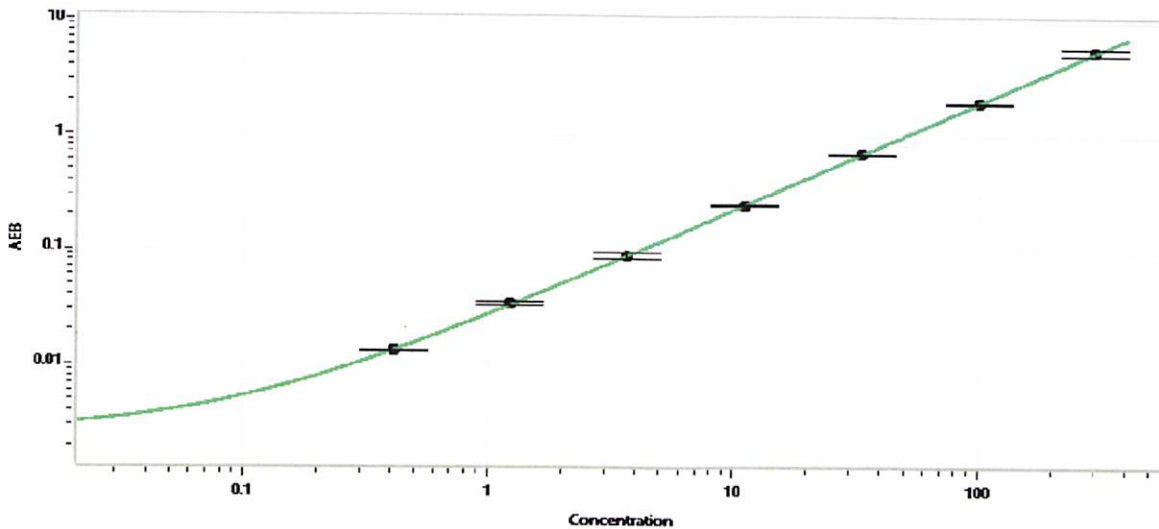
Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number
Bead Stock	101929	228003
Detector Stock	101930	228302
Calibrator Concentrate	101819	127003
Calibrator Diluent	101943	227714
Sample Diluent	101944	227231

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 (pg/mL)	6000 pg/mL
---	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.412
C	1.23
D	3.70
E	11.1
F	33.3
G	100
H	300

Review/Approval

*Sonia Khillan*  
Name

QC  
Title

*Sonia Khillan 1/15 NOV 2022*  
Signature/Date