

Product Number:	102475
Lot Number:	502859
Expiration:	26-Jan-2022
Platform(s):	HD-1/HD-X



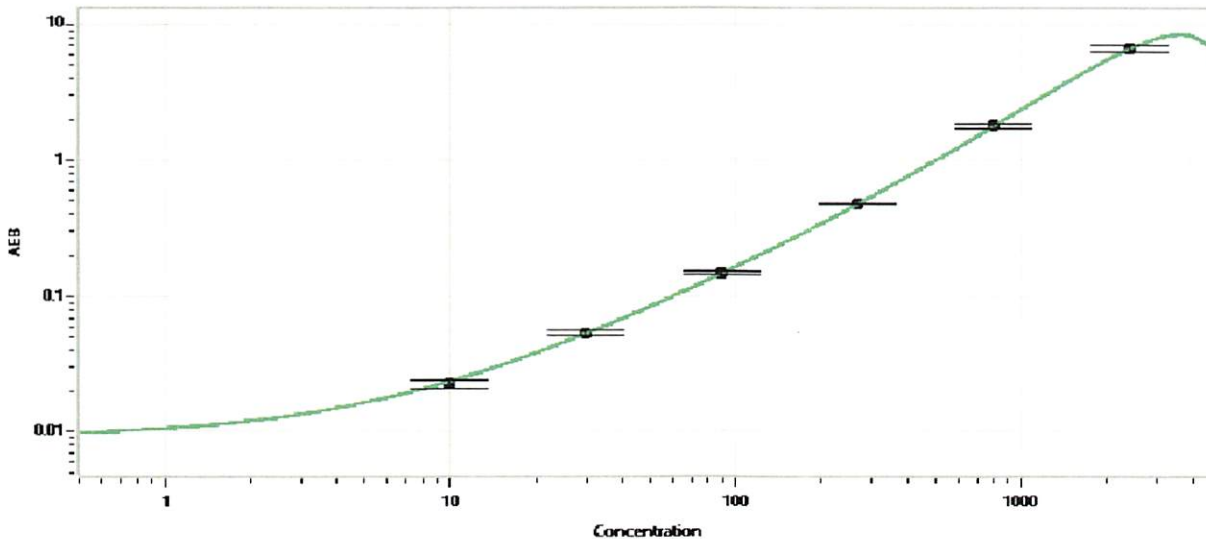
Data below represents results generated on the Simoa™ HD-X Analyzer.

Component	Part Number	Lot Number
Bead Stock	102468	109817
Detector Stock	102469	110004
Calibrator Concentrate	102472	004511
Calibrator Diluent	102470	108128
Sample Diluent	102471	109021

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	9.88
C	29.6
D	88.9
E	267
F	800
G	2400

Review/Approval



 23 Apr 2021
 Name Title Signature/Date