

Product Number:	102700
Lot Number:	503728
Expiration:	10-Aug-2023
Platform (s):	HD-1/HD-X



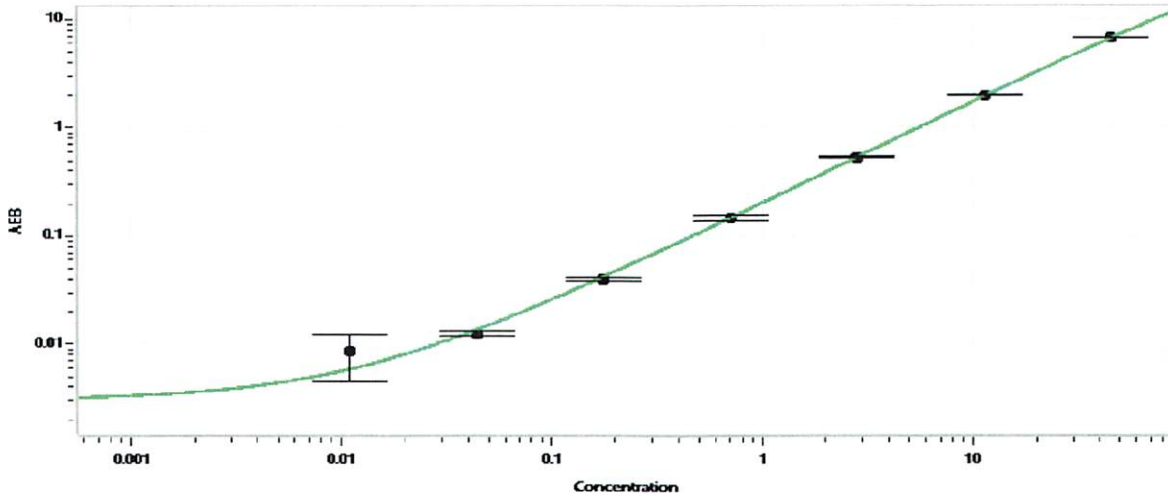
Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number
Bead Stock	102693	229905
Detector Stock	102694	230105
Calibrator Concentrate	102697	206403
Calibrator Diluent	102695	230107
Sample Diluent	102696	230108

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)	1000
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.011
C	0.044
D	0.176
E	0.703
F	2.81
G	11.3
H	45.0

Review/Approval

*Matt Koubouris*  
Name

QC Supervisor  
Title

*[Signature]* 06Feb2023  
Signature/Date