

Product Number:	102209
Lot Number:	503279
Expiration:	18-Aug-2022
Platform(s):	HD-1/HD-X & SR-X



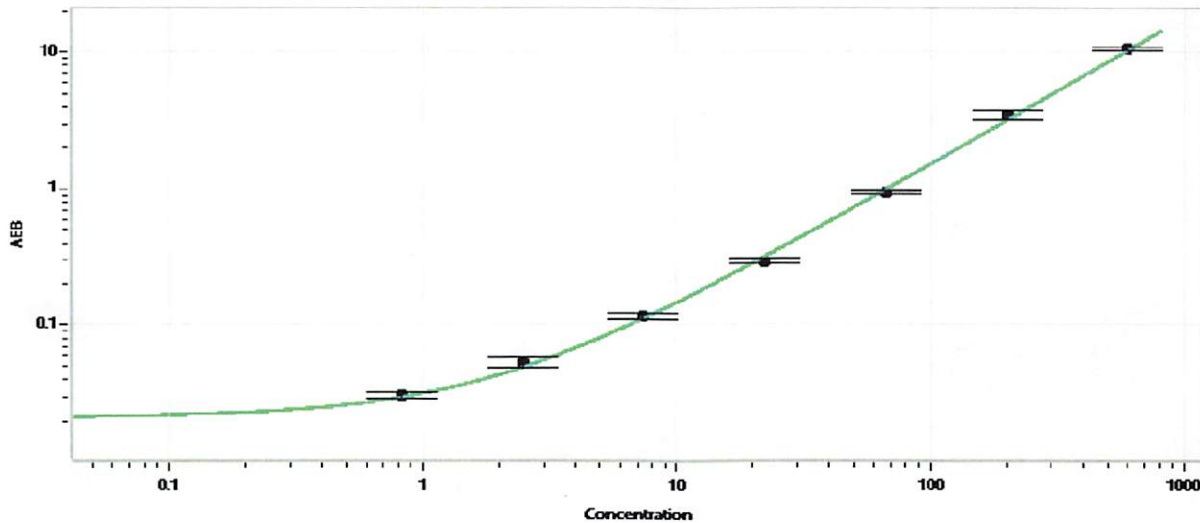
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102201	135303
Detector Stock	102202	135405
Calibrator Concentrate	102205	023103
Calibrator Diluent	102203	135309
Sample Diluent	102204	135308

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.823
C	2.47
D	7.41
E	22.2
F	66.7
G	200
H	600

Review/Approval

*Wenkin, Tu*

Name

QC

Title

*[Signature]* 14 Jun 2022

Signature/Date