

Product Number:	102570
Lot Number:	503368
Expiration:	28-Feb-2023
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	102563	225114
Detector Stock	102564	206006
Detector Diluent	102590	205908
Calibrator Concentrate	102567	133306
Calibrator Diluent	102565	217220
Sample Diluent	102566	217219

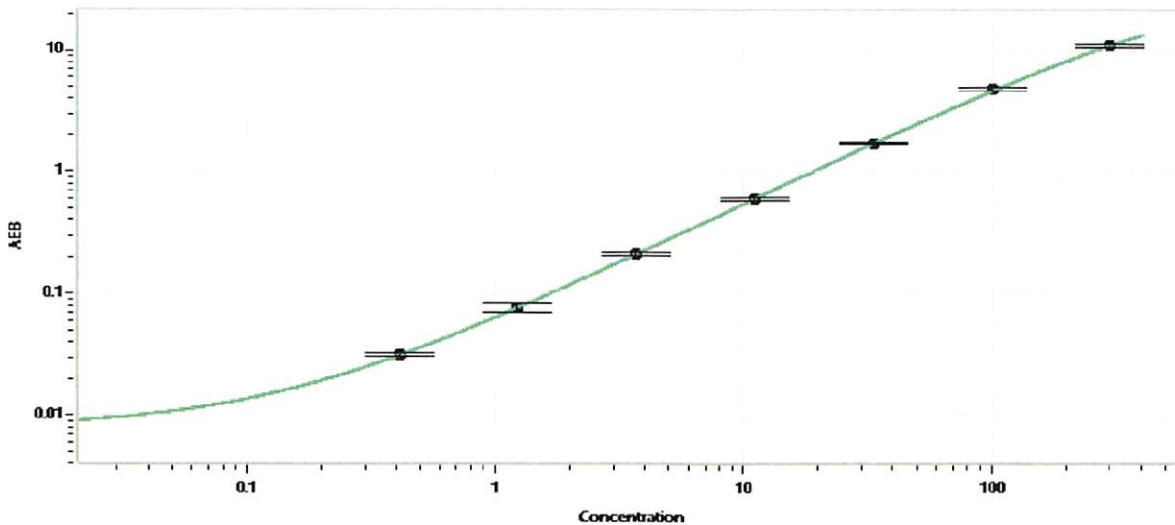


Data below represents results generated on the Simoa® HD-X Analyzer.

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159

Calibrator Concentrate Target Value (ng/mL)	6.0 ng/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 Manager Title, Sonia 12/16/12012022 Signature/Date