

Product Number:	102336
Lot Number:	503879
Expiration:	4-Apr-2024
Platform(s):	HD-1 / HD-X & SR-X

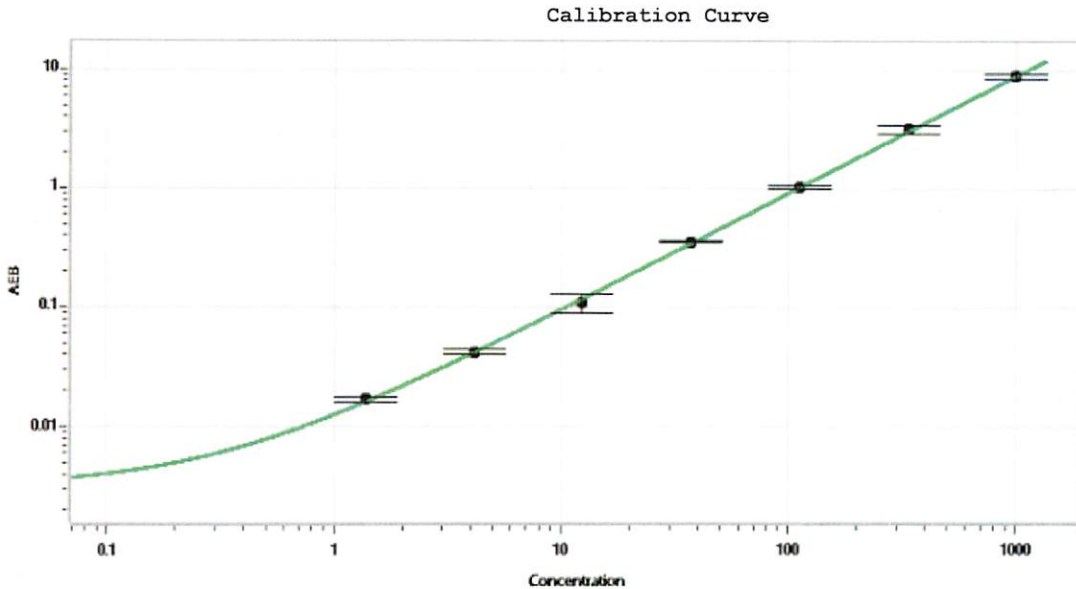


Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number	Control Lot Number
Bead Stock	102330	312103	312103A
Detector Stock	102331	312303	312303A
Calibrator Concentrate	102335	312406	312406A
Calibrator Diluent	102332	312316	312316A
Sample Diluent	102333	312310	312310A

Generic Components Included With Kit But Not Lot Number Specific		
Discovery Bead Diluent	101954	N/A
Discovery Detector Diluent	101955	N/A
Discovery SBG Diluent	101956	N/A
Discovery SBG Concentrate	103297	N/A
RGP Reagent	103159	N/A

Calibrator Concentrate Target Value (ng/mL)	20
--	----



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.11
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia K. P. Khillan 16 MAY 2023
Signature/Date