

Product Number:	102336
Lot Number:	503710
Expiration:	30-Sep-2023
Platform(s):	HD-1 & HD-X



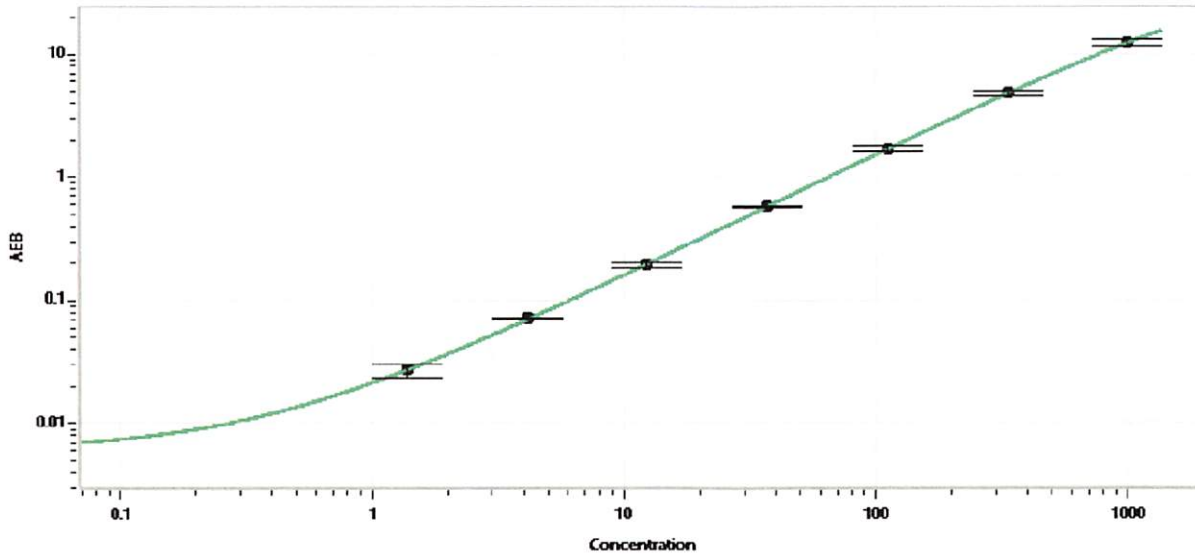
Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number
Bead Stock	102330	227310
Detector Stock	102331	227610
Calibrator Concentrate	102335	227614
Calibrator Diluent	102332	227613
Sample Diluent	102333	227611

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
---	----

Calibration Curve



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 Khillan / 10 NOV 2022* Signature/Date