

Product Number:	102039
Lot Number:	503450
Expiration:	14-Apr-2023
Platform(s):	HD-1/HD-X & SR-X



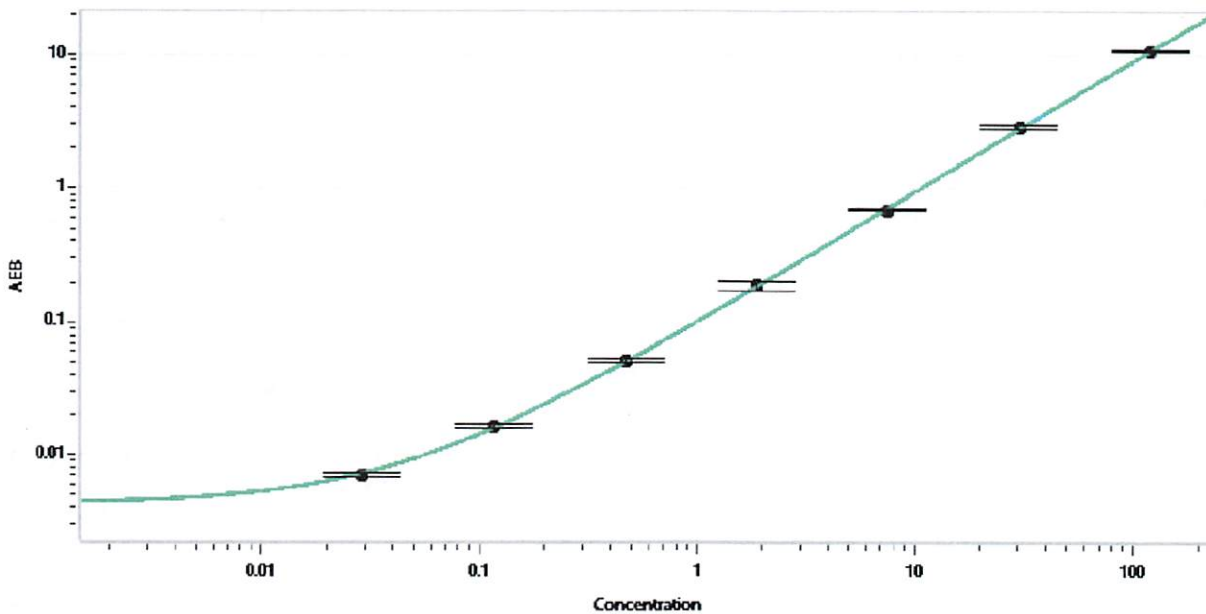
Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number
Bead Stock	102028	211311
Detector Stock	102029	211401
Calibrator Concentrate	102032	119607
Calibrator Diluent	102030	210410
Sample Diluent	103402	211312

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 (ng/mL)	2
---	---

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.029
C	0.117
D	0.469
E	1.88
F	7.50
G	30.0
H	120

Review/Approval: Sonia Khillan Name, AC Manager Title, Sonia Khillan 08 JUL 2022 Signature/Date