

Product Number:	102039
Lot Number:	503934
Expiration:	14-Jun-2024
Platform(s):	HD-1/HD-X & SR-X



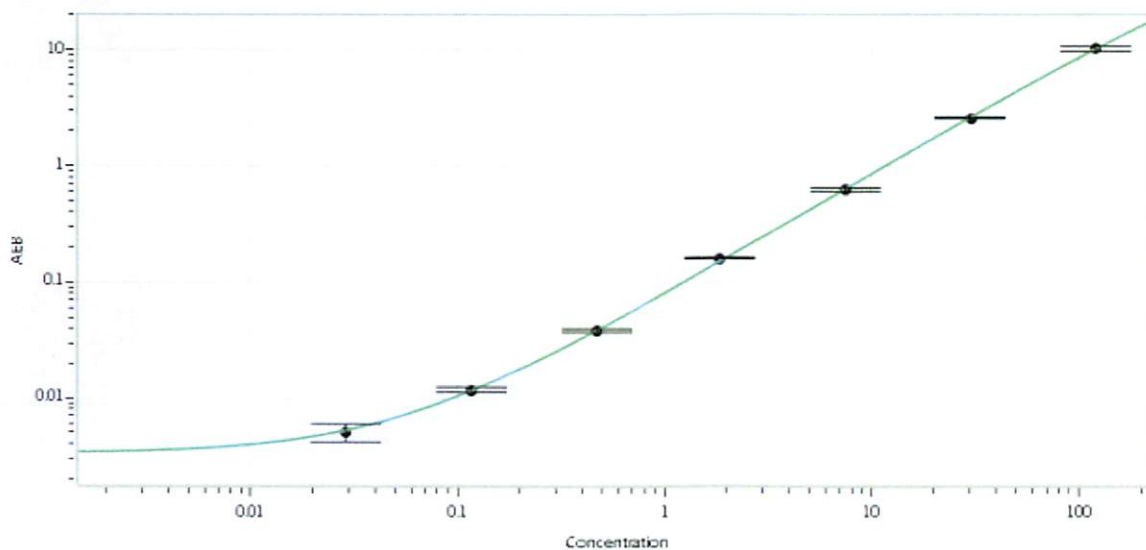
Data below represents results generated on the Simoa® HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102028	323311
Detector Stock	102029	323307
Calibrator Concentrate	102032	323701
Calibrator Diluent	102030	322622
Sample Diluent	103402	323414

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, QC Manager Title, [Signature] 26 SEP 2023 Signature/Date