

Product Number:	102570
Lot Number:	503558
Expiration:	23-May-2023
Platform(s):	HD-1/HD-X



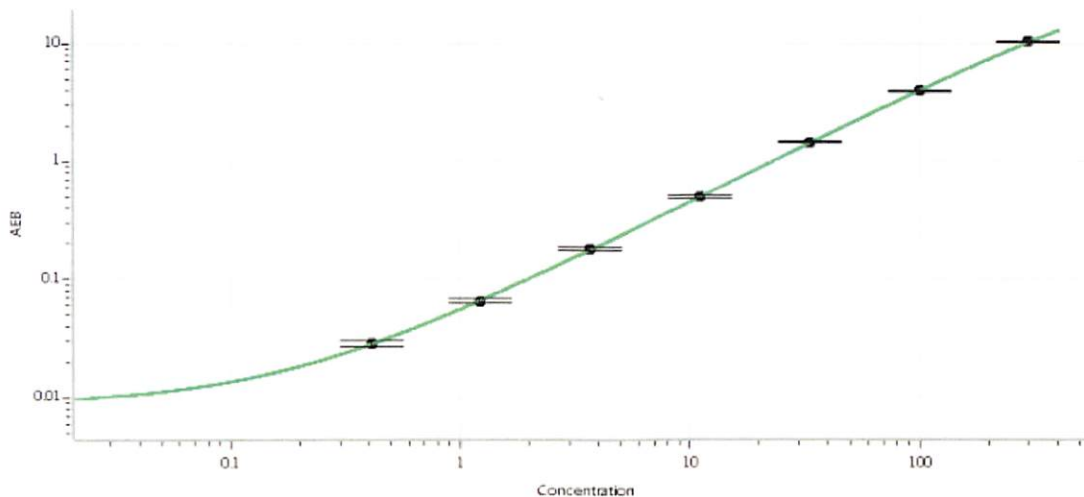
Data below represents results generated on the Simoa® HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102563	221008
Detector Stock	102564	221309
Detector Diluent	102590	221310
Calibrator Concentrate	102567	133306
Calibrator Diluent	102565	220719
Sample Diluent	102566	220718

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)	6.00 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 QC manager Sonia Khillan 1/35EP2022
 Name Title Signature/Date