

Product Number:	102336
Lot Number:	502587
Expiration:	14-Jul-2021
Platform(s):	HD-1/HD-X



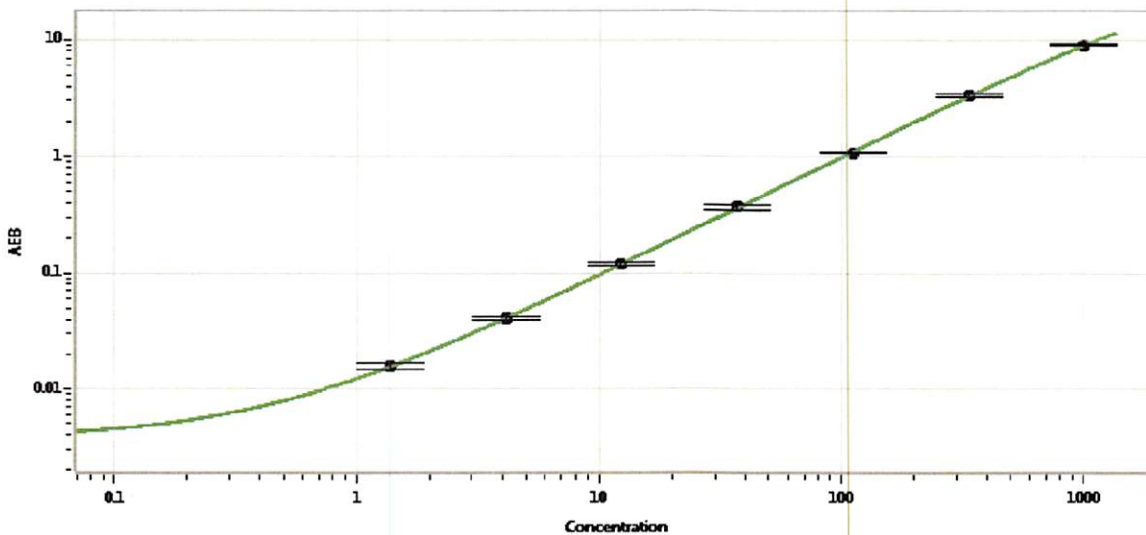
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102330	029008
Detector Stock	102331	029306
Calibrator Concentrate	102335	006004
Calibrator Diluent	102332	024414
Sample Diluent	102333	019704

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	20 ng/mL
-------------------------------------	----------

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

Shivani Goel
Name

QC Manager
Title

Shivani Goel 02 Nov 2020
Signature/Date