

Product Number:	102336
Lot Number:	502477
Expiration:	18-Feb-2021
Platform(s):	HD-1/HD-X



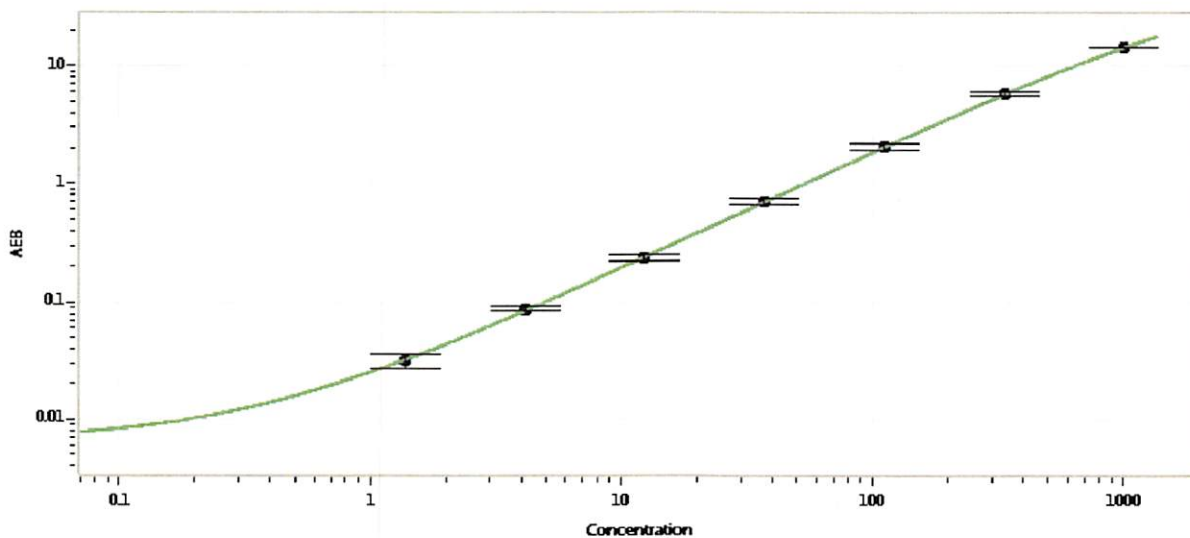
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102330	006003
Detector Stock	102331	006001
Calibrator Concentrate	102335	006004
Calibrator Diluent	102332	004912
Sample Diluent	102333	006202

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 28 Jul 2020
Signature/Date