

Product Number:	101419
Lot Number:	502416
Expiration:	18-Feb-2021
Platform(s):	HD-1/HD-X



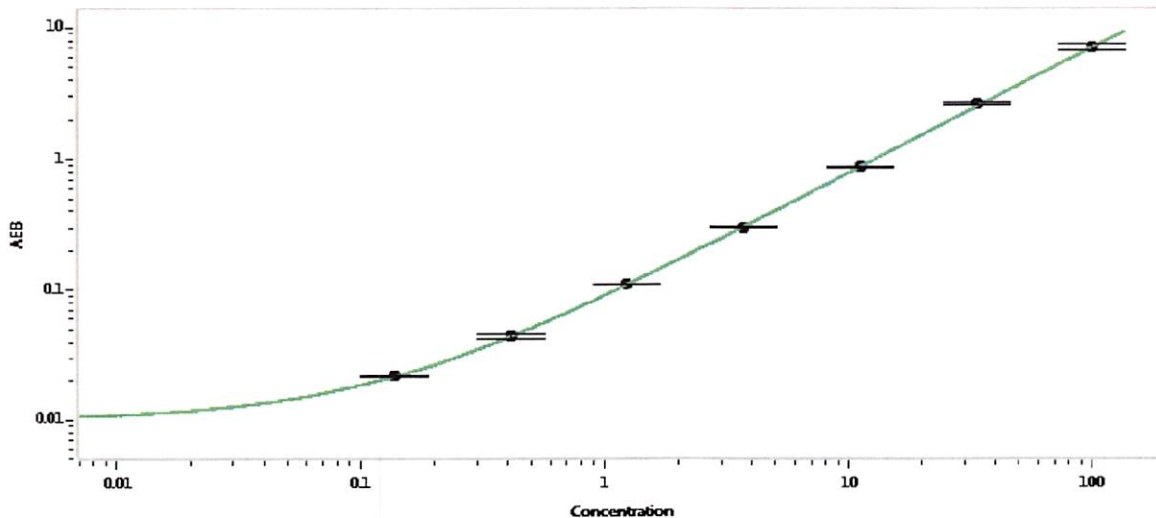
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101899	016009
Detector Stock	101900	015507
Calibrator Concentrate	101420	016014
Calibrator Diluent	101418	015104
Sample Diluent	101416	004919

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	2 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.24
E	3.70
F	11.1
G	33.3
H	100

Review/Approval: Shivani Goel Quality Control Manager Shivani Goel 07/11/2020
 Name Title Signature/Date