

Product Number:	101419
Lot Number:	502969
Expiration:	4-Mar-2022
Platform(s):	HD-1/HD-X



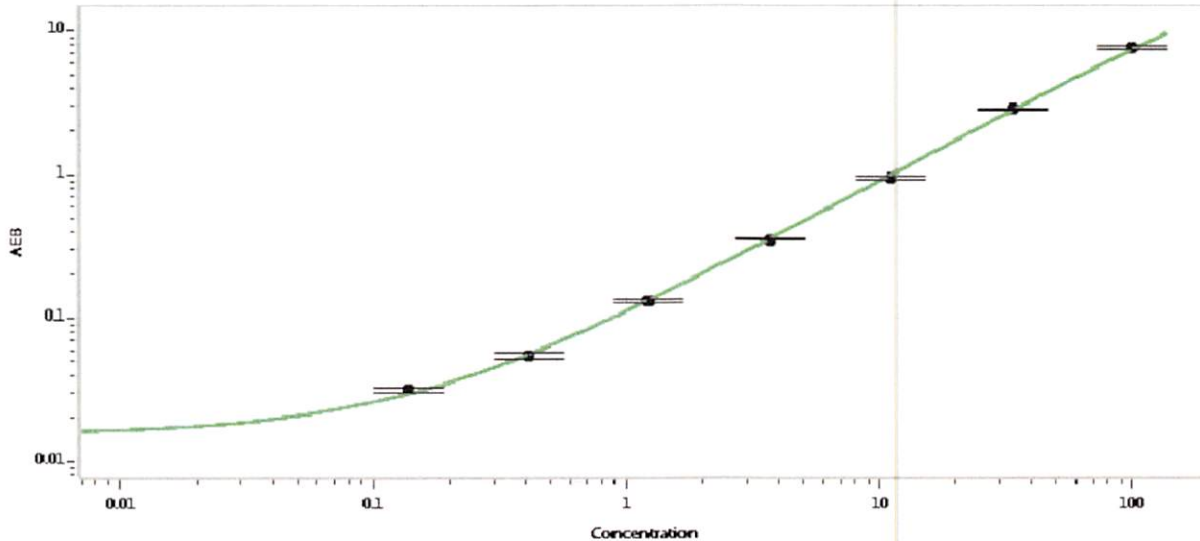
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101899	115002
Detector Stock	101900	115003
Calibrator Concentrate	101420	016014
Calibrator Diluent	101418	111727
Sample Diluent	101416	115219

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

Junpil Park
Name

QC Supervisor
Title

[Signature] 09 JUN 2021
Signature/Date