

Product Number:	102336
Lot Number:	502943
Expiration:	17-Mar-2022
Platform(s):	HD-1 / HD-X



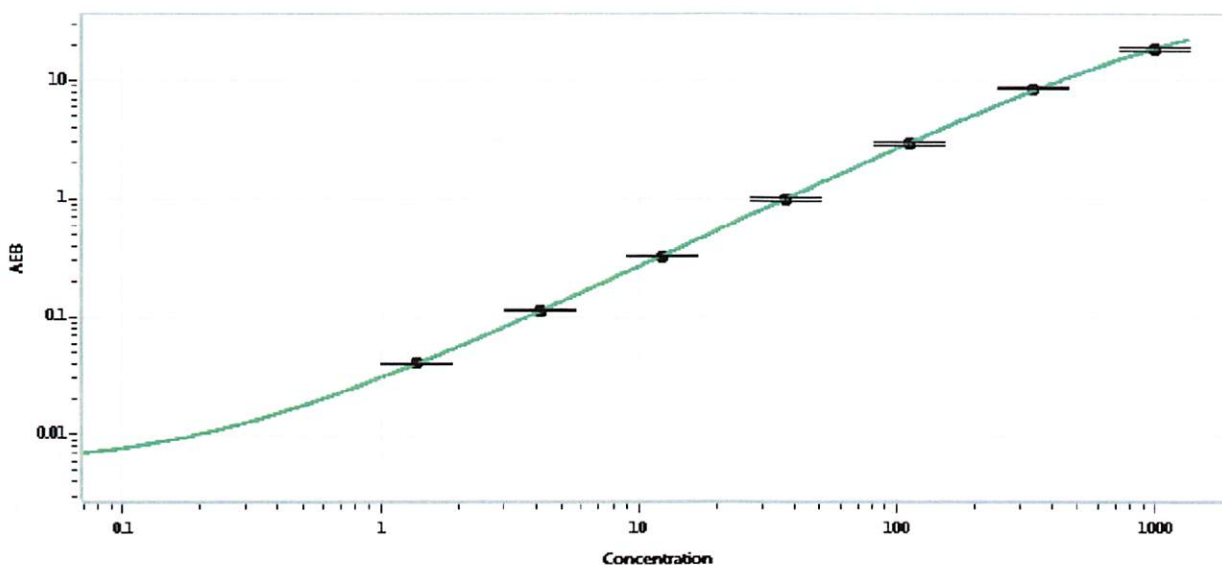
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102330	114002
Detector Stock	102331	114014
Calibrator Concentrate	102335	114016
Calibrator Diluent	102332	111410
Sample Diluent	102333	114015

Generic Components - Included With Kit But Not Lot Number Specified	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Concentrate	101957
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

Junpil Park
Name

QC Supervisor
Title

02 Jun 2021
Signature/Date