

Product Number:	102700
Lot Number:	503069
Expiration:	1-Jun-2022
Platform(s):	HD-1/HD-X

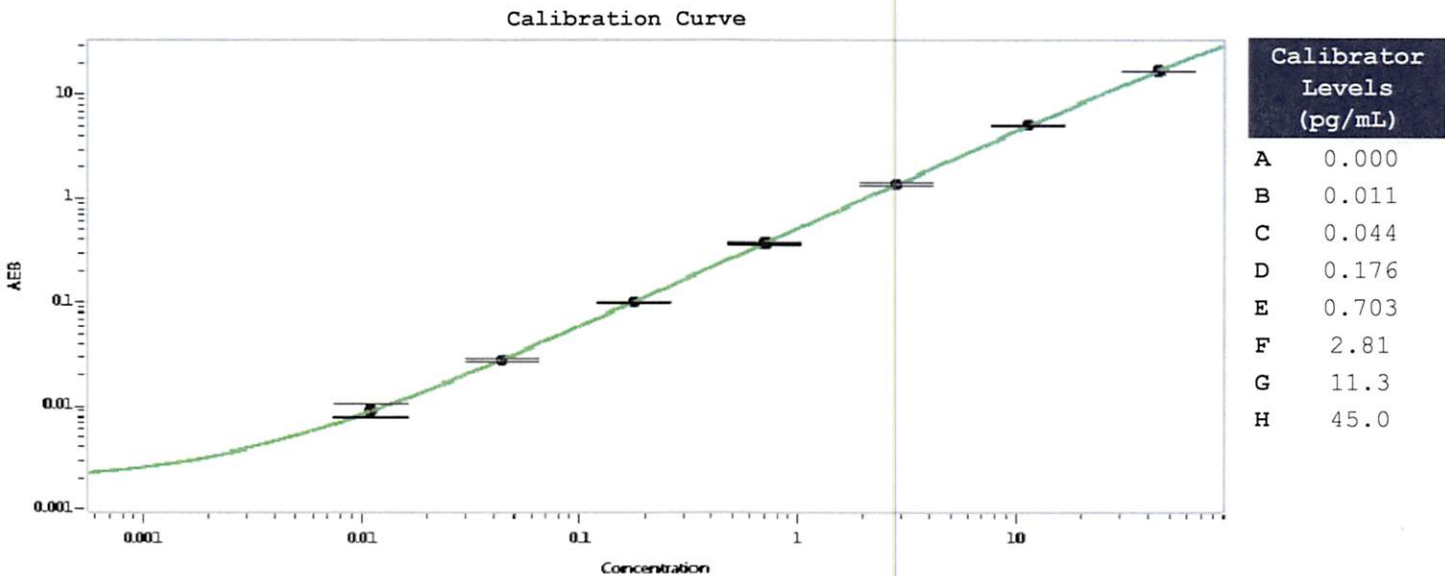


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102693	120501
Detector Stock	102694	120603
Calibrator Concentrate	102697	016302
Calibrator Diluent	102695	115226
Sample Diluent	102696	115225

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	1,000 pg/mL
-------------------------------------	-------------



Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 03 Aug 2021
Signature/Date