

Product Number:	102700
Lot Number:	502991
Expiration:	10-Apr-2022
Platform(s):	HD-1/HD-X



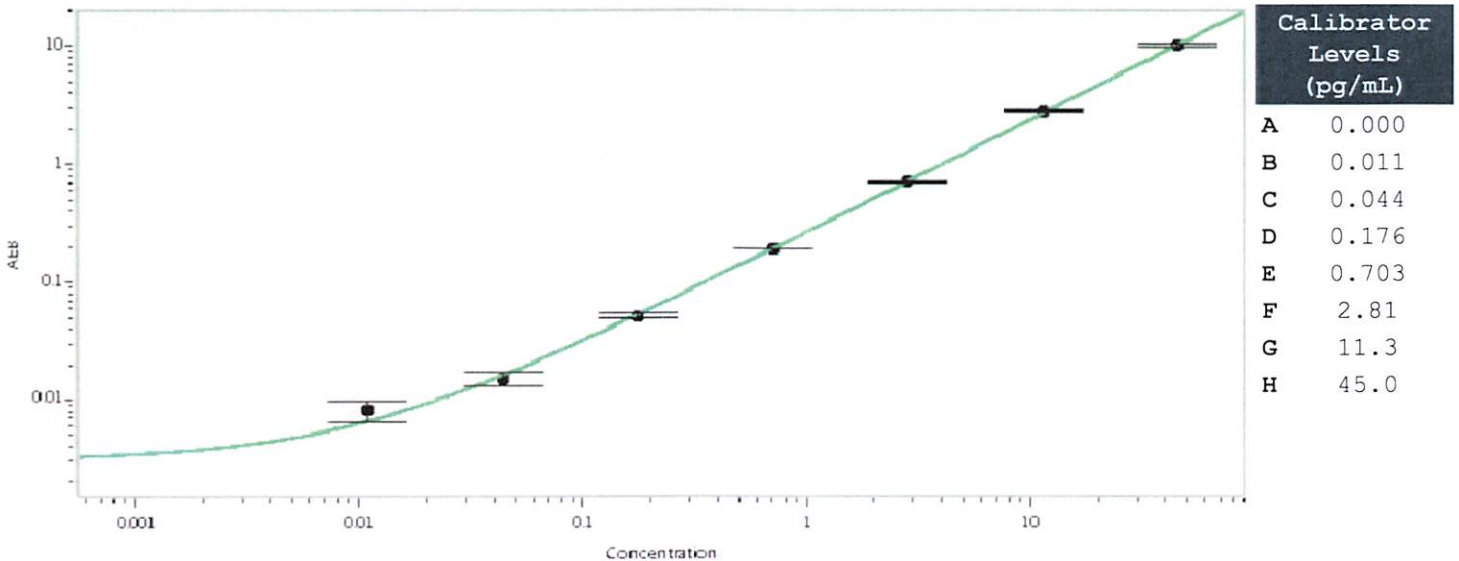
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102693	116504
Detector Stock	102694	116512
Calibrator Concentrate	102697	016302
Calibrator Diluent	102695	111914
Sample Diluent	102696	111915

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
-------------------------------------	-------------

Calibration Curve



Review/Approval

Junpil Park
Name

QC Supervisor
Title

Signature/Date 28 Jun 2021