

Product Number:	102583
Lot Number:	502446
Expiration:	25-Nov-2020
Platform(s):	HD-1/HD-X



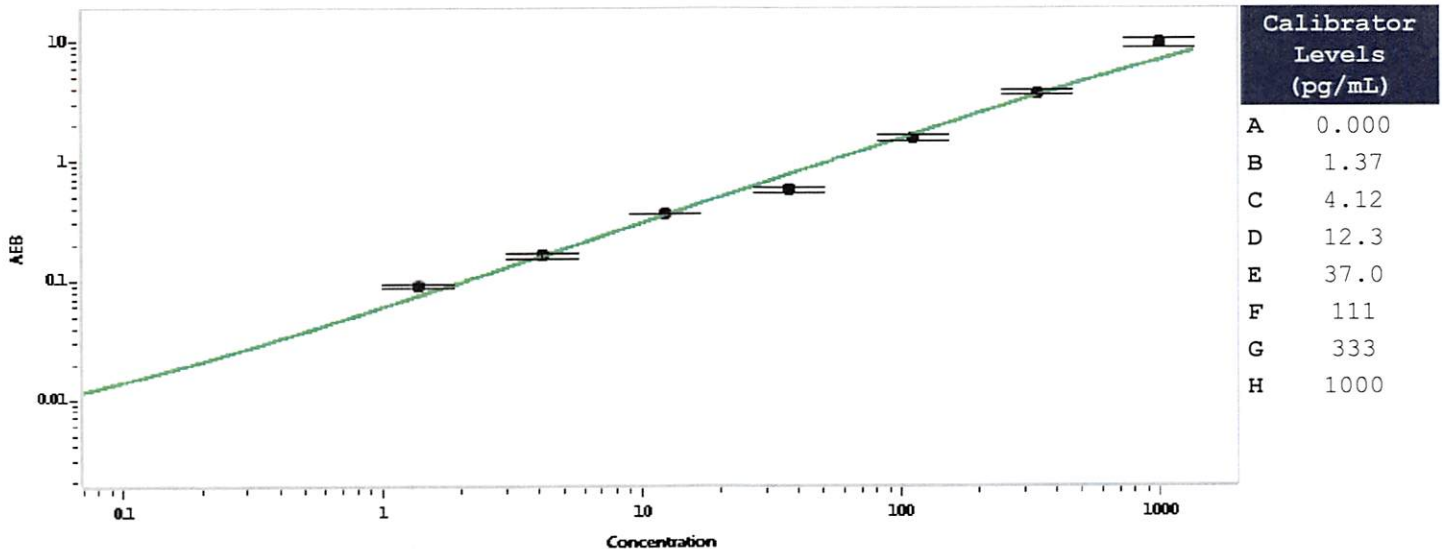
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102576	002209
Detector Stock	102577	002205
Calibrator Concentrate	102580	912804
Calibrator Diluent	102578	932925
Sample Diluent	102579	933813

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	10 ng/mL
-------------------------------------	----------

Calibration Curve



Review/Approval

Shivani Goel

QC Manager

Shivani Goel 01/17/2020

Name

Title

Signature/Date