

Product Number:	102233
Lot Number:	501913
Expiration:	10-Jan-2020
Platform(s):	HD-1/HD-X



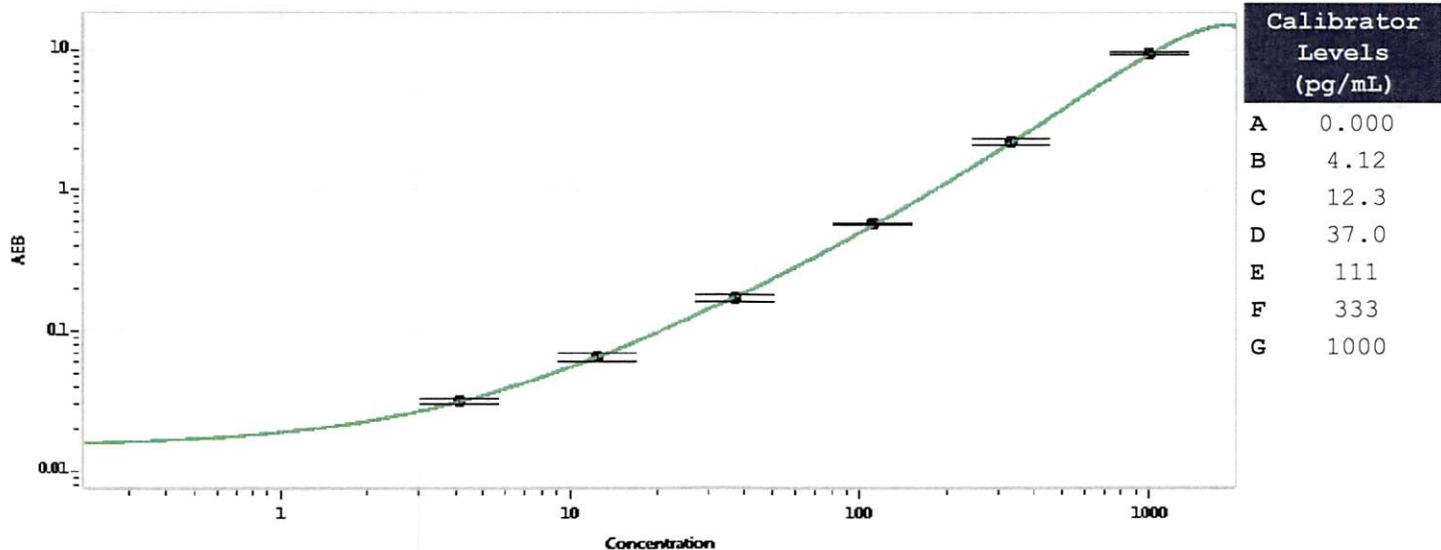
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102223	919302
Detector Stock	102224	919204
Calibrator Concentrate	102229	835002
Calibrator Diluent	102225	919106
Sample Diluent	102226	919105

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	20 ng/mL
-------------------------------------	----------

Calibration Curve



Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 17 Jul 2019
Signature/Date