

Product Number:	102635
Lot Number:	502615
Expiration:	29-Jul-2021
Platform(s):	HD-1/ HD-X

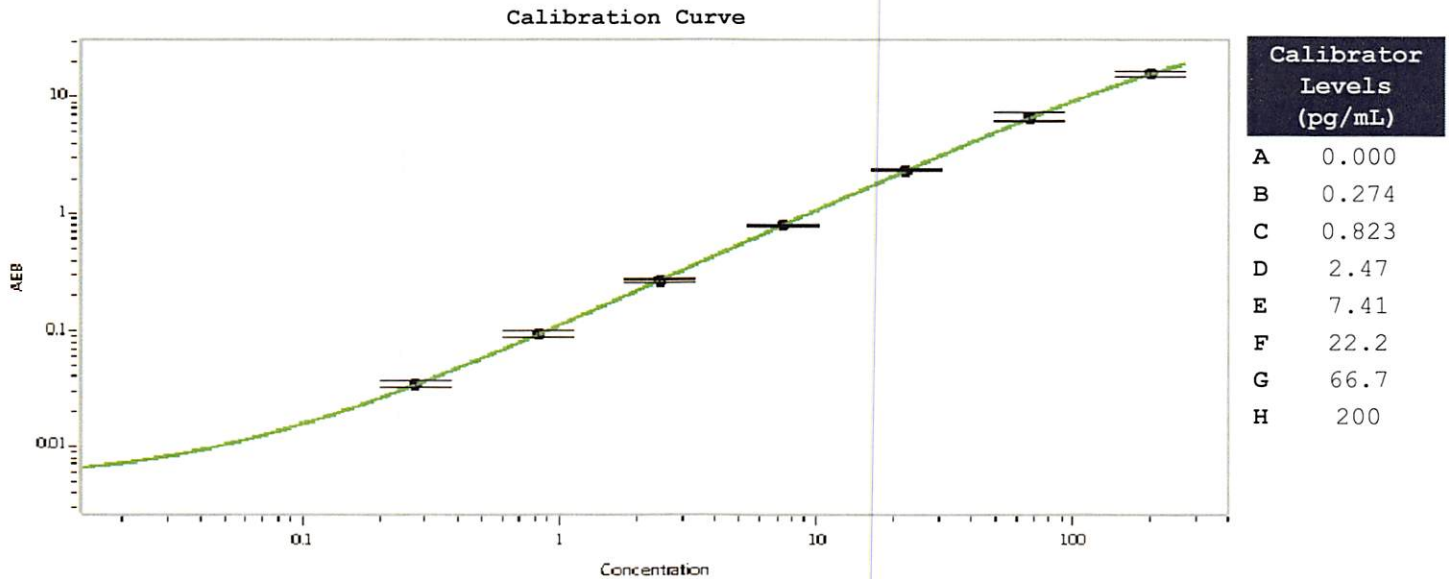


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102628	031701
Detector Stock	102629	021104
Calibrator Concentrate	102632	013308
Calibrator Diluent	102630	031413
Sample Diluent	102631	032507

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	4 ng/mL
-------------------------------------	---------



Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 25Nov2020
Signature/Date