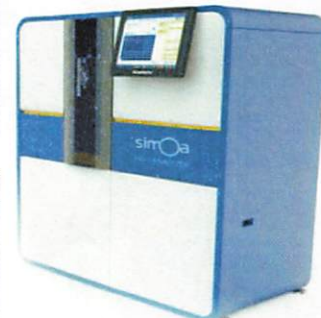


Product Number:	102064
Lot Number:	503280
Expiration:	3-Sep-2022
Platform(s):	HD-1/HD-X

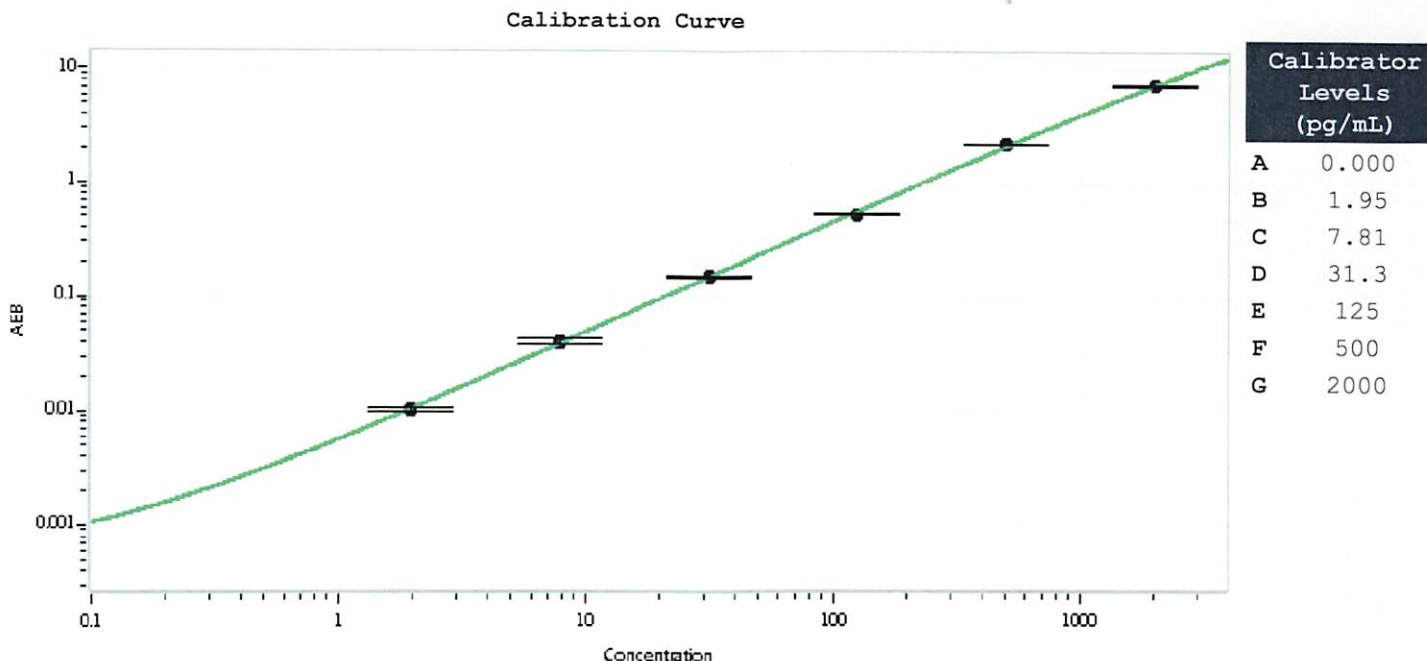
Component	Part Number	Lot Number
Bead Stock	102056	135306
Detector Stock	102057	136210
Calibrator Concentrate	102060	110408
Calibrator Diluent	102058	135419
Sample Diluent	102059	135418



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

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	40 ng/mL
-------------------------------------	----------



Review/Approval: Wenku Tsui Name: Wenku Tsui Title: QC Signature/Date: [Signature] 14 Jun 2022