

Product Number:	102669
Lot Number:	502562
Expiration:	9-Jun-2021
Platform(s):	HD-1/HD-X

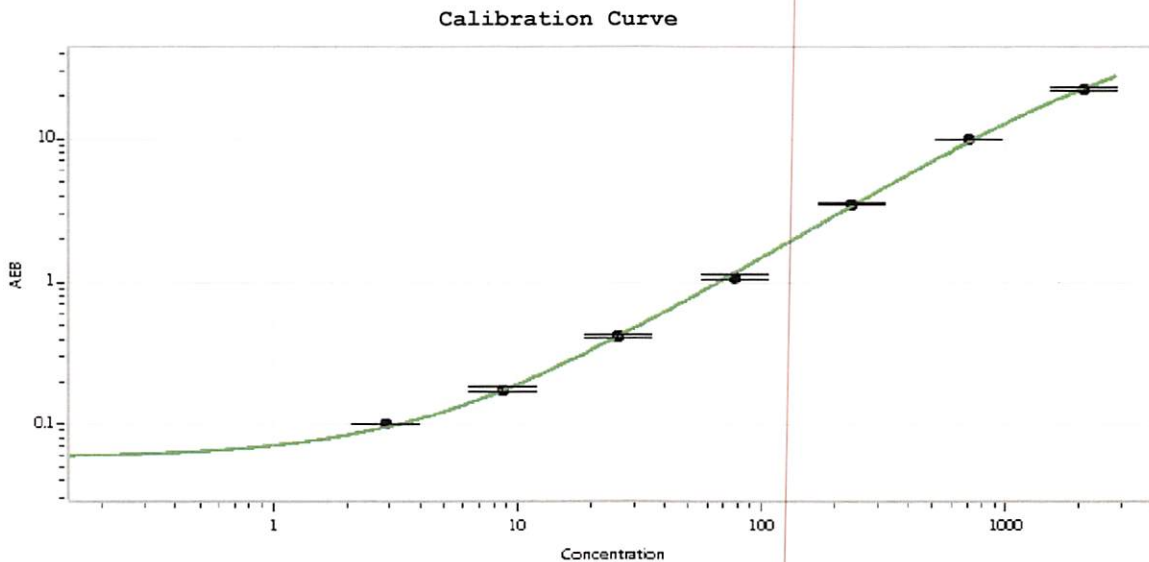


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102662	016705
Detector Stock	102663	017702
Calibrator Concentrate	102666	934501
Calibrator Diluent	102664	016110
Sample Diluent	102665	016701

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	60 ng/mL
-------------------------------------	----------



Review/Approval	Shivani Goel	QC Manager	Shivani Goel 06Oct2020
	Patrizia Stadler	QC Assoc. Director	
	Name	Title	Signature/Date