

Product Number:	101808
Lot Number:	502875
Expiration:	26-Jan-2022
Platform(s):	HD-1/HD-X

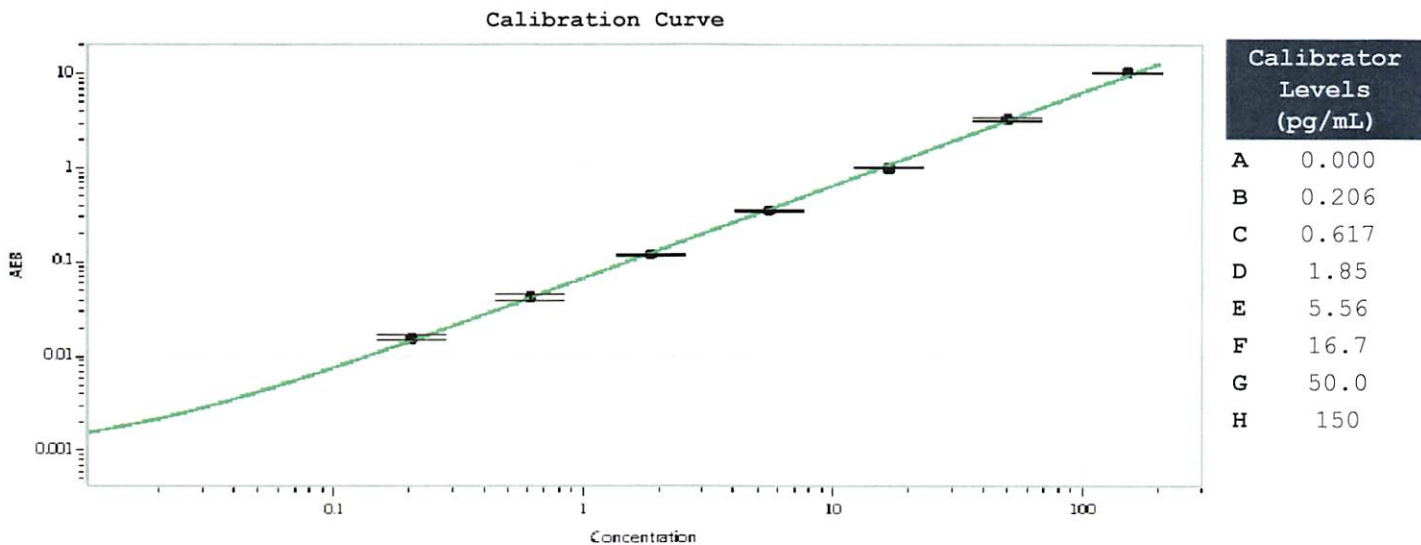


Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	101922	102612
Detector Stock	101923	103002
Calibrator Concentrate	101803	110512
Calibrator Diluent	101925	103805
Sample Diluent	101926	103001

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	3 ng/mL
-------------------------------------	---------



Review/Approval: Junpil Park Name, QC Supervisor Title, [Signature] 21 Apr 2021 Signature/Date