

Product Number:	102504
Lot Number:	501830
Expiration:	10-May-2020
Platform(s):	HD-1



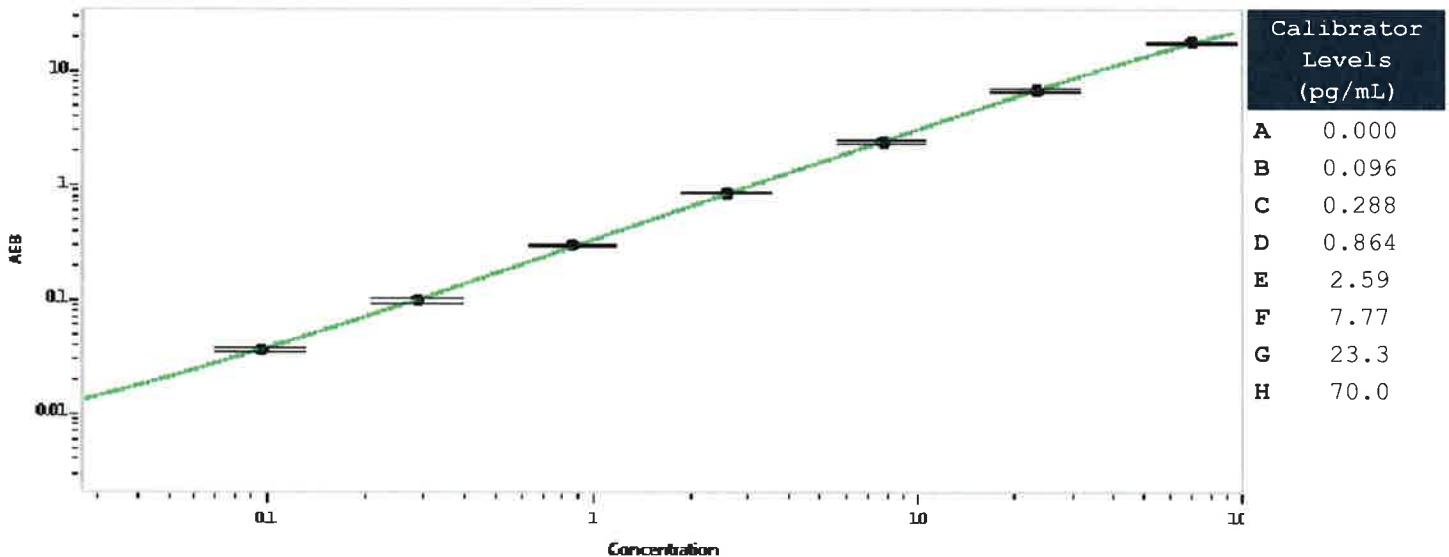
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102497	913008
Detector Stock	102498	913304
Calibrator Concentrate	102501	913702
Calibrator Diluent	102499	913305
Sample Diluent	102500	913306

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	1.4 ng/mL
-------------------------------------	-----------

Calibration Curve



Review/Approval

Patrizia Stadler

QC Assoc. Director

P. Stadler 11 Jun 2019

Name

Title

Signature/Date