

Product Number:	102504
Lot Number:	501830
Expiration:	30-Mar-2020
Platform(s):	HD-1



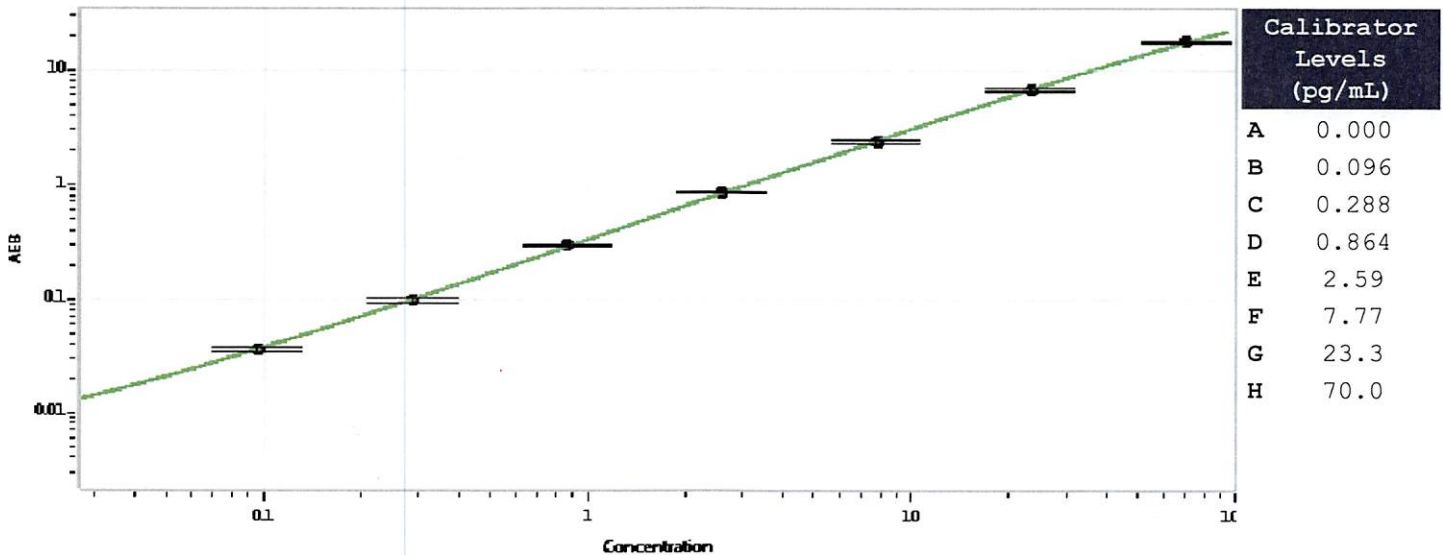
Data below represents results generated on the

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
Name

QC Assoc. Director
Title

P. Stadler 12 Jun 2019
Signature/Date