

Product Number:	101847
Lot Number:	501761
Expiration:	11-Feb-2020
Platform(s):	HD-1



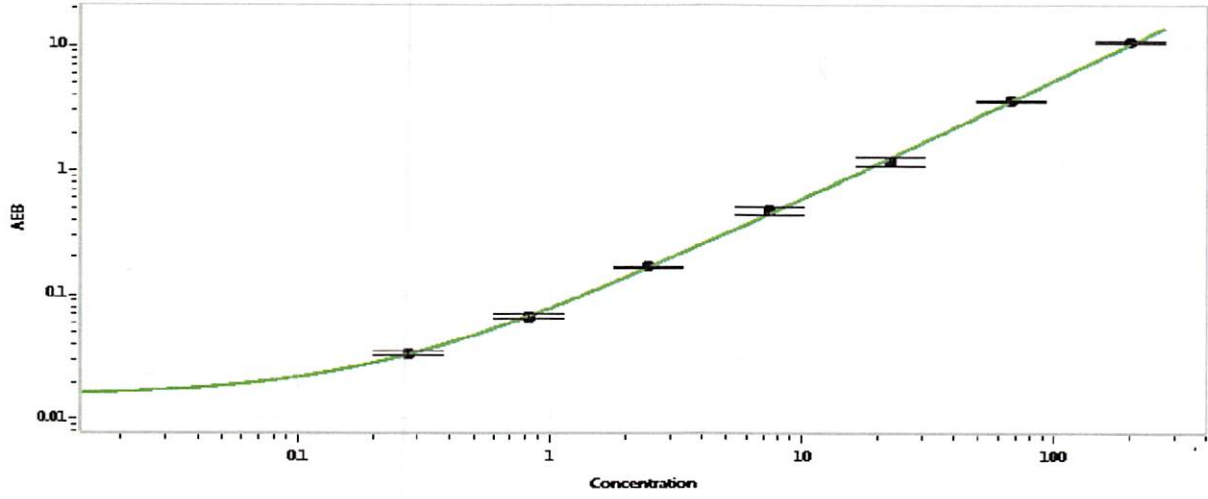
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	101935	908802
Detector Stock	101936	908806
Calibrator Concentrate	101842	909303
Calibrator Diluent	101951	908205
Sample Diluent	101952	907311

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	4000 pg/mL
-------------------------------------	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 05 April 2019
Signature/Date