

Product Number:	101839
Lot Number:	501874
Expiration:	30-Mar-2020
Platform(s):	HD-1/HD-X



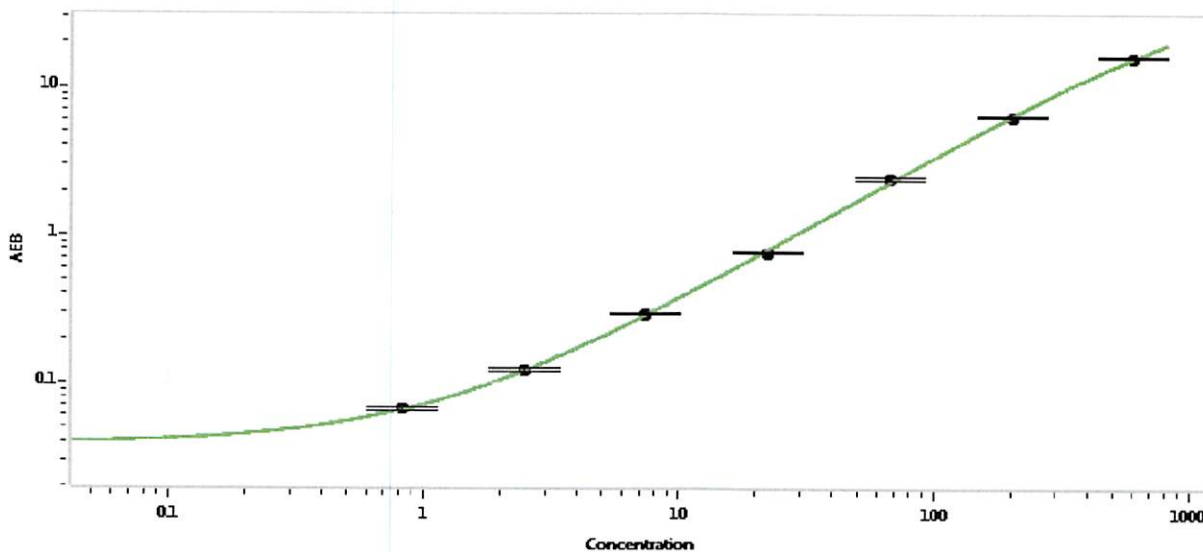
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	101931	916407
Detector Stock	101932	916505
Calibrator Concentrate	101835	916807
Calibrator Diluent	101947	916310
Sample Diluent	101948	908912

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	12,000 pg/mL
-------------------------------------	--------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.823
C	2.47
D	7.41
E	22.2
F	66.7
G	200
H	600

Review/Approval

Patrizia Stadler  
Name

Assoc. Director, QC  
Title

*P. Stadler* 25 JUN 2019  
Signature/Date