

Product Number:	101968
Lot Number:	501816
Expiration:	8-Nov-2019
Platform(s):	HD-1

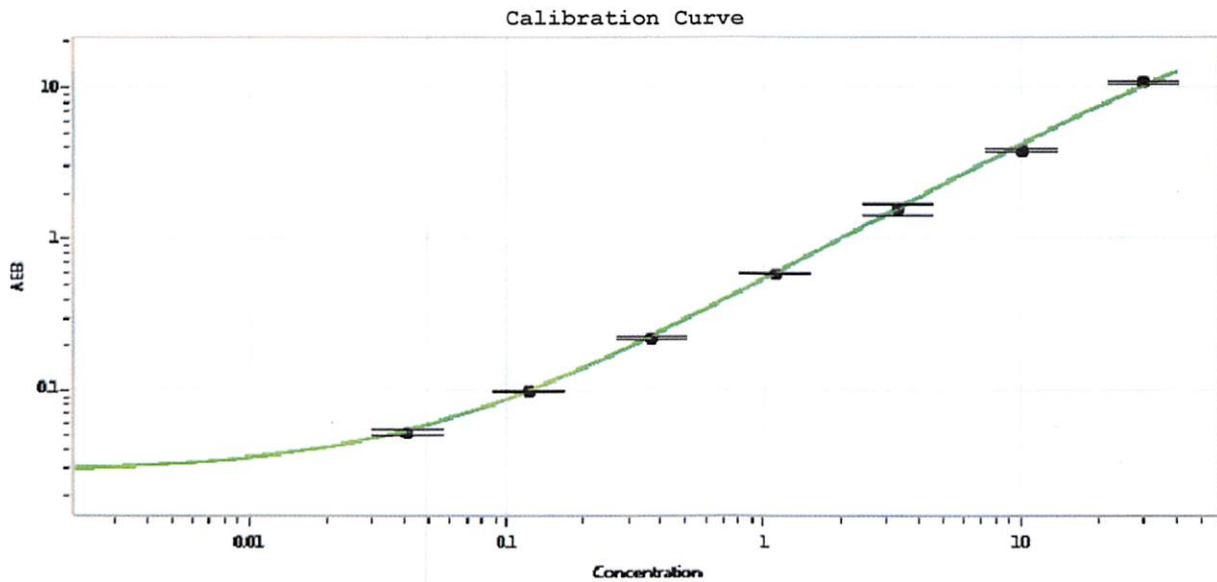


Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	101966	912211
Detector Stock	101967	912210
Calibrator Concentrate	101965	MF-110817-03
Calibrator Diluent	101963	912013
Sample Diluent	101964	835511

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	600 pg/mL
-------------------------------------	-----------



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 13 May 2019
Signature/Date