

Product Number:	102700
Lot Number:	501771
Expiration:	11-Feb-2020
Platform(s):	HD-1



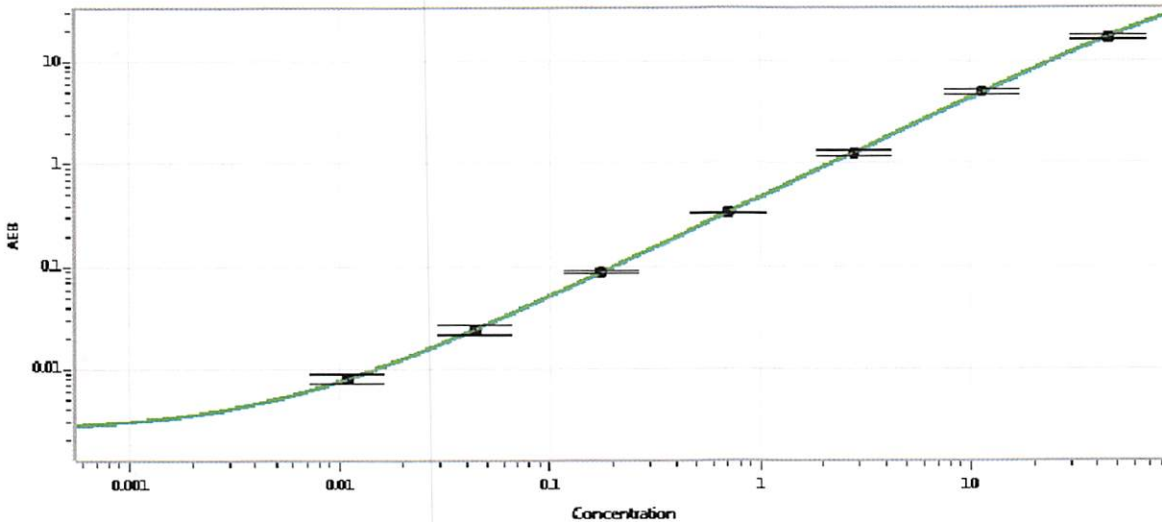
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102693	908606
Detector Stock	102694	910109
Calibrator Concentrate	102697	MF-060418-01
Calibrator Diluent	102695	907308
Sample Diluent	102696	907307

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	1000 pg/mL
-------------------------------------	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.011
C	0.044
D	0.176
E	0.70
F	2.81
G	11.3
H	45.0

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 16 Apr 2019
Signature/Date