

Product Number:	101984
Lot Number:	501634
Expiration:	14-Dec-2019
Platform(s):	HD-1



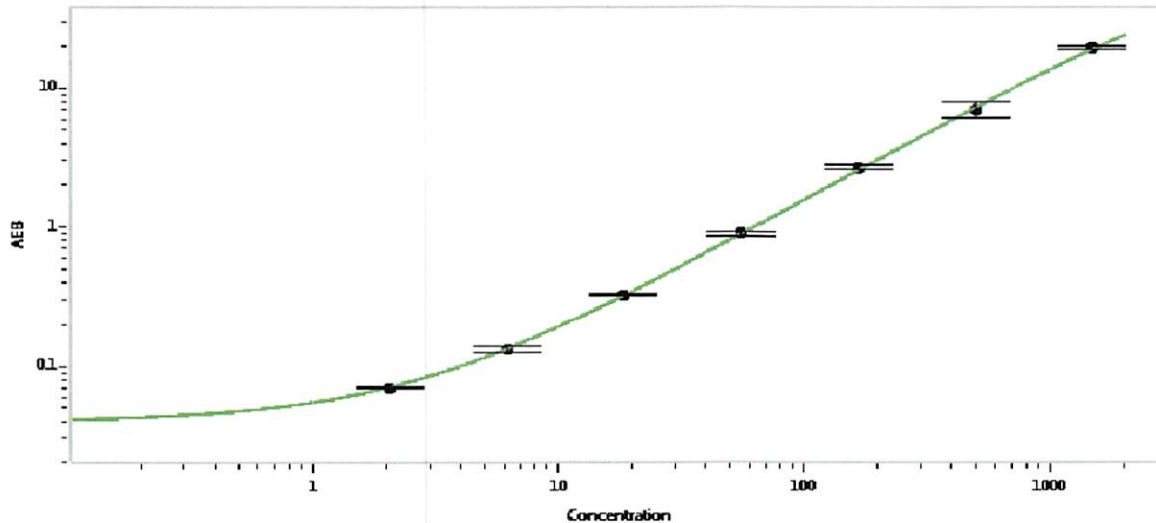
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	101910	902504
Detector Stock	101911	903303
Calibrator Concentrate	101983	MF-121417-01
Calibrator Diluent	101908	902603
Sample Diluent	101909	902602

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	30 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.06
C	6.17
D	18.5
E	55.6
F	167
G	500
H	1500

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 14Feb2019
Signature/Date