

Product Number:	101824
Lot Number:	502236
Expiration:	7-Oct-2020
Platform(s):	HD-1/ HD-X



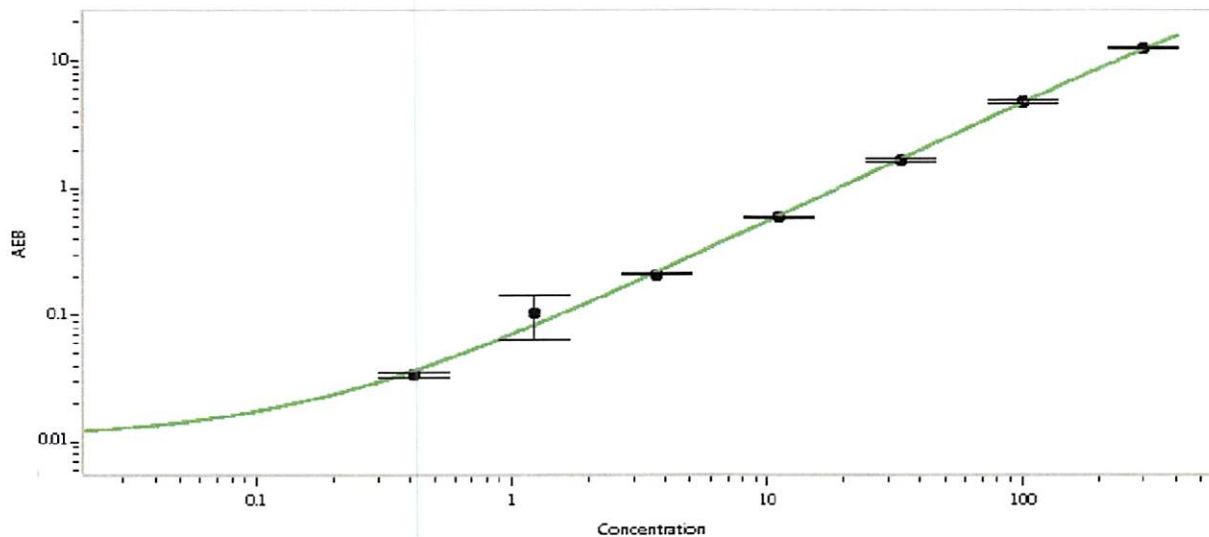
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	101929	004106
Detector Stock	101930	004107
Calibrator Concentrate	101819	907806
Calibrator Diluent	101943	003511
Sample Diluent	101944	928027

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	6000 pg/mL
-------------------------------------	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.412
C	1.23
D	3.70
E	11.1
F	33.3
G	100
H	300

Review/Approval

Patrizia Stadler

Assoc. Director QC

P. Stadler 14 Feb 2020

Name

Title

Signature/Date