

Product Number:	101808
Lot Number:	501970
Expiration:	19-Jun-2020
Platform(s):	HD-1/HD-X



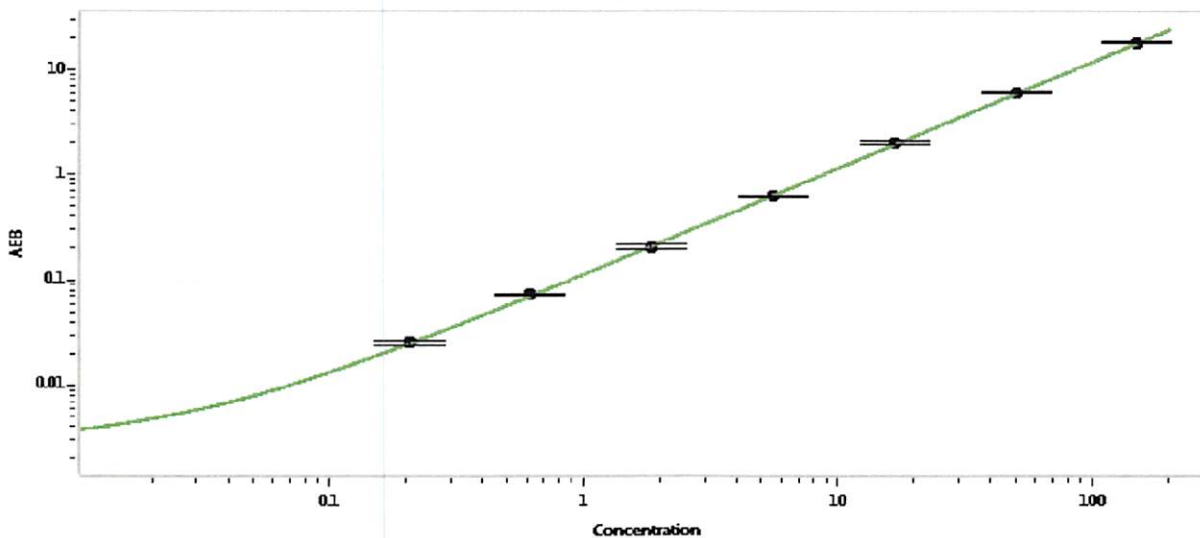
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	101922	921404
Detector Stock	101923	921308
Calibrator Concentrate	101803	831707
Calibrator Diluent	101925	921304
Sample Diluent	101926	921305

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	3 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.206
C	0.617
D	1.85
E	5.56
F	16.7
G	50.0
H	150

Review/Approval	Patrizia Stadler	Associate Director, QC	<i>P. Stadler</i>	14 Aug 2019
	Name	Title	Signature/Date	