

Product Number:	101968
Lot Number:	502173
Expiration:	21-Nov-2020
Platform(s):	HD-1/HD-X



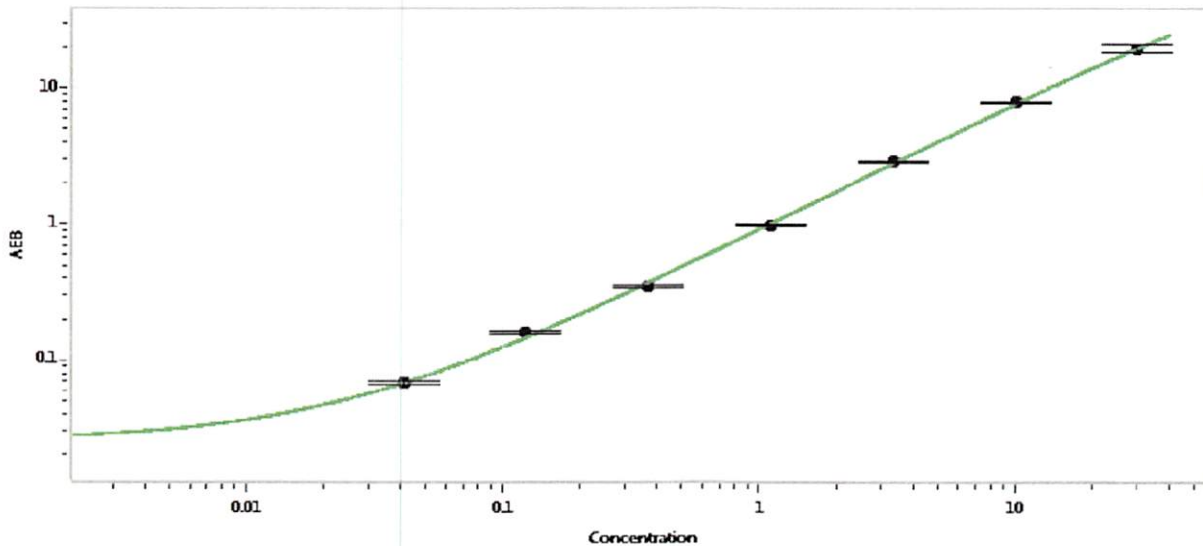
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	101966	002304
Detector Stock	101967	002109
Calibrator Concentrate	101965	000719
Calibrator Diluent	101963	932922
Sample Diluent	101964	932622

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
-------------------------------------	-----------

Calibration Curve



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	QC Asso. Director	<i>P. Stadler</i> 29 Jan 2020
	Name	Title	Signature/Date