

Product Number:	102039
Lot Number:	501774
Expiration:	14-Mar-2020
Platform(s):	HD-1 & SR-X



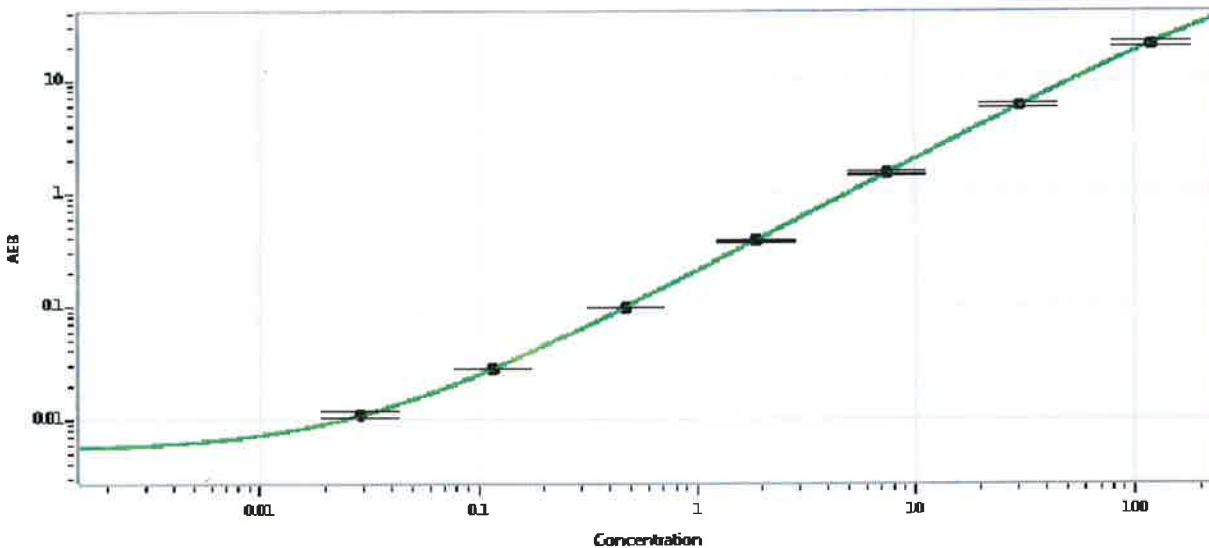
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102028	910809
Detector Stock	102029	910506
Calibrator Concentrate	102032	910804
Calibrator Diluent	102030	910306
Sample Diluent	103402	907314

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	2 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.029
C	0.117
D	0.469
E	1.88
F	7.50
G	30.0
H	120

Review/Approval

Patrizia Stadler

QC Assoc. Director

p. stodler 24 Apr 2019

Name

Title

Signature/Date

