

Product Number:	102648
Lot Number:	501778
Expiration:	11-Feb-2020
Platform(s):	HD-1 & SR-X



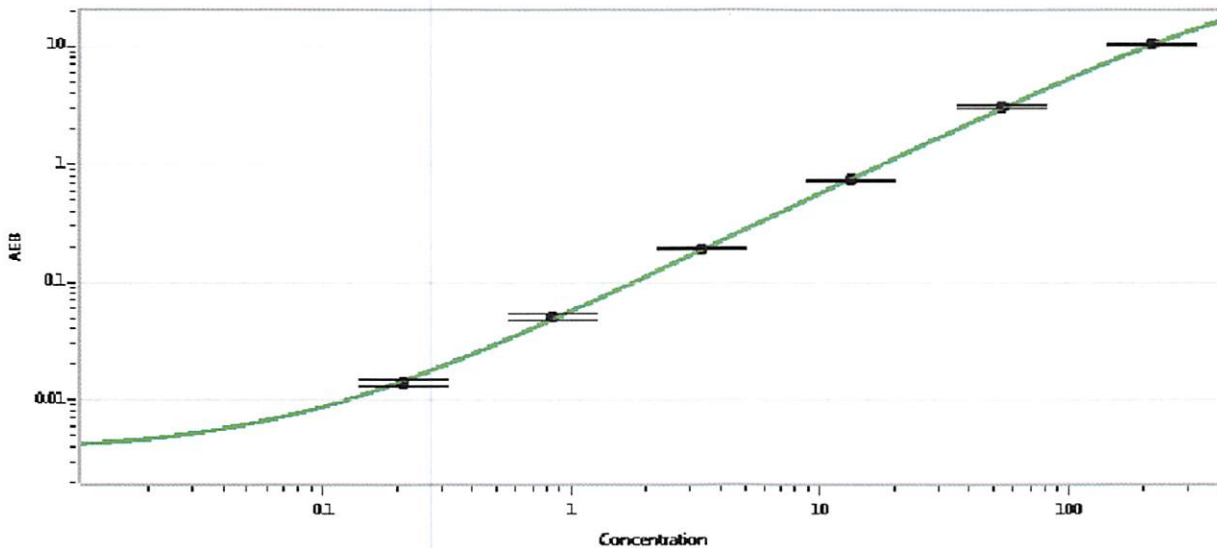
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102641	909410
Detector Stock	102642	910305
Detector Diluent	102734	910212
Calibrator Concentrate	102645	902915
Calibrator Diluent	102643	908819
Sample Diluent	102644	907313

Generic Components - Included With Kit But Not Lot Number Specific		
Discovery Bead Diluent		101954
Discovery SBG Diluent		101956
Discovery SBG Stock		103297
RGP Reagent		103159

Calibrator Concentrate Target Value	3000 pg/mL
-------------------------------------	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.210
C	0.840
D	3.36
E	13.4
F	53.8
G	215

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 22 Apr 2019
Signature/Date