

Product Number:	102039
Lot Number:	501774
Expiration:	11-Feb-2020
Platform(s):	HD-1 & SR-X



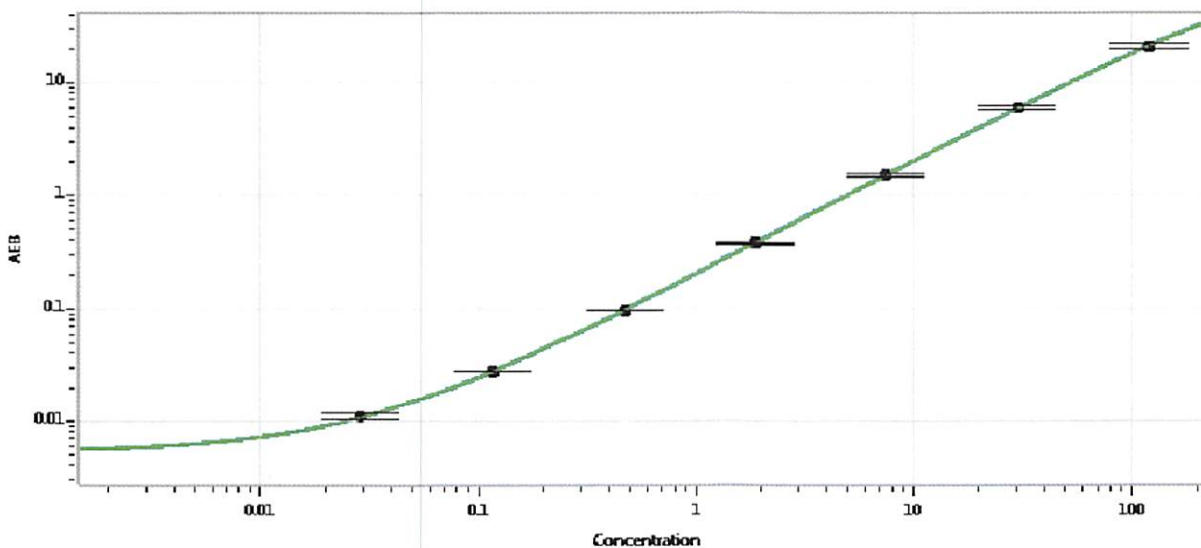
Data below represents results generated on the

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  
Name

QC Assoc. Director  
Title

*P. Stadler* 24 Apr 2019  
Signature/Date