

Product Number:	102517
Lot Number:	501623
Expiration:	3-Jul-2019
Platform(s):	HD-1 & SR-X



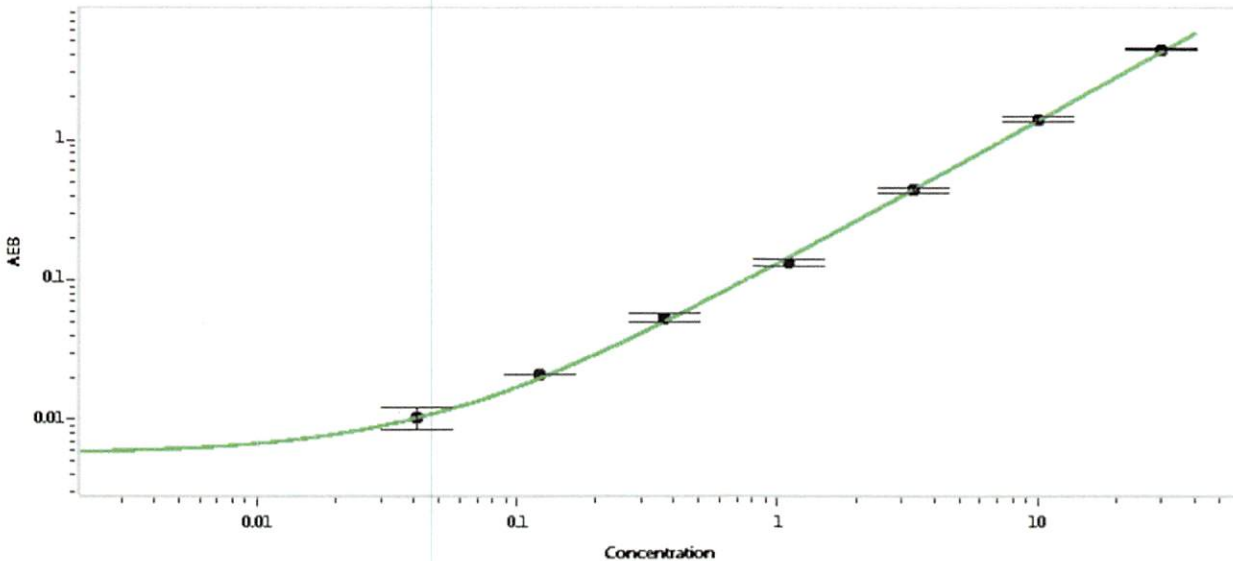
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102510	822608
Detector Stock	102511	822609
Calibrator Concentrate	102514	MF-022818-09
Calibrator Diluent	102512	818420
Sample Diluent	102513	818419

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
Name

QC Assoc. Director
Title

P. Stadler 29 Jan 2019
Signature/Date