

Product Number: 102700
 Lot Number: 501631
 Expiration: 10-Jan-2019
 Platform(s): HD-1



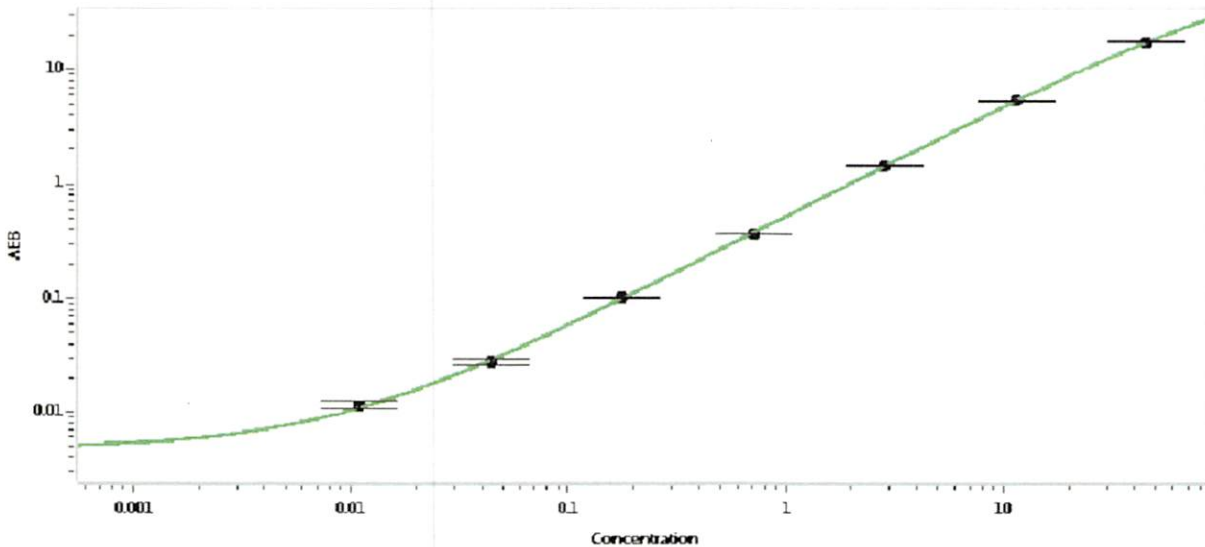
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102693	903205
Detector Stock	102694	903107
Calibrator Concentrate	102697	MF-060418-01
Calibrator Diluent	102695	902922
Sample Diluent	102696	902921

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: 1 ng/mL

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.011
C	0.044
D	0.176
E	0.703
F	2.81
G	11.3
H	45.0

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Scollin 13 Feb 2019
Signature/Date