

Product Number:	102556
Lot Number:	501564
Expiration:	19-Sep-2019
Platform(s):	HD-1



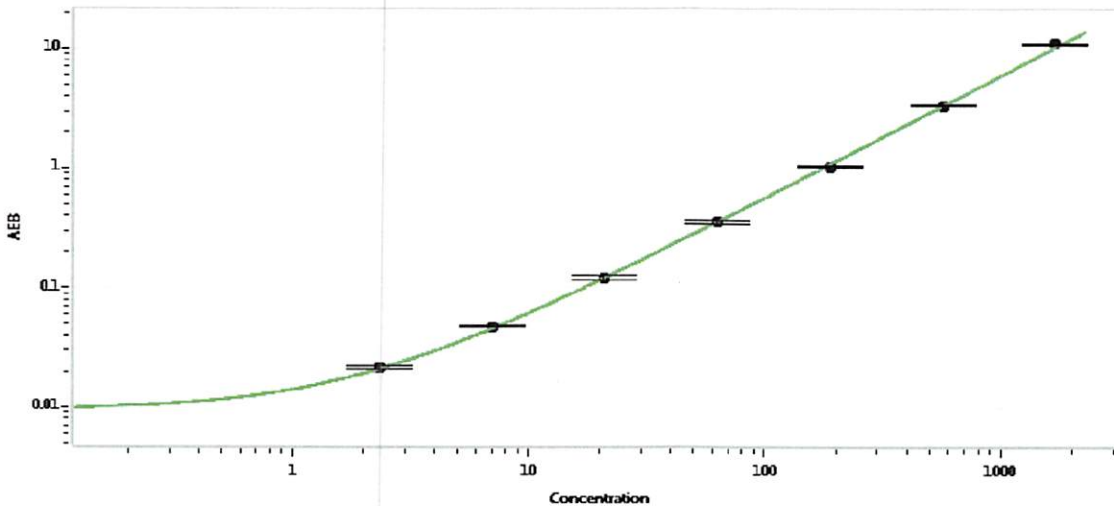
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102549	835107
Detector Stock	102550	834901
Calibrator Concentrate	102553	900707
Calibrator Diluent	102551	834710
Sample Diluent	102552	826224

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	40 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	2.33
C	7.00
D	21.0
E	63.0
F	189
G	567
H	1700

Review/Approval

Patrizia Stadler  
Name

QC Assoc. Director  
Title

*P. Stadler* 30 Jun 2019  
Signature/Date