

Product Number:	102336
Lot Number:	501887
Expiration:	30-Mar-2020
Platform(s):	HD-1 / HD-X



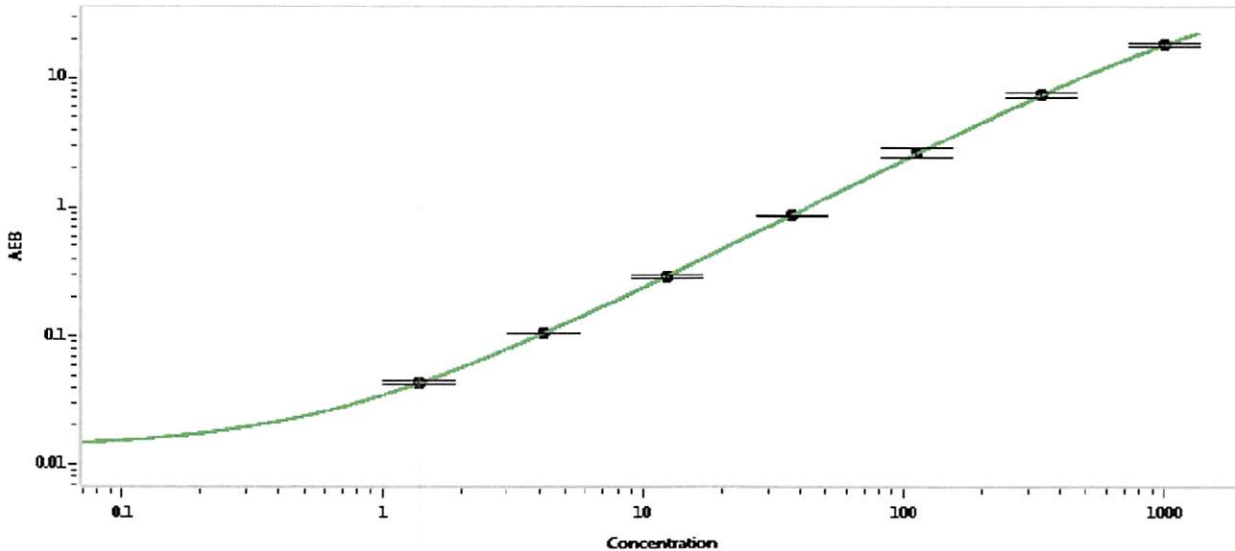
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102330	917901
Detector Stock	102331	917902
Calibrator Concentrate	102335	912103
Calibrator Diluent	102332	908915
Sample Diluent	102333	915702

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	20 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.12
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Patrizia Stadler
Name

QC Assoc. Director
Title

P. Stadler 03 JUL 2019
Signature/Date