

Product Number:	102462
Lot Number:	502341
Expiration:	15-Jan-2021
Platform(s):	HD-1/ HD-X

Component	Part Number	Lot Number
Bead Stock	102455	011106
Detector Stock	102456	011107
Calibrator Concentrate	102459	911304
Calibrator Diluent	102457	011004
Sample Diluent	102458	011005

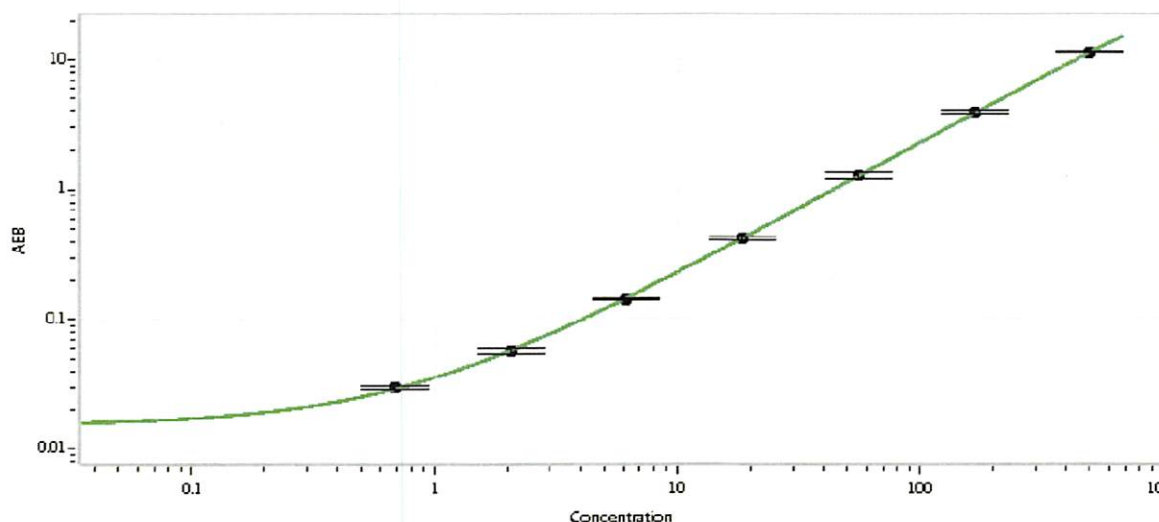


Data below represents results generated on the

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	10 ng/mL
--	----------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.69
C	2.1
D	6.2
E	19
F	56
G	167
H	500

Review/Approval

Patrizia Stadler
Name

QC Assoc. Director
Title

P. Stadler 23 Apr 2020
Signature/Date