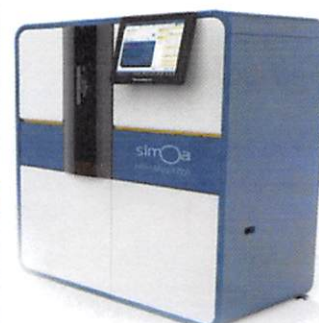


Product Number:	102462
Lot Number:	501795
Expiration:	11-Feb-2020
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	102455	911208
Detector Stock	102456	910904
Calibrator Concentrate	102459	911304
Calibrator Diluent	102457	910903
Sample Diluent	102458	910902

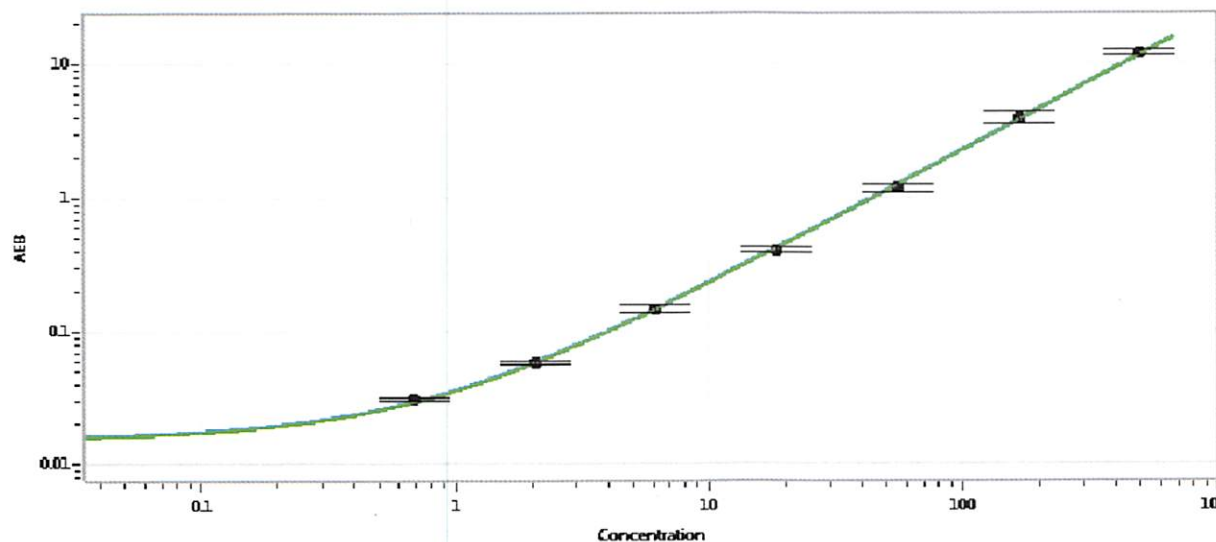


Data below represents results generated on the

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	0.686
C	2.06
D	6.17
E	18.5
F	55.6
G	167
H	500

Review/Approval

Patrizia Stadler
Name

QC Assoc. Director
Title

P. Stadler 29 Apr 2019
Signature/Date