

Product Number:	102136
Lot Number:	501637
Expiration:	10-Jan-2020
Platform(s):	HD-1



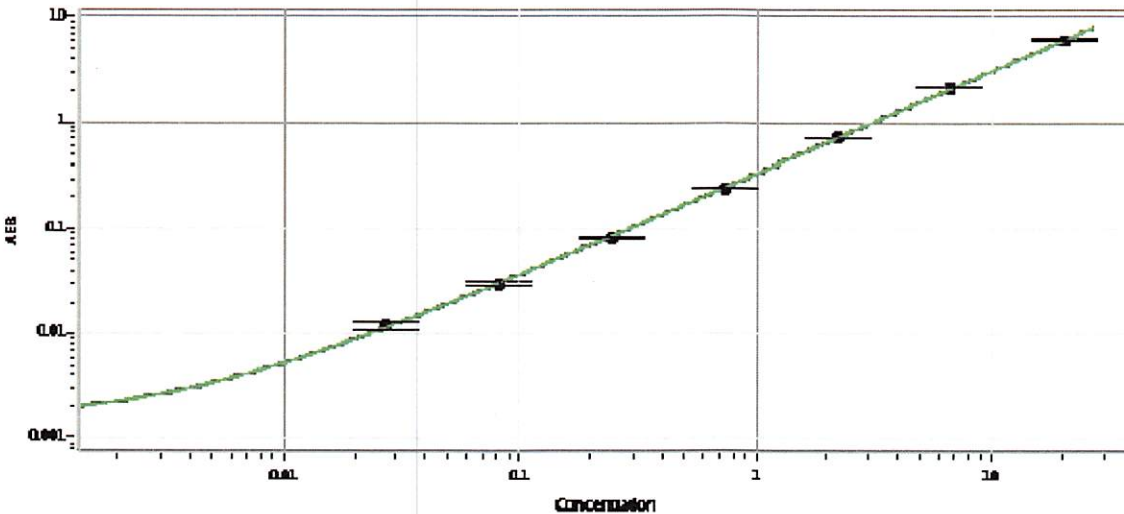
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102128	903504
Detector Stock	102129	903611
Calibrator Concentrate	102132	904303
Calibrator Diluent	102130	901017
Sample Diluent	102131	902924

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	1000 U/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.027
C	0.082
D	0.247
E	0.741
F	2.22
G	6.67
H	20.0

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i> 22 Feb 2019
	Name	Title	Signature/Date