

Product Number:	102543
Lot Number:	501674
Expiration:	10-Jan-2020
Platform(s):	HD-1



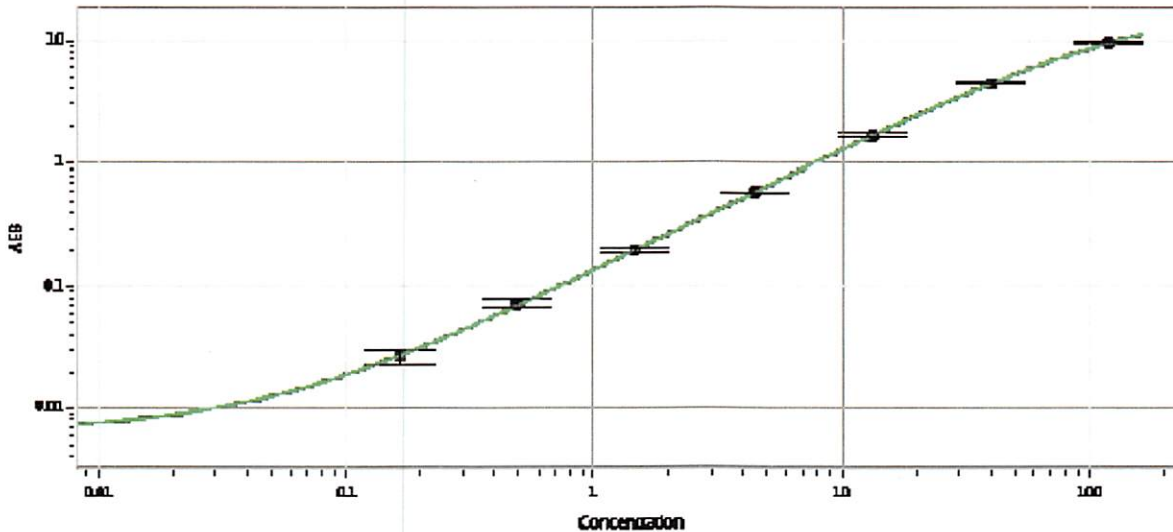
Data below represents results generated on the

Component	Part Number	Lot Number
Bead Stock	102536	905104
Detector Stock	102537	903612
Calibrator Concentrate	102540	904304
Calibrator Diluent	102538	901018
Sample Diluent	102539	902926

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	2400 U/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.165
C	0.494
D	1.48
E	4.44
F	13.3
G	40.0
H	120

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