

Product Number:	101824
Lot Number:	501724
Expiration:	29-Jan-2020
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	101929	907216
Detector Stock	101930	906607
Calibrator Concentrate	101819	907806
Calibrator Diluent	101943	905822
Sample Diluent	101944	902934

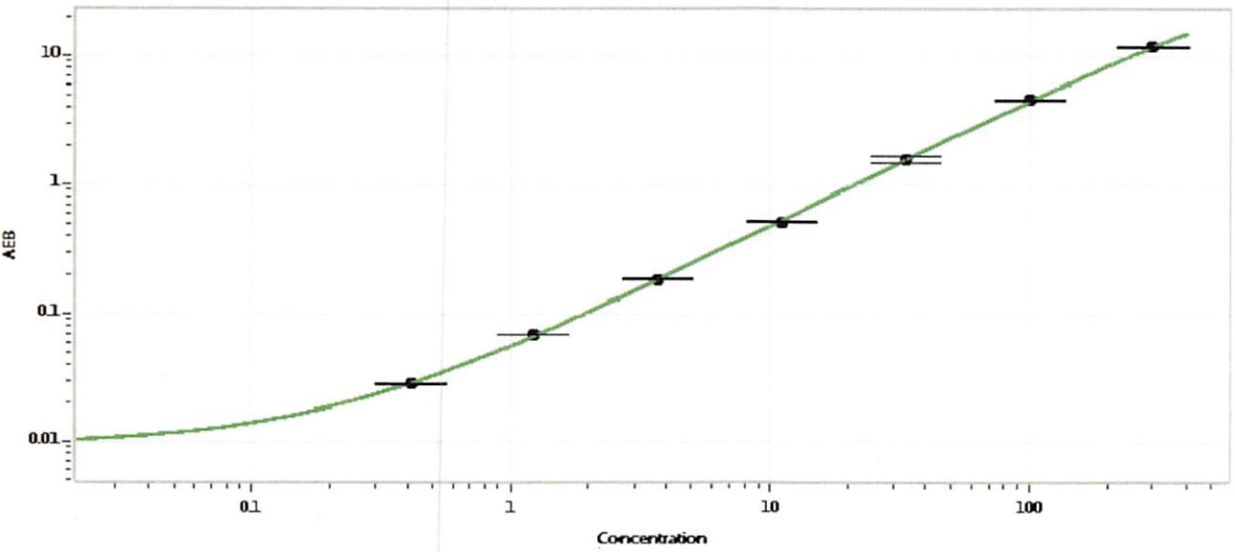
Generic Components - Included With Kit But Not Lot Number Specific		
Discovery Bead Diluent		101954
Discovery Detector Diluent		101955
Discovery SBG Diluent		101956
Discovery SBG Stock		103297
RGP Reagent		103159



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Calibrator Concentrate Target Value	6000 pg/mL
-------------------------------------	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.412
C	1.23
D	3.70
E	11.1
F	33.3
G	100
H	300

Review/Approval	Patrizia Stadler	Assoc. Director QC	P. Stadler	27 MAR 2019
	Name	Title	Signature/Date	