

Product Number:	101863
Lot Number:	501956
Expiration:	15-Mar-2020
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	101904	921009
Detector Stock	101905	921113
Calibrator Concentrate	101858	MF-031518-02
Calibrator Diluent	101990	917621
Sample Diluent	101991	921004

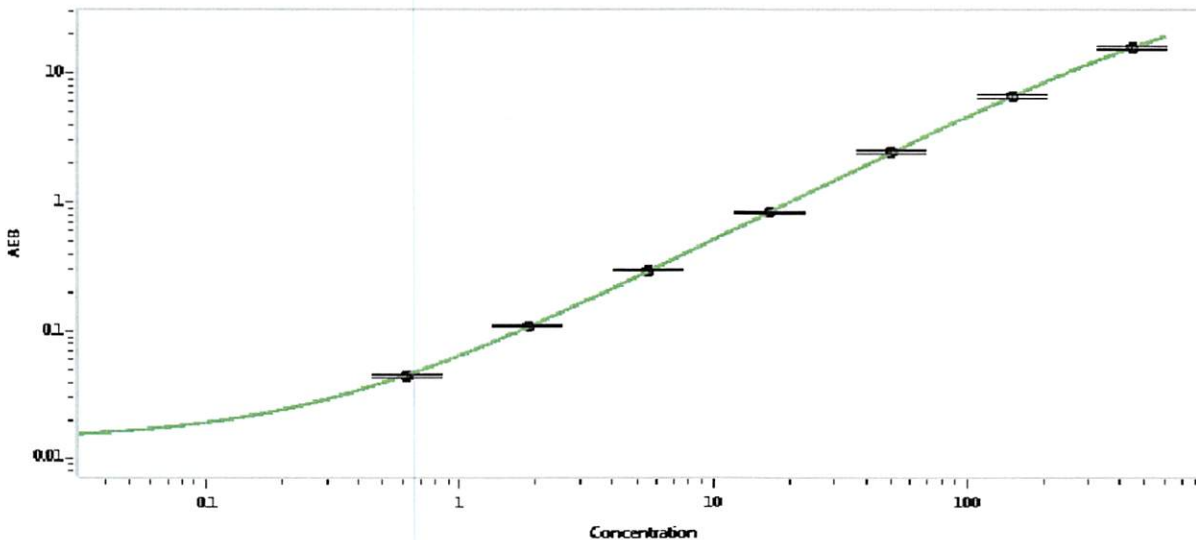


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	9 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.617
C	1.85
D	5.56
E	16.7
F	50.0
G	150
H	450

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i> 07 Aug 2019
	Name	Title	Signature/Date