

Product Number:	102929
Lot Number:	501602
Expiration:	18-Aug-2019
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	102922	902210
Detector Stock	102923	901806
Calibrator Concentrate	102926	MF-081817-01
Calibrator Diluent	102924	901805
Sample Diluent	102925	901804

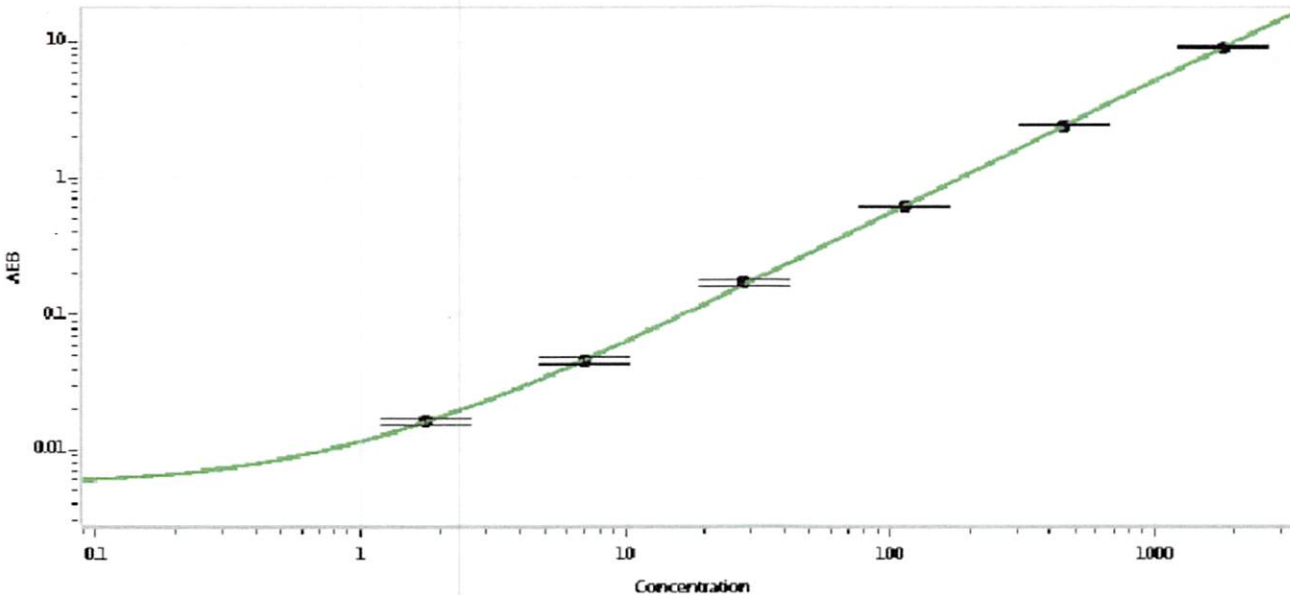
Generic Components - Included With Kit But Not Lot Number Specific		
Discovery Bead Diluent		101954
Discovery Detector Diluent		101955
Discovery SBG Diluent		101956
Discovery SBG Stock		103297
RGP Reagent		103159

Calibrator Concentrate Target Value	36 ng/mL
-------------------------------------	----------



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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.76
C	7.03
D	28.1
E	113
F	450
G	1800

Review/Approval	Patrizia Stadler	Associate Director, QC	<i>p. Stadler</i> 28 Jan 2019
	Name	Title	Signature/Date