

<b>Product Number:</b>	102794
<b>Lot Number:</b>	501632
<b>Expiration:</b>	10-Jan-2020
<b>Platform(s):</b>	HD-1

Component	Part Number	Lot Number
Bead Stock	102787	903204
Detector Stock	102788	903108
Calibrator Concentrate	102791	MF-021018-04
Calibrator Diluent	102789	901914
Sample Diluent	102790	903203

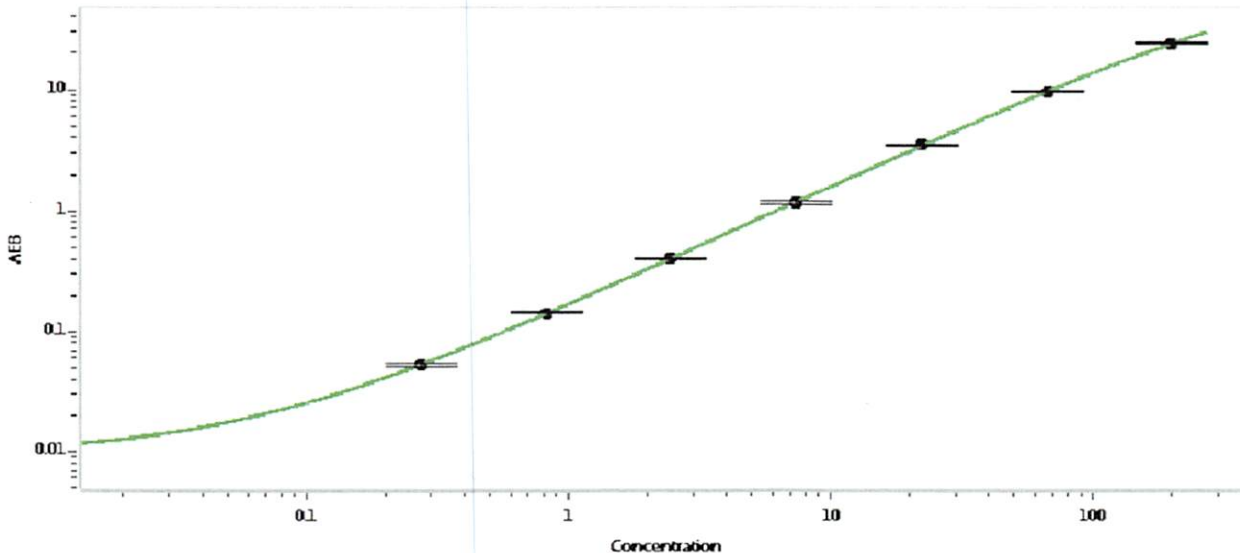
Generic Components - Included With Kit But Not Lot Number Specific		
Discovery Bead Diluent		101954
Discovery Detector Diluent		101955
Discovery SBG Diluent		101956
Discovery SBG Stock		103297
RGP Reagent		103159

<b>Calibrator Concentrate Target Value</b>	4 ng/mL
--	---------



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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	15 Feb 2019
	Name	Title	Signature/Date	