

Product Number:	102635
Lot Number:	502401
Expiration:	17-Mar-2021
Platform(s):	HD-1/ HD-X

Component	Part Number	Lot Number
Bead Stock	102628	013504
Detector Stock	102629	014804
Calibrator Concentrate	102632	013308
Calibrator Diluent	102630	007814
Sample Diluent	102631	013601

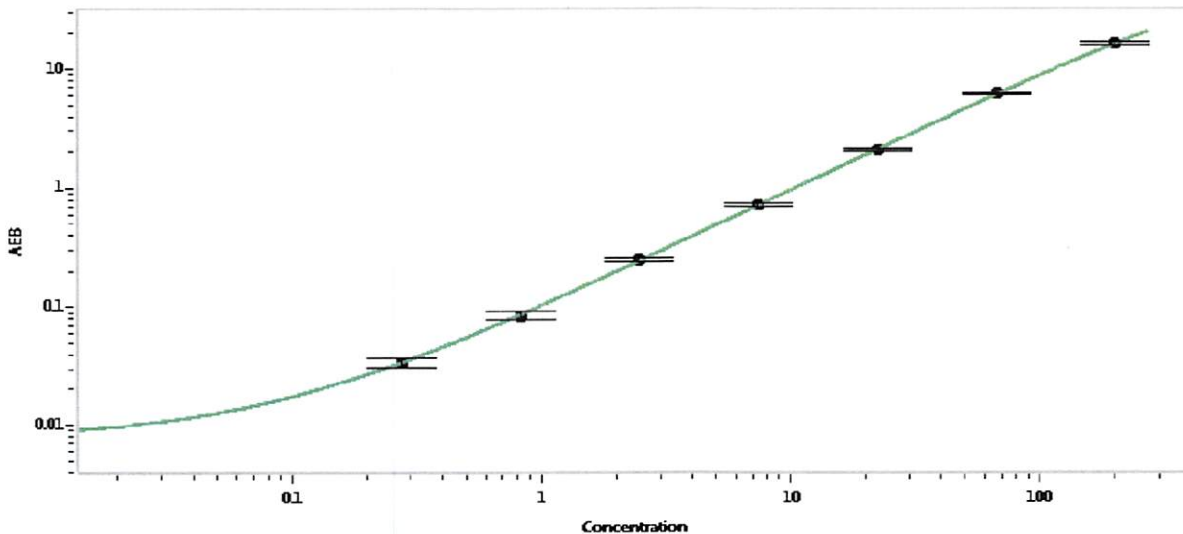


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	4 ng/mL
-------------------------------------	---------

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	Shivani Goel	QC Manager	Shivani Goel 02Jun2020
	Name	Title	Signature/Date