

Product Number:	102336
Lot Number:	503164
Expiration:	13-Jun-2022
Platform(s):	HD-1 / HD-X

Component	Part Number	Lot Number
Bead Stock	102330	127202
Detector Stock	102331	127505
Calibrator Concentrate	102335	127205
Calibrator Diluent	102332	124009
Sample Diluent	102333	127302

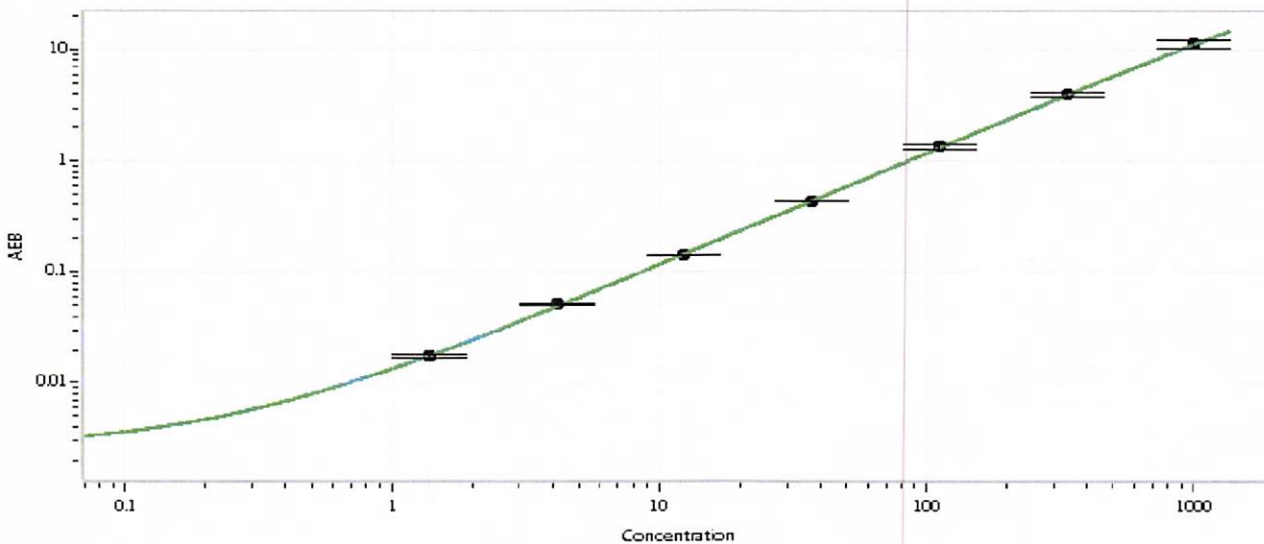


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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

Calibrator Concentrate Target Value	20 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.11
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval: Shivani Goel Name: Shivani Goel Title: QC Signature/Date: Shivani Goel 20oct2021