

Product Number:	101984
Lot Number:	502388
Expiration:	20-Mar-2021
Platform(s):	HD-1/HD-X



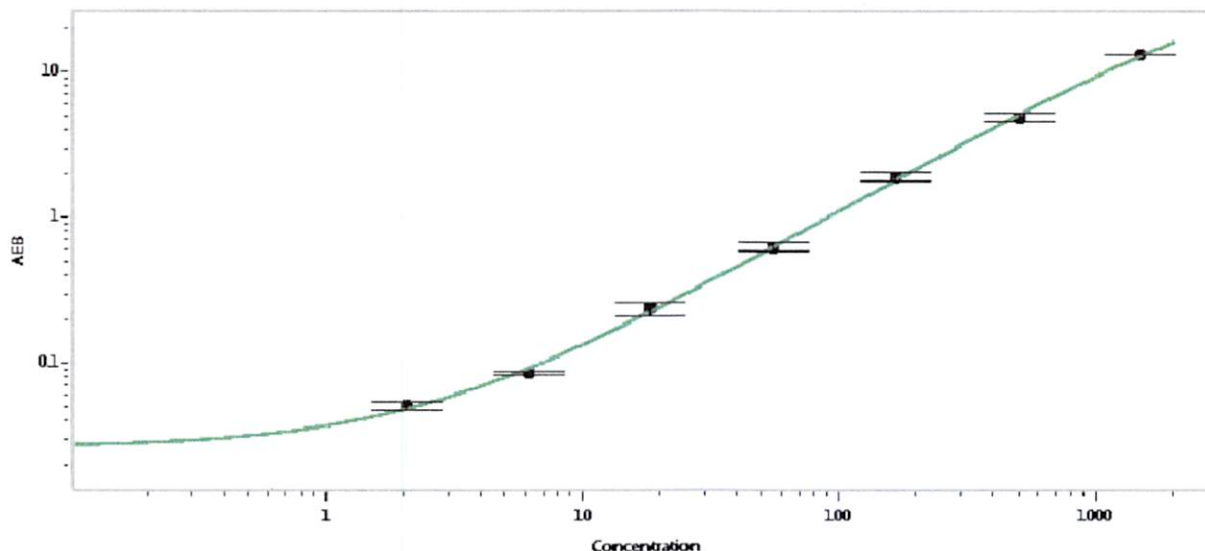
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101910	013802
Detector Stock	101911	013804
Calibrator Concentrate	101983	910010
Calibrator Diluent	101908	014501
Sample Diluent	101909	014502

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	30 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.06
C	6.17
D	18.5
E	55.6
F	167
G	500
H	1500

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 08/02/2020  
Signature/Date