

Product Number:	101984
Lot Number:	502550
Expiration:	10-Apr-2021
Platform(s):	HD-1/HD-X



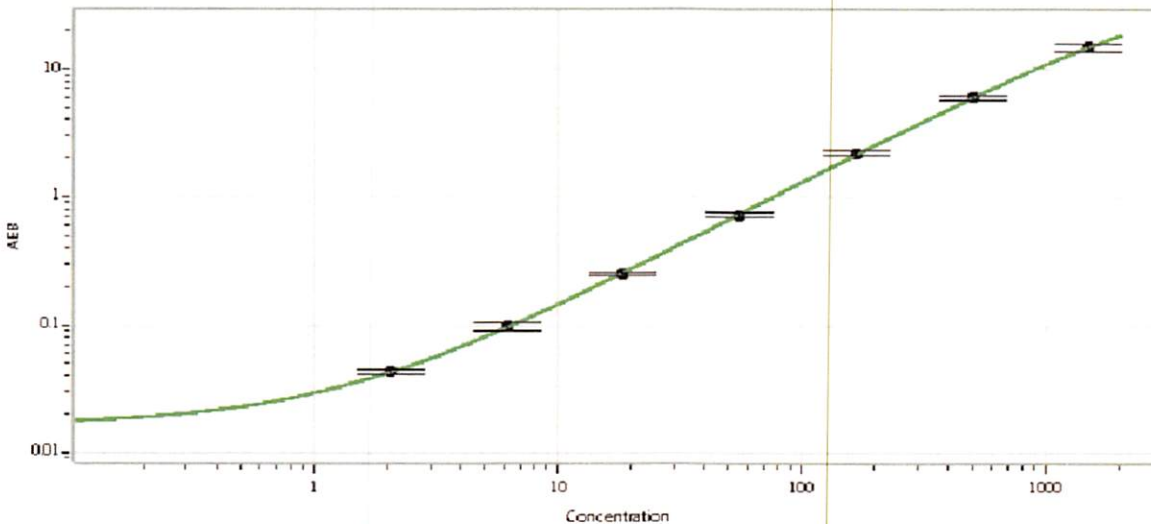
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101910	027205
Detector Stock	101911	027303
Calibrator Concentrate	101983	910010
Calibrator Diluent	101908	027305
Sample Diluent	101909	027304

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 23 Oct 2020
Signature/Date