

Product Number:	102491
Lot Number:	502498
Expiration:	13-Feb-2021
Platform(s):	HD-1/HD-X



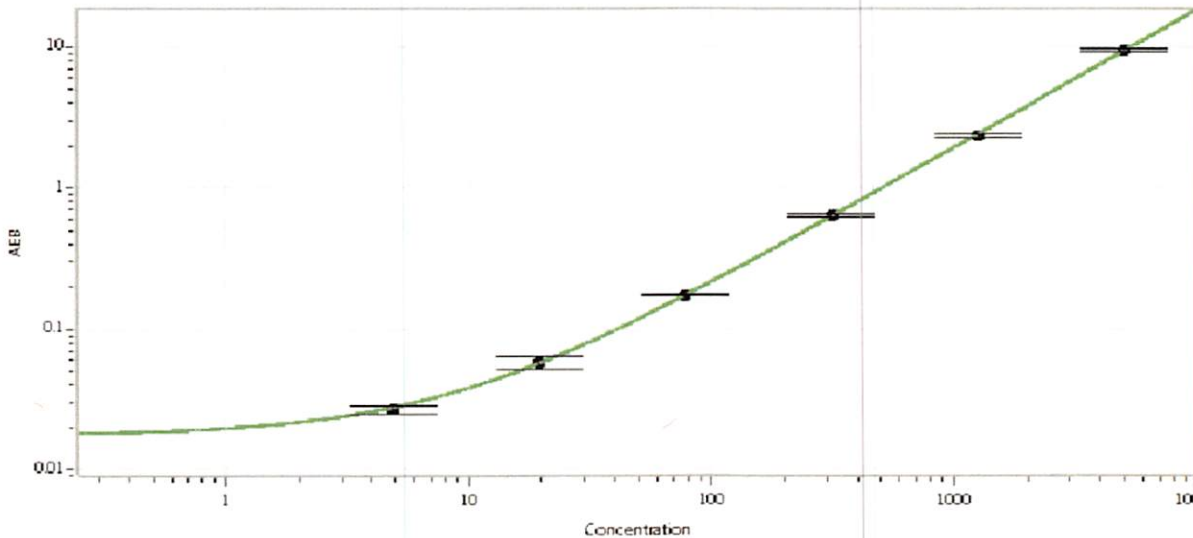
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102484	022601
Detector Stock	102485	023001
Detector Diluent	103084	023006
Calibrator Concentrate	102488	907004
Calibrator Diluent	102486	023003
Sample Diluent	102487	023002

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

Calibrator Concentrate Target Value	100 ng/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	4.88
C	19.5
D	78.1
E	313
F	1250
G	5000

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 24 Aug 2020
	Name	Title	Signature/Date