

Product Number:	103059
Lot Number:	502363
Expiration:	21-Feb-2021
Platform(s):	HD-1 / HD-X



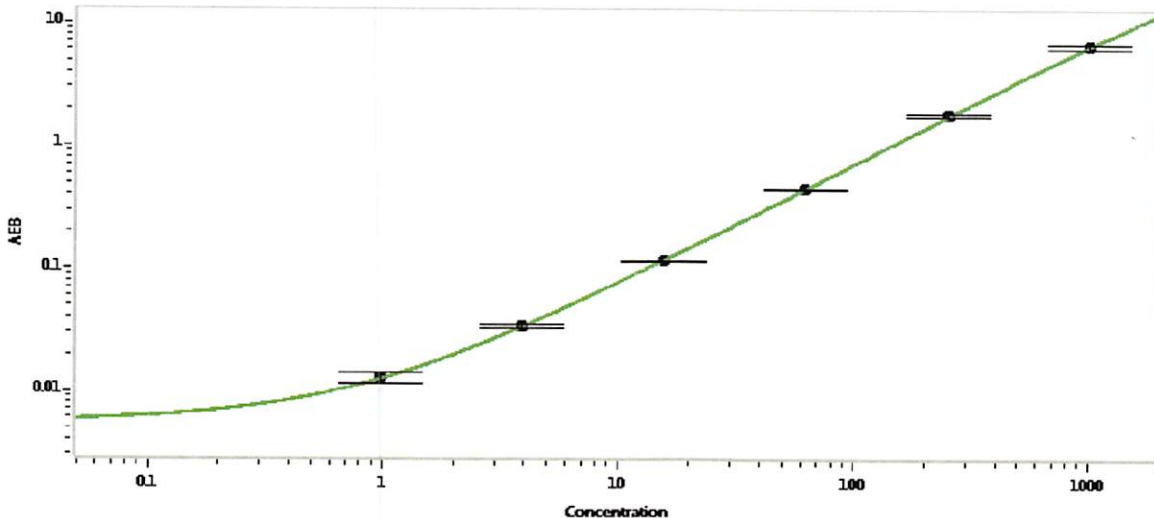
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	103052	012607
Detector Stock	103053	012502
Calibrator Concentrate	103054	012501
Calibrator Diluent	103056	010811
Sample Diluent	103055	005229
Detector Diluent	103142	010812

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	20,000 pg/mL
-------------------------------------	--------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.977
C	3.91
D	15.6
E	62.5
F	250
G	1000

Review/Approval	Shivani Goel	QC Manager	Shivani Goel (15 May 2020)
	Name	Title	Signature/Date