

Product Number:	102318
Lot Number:	502391
Expiration:	7-Jan-2021
Platform(s):	HD-1/HDX



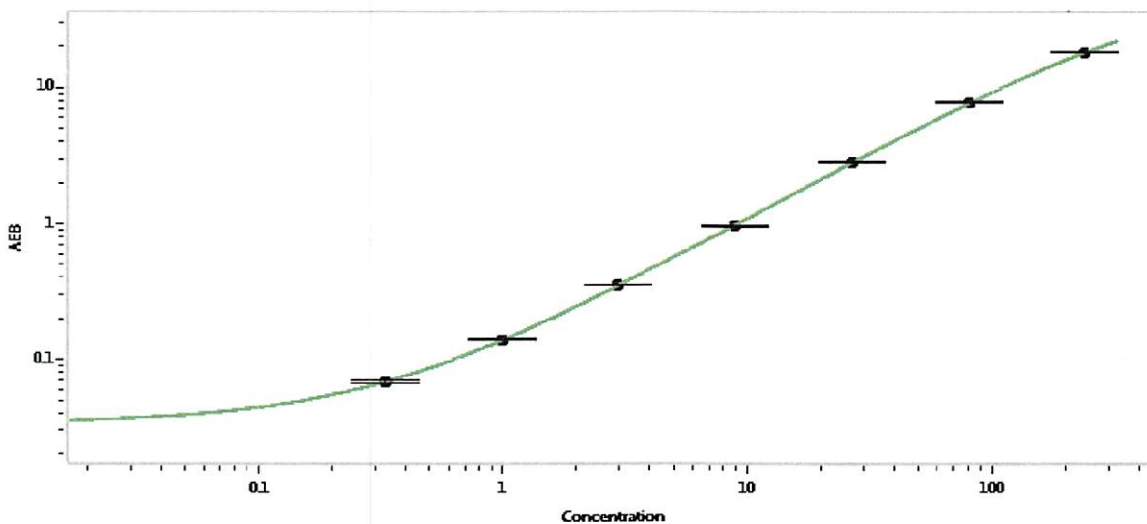
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102309	014008
Detector Stock	102310	000722
Calibrator Concentrate	102314	929702
Calibrator Diluent	102311	014001
Sample Diluent	102312	014002

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	4,800 pg/mL
-------------------------------------	-------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.329
C	0.988
D	2.96
E	8.89
F	26.7
G	80.0
H	240

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 02Jun2020
	Name	Title	Signature/Date