

Product Number:	101968
Lot Number:	503317
Expiration:	3-Sep-2022
Platform(s):	HD-1/HD-X



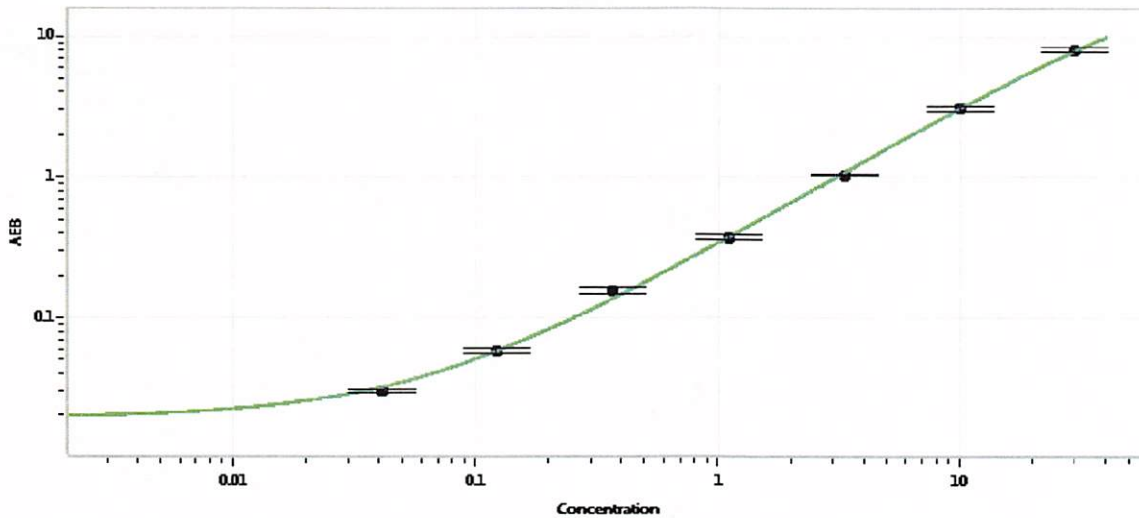
Data below represents results generated on the Simoa® HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101966	202104
Detector Stock	101967	202203
Calibrator Concentrate	101965	206206
Calibrator Diluent	101963	129130
Sample Diluent	101964	133917

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	600 pg/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval: Shivani Goel Name, QC Manager Title, Shivani Goel 05Apr2022 Signature/Date