

Product Number:	102039
Lot Number:	503297
Expiration:	30-Jul-2022
Platform(s):	HD-1/HD-X & SR-X



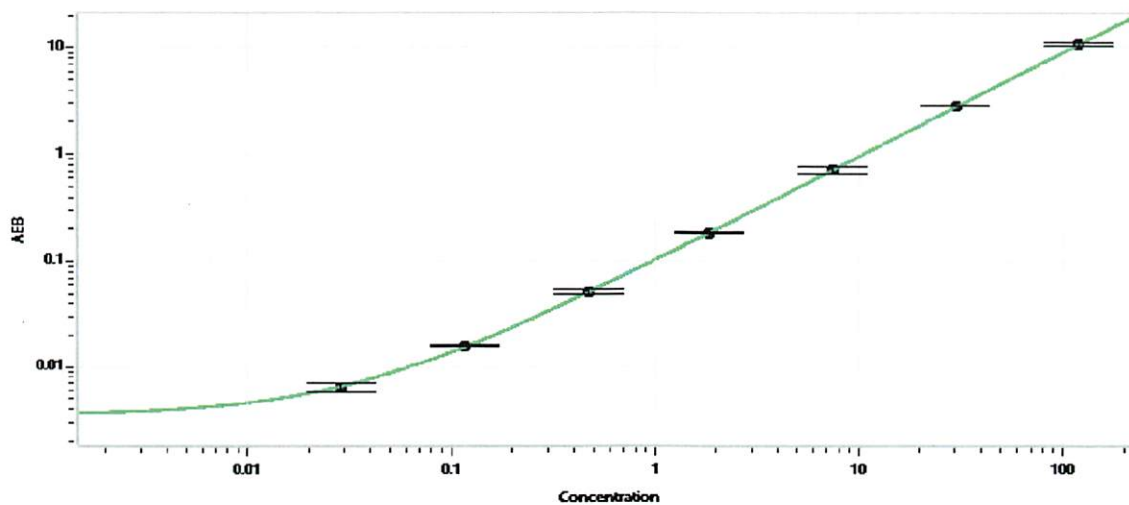
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102028	201214
Detector Stock	102029	136403
Calibrator Concentrate	102032	119607
Calibrator Diluent	102030	128017
Sample Diluent	103402	121124

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	2 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.029
C	0.117
D	0.469
E	1.88
F	7.50
G	30.0
H	120

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 20Jan2022

Name

Title

Signature/Date