

Product Number:	102039
Lot Number:	502786
Expiration:	30-Oct-2021
Platform(s):	HD-1/HD-X & SR-X



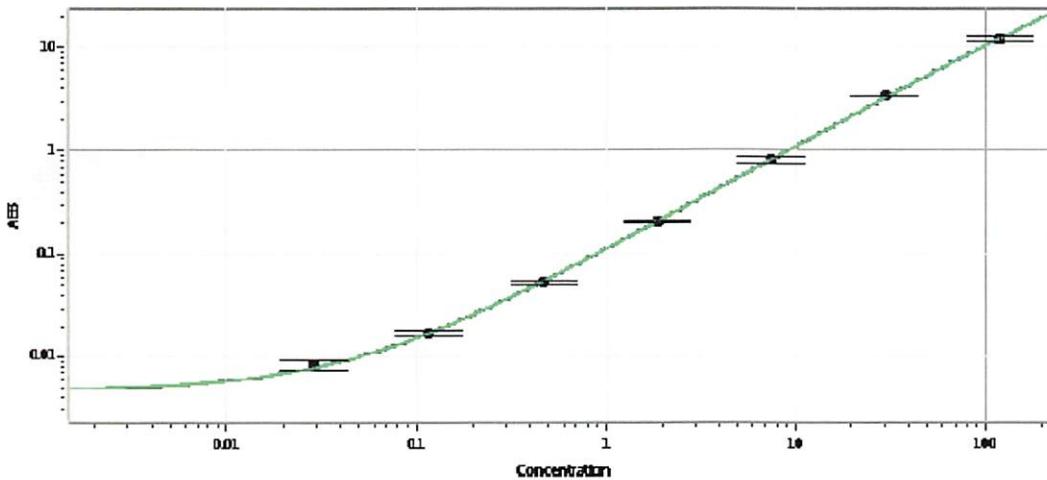
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102028	032207
Detector Stock	102029	032407
Calibrator Concentrate	102032	032408
Calibrator Diluent	102030	031412
Sample Diluent	103402	030405

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

Junpil Park
Name

QC Supervisor
Title

 04Mar2021
Signature/Date