

Product Number:	102394
Lot Number:	503009
Expiration:	4-Mar-2022
Platform(s):	HD-1/ HDX



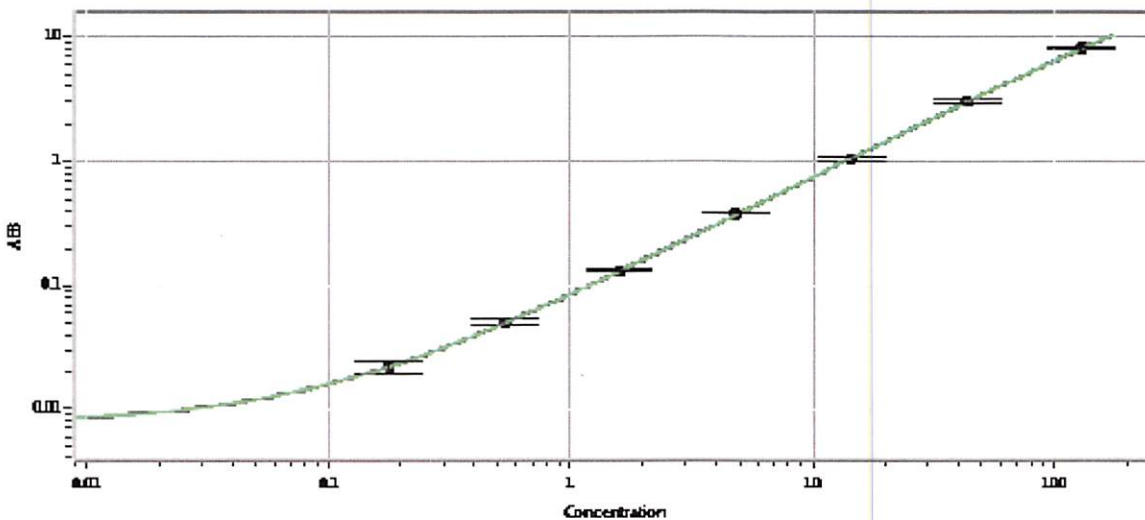
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102387	116706
Detector Stock	102388	109302
Calibrator Concentrate	102391	015401
Calibrator Diluent	102389	106320
Sample Diluent	102390	109020

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	4,000 pg/mL
-------------------------------------	-------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.178
C	0.535
D	1.60
E	4.81
F	14.4
G	43.3
H	130

Review/Approval

Junpil Park  
Name

QC Supervisor  
Title

*[Signature]* 24 Jun 2021  
Signature/Date