

Product Number:	102648
Lot Number:	502942
Expiration:	17-Mar-2022
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Stock	102641	113807
Detector Stock	102642	113913
Calibrator Concentrate	102645	114006
Calibrator Diluent	102643	112910
Sample Diluent	102644	111912
Detector Diluent	102734	113912

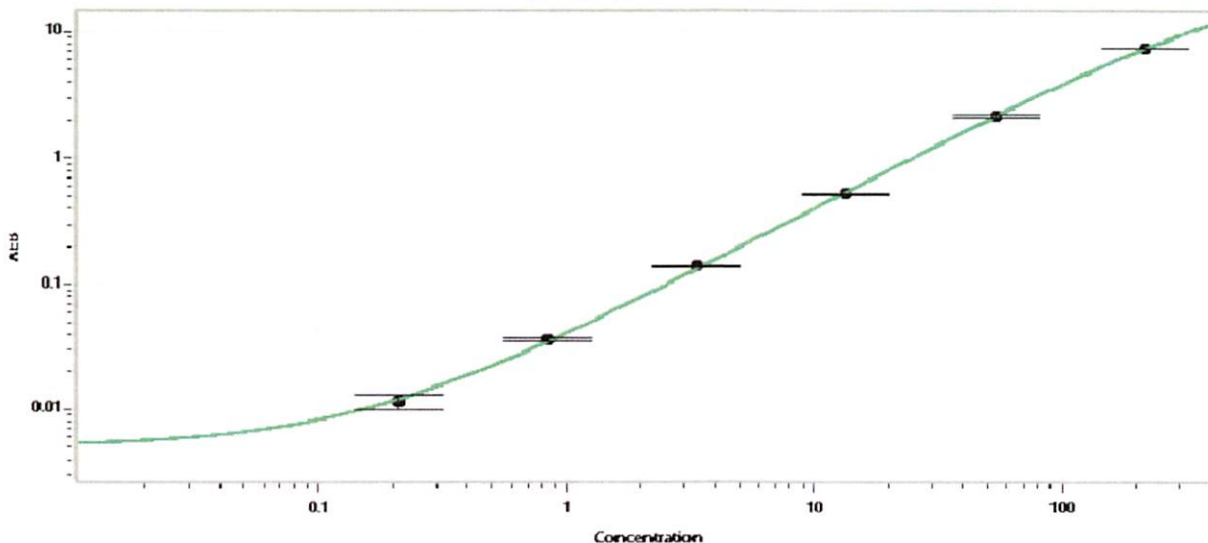


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159

Calibrator Concentrate Target Value	3,000 pg/mL
-------------------------------------	-------------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.210
C	0.840
D	3.36
E	13.4
F	53.8
G	215

Review/Approval

Junpil Park  
Name

QC Supervisor  
Title

*[Signature]* 07 Jun 2021  
Signature/Date