

Product Number:	103071
Lot Number:	502660
Expiration:	28-Sep-2021
Platform(s):	HD-1/ HD-X & SR-X

Component	Part Number	Lot Number
Bead Stock	103064	034901
Detector Stock	103065	035002
Calibrator Concentrate	103068	934309
Calibrator Diluent	103066	031907
Sample Diluent	103067	033809

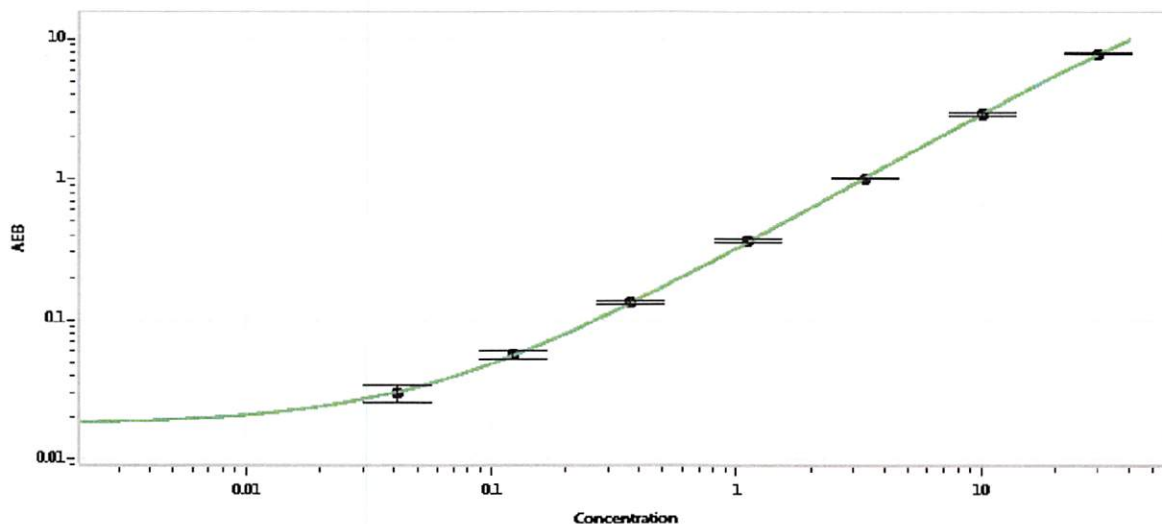
Generic Components - Included With Kit But Not Lot Number Specific		
Discovery Bead Diluent		101954
Discovery Detector Diluent		101955
Discovery SBG Diluent		101956
Discovery SBG Stock		103297
RGP Reagent		103159



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

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

Junpil Park

Name

QC Supervisor

Title

30 Dec 2020

Signature/Date