

Product Number:	103071
Lot Number:	503109
Expiration:	17-Mar-2022
Platform(s) :	HD-1/ HD-X & SR-X

Component	Part Number	Lot Number
Bead Stock	103064	122601
Detector Stock	103065	114001
Calibrator Concentrate	103068	114717
Calibrator Diluent	103066	112911
Sample Diluent	103067	111913

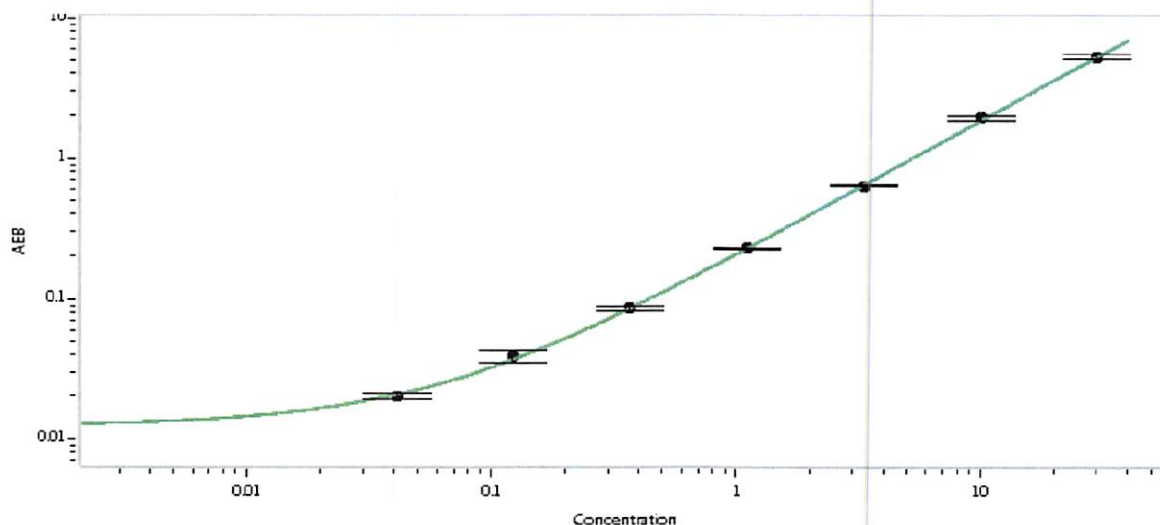
Generic Components - Included With Kit But Not Lot Number Specific		
Discovery Bead Diluent		101954
Discovery Detector Diluent		101955
Discovery SBG Diluent		101956
Discovery SBG Stock		103297
RGP Reagent		103159



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

[Signature] 25 Aug 2021
Signature/Date