

Product Number:	103071
Lot Number:	503133
Expiration:	20-May-2022
Platform(s):	HD-1/ HD-X & SR-X

Component	Part Number	Lot Number
Bead Stock	103064	125308
Detector Stock	103065	125806
Calibrator Concentrate	103068	114717
Calibrator Diluent	103066	121311
Sample Diluent	103067	121113

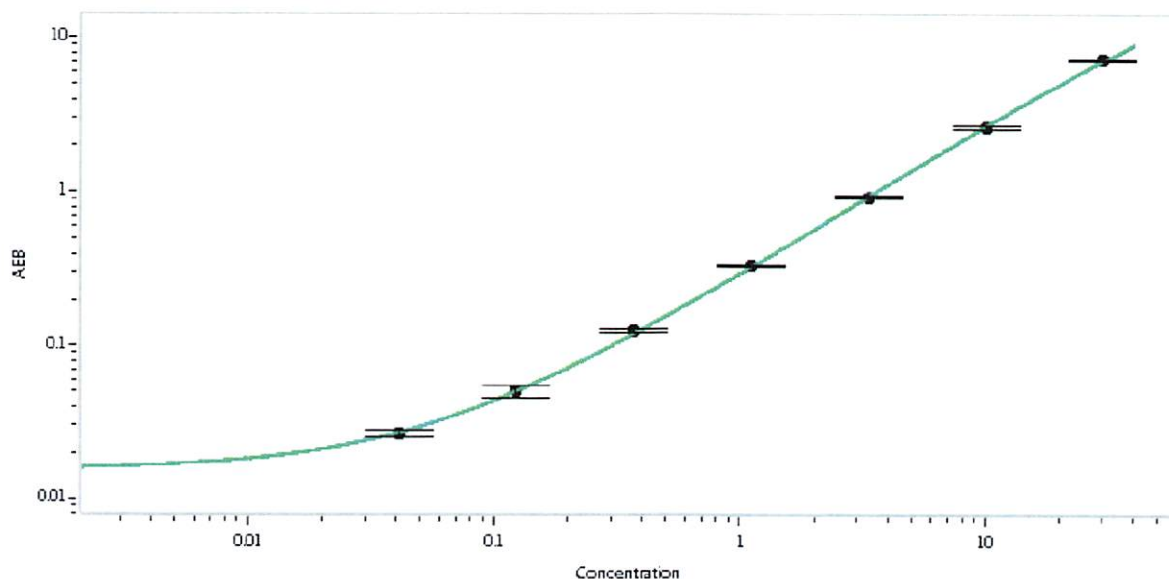
Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159



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/Date