

Product Number:	101968
Lot Number:	502491
Expiration:	18-Feb-2021
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	101966	022501
Detector Stock	101967	022602
Calibrator Concentrate	101965	000719
Calibrator Diluent	101963	020423
Sample Diluent	101964	004922

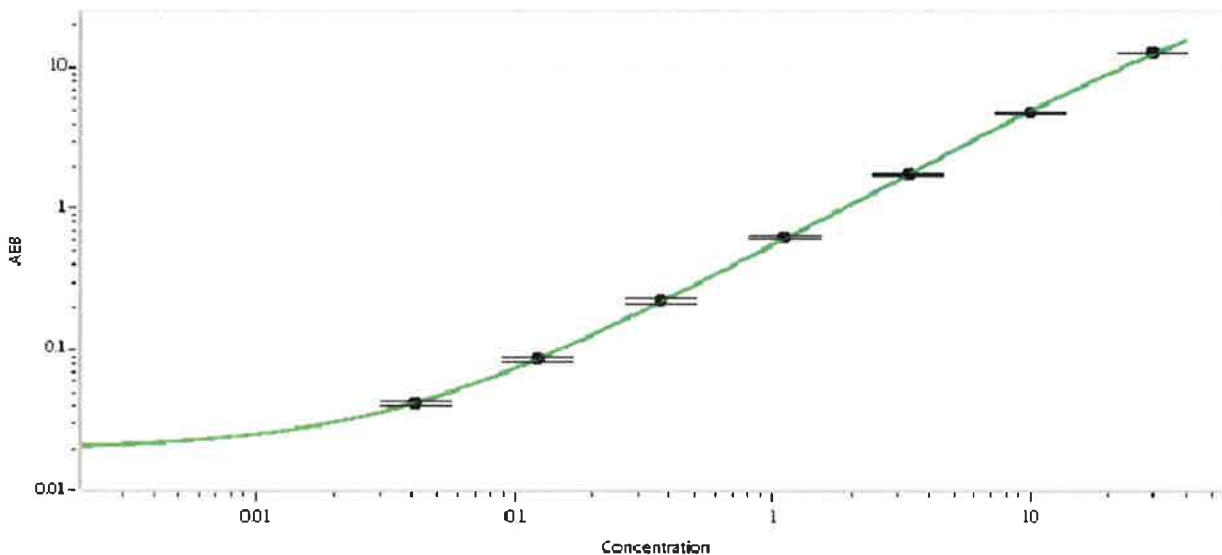
Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159



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

Shivani Goel
Name

QC Manager
Title

Shivani Goel 20 Aug 2020
Signature/Date