

Product Number:	102517
Lot Number:	503701
Expiration:	14-Jul-2023
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Stock	102510	227305
Detector Stock	102511	227308
Calibrator Concentrate	102514	209801
Calibrator Diluent	102512	219514
Sample Diluent	102513	219513

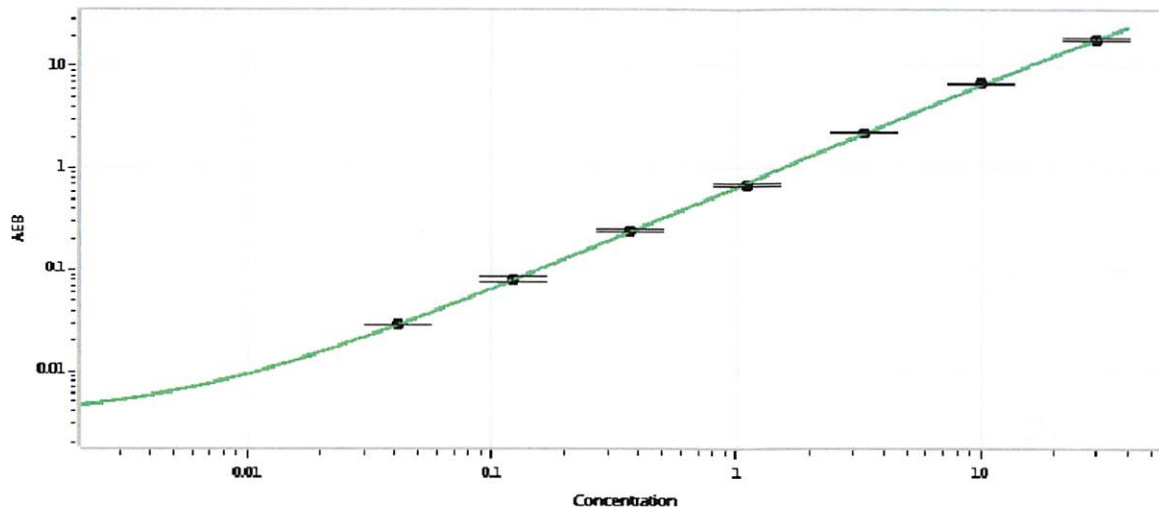
Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159



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

Calibrator Concentrate Target Value (pg/mL)	600
---	-----

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

*Sonia Khillan*  
Name

QC  
Title

*Sonia Khillan / 14/10/2022*  
Signature/Date