

Product Number:	102517
Lot Number:	503274
Expiration:	24-Oct-2022
Platform(s):	HD-1/HD-X & SR-X



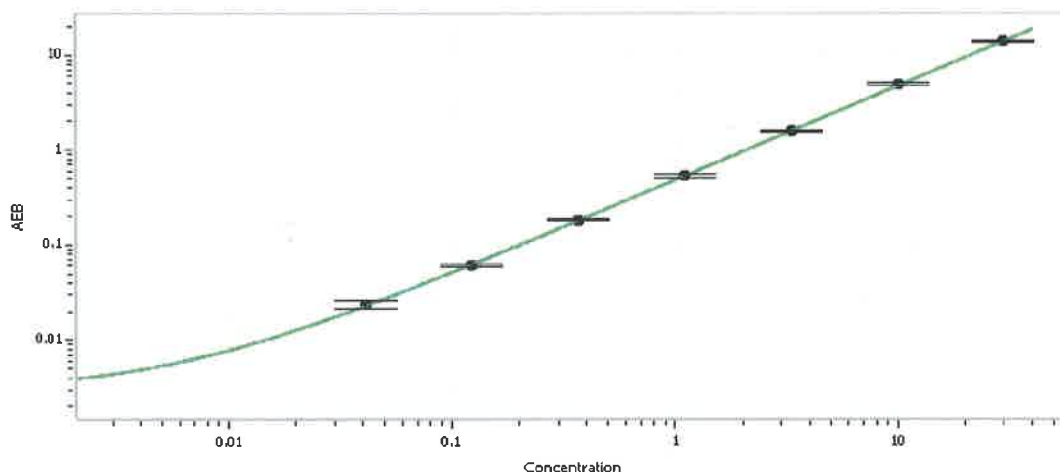
Component	Part Number	Lot Number
Bead Stock	102510	134914
Detector Stock	102511	135001
Calibrator Concentrate	102514	209801
Calibrator Diluent	102512	129708
Sample Diluent	102513	129707

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

Richa Shrestha

QC Supervisor

Name

Title

Signature/Date

Richa Shrestha 22 Apr 2022