

Product Number:	102700
Lot Number:	503354
Expiration:	10-Nov-2022
Platform(s):	HD-1/HD-X



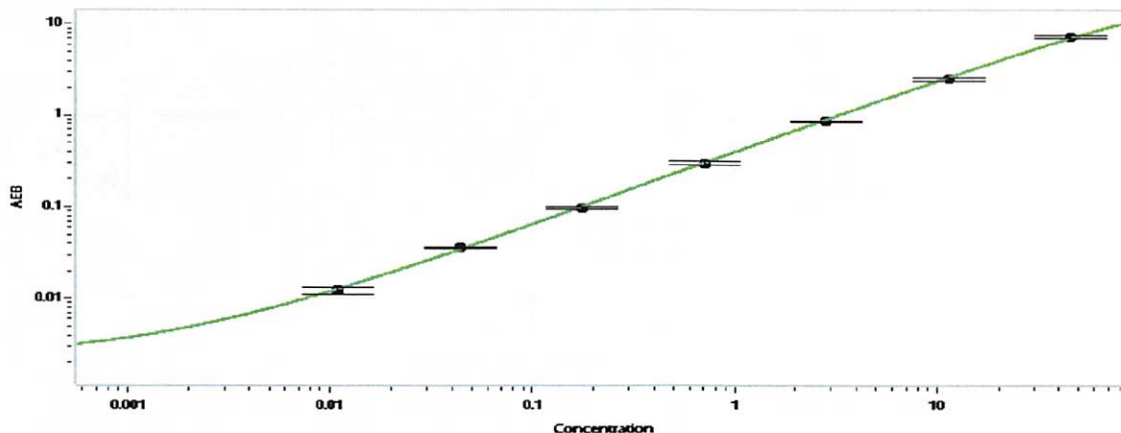
Component	Part Number	Lot Number
Bead Stock	102693	206001
Detector Stock	102694	206008
Calibrator Concentrate	102697	206403
Calibrator Diluent	102695	133935
Sample Diluent	102696	133936

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

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

Calibrator Concentrate Target Value	1000 pg/mL
-------------------------------------	------------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.011
C	0.044
D	0.176
E	0.703
F	2.81
G	11.3
H	45.0

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 31 Mar 2022

Name

Title

Signature/Date