

Product Number:	102336
Lot Number:	503383
Expiration:	15-Feb-2023
Platform(s):	HD-1 & HD-X



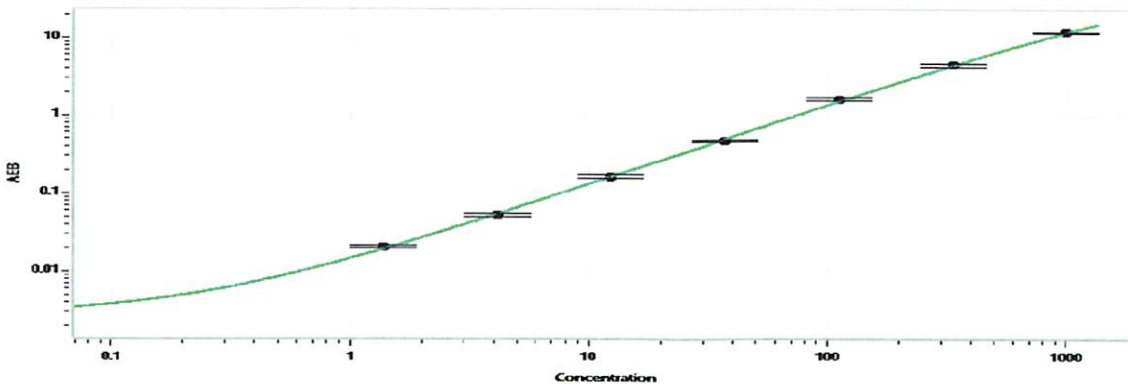
Component	Part Number	Lot Number
Bead Stock	102330	206404
Detector Stock	102331	206503
Calibrator Concentrate	102335	203112
Calibrator Diluent	102332	206505
Sample Diluent	102333	206506

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	N/A
Discovery Detector Diluent	101955	N/A
Discovery SBG Diluent	101956	N/A
Discovery SBG Concentrate	103297	N/A
RGP Reagent	103159	N/A

Calibrator Concentrate Target Value	20 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.12
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 30 Mar 2022
Signature/Date