

Product Number:	102336
Lot Number:	503309
Expiration:	10-Nov-2022
Platform(s):	HD-1 & HD-X



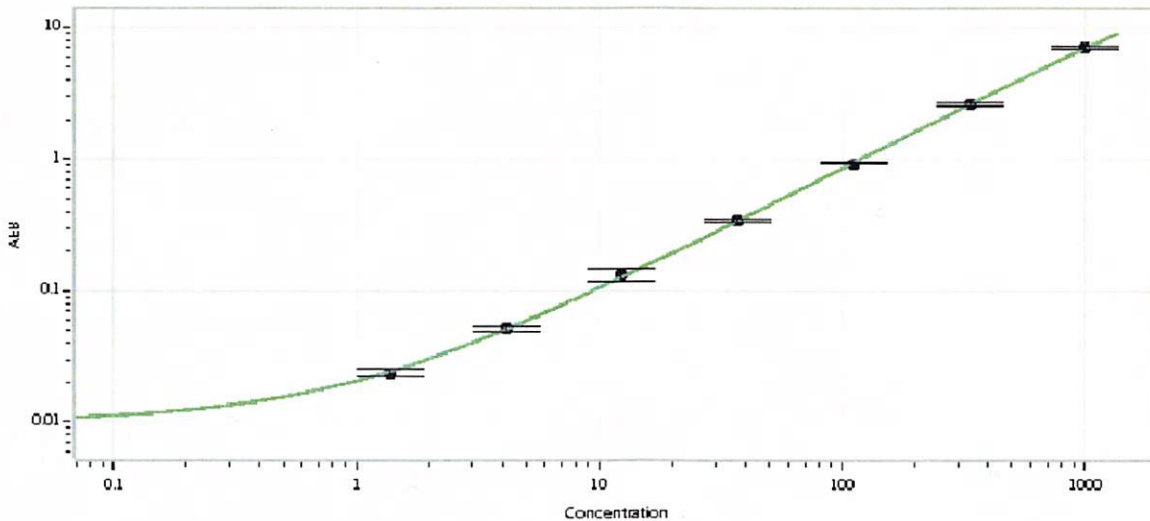
Data below represents results generated on the Simoa® HD-1 & HD-X Analyzer.

Component	Part Number	Lot Number
Bead Stock	102330	202807
Detector Stock	102331	203003
Calibrator Concentrate	102335	203112
Calibrator Diluent	102332	133920
Sample Diluent	102333	203010

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	4.11
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 08 Mar 2022  
Signature/Date