

Product Number:	101984
Lot Number:	504389
Expiration:	13-Dec-2025
Platform(s):	HD-1/HD-X



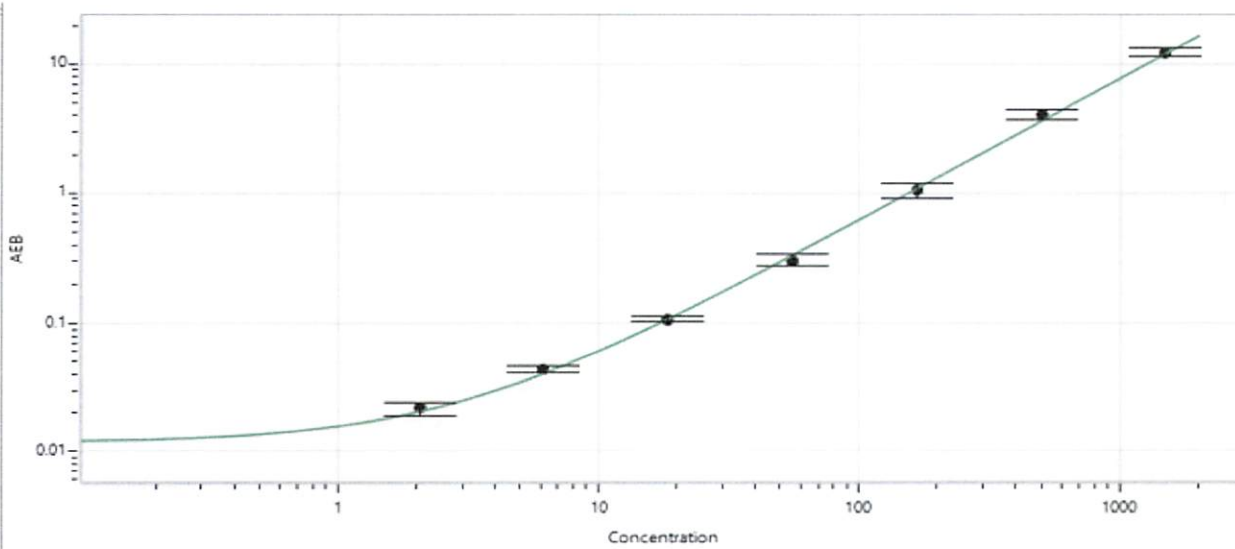
Component	Part Number	Lot Number
Bead Stock	101910	425305
Detector Stock	101911	425505
Calibrator Concentrate	101983	427002
Calibrator Diluent	101908	425308
Sample Diluent	101909	425307

Data below represents results generated on the Simoa® 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 (ng/mL)	30.0 ng/mL
----------------------------------------------------	------------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	2.06
C	6.17
D	18.5
E	55.6
F	167
G	500
H	1500

Review/Approval

Dipesh Patel *Documentation II QA* *et 18 OCT 2024*
 Name Title Signature/Date