

Product Number:	102318
Lot Number:	504328
Expiration:	2-Nov-2025
Platform(s):	HD-1/HD-X



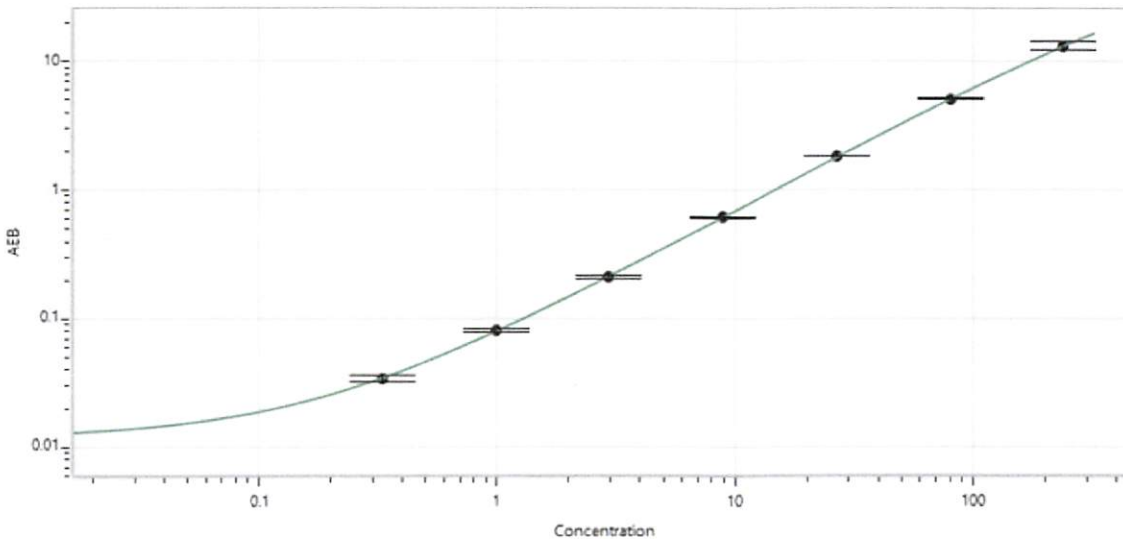
Component	Part Number	Lot Number
Bead Stock	102309	421513
Detector Stock	102310	421404
Calibrator Concentrate	102314	422202
Calibrator Diluent	102311	421208
Sample Diluent	102312	421202

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 Stock	103297	N/A
RGP Reagent	103159	N/A

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

Calibrator Concentrate Target Value (ng/mL)	4.8 ng/mL
--	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.329
C	0.988
D	2.96
E	8.89
F	26.7
G	80.0
H	240

Review/Approval

Robert Somogyi

QC Manager

[Handwritten Signature] 19 AUG 2024

Name

Title

Signature/Date