

Product Number:	102318
Lot Number:	504271
Expiration:	10-Nov-2025
Platform(s):	HD-1/HD-X



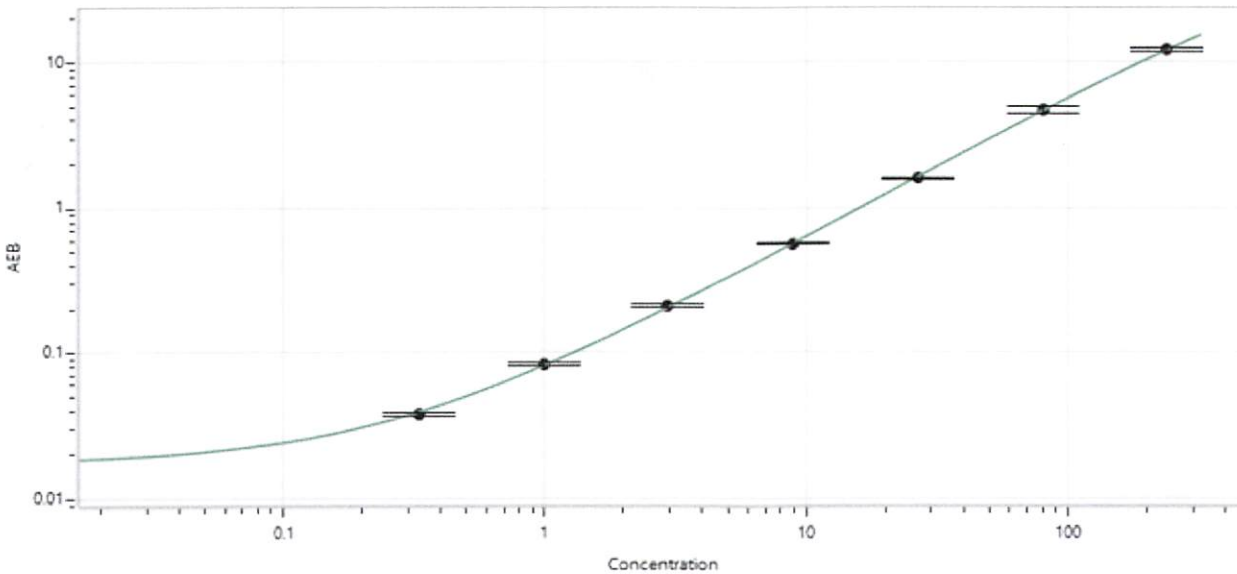
Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number
Bead Stock	102309	413704
Detector Stock	102310	415006
Calibrator Concentrate	102314	417208
Calibrator Diluent	102311	413804
Sample Diluent	102312	414110

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	4,800 pg/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
Name

QC Manager
Title

[Handwritten Signature]
Signature/Date
24 JUN 2024