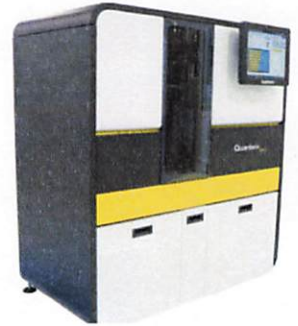


Product Number:	102109
Lot Number:	503707
Expiration:	29-Aug-2023
Platform(s):	HD-1 / HD-X & SR-X



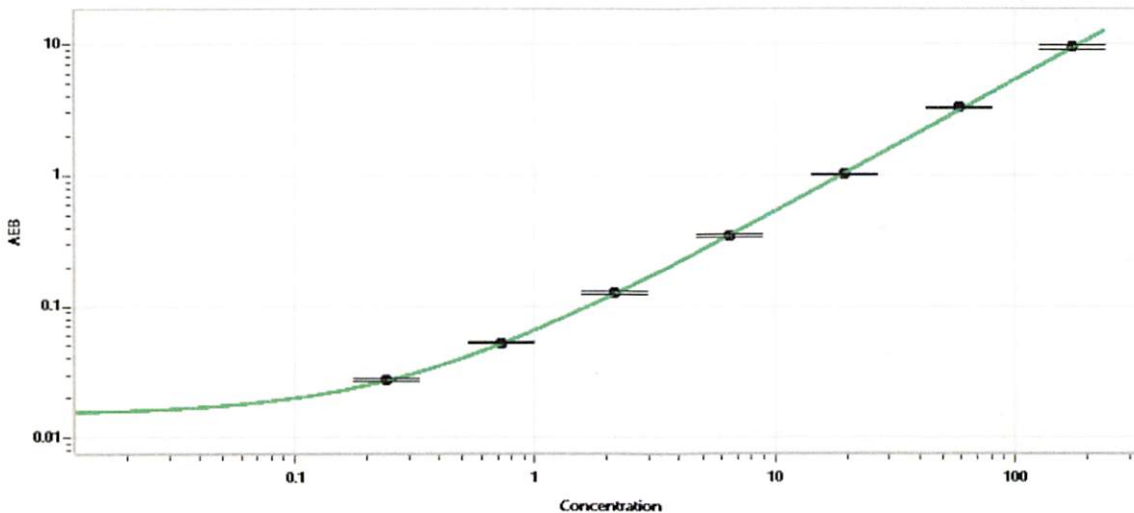
Component	Part Number	Lot Number
Bead Stock	102101	228007
Detector Stock	102102	230610
Calibrator Concentrate	102105	228608
Calibrator Diluent	102103	227715
Sample Diluent	102104	227232

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 (pg/mL)	3,500 pg/ml
---	-------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.240
C	0.720
D	2.16
E	6.48
F	19.4
G	58.3
H	175

Review/Approval

*Sonia Khillan*  
Name

QC

Title

*Sonia Khillan / 18 Nov 2022*  
Signature/Date