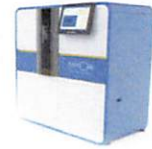


Product Number:	102475
Lot Number:	503355
Expiration:	10-Nov-2022
Platform(s):	HD-1/HD-X



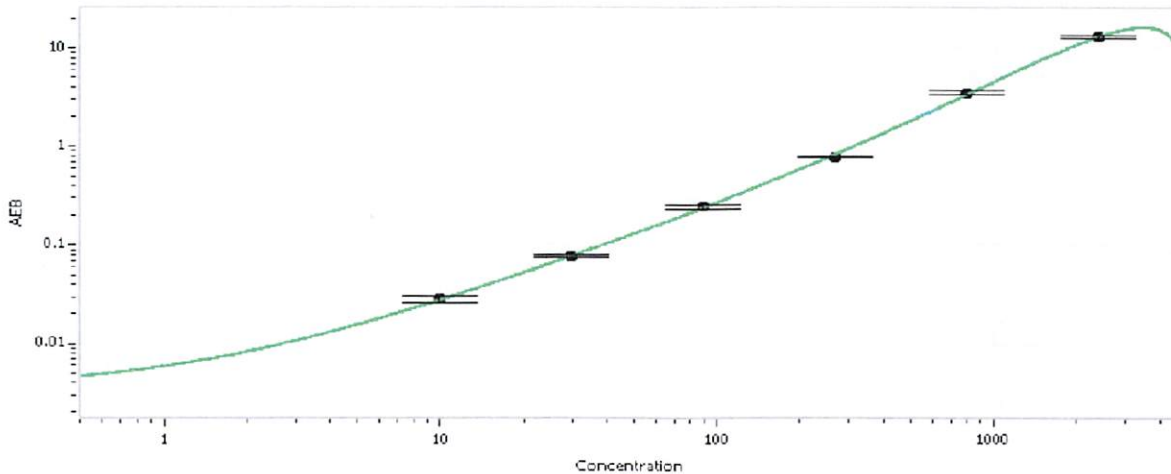
Component	Part Number	Lot Number
Bead Stock	102468	205306
Detector Stock	102469	205401
Calibrator Concentrate	102472	214307
Calibrator Diluent	102470	204612
Sample Diluent	102471	133930

Data below represents results generated on the Simoa® HD-1 / 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	60 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	9.88
C	29.6
D	88.9
E	267
F	800
G	2400

Review/Approval

Matt Smart
Name

Director, QC
Title

MRA 13 Jun 2022
Signature/Date