

Product Number:	102336
Lot Number:	503548
Expiration:	23-May-2023
Platform(s):	HD-1 & HD-X

Component	Part Number	Lot Number
Bead Stock	102330	220712
Detector Stock	102331	220801
Calibrator Concentrate	102335	220811
Calibrator Diluent	102332	220207
Sample Diluent	102333	220802

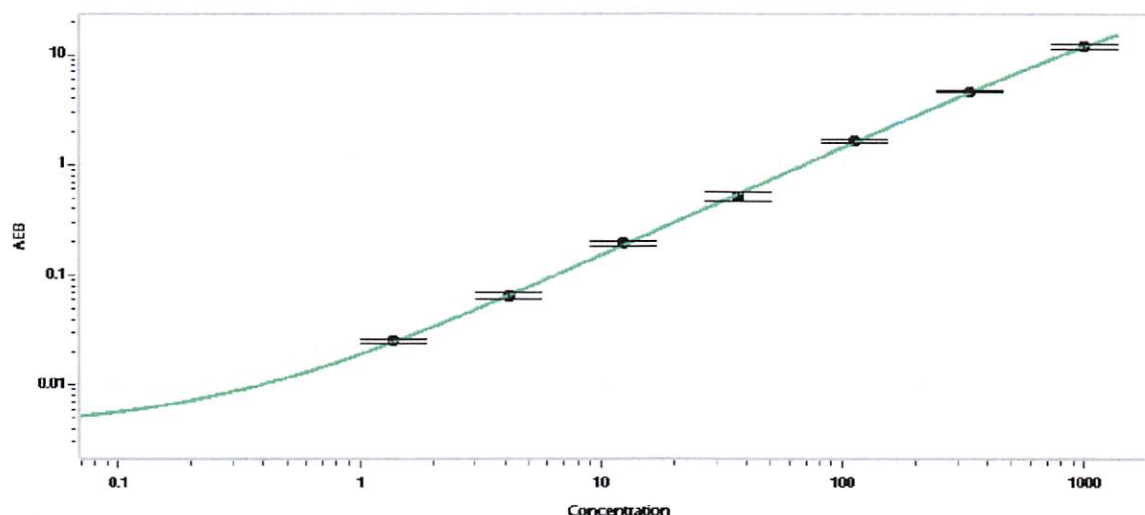


Data below represents results generated on the Simoa® HD-X Analyzer.

Generic Components - Included With Kit But Not Lot Number Specific		
Discovery Bead Diluent	101954	N/A
Discovery Detector Diluent	101955	N/A
Discovery SBG Diluent	101956	N/A
Discovery SBG Concentrate	103297	N/A
RGP Reagent	103159	N/A

Calibrator Concentrate Target Value (ng/mL)	20
--	----

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	1.37
C	4.11
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Matt Smart

Name

QC Director

Title

MDA 15 Aug 2022
Signature/Date