

Product Number:	103071
Lot Number:	503713
Expiration:	22-May-2023
Platform(s):	HD-1/ HD-X & SR-X



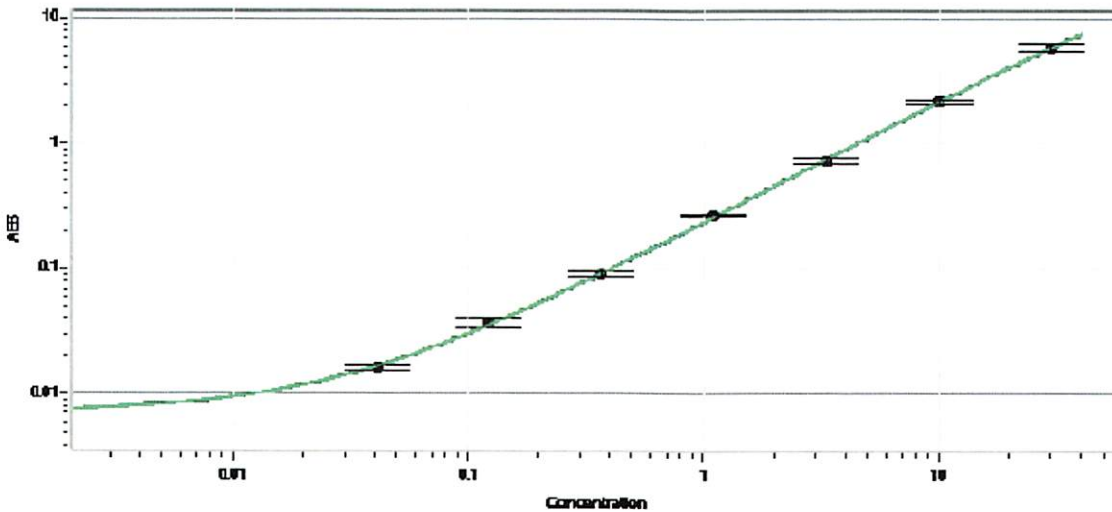
Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number
Bead Stock	103064	228505
Detector Stock	103065	228511
Calibrator Concentrate	103068	212505
Calibrator Diluent	103066	227717
Sample Diluent	103067	227234

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 (pg/mL)	600
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Matt Koulouris
Name

QC Supervisor
Title

[Signature] 03NOV2022
Signature/Date