

Product Number:	101968
Lot Number:	503520
Expiration:	15-Feb-2023
Platform(s):	HD-1/HD-X



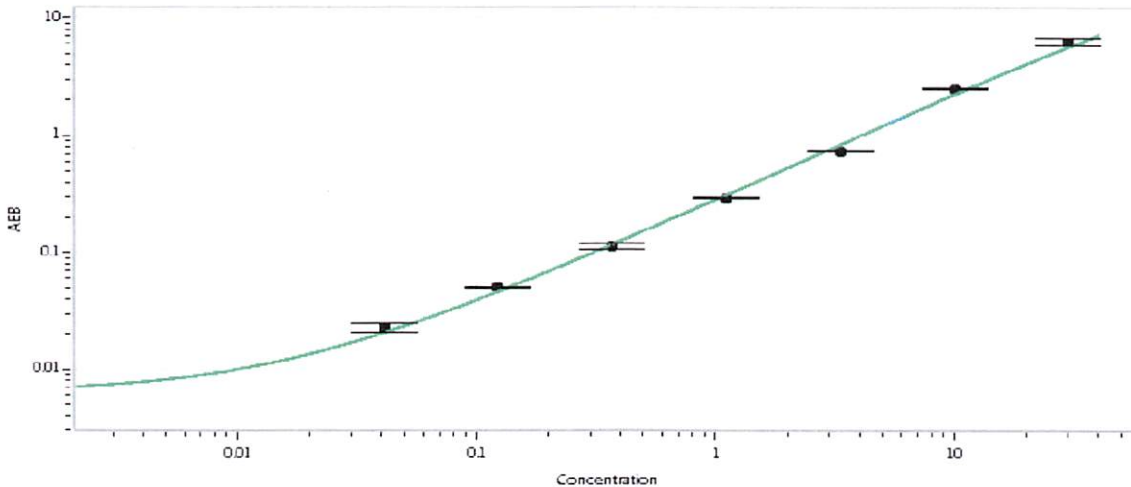
Component	Part Number	Lot Number
Bead Stock	101966	218814
Detector Stock	101967	218605
Calibrator Concentrate	101965	206206
Calibrator Diluent	101963	210415
Sample Diluent	101964	204619

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

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 Smart
Name

QC Director
Title

M. Smart 15 Aug 2022
Signature/Date