

Product Number:	102039
Lot Number:	503562
Expiration:	14-Apr-2023
Platform(s):	HD-1/HD-X & SR-X



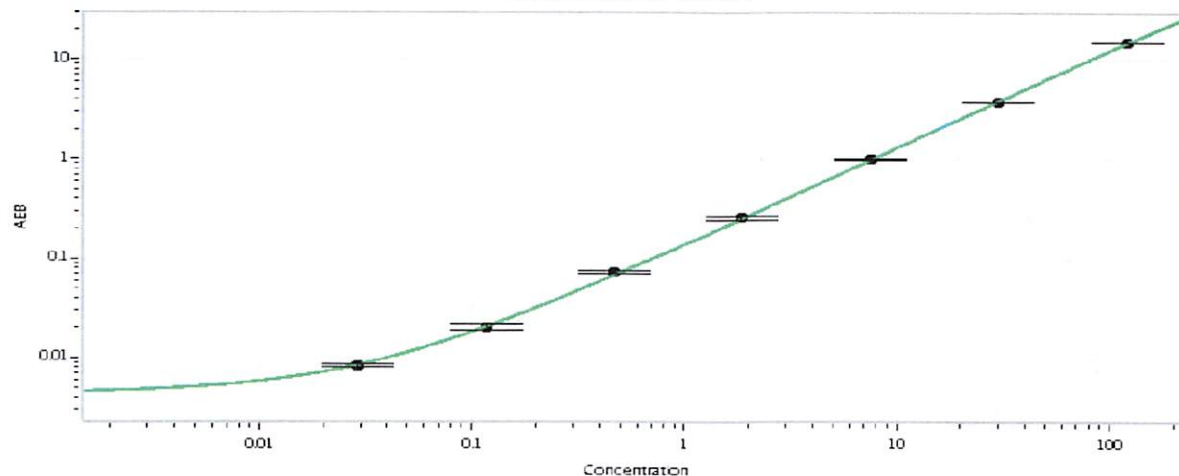
Component	Part Number	Lot Number
Bead Stock	102028	211311
Detector Stock	102029	211401
Calibrator Concentrate	102032	220308
Calibrator Diluent	102030	210410
Sample Diluent	103402	211312

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

Data below represents results generated on the Simoa® HD-1 Analyzer.

Calibrator Concentrate Target Value (ng/mL)	2.00
--	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.029
C	0.117
D	0.469
E	1.88
F	7.50
G	30.0
H	120

Review/Approval

<i>Sonia Khillan</i>	<i>QC Manager</i>	<i>Sonia Khillan / 03AUG2022</i>
Name	Title	Signature/Date