

Product Number:	102109
Lot Number:	503315
Expiration:	24-Jul-2022
Platform(s):	HD-1 / HD-X & SR-X



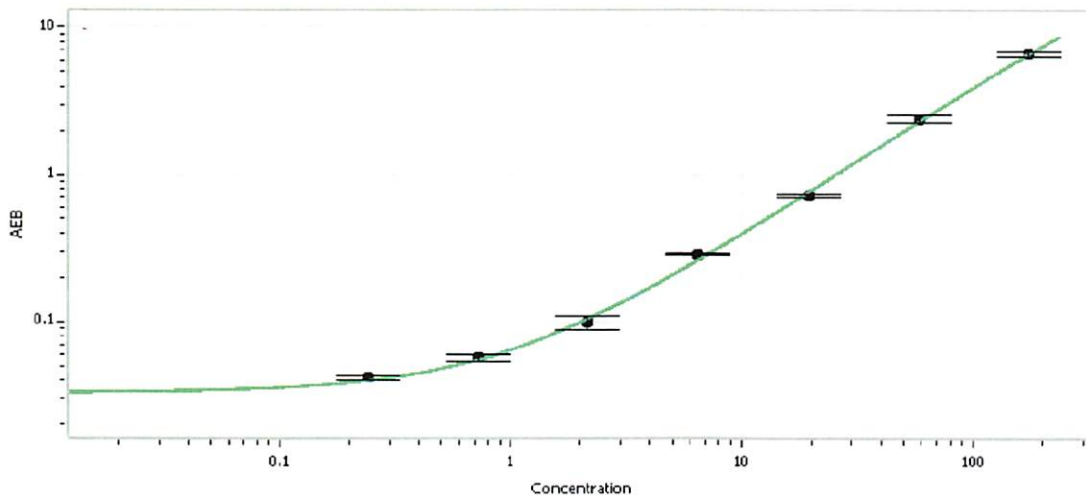
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102101	201808
Detector Stock	102102	202202
Calibrator Concentrate	102105	025403
Calibrator Diluent	102103	129129
Sample Diluent	102104	121126

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 Stock	103297	N/A
RGP Reagent	103159	N/A

Calibrator Concentrate Target Value	3500 pg/mL
-------------------------------------	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.240
C	0.720
D	2.16
E	6.48
F	19.4
G	58.3
H	175

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 08Feb2022
Signature/Date