

Product Number:	102109
Lot Number:	502577
Expiration:	2-Jun-2021
Platform(s):	HD-1 / HD-X & SR-X



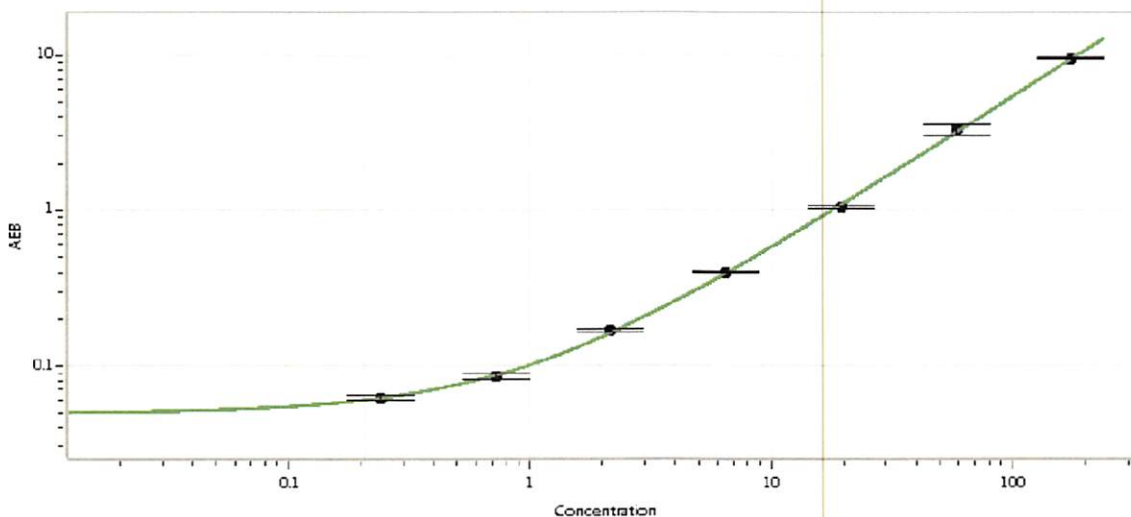
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102101	028709
Detector Stock	102102	029006
Calibrator Concentrate	102105	025403
Calibrator Diluent	102103	024717
Sample Diluent	102104	015433

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	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 Quality Control Manager Shivani Goel 28oct2020

Name: Shivani Goel Title: Quality Control Manager Signature/Date: Shivani Goel 28oct2020