

Product Number:	102700
Lot Number:	502540
Expiration:	2-Jun-2021
Platform(s):	HD-1/HD-X



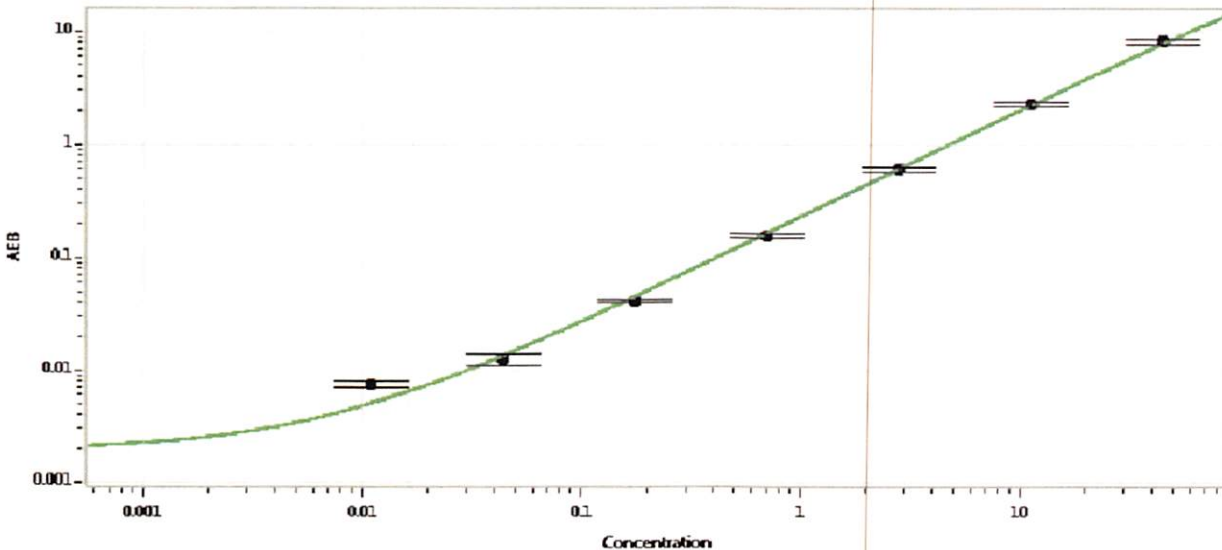
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102693	026203
Detector Stock	102694	026607
Calibrator Concentrate	102697	016302
Calibrator Diluent	102695	015429
Sample Diluent	102696	015428

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	1,000 pg/mL
-------------------------------------	-------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.011
C	0.044
D	0.176
E	0.703
F	2.81
G	11.3
H	45.0

Review/Approval	<u>Shivani Goel</u> <u>Patrizia Stadler</u>	<u>QC Manager</u> <u>Assoc. Director, QC</u>	<u>Shivani Goel 06Oct2020</u>
	Name	Title	Signature/Date