

Product Number:	102700
Lot Number:	502425
Expiration:	4-Dec-2020
Platform(s):	HD-1/HD-X



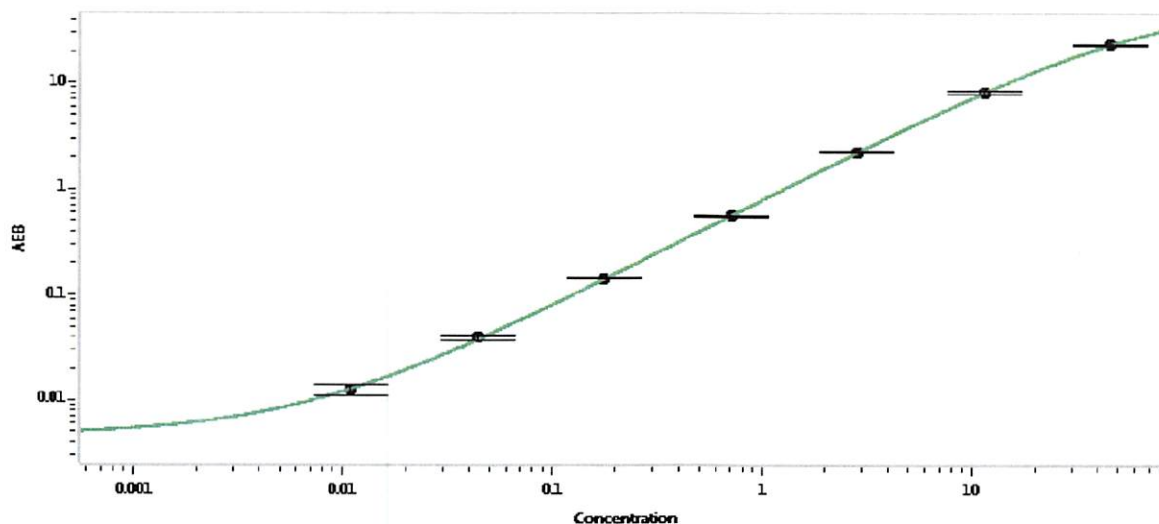
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102693	001306
Detector Stock	102694	001310
Calibrator Concentrate	102697	016302
Calibrator Diluent	102695	933811
Sample Diluent	102696	933812

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

Shivani Goel

QC Manager

Shivani Goel 22Jun2020

Name

Title

Signature/Date