

Product Number:	102648
Lot Number:	502489
Expiration:	21-Feb-2021
Platform(s):	HD-1/HD-X & SR-X



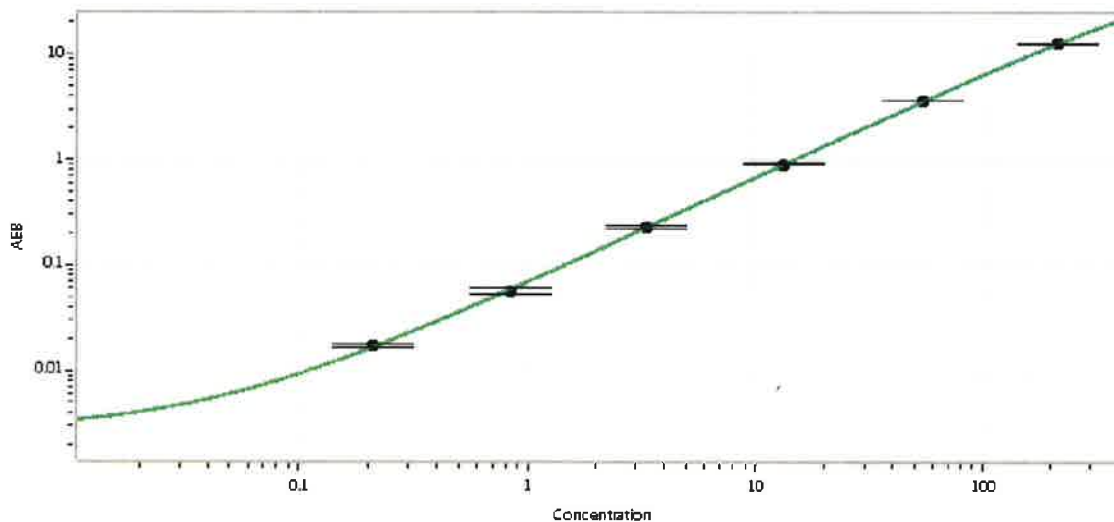
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102641	021303
Detector Stock	102642	022304
Calibrator Concentrate	102645	924602
Calibrator Diluent	102643	020418
Sample Diluent	102644	005231
Detector Diluent	102734	022004

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	3,000 pg/mL
-------------------------------------	-------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.210
C	0.840
D	3.36
E	13.4
F	53.8
G	215

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 18 Aug 2020
Signature/Date