

Product Number:	102648
Lot Number:	502718
Expiration:	3-Sep-2021
Platform(s):	HD-1/HD-X & SR-X



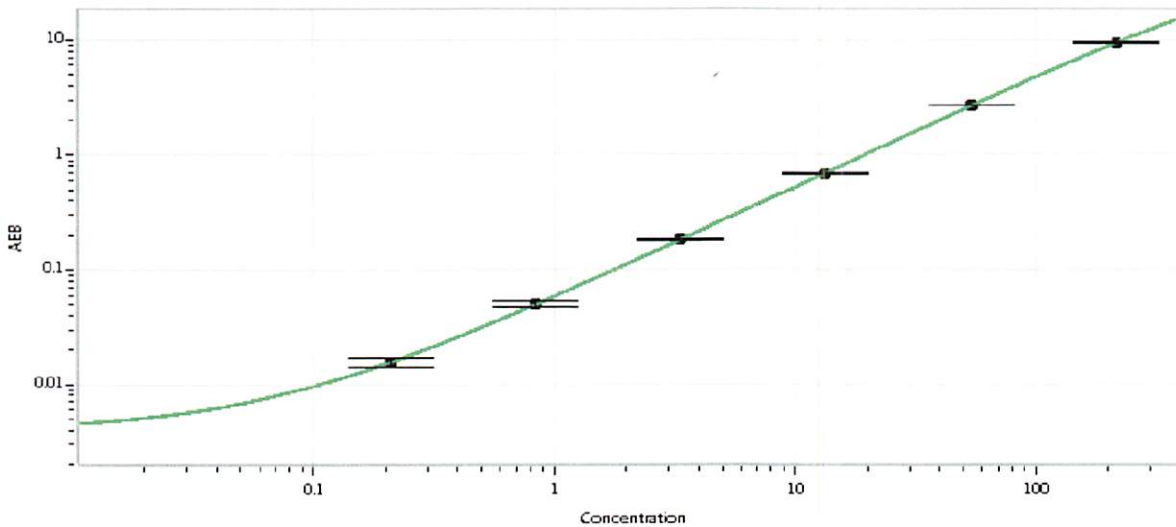
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102641	103210
Detector Stock	102642	106307
Calibrator Concentrate	102645	924602
Calibrator Diluent	102643	102709
Sample Diluent	102644	033811
Detector Diluent	102734	103506

Generic Components - Included With Kit But Not Lot	Number Specific
Discovery Bead Diluent	101954
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

Jungil Park
Name

QC supervisor
Title

[Signature] 23 Mar 2021
Signature/Date