

Product Number:	102039
Lot Number:	502910
Expiration:	4-Mar-2022
Platform(s):	HD-1/HD-X & SR-X



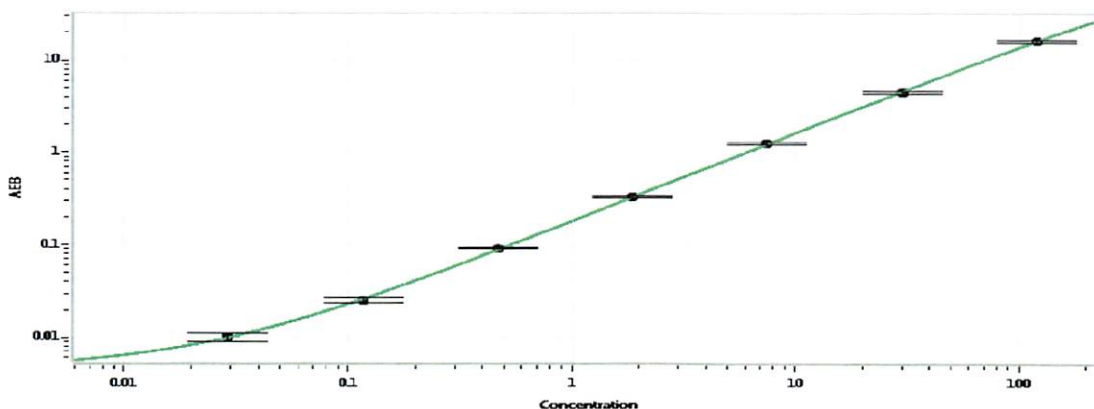
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102028	111604
Detector Stock	102029	111606
Calibrator Concentrate	102032	110318
Calibrator Diluent	102030	111313
Sample Diluent	103402	110013

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	2 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.029
C	0.117
D	0.469
E	1.88
F	7.50
G	30.0
H	120

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 12 May 2021
Signature/Date