

Product Number:	102209
Lot Number:	502502
Expiration:	24-Mar-2021
Platform(s):	HD-1/HD-X & SR-X



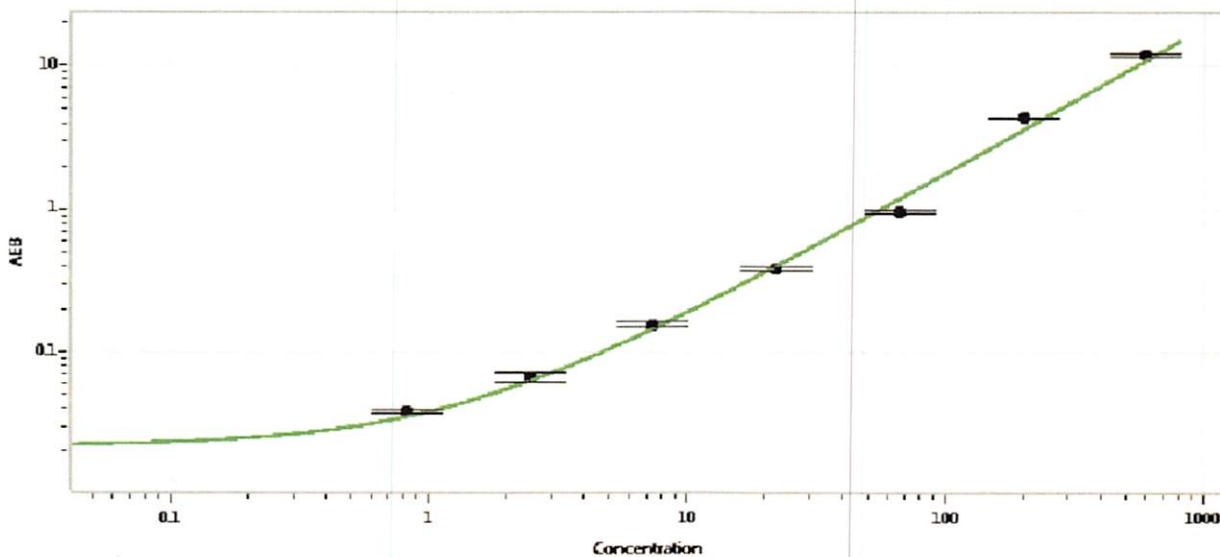
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102201	023004
Detector Stock	102202	023104
Calibrator Concentrate	102205	023103
Calibrator Diluent	102203	023101
Sample Diluent	102204	023102

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	12 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.823
C	2.47
D	7.41
E	22.2
F	66.7
G	200
H	600

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 02Sep2020

Name

Title

Signature/Date