

Product Number:	102475
Lot Number:	502438
Expiration:	4-Dec-2020
Platform(s):	HD-1/HD-X



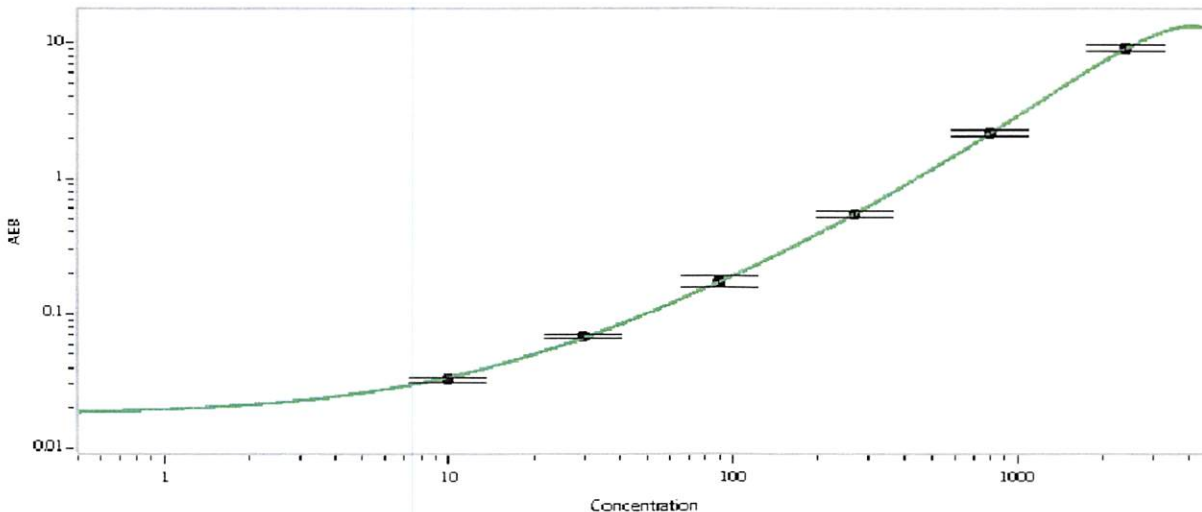
Data below represents results generated on the Simoa™ HD-X Analyzer.

Component	Part Number	Lot Number
Bead Stock	102468	017409
Detector Stock	102469	016407
Calibrator Concentrate	102472	004511
Calibrator Diluent	102470	934510
Sample Diluent	102471	933809

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	60 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	9.88
C	29.6
D	88.9
E	267
F	800
G	2400

Review/Approval

Shivani Goel
Name

Manager, Quality Control
Title

Shivani Goel 26 Jun 2020
Signature/Date