

Product Number:	101855
Lot Number:	502319
Expiration:	18-Feb-2021
Platform(s):	HD-1/HD-X



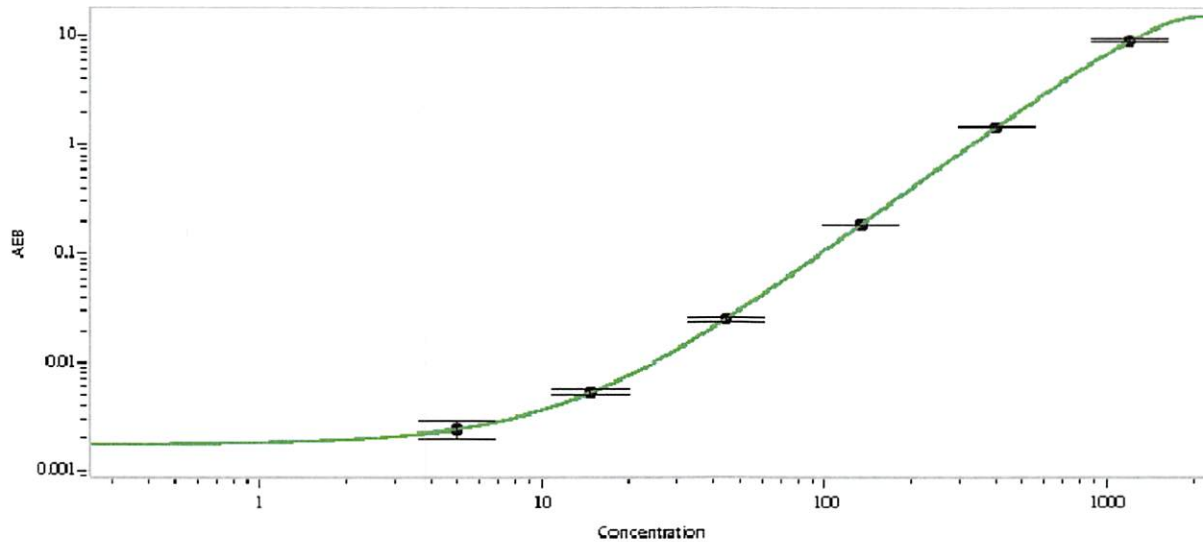
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101902	009806
Detector Stock	101903	009906
Calibrator Concentrate	101850	010102
Calibrator Diluent	101988	009506
Sample Diluent	101989	004914

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	30 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	4.94
C	14.8
D	44.4
E	133
F	400
G	1200

Review/Approval: Shivani Goel Name: Shivani Goel Title: QC Manager Signature/Date: Shivani Goel 16Apr2020