

Product Number:	102109
Lot Number:	502533
Expiration:	2-May-2021
Platform(s):	HD-1 / HD-X & SR-X



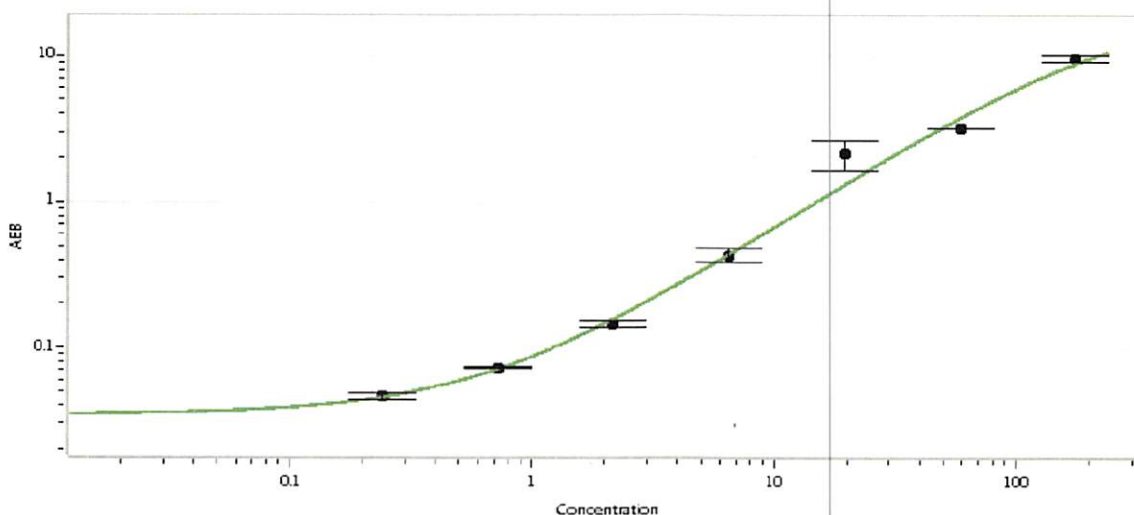
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102101	025801
Detector Stock	102102	025404
Calibrator Concentrate	102105	025403
Calibrator Diluent	102103	024707
Sample Diluent	102104	015427

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	3500 pg/mL
-------------------------------------	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.240
C	0.720
D	2.16
E	6.48
F	19.4
G	58.3
H	175

Review/Approval

Shivani Goel  
Name

Quality Control Manager  
Title

Shivani Goel 24 Sep 2020  
Signature/Date