

Product Number:	103093
Lot Number:	502534
Expiration:	2-Mar-2021
Platform(s):	HD-1/ HD-X & SR-X

Component	Part Number	Lot Number
Bead Stock	102021	025308
Detector Stock	103043	025505
Calibrator Concentrate	103047	935308
Calibrator Diluent	103044	025507
Sample Diluent	103046	025506

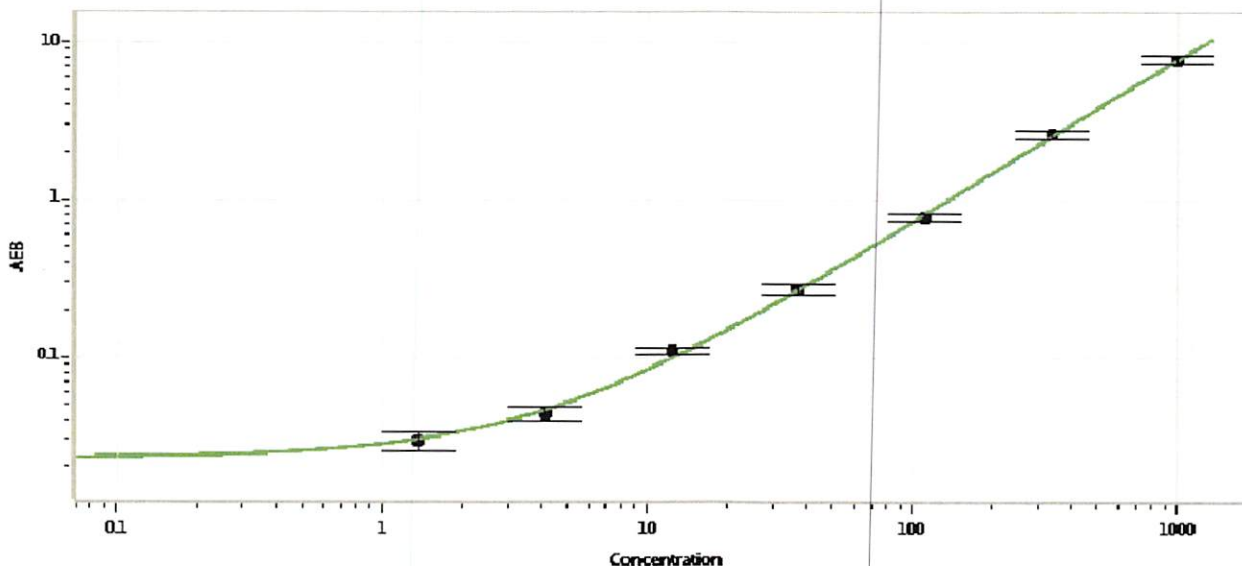


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	1.37
C	4.12
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 24Sep
Signature/Date 2020