

Product Number:	102184
Lot Number:	502569
Expiration:	2-Jun-2021
Platform(s):	HD-1/ HD-X

Component	Part Number	Lot Number
Bead Stock	102176	022408
Detector Stock	102177	028003
Calibrator Concentrate	102180	005605
Calibrator Diluent	102178	024715
Sample Diluent	102179	015431

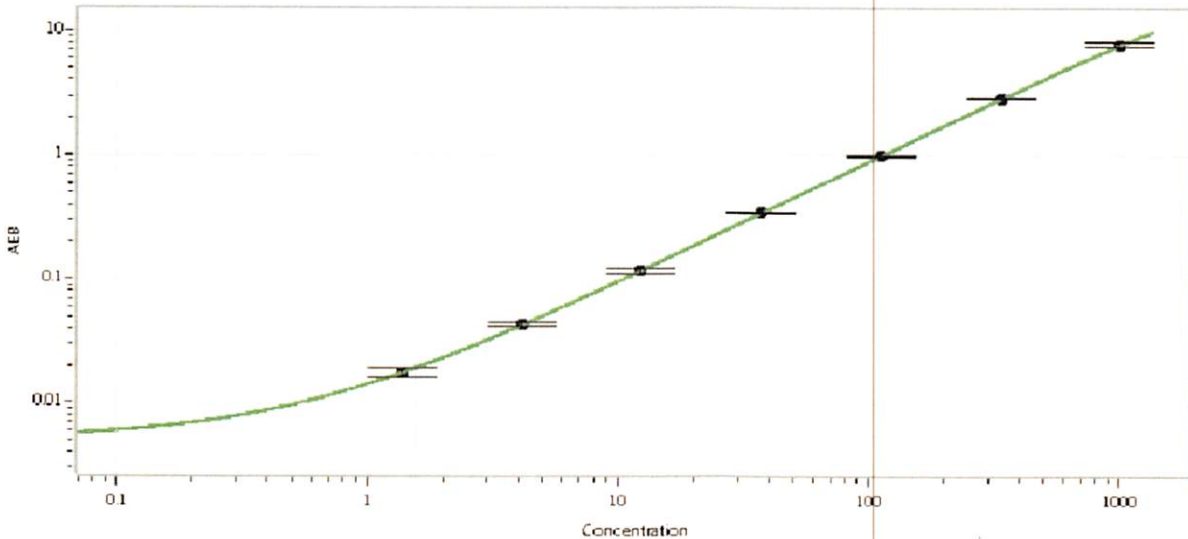


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	40 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

QC Manager

Shivani Goel 02NOV 2020

Name

Title

Signature/Date