

Product Number:	102233
Lot Number:	502501
Expiration:	5-Nov-2020
Platform(s):	HD-1/HD-X, SRX



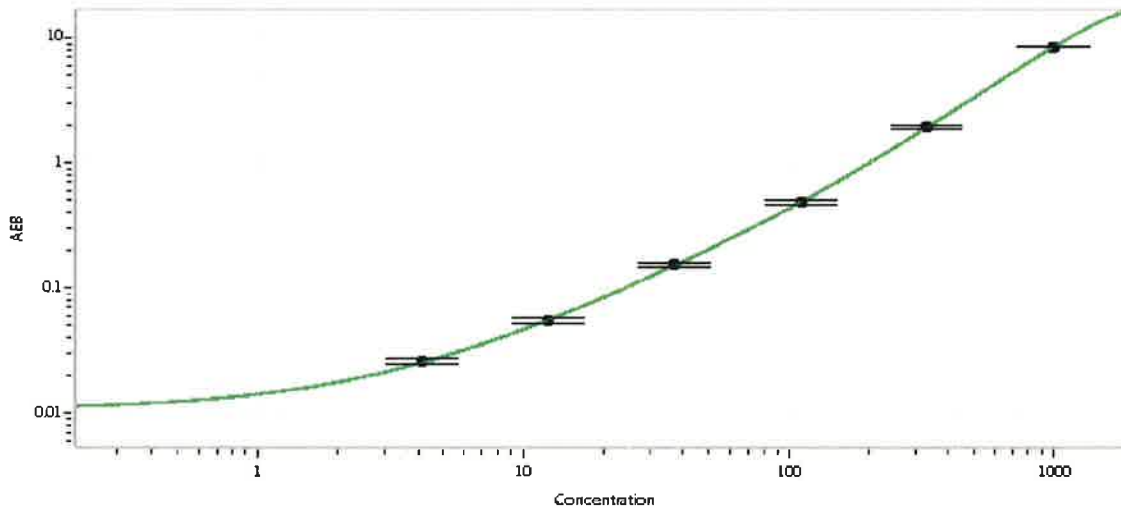
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102223	022707
Detector Stock	102224	020206
Calibrator Concentrate	102229	835303
Calibrator Diluent	102225	012605
Sample Diluent	102226	012604

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	20 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	4.12
C	12.3
D	37.0
E	111
F	333
G	1000

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 20 Aug 2020

Name

Title

Signature/Date