

Product Number:	102475
Lot Number:	503454
Expiration:	10-Nov-2022
Platform(s):	HD-1/HD-X



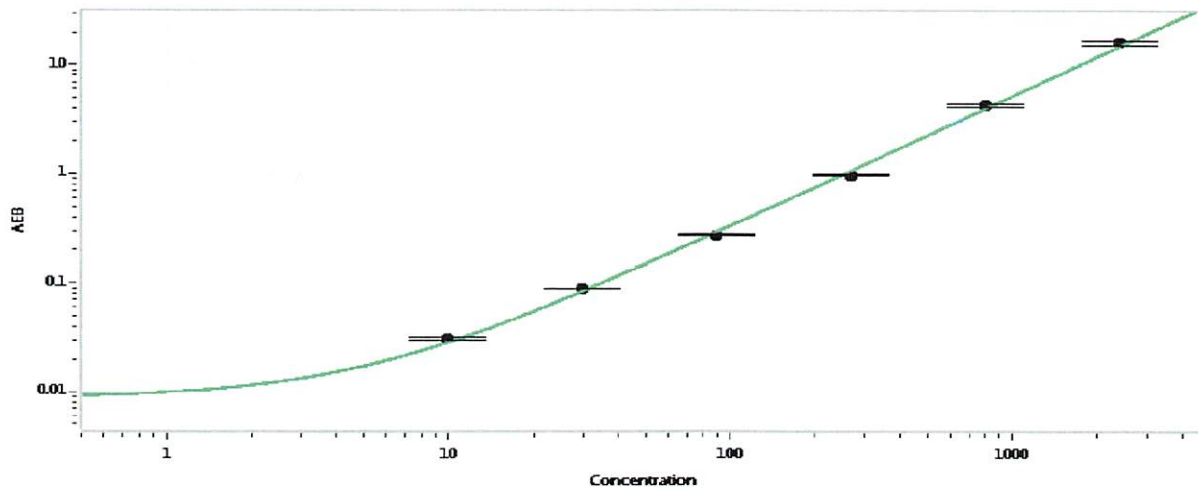
Data below represents results generated on the Simoa® HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102468	215205
Detector Stock	102469	215306
Calibrator Concentrate	102472	214307
Calibrator Diluent	102470	212706
Sample Diluent	102471	211314

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	60ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	9.88
C	29.6
D	88.9
E	267
F	800
G	~2400

Review/Approval

*Matt Smart*  
Name

*Director, QC*  
Title

*MAT 16 Jun 2022*  
Signature/Date