

Product Number:	102382
Lot Number:	503429
Expiration:	12-Dec-2022
Platform(s):	HD-1



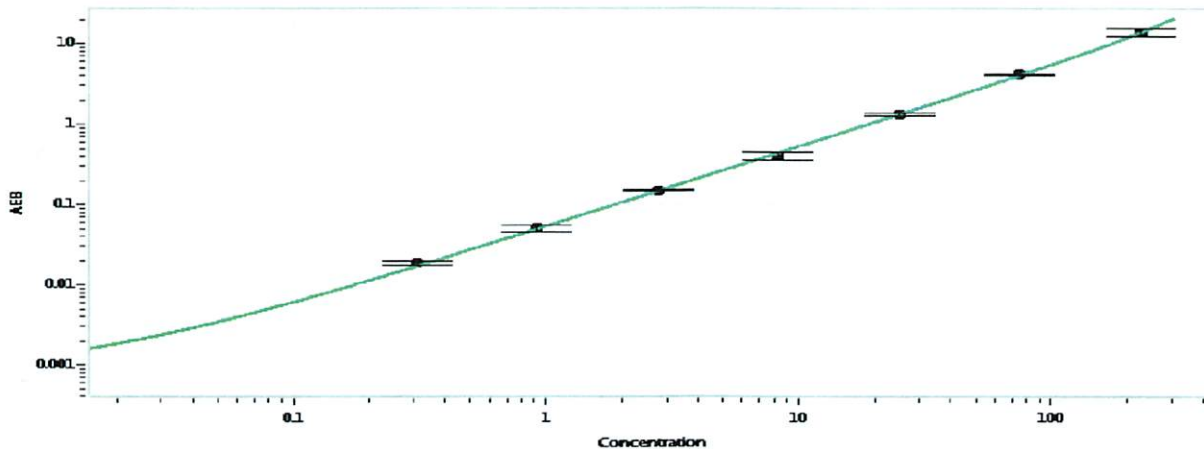
Data below represents results generated on the Simoa® HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102374	210501
Detector Stock	102375	210601
Calibrator Concentrate	102379	210402
Calibrator Diluent	102376	210603
Sample Diluent	102378	210602

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	4500 pg/mL
-------------------------------------	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.309
C	0.926
D	2.78
E	8.33
F	25.0
G	75.0
H	225

Review/Approval

Matt Smart

Director, Quality Control

M. A. Smart 20 May 2022

Name

Title

Signature/Date