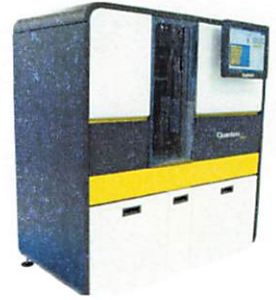


Product Number:	102794
Lot Number:	503366
Expiration:	10-Nov-2022
Platform(s):	HD-1/HD-X



Component	Part Number	Lot Number
Bead Stock	102787	205806
Detector Stock	102788	206103
Calibrator Concentrate	102791	202306
Calibrator Diluent	102789	133932
Sample Diluent	102790	205805

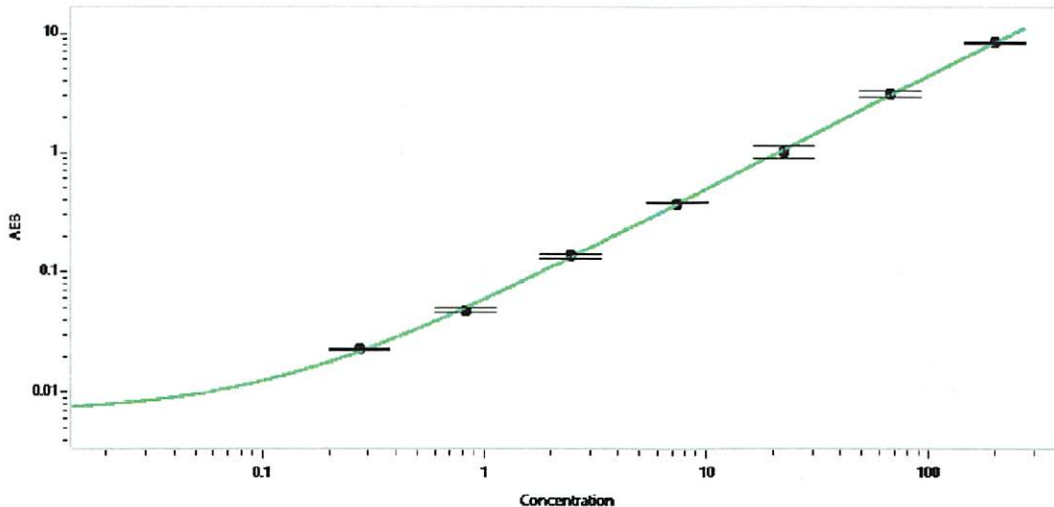
Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159

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

Calibrator Concentrate Target Value	4 ng/mL
-------------------------------------	---------

Calibration Curve

2022-03-20 VegF Discovery Kit lot 503366 Release QC (HDX-83) PR Curve 1



Calibrator Levels (pg/mL)

A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 31 Mar 2022  
Signature/Date