

Product Number:	102039
Lot Number:	502409
Expiration:	21-Feb-2021
Platform(s):	HD-1/HD-X & SR-X



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

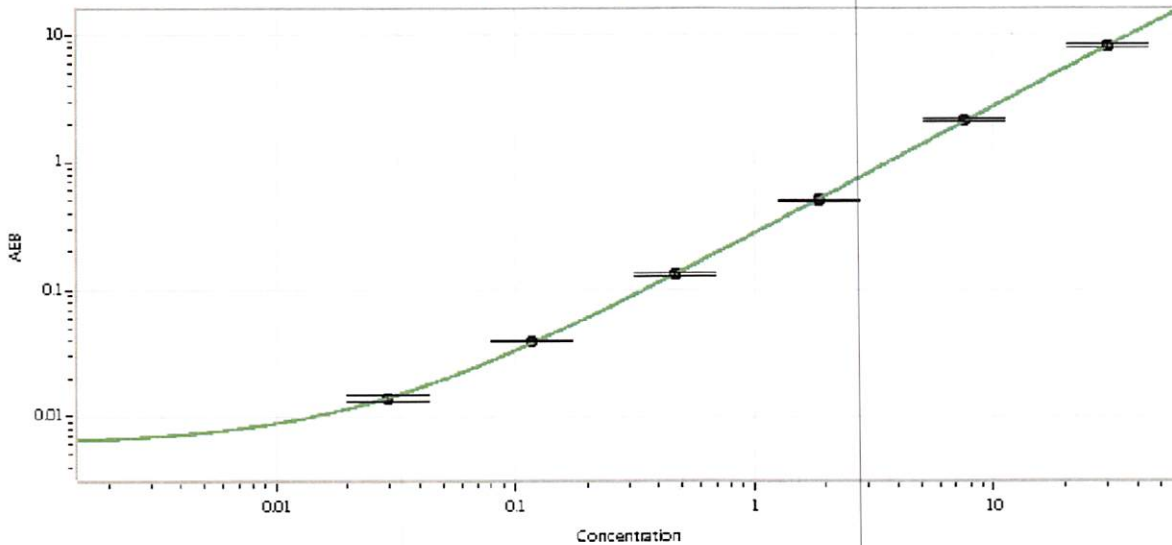
Component	Part Number	Lot Number
Bead Stock	102028	026205
Detector Stock	102029	015904
Calibrator Concentrate	102032	910804
Calibrator Diluent	102030	015112
Sample Diluent	103402	005236

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	2 ng/mL
-------------------------------------	---------

Calibration Curve

2020-09-21 BDNF D Kit (Lot 502409) Release QC (CU-122) Curve 1



Calibrator Levels (pg/mL)

A	0.000
B	0.029
C	0.117
D	0.469
E	1.88
F	7.50
G	30.0
H	120

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 25 Sep 2020

Name

Title

Signature/Date