

Product Number:	102209
Lot Number:	502266
Expiration:	2-Oct-2020
Platform(s):	HD-1/HD-X & SR-X



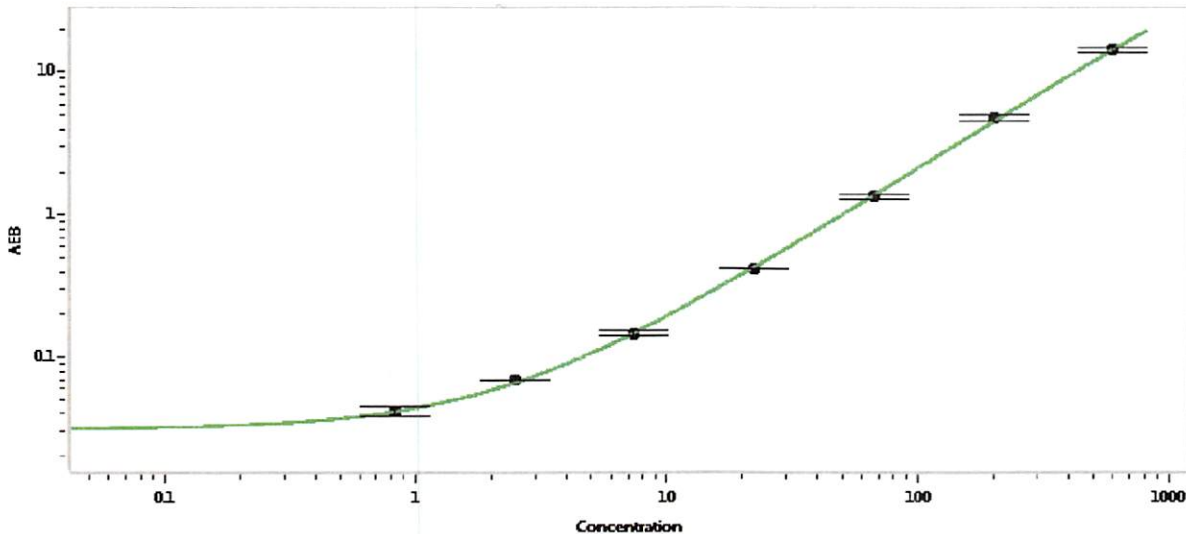
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102201	006409
Detector Stock	102202	006301
Calibrator Concentrate	102205	827108
Calibrator Diluent	102203	006504
Sample Diluent	102204	006505

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	12 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.823
C	2.47
D	7.41
E	22.2
F	66.7
G	200
H	600

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	11/10/2020
	Name	Title	Signature/Date	