

Product Number:	102064
Lot Number:	501684
Expiration:	27-Sep-2019
Platform(s):	HD-1



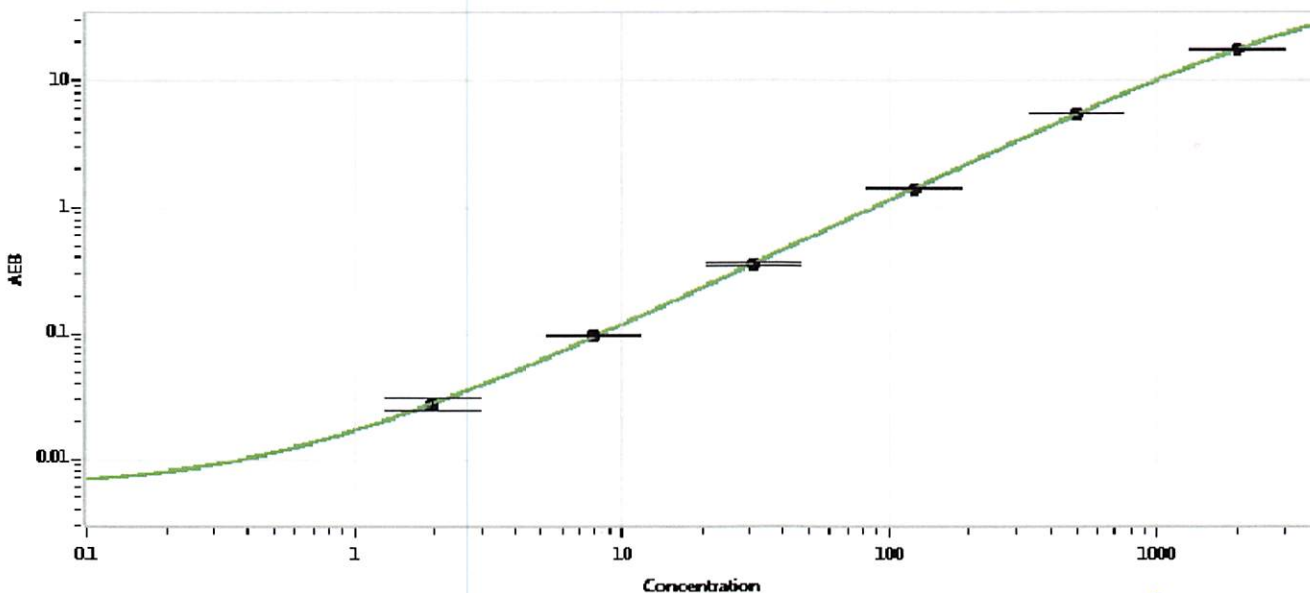
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102056	831817
Detector Stock	102057	827403
Calibrator Concentrate	102060	833709
Calibrator Diluent	102058	827014
Sample Diluent	102059	827015

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	40 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.95
C	7.81
D	31.3
E	125
F	500
G	2000

Review/Approval _____ *Junpil Park QC Analyst II  05Mar2019

Name Title Signature/Date

*On behalf of Patrizia Stadler Assoc. Director, QC