

Product Number:	102929
Lot Number:	502022
Expiration:	01-AUG-202
Platform(s):	HD-1/HD-X



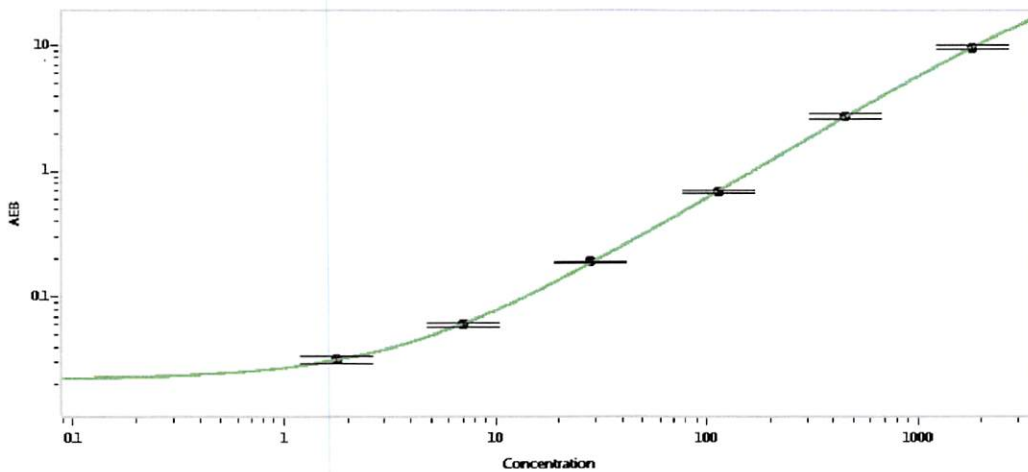
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102922	926312
Detector Stock	102923	926311
Calibrator Concentrate	102926	926313
Calibrator Diluent	102924	926310
Sample Diluent	102925	926210

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	36 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.76
C	7.03
D	28.1
E	113
F	450
G	180

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i> 26 Sep 2019
	Name	Title	Signature/Date