

Product Number:	103093
Lot Number:	502611
Expiration:	16-Oct-2021
Platform(s):	HD-1/ HD-X & SR-X



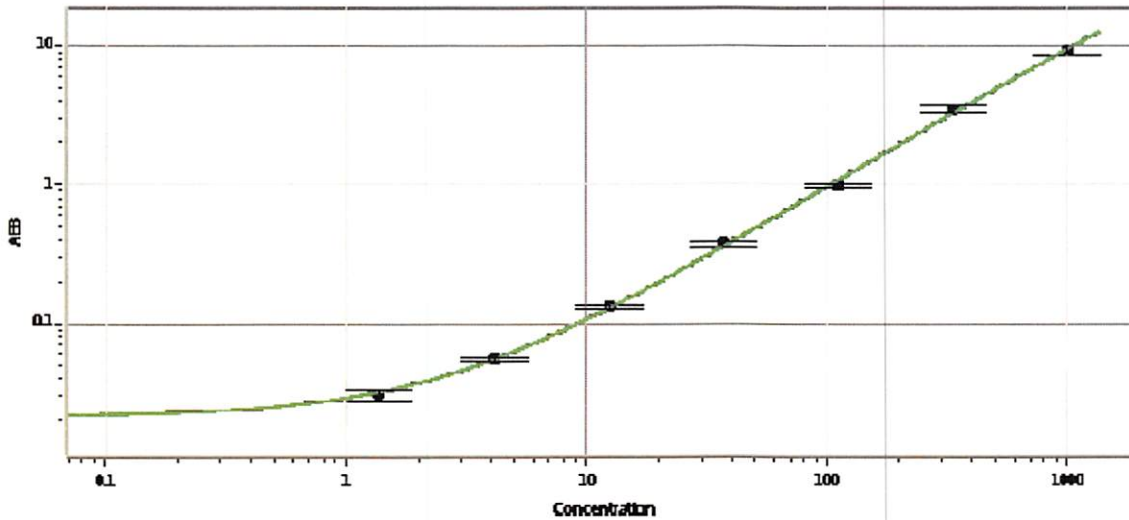
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102021	031011
Detector Stock	103043	031405
Calibrator Concentrate	103047	935308
Calibrator Diluent	103044	032111
Sample Diluent	103046	031110

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	1.37
C	4.12
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 25 Nov 2020
	Name	Title	Signature/Date