

Product Number:	102394
Lot Number:	502407
Expiration:	21-Feb-2021
Platform(s):	HD-1/ HDX



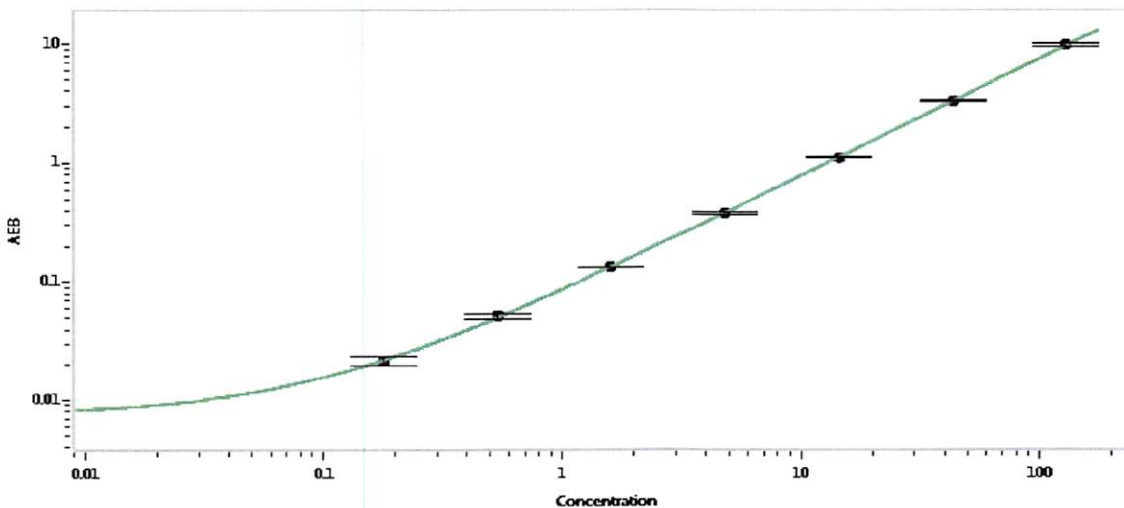
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102387	015101
Detector Stock	102388	015008
Calibrator Concentrate	102391	015401
Calibrator Diluent	102389	012513
Sample Diluent	102390	005234

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	4,000 pg/mL
-------------------------------------	-------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.178
C	0.535
D	1.60
E	4.81
F	14.4
G	43.3
H	130

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 08 Jun 2020
	Name	Title	Signature/Date