

Product Number:	102669
Lot Number:	502536
Expiration:	3-Sep-2021
Platform(s):	HD-1/HD-X



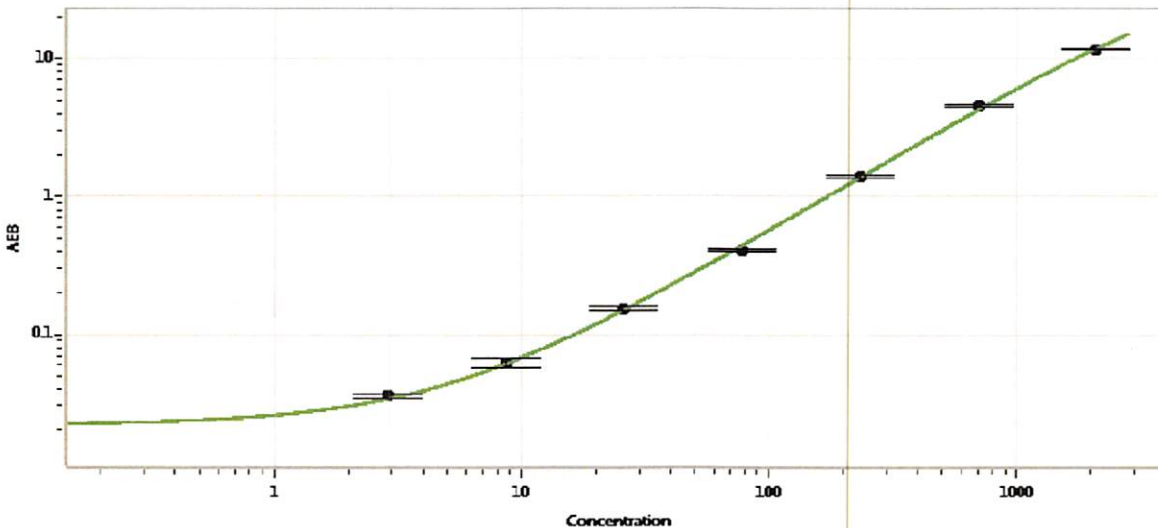
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102662	024704
Detector Stock	102663	030804
Calibrator Concentrate	102666	031803
Calibrator Diluent	102664	025214
Sample Diluent	102665	025206

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	60 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.88
C	8.64
D	25.9
E	77.8
F	233
G	700
H	2100

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 20 Nov 2020
	Name	Title	Signature/Date