

Product Number:	102336
Lot Number:	502365
Expiration:	18-Feb-2021
Platform(s):	HD-1/HD-X



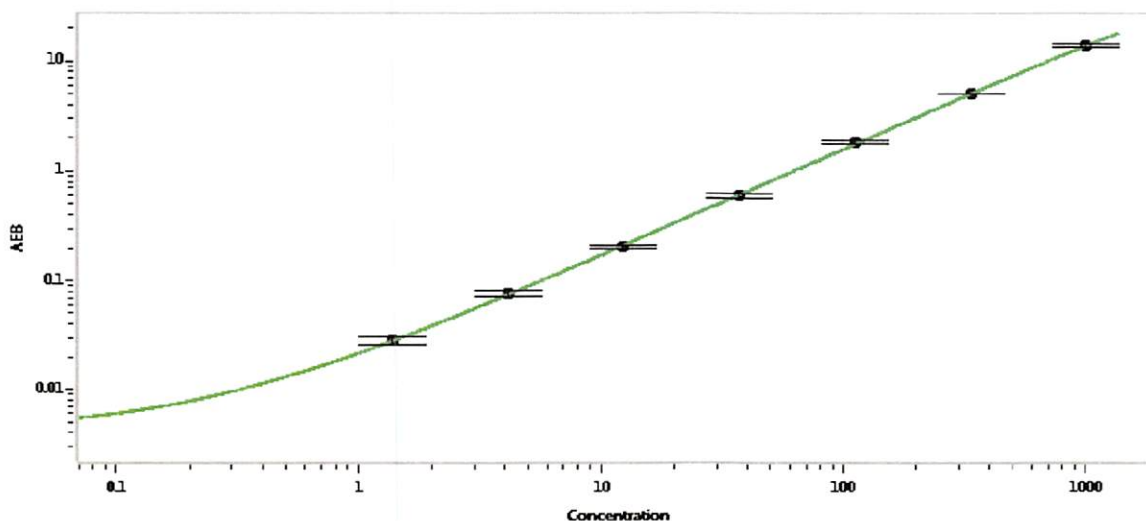
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102330	012401
Detector Stock	102331	012402
Calibrator Concentrate	102335	012106
Calibrator Diluent	102332	004916
Sample Diluent	102333	012103

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	20 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.11
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 13 May 2020
	Name	Title	Signature/Date