

Product Number:	102073
Lot Number:	502485
Expiration:	27-May-2021
Platform(s):	HD-1/HD-X

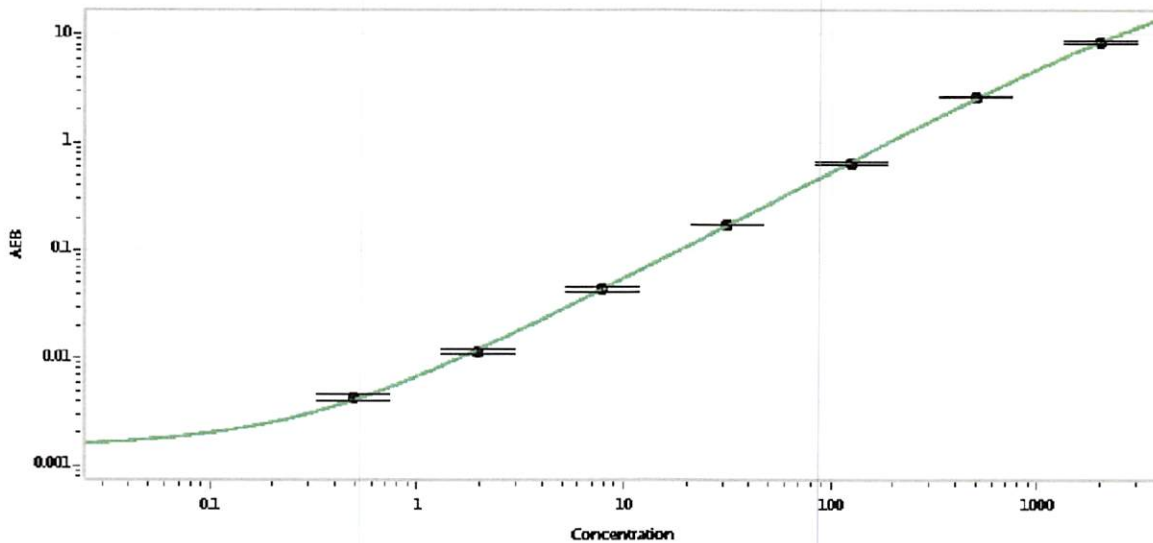


Component	Part Number	Lot Number
Bead Stock	102065	018906
Detector Stock	102066	021103
Calibrator Concentrate	102069	019101
Calibrator Diluent	102067	019006
Sample Diluent	102068	019005

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

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

Calibrator Concentrate Target Value	40 ng/mL
-------------------------------------	----------



Calibrator Levels (pg/mL)	
A	0.000
B	0.488
C	1.95
D	7.81
E	31.3
F	125
G	500
H	2000

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 06 Aug 2020
	Name	Title	Signature/Date