

Product Number:	102233
Lot Number:	502144
Expiration:	23-May-2020
Platform(s):	HD-1/HD-X



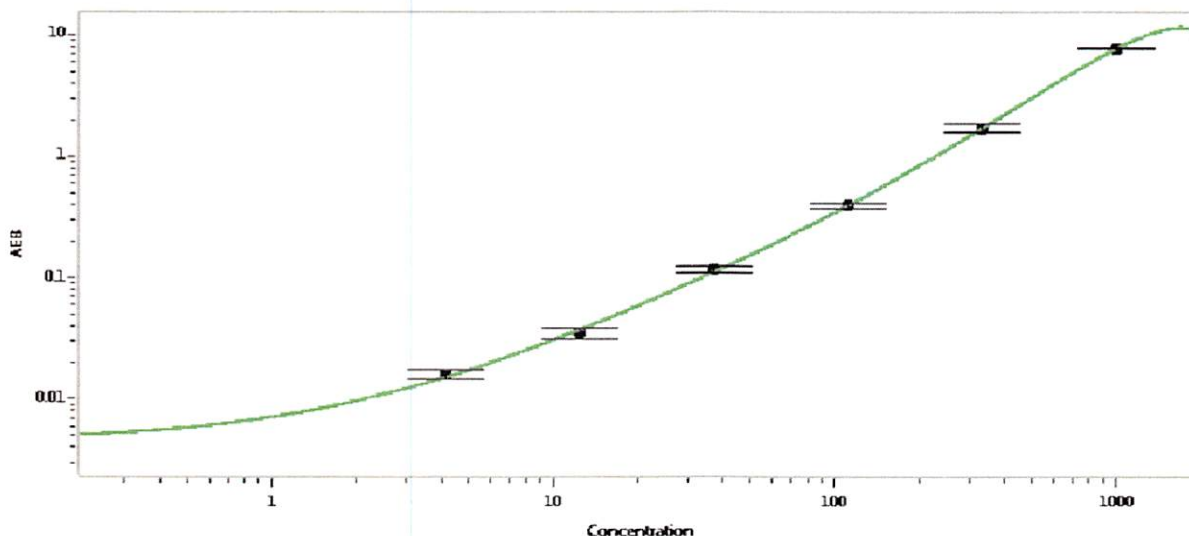
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102223	933608
Detector Stock	102224	932705
Calibrator Concentrate	102229	835303
Calibrator Diluent	102225	932902
Sample Diluent	102226	932905

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	4.12
C	12.3
D	37.0
E	111
F	333
G	1000

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>p.stadler</i> 19 Dec 2019
	Name	Title	Signature/Date