

Product Number:	102233
Lot Number:	502566
Expiration:	20-Jan-2021
Platform(s):	HD-1/HD-X, SRX

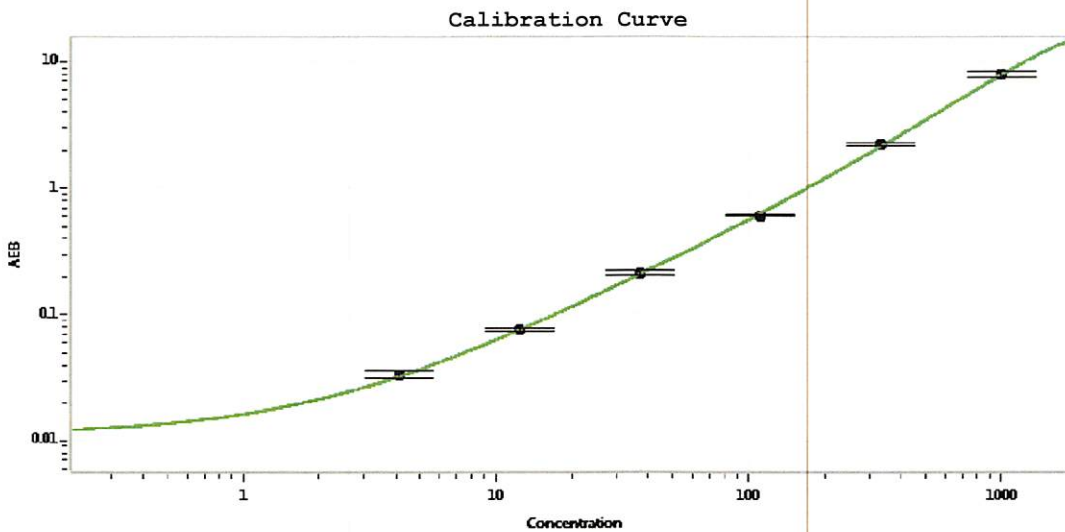


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102223	022707
Detector Stock	102224	020206
Calibrator Concentrate	102229	027601
Calibrator Diluent	102225	025410
Sample Diluent	102226	025409

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
-------------------------------------	----------



Calibrator Levels (pg/mL)	
A	0.000
B	4.12
C	12.3
D	37.0
E	111
F	333
G	1000

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 09 Nov 2020
	Name	Title	Signature/Date