

Product Number:	102318
Lot Number:	502588
Expiration:	24-Aug-2021
Platform(s):	HD-1/HDX



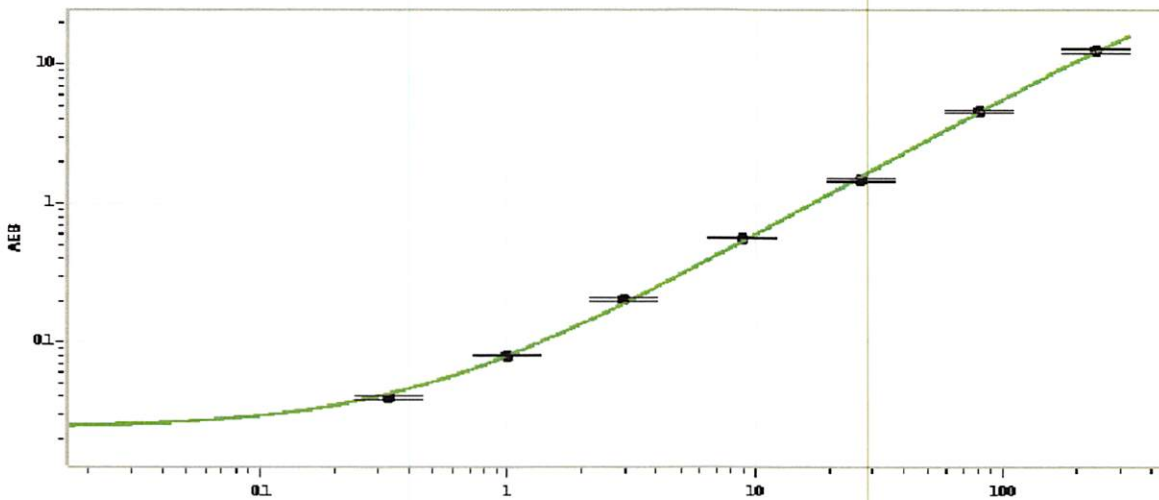
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102309	029312
Detector Stock	102310	029406
Calibrator Concentrate	102314	030204
Calibrator Diluent	102311	029402
Sample Diluent	102312	029404

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	4,800 pg/mL
--	-------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.329
C	0.988
D	2.96
E	8.89
F	26.7
G	80.0
H	240

Review/Approval

Shivani Goel
Name

Manager, QC
Title

Shivani Goel 09 Nov 2020
Signature/Date