

Product Number:	101871
Lot Number:	502201
Expiration:	21-Nov-2020
Platform(s):	HD-1/HD-X



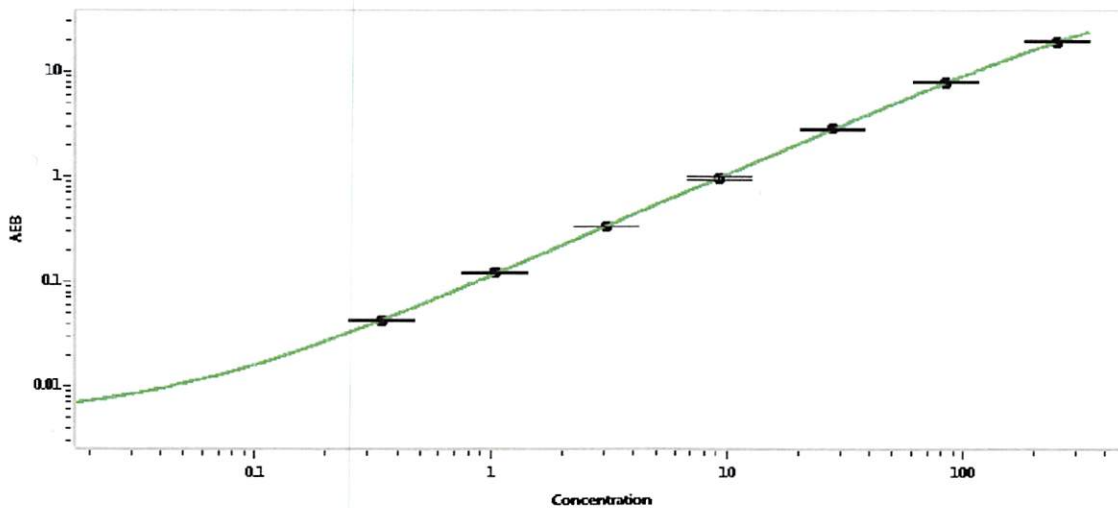
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101893	001605
Detector Stock	101894	001601
Calibrator Concentrate	101866	926009
Calibrator Diluent	101959	001316
Sample Diluent	101958	932625

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	5 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.343
C	1.03
D	3.09
E	9.26
F	27.8
G	83.3
H	250

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler 27 Jan 2020

Name

Title

Signature/Date