

Product Number:	102794
Lot Number:	502176
Expiration:	21-Nov-2020
Platform(s):	HD-1/HD-X



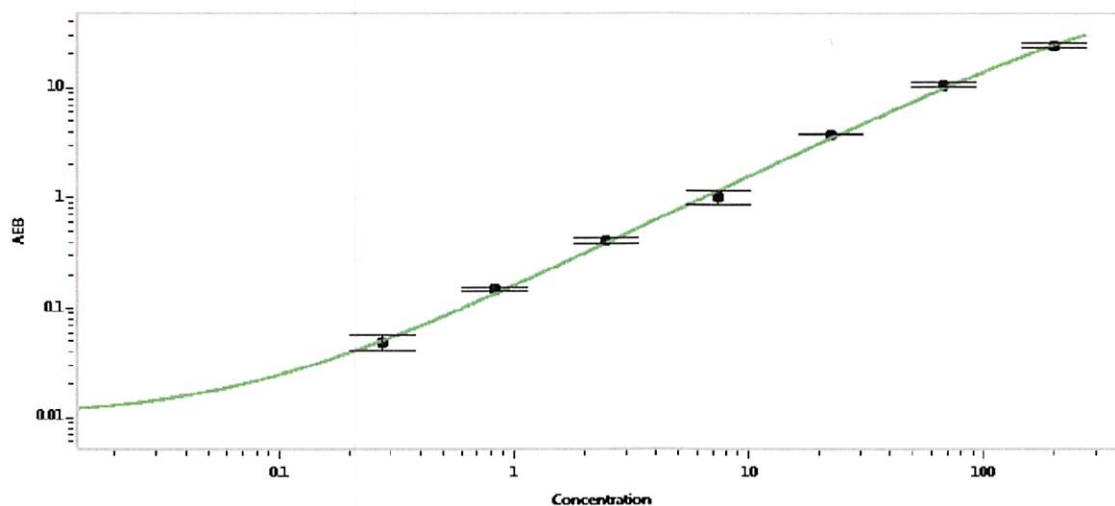
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102787	935111
Detector Stock	102788	935305
Calibrator Concentrate	102791	936007
Calibrator Diluent	102789	935328
Sample Diluent	102790	935320

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Patrizia Stadler

Assoc. Director, QC

*P. Stadler* 14 Jan 2020

Name

Title

Signature/Date