

Product Number:	102669
Lot Number:	502078
Expiration:	11-Sep-2020
Platform(s):	HD-1/HD-X



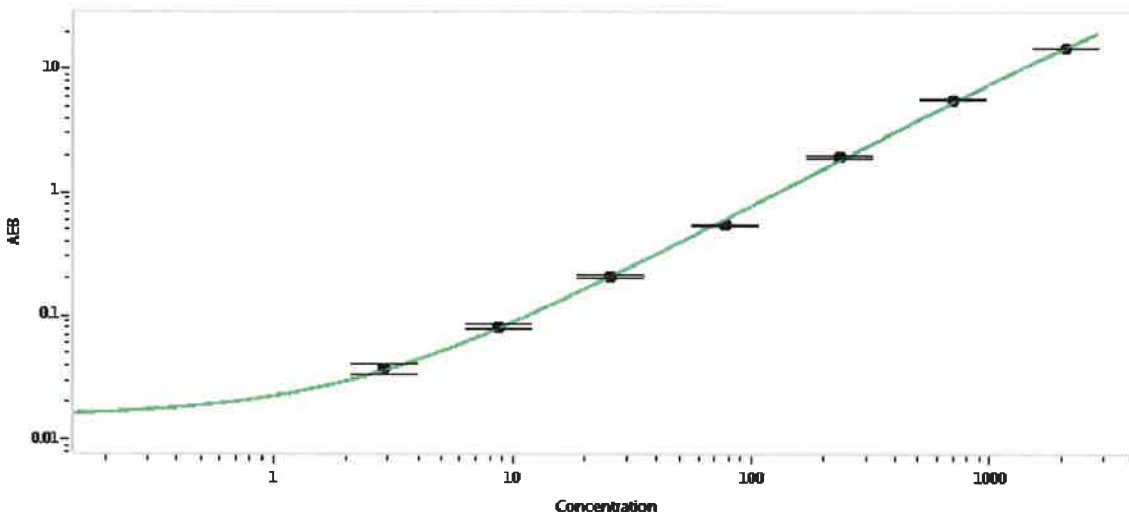
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102662	925508
Detector Stock	102663	928209
Calibrator Concentrate	102666	913401
Calibrator Diluent	102664	926209
Sample Diluent	102665	925409

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.88
C	8.64
D	25.9
E	77.8
F	233
G	700
H	2100

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler 18 Oct 2019

Name

Title

Signature/Date