

Product Number:	102669
Lot Number:	502086
Expiration:	17-Sep-2020
Platform(s):	HD-1/HD-X



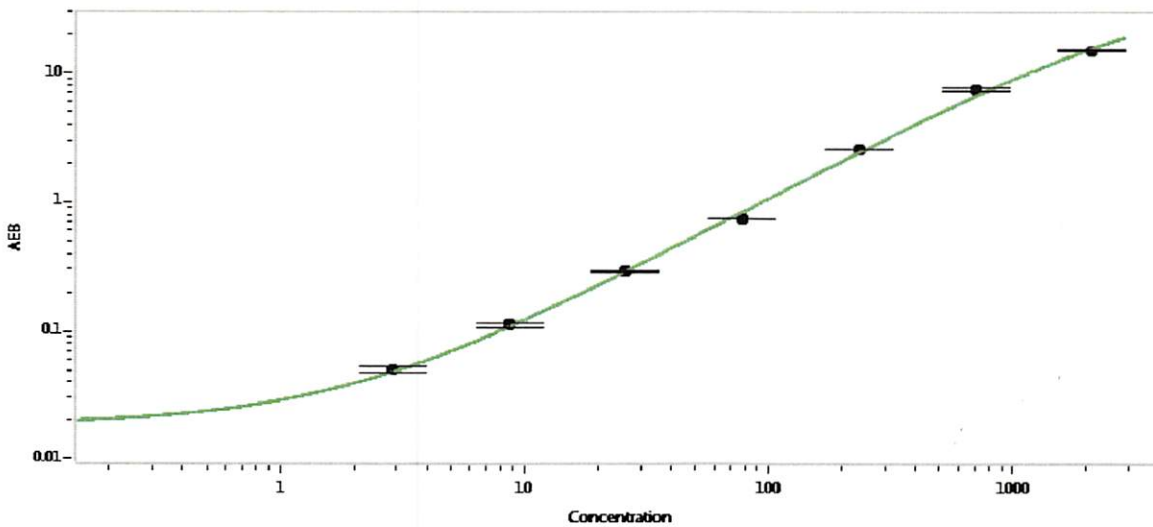
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102662	929111
Detector Stock	102663	928213
Calibrator Concentrate	102666	913401
Calibrator Diluent	102664	929112
Sample Diluent	102665	929105

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
Name

Assoc. Director, QC
Title

p. Stoollev 01 NOV 2019
Signature/Date