

Product Number:	102570
Lot Number:	502116
Expiration:	1-Aug-2020
Platform(s):	HD-1/HD-X



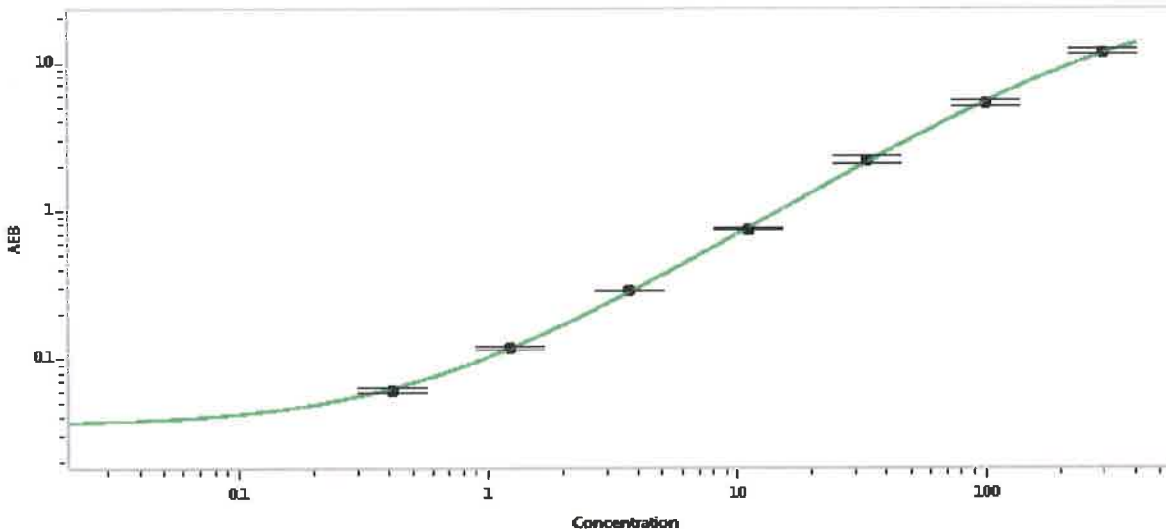
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102563	931907
Detector Stock	102564	931812
Detector Diluent	102590	931814
Calibrator Concentrate	102567	829005
Calibrator Diluent	102565	927718
Sample Diluent	102566	927717

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159

Calibrator Concentrate Target Value	6 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.412
C	1.23
D	3.70
E	11.1
F	33.3
G	100
H	300

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 06 Dec 2019
Signature/Date