

Product Number:	102336
Lot Number:	502263
Expiration:	21-Nov-2020
Platform(s):	HD-1/HD-X



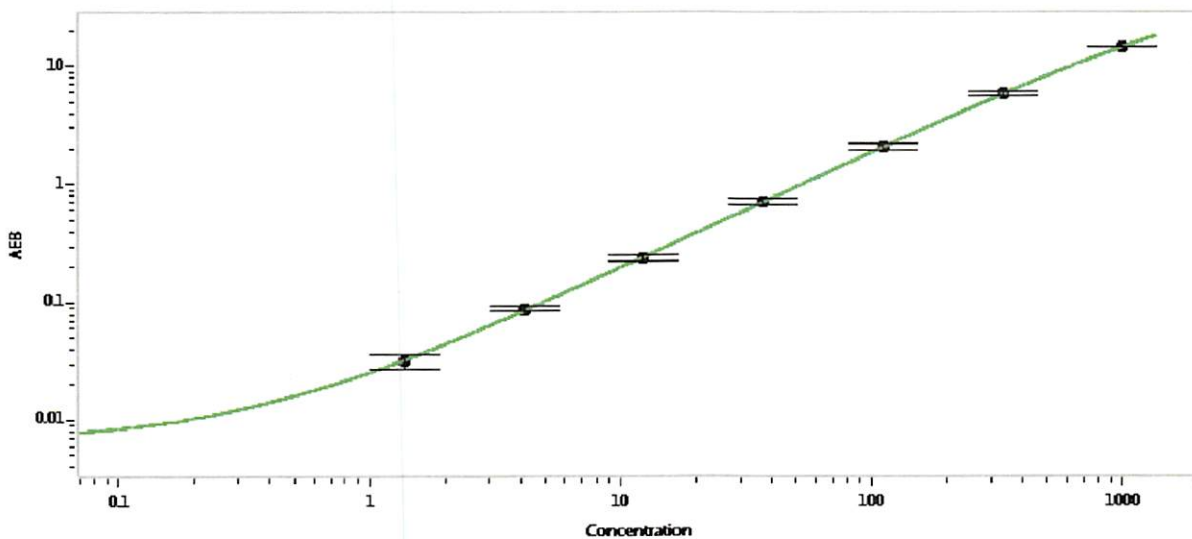
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102330	006003
Detector Stock	102331	006001
Calibrator Concentrate	102335	006004
Calibrator Diluent	102332	004912
Sample Diluent	102333	006202

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.11
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler 06 MAR 2020

Name

Title

Signature/Date